thermoscientific

Maps 3.32.1 User Guide

PN 1551059

Revision A • March 2025 • Original Instructions Limited Rights



Copyright and Trademarks

Technical Publications

Technical Publications Team - Hillsboro

Copyright © 2025 by FEI Company, a part of Thermo Fisher Scientific. The information and materials contained herein are confidential and proprietary to Thermo Fisher and are provided for your organization's internal use on a need-to-know basis. They cannot be duplicated, published or disseminated to any third party without the express written consent of Thermo Fisher.

Trademark Acknowledgments

All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. FEI and the FEI logo are registered trademarks of FEI Company (part of Thermo Fisher Scientific and its affiliates). All other trademarks belong to their respective owners.

Excel and Microsoft are registered trademarks of Microsoft Corporation.

Limited Rights

The following notice applies to the U.S. Government and other purchases with federal funds:

Contractor Name: Thermo Fisher Scientific

Contractor Address: 5350 NE Dawson Creek Drive, Hillsboro, OR 97124. The Government's rights to use, modify, reproduce, release, perform, display, or disclose these technical data are restricted to those rights specified in DFARS 252.227-7015(b)(2), FAR 52.227-14(g)(2)(Alternate II) and FAR 12.211. Any reproduction of technical data or portions thereof marked with this legend must also reproduce the markings. Any person, other than the Government, who has been provided access to such data, must promptly notify the above named Contractor.

Document History

Rev A: March 2025

Maps 3.32.1 Release Notes

This document specifies changes made to the Thermo Scientific[™] Maps 3.32.1 software application.

New Features

Phenom: Manual selection of colors for EDS

Maps can now assign predefined colors with EDS elements.

See Phenom Tile Set Tab for more information.

Phenom: EDS parameters in tile set templates

Maps now includes EDS parameters in tile set templates.

See Tile Set Templates for more information.

TEM: Support for TEM 3.23 and TEM 2.23

Maps now supports TEM server versions 3.18-3.23 for Titan and 2.18-2.23 for Talos.

Known Issues

See Known issues for known issues introduced with this version of Maps software.

Bug Fixes

The following bugs were fixed in this release of the Maps software.

Bug ID	Description
MAPS-10481	Maps EDS colors for Phenom are now being taken from the same registry in which Phenom UI stores them. Therefore, colors in both MAPS and Phenom match one another.
MAPS-10655	The Phenom metadata is now retrieved and saved in Maps. This information is available under the Show Metadata button on snapshots and stitched tile sets.

Bug ID	Description
MAPS-11408	Now when you export a stitched tile set, the scale bar option is available in the context menu.
MAPS-11434	If you import an image into Maps and the file name does not include the image resolution, an error message now appears to indicate that the file name must include the image resolution.
MAPS-12291	All EDS fields are now displayed regardless of their checkbox status (that is, activated/deactivated). To change a checkbox status, click the checkbox of the desired EDS field.
MAPS-12654	Temporary files are now deleted correctly during time-intensive runs, which prevents crashes in the Maps software application.
MAPS-12980	The naming convention has been updated to align with the labeling of tiles in Maps. Tiles on Phenom are now labeled in the same manner.
MAPS-12995	The interface between Maps and ChemiSEM has been extended, so that both Regular and Object-based segmentation types are properly recognized. Selecting Regular segmentation in Maps now works as expected.

Overview

This document describes the Maps 3.32.1 application software used on SEM/SDB and TEM systems.

Topics in this section include:

Introduction to Maps

Terminology

System Requirements

Data Storage Requirements

Project Directory

Package options

Introduction to Maps

The Maps software application provides users with the ability to focus on areas of interest within a sample region by organizing the sample into tiles and then navigating, selecting, and overlapping tiles so they can be viewed and analyzed in a higher level of detail than previously available.

The application will:

- Create a grid of a sample to enable efficient navigation and previews of selected tiles.
- Acquire tiled images in an automated fashion from multiple sample locations, with correlation between multiple layers.
- Find stitching information between tiles in an automated or semi-automated fashion using cross-correlation.
- Stitch and blend tiles into larger images with the capability to export them into common industry formats for viewing.
- Provide the capability to interactively display and investigate images constructed from multiple locations and layers of the sample.
- Process powerful workflows to do alignments or array tomography.
- Use the mineralogy workflow to collect and analyze EDS and mineralogy data from geological samples.

Maps complements the Correlative Workflow features that enable you to move samples between light microscopes and electron microscopes with the help of automated alignment on supported sample holders and a suite of powerful correlation tools.

Maps provides a solution for correlative microscopy by facilitating workflows that involve optical microscopy, SEM, and TEM. Both automated and manual alignments are used for correlation of the data to truly understand the sample

i Note: When upgrading from version 1 or 2 of Maps to version 3, a new license key is required.

Terminology

Common terms used in Maps and Array Tomography are defined below.

Maps Terms

Tomography Terms

Maps Terms

Common Maps terms are listed below.

Alignment

In Maps, adjusts project data to the current stage position.

Blending

The Maps software will create a smooth transition between images during stitching.

Correlation

In Maps context, correlation is used for two things: 1. finding back areas of interest that show up in different modalities, and 2. for visually overlaying data sets that correspond to the same area on the sample.

EM

Electron microscopy

FIB

Focused Ion Beam

Holder Calibration

This procedure is used to adjust the expected fiducial positions for automated holder alignment on individual systems. It must be performed the first time alignment is run on a supported holder. See the Automated Holder Alignment section on Calibration.

Image Registration

The computational process Maps uses to determine the position of a single tile within the greater tile set. The Image Registration process is run during tile set acquisition.

LM

Low Magnification mode on TEM.

Maps Offline Viewer

A version of Maps that has read-only capabilities for Maps projects.

MIP

Maximum intensity projection. Collapses a Z-Stack of images into a single image using the highest intensity pixels from the stack.

Nav-Cam

Navigation Camera

SDB

Small DualBeam; a family of microscopes that have both scanning electron and ion columns.

Segmentation

In the context of images, segmentation is the process of assigning a label to specified pixels in an image such that pixels with the same label share certain visual characteristics, such as color.

SEM

Scanning Electron Microscope

Stitching

The computational task of combining a set of individual images to form a single image.

STEM

Scanning Transmission Electron Microscopy. STEM mode exists on SEMs, SDBs and TEMs (depending on the exact configuration).

TEM

Transmission Electron Microscope

TIA

On TEM, the instrument user interface consists of two parts: TEM User Interface for instrument control and TEM Imaging and Analysis for acquisition and analysis.

Tile Acquisition

The automated task of acquiring, aligning, and registering tiled images, as defined by the user.

Tile Grid

The empty grid that defines where and how a set of tiles will be acquired.

Tile Set

The output of a tile grid once acquisition is run, or the set of actual tiled images.

xT Software

Platform software for the SEM/SDB microscopes.

xT UI

xT software user interface

Z-Stack

A sequence of images through the z-axis, used to build a 3D volume.

Z-Stack Browser

The Maps UI component that allows you to view and select image planes from a z-stack.

Tomography Terms

Common Array Tomography terms are listed below.

Classic Ribbons

(Array Tomography) Sample sections are attached to each other and the ribbon is always slightly bent. The angle between each section is constant, for the most part.

Serial Section Array

The layer group that contains tile set arrays and section outlines. The purpose of the group is to easily contain most of the array tomography data about a sample in one section of the Layer tree.

Tape Ribbons

Sample sections are regularly spaced and the angle between each slice varies slightly.

Tile Set Array

A collection of tile sets that all share the same acquisition parameters, but can differ in location and rotation.

Workflow

A series of steps that are necessary to complete a task.

System Requirements

- 3 GB of physical RAM
- 4.5 GB of memory paging allocated in Windows®
- Dual-core processor
- For Microscope Operation, microscope system software (xT, TEM Server, or LA) must be running before launching Maps

Data Storage Requirements

6

You can select a default location for project data. We suggest that you place the Project Data Directory on the D drive or a supplemental data storage device such as a network drive, external USB drive, or additional internal hard drive.

- You may need additional internal or external storage capacity to acquire large tile sets.
- Acquisition speed might be reduced when saving project data to USB or network locations.

The table below indicates the number of tile sets you can acquire at various image resolutions and dimensions, provided that the Project Data Directory is configured to save to a standard empty 350 GB drive partition.

Note: For multiple channel acquisitions, the storage requirements increase incrementally with the number of channels.

Data Storage Requirements for Single Channel Image Acquisition				
Resolution	Tile Set Size			
	5x5	10x10	50x50	100x100
512	15 MB	100 MB	1.25 GB	5 GB
1 K	55 MB	200 MB	5 GB	20 GB
2 K	200 MB	700 MB	20 GB	75 GB
4 K	750 MB	2 GB	75 GB	250 GB
8 K	4 GB	10 GB	250 GB	1.00 TB
Color Key:				
Red	< 1 Tile Set			
Orange	1 to 2 Tile Sets			
Yellow	2 to 5 Tile Sets			

Data Storage Requirements for Single Channel Image Acquisition(Continued)			
Resolution	Tile Set Size		
Green	> 4 Tile Sets		

Project Directory

At the root directory, you will see the following:

퉬 LayersData	4/12/2016 6:58
鷆 MetaData	4/13/2016 7:08
MapsProject.xml	4/13/2016 7:08
MapsProject.xml.backup	4/13/2016 7:07
Project.mapsxml	4/12/2016 6:58
🗎 ProjectLog.log	4/13/2016 7:08
)

The LayersData folder contains a directory structure that mimics the Layer tree in Maps, so you can manually locate the image files.

Launch Maps from a New File in the Project Directory

Double-click **Project.mapsxml** to launch Maps and automatically open the project from the directory in which this mapsxml file exists.

Package options

The table below lists all the package options available to you with the Maps software. To view your installed licenses and their expiration date(s), click the **Help** menu, and then click **About**. The expiration date(s) are determined by your product license key(s).

Contact your Thermo Fisher representative for licensing information.



Maps Package Options			
Option	Description		
Analytics for TEM	Licensed to run the Analytics plug-in and acquire EDS data through a first- party device interface (FDI). See TEM Analytics for more information.		
Analytics for SEM	Licensed to run the Analytics plug-in and acquire EDS data through a software application interface on a SEM/SDB microscope. See SEM Analytics for more information.		
Annotations	Licensed to create annotations in the project.		
Array Tomography	Licensed to enable the Array Tomography workflow. See Array Tomography Workflow for more information.		
Big Snapshot	Licensed to acquire high resolution square images (4k and 8k) on supported xT microscopes.		
Color Correlation	Licensed to allow correlation features including segmentation and color maps.		
External Image Import	Licensed to allow importing of images from any source such as other microscopes, cameras, CAD exports, etc.		
Maps Offline Viewer	Installations of Maps Offline Viewer are licensed only for viewing datasets that have been generated by the full version of Maps.		
Maps Min	Licensed to set up and acquire mineral EDS data with automated mineral classification post processing. See Maps Min for more information.		
Phenom	Licensed to use the Maps application software on Phenom desktop SEM systems.		
Project Editing	Licensed to edit projects. Without this projects are opened read-only.		
Stage Correlation	Licensed to allow 1-3 point alignment.		
Stitching	Licensed to allow stitching of acquired tile sets.		
TEM Online	Licensed to run on TEM family of microscopes, such as Titan, Talos, etc.		

	Maps Package Options (Continued)
Option	Description
xT Online	Licensed to run on the xT family of SEM and SDB microscopes, such as Helios NanoLab™, Quanta™, Versa, and so on.

Maps User Interface

This section describes the Maps software user interface.

Topics include:

Default Shortcut Keys Viewer Shortcut Keys Main Screen Menus Toolbar Layer/Holder Definition Control Layer Tree Tabs Layer Control Job Control Viewer Rotation Control in the Viewer

Default Shortcut Keys

This topic lists default shortcut keys in Maps.

Shortcut Keys			
Keys	Description		
F1	Opens the current Maps help system.		
F2	Saves the screenshot with an automatically incremented file name. Same as selecting File > Save Screenshot.		
Space	View all objects (zoom out).		
Ctrl+Shift+C	Centers the view on the current layer, same as Center View on Selection and Center View .		
Ctrl+G	Shows/hides grid lines for the current layer, same as Show Grid Lines (Tile Set right-click menu selection).		
Ctrl+H	Hides all graphic elements in the viewer, leaving only the images. Same as selecting Options > Hide Annotations .		
Ctrl+L	Opens the Log Viewer. Same as selecting Help > Open Logs .		
Ctrl+O	Opens the project folder. Same as selecting File > Open Project Folder.		
Arrows	Moves selected tile set in small increments.		
Shift + arrow	Moves selected tile set in finer increments.		
Ctrl + click or drag	Selects tiles.		
Shift + click	Selects tiles by row.		

Viewer Shortcut Keys

There are three profiles defined for Viewer's shortcut keys (Application settings).

- **Default Profile**
- Classic Maps Profile
- Avizo Profile

Default Profile

Viewer Default Profile Shortcut Keys		
Keys	Description	
Ctrl + mouse wheel +/-	Zooms in and out.	
Mouse wheel	Zooms in and out slowly.	
Shift + Ctrl + left-click and drag	Zooms the image to the selected region.	
Left-click and drag	Pan (move)	
Shift + left-click and drag	Rotates the viewer without moving the stage.	
Alt + left-click and drag	Draw a feature of interest rectangle with a number of available options.	
Left-click	Selects an object.	
Ctrl + left-click	Selects multi objects.	
Escape	Deselects all objects.	
Shift + left-click	Selects tiles by row.	
L + left-click and drag	Draw a line measurement tool.	
A + left-click and drag	Draw an angle measurement tool.	

Classic Maps Profile

Viewer Classic Maps Profile Shortcut Keys		
Keys	Description	
Ctrl + mouse wheel	Zooms in and out.	
Shift + mouse wheel	Zooms in and out slowly.	
Shift + Ctrl + left-click and drag	Zooms the image to the selected region.	
Shift + left-click and drag	Pan (move)	
Shift + right-click and drag	Rotates the viewer without moving the stage.	
Left-click and drag	Draw a feature of interest rectangle with a number of available options.	
Left-click	Selects an object.	
Ctrl + left-click	Selects multi objects.	
L + left-click and drag	Draw a line measurement tool.	
A + left-click and drag	Draw an angle measurement tool.	

Avizo Profile

This profile defines the input as in Avizo® software.

Viewer Classic Maps Profile Shortcut Keys		
Keys	Description	
Ctrl + middle-click and drag	Zooms in and out.	
Mouse wheel	Zooms in and out slowly.	

Viewer Classic Maps Profile Shortcut Keys(Continued)		
Keys	Description	
Shift + Ctrl + left-click and drag	Zooms the image to the selected region.	
Middle-click and drag	Pan (move)	
Left-click and drag	Rotates the viewer without moving the stage.	
Ctrl + Left-click and drag	Draw a feature of interest rectangle with a number of available options.	
P + Left-click	Selects an object.	
Ctrl + left-click	Selects multiple objects.	
L + left-click and drag	Draw a line measurement tool.	
A + left-click and drag	Draw an angle measurement tool.	

Main Screen

This topic shows and explains the Main screen of the user interface.

Maps User Interface, Control Panel



- 1. Menus
- 2. Toolbar
- 3. Layer Control
- 4. Overview Viewer
- 5. Tile Set Tab
- 6. Visualization Tab
- 7. Viewer
- 8. Micron Bar

Show or hide main controls

Click the last icon on the left side of the tool bar to hide the main controls. Click the icon again to show the main controls.



Microscope connection states

The bottom of the main screen contains two status icons that indicate the connection states between the Maps application and the microscope.

Green connection states include:



- Solid Green: If the status icon is solid green, then the connection is active and you can use the Maps application online.
- **Spinning Green**: If the status icon is green with a spinning wheel, then the connection is currently occupied with a task. To use the Maps application online, wait until the status icon becomes solid green.

Red connection states include:



- **Spinning Red**: If the status icon is red with a spinning wheel, then the connection is currently unavailable and the Maps application is attempting to secure the online connection.
- **Solid Red**: If the status icon is solid red, then the connection could not be established and the Maps application is no longer attempting to connect to the microscope. Click the status icon to initiate a new online connection.

Menus

This section describes options listed in Maps software menus.

Topics include:

File menu

View menu

Phenom menu

Options menu

Help menu

Tomography menu

File menu

This section describes options listed in the File menu.

File		
	New Project	
	Open Project	
	Import Project	
	Project Properties	
	Open Project Folder	Ctrl+O
	Import Grid	
	Import Images	
	Save Screenshot As	
	Save Screenshot	F2
	Quit	

	File Menu Overview
Menu Selection	Description
New Project	Displays the New Project window. See New Project for full instructions.

	File Menu Overview(Continued)		
Menu Selection	Description		
Open Project	Displays the Project History window showing all previous projects. See Open a Project for more information.		
Import Project	Displays the Import Project window. Browse to a Source directory and select a project, rename it if appropriate, and then click Import . The imported project name will appear in the Project History list. This can be used for passing projects between computers.		
	Ma Import Project		
	Source directory:		
	C:\Data\Project Demo 2000		
New project name:			
	Project Demo 2000		
	Import Cancel		
	See Import a Project for full instructions.		
Project Properties	Displays the Project Properties window for selecting the Project Path, Sample Holder, and a project description. See Project properties for more information.		
Open Project Folder	Displays the project folder. This can also be accessed by selecting Open Folder in the Project Properties window.		
(001.0)			
Import Grid	Creates a blank tile set that matches the settings of a previously exported tile set. An exported tile set is saved as a .gridx file.		

File Menu Overview(Continued)				
Menu Selection	Description			
Import Images	Opens a standard open file dialog that enables the user to import a image file into the current Maps project.			
Save Screenshot As	Displays a standard Windows Save As dialog box for saving a screen capture of the Maps viewer. Used for selecting a file name and a file type for saved viewer screenshots.			
Save Screenshot	Saves the screenshot with an automatically incremented file name. Briefly displays a dialog box with the saved (automatically incremented) file name.			
(F2)	Screenshot saved X Image: Screenshot saved to C:\Users\User\AppData\Loca\FEI\Maps\Demo project\screenshot 001.tif Image: Screenshot saved to C:\Users\User\AppData\Loca\FEI\Maps\Demo project\screenshot 001.tif			
Quit	Closes the Maps software.			

Project properties

View the attributes of the current project and modify the selected sample holder and project description here.

Ma Project properties			×
Project Name:			
Demo project			
Project Path:			
C:\Maps Projects\Demo project			
Disk Space Usage:			
279.56 MB (293,141,898 bytes)			
Total Number of Images:			
25 tiles, 2 imported images			
Project Description:			
Sample Holder:			
3mm Grid holder			~
ក្រសួសស្នា	Holder:	3mm Grid holder	
_	Version:	1.1	
	Manufacturer:	FEI	
	Part number:	1046764	
	Ok	Apply	Cancel

Project Properties Overview			
Interface Item	Description		
Project Name	Displays the name given to the project.		
Project Path	Displays the folder and location of the selected project.		
Disk Space Usage	Displays the amount of disk space used by the project.		
Total Number of Images	Displays the number of tiles and imported images for the project.		
Project Description	Displays the description of the current project.		
Sample Holder	Displays a dropdown list of possible sample holders that support automated alignment on SEM/SDB microscopes.		
OK	Applies the selections and closes the window.		
Apply	Applies the selection only.		

	Project Properties Overview(Continued)
Interface Item	Description
Cancel	Cancels any changes to the Project Properties and closes the window.

View menu

This section describes options listed in the View menu.

All selections

	View	
	Hide All Annotations	Ctrl+H
~	Show Layer Name Annotations	
	Show Annotations on Top	
~	Animate Pan & Zoom	
	Presentation Mode	

View Menu, All Selections Overview				
Menu Selection	Description			
Hide All Annotations (Ctrl+H)	Hides all graphic elements in the viewer, leaving only the images.			
Show Layer Name Annotations	Displays an additional annotation of the tile set names to help identify data during viewing. This option is activated by default and persists when you restart Maps. To deactivate this feature, click the Show Layer Name Annotations menu item to clear the check mark associated with it.			

View Menu, All Selections Overview(Continued)				
Menu Selection	Description			
Show Annotations on Top	When activated, this menu option displays annotations on top of the image data. Annotations include areas of interest; sites of interest; and line and angle measurements.			
	Note: This option persists when you restart Maps.			
Animate Pan & Zoom	Allows a smoother transition of panning and zooming. Note that this is a slower process and takes more memory. This option is activated by default and persists when you restart Maps. To deactivate this feature, click the Show Layer Name Annotations menu item to clear the check mark associated with it.			
	• Pan : To make a sweeping movement			
	 Zoom: To simulate movement rapidly away from or toward a subject using a zoom lens 			
	The "Center View on Sample," "Center View on Selection," and "Go to Microscope FOV" viewer functions animate pan and zoom by default.			
Presentation Mode	This option is available when you have two monitors. When activated, a second full screen view of the project data displays on the other screen. This is useful for demos and presentations.			

Phenom menu

The **Phenom** menu is only available for Maps software on Phenom desktop SEM systems.



Phenom Menu Overview			
Menu Selection	Description		
Phenom instrument	Opens the Phenom instrument dialog. The dialog shows details of the connected Phenom microscope. You can also use the dialog to disconnect from the microscope. See Phenom Connectivity for more information.		

Options menu

This section describes options listed in the **Options** menu.

System views

The common selections are described in the table below. For system-specific menu selections outlined in purple, click the following links.

- SEM/SDB Options Menu
- TEM Options Menu

Options Menu, Specific Selections by System

SEM/SDB View



TEM View



Ontions Menu All Selections Overview				
Menu Selection	Description			
Application Settings	Displays the projects.	Application Settings window f	or changing de	efaults for all new
	Application	Default project data directory:		
	Application	D:\Data Sets\Projects		
	XPS	Appearance		
	Mineralogy	Stage cursor color: Image: Color: Measurements default color: Image: Color:	Site/area of interest defaul	lt color:
	Velox	Viewer		
	Сгуо	 Display 360° Scan Rotation Viewer Fonts 	Viewer Profile:	Default 🗸
	Stitching	Times New Roman 🗸	۵aF	3677123
	Internal	14 v B <i>I</i>		
		Correlation Default Holder: none Histogram white tail: 1% Histogram black tail: 1% Screen Saver 2 ✓ Wait: 2 Improvement Program ClientID: ? Allow usage tracking	is busy and there is no user ir g the remaining time and late 1A2B3C4D5E6F7G89 Apply	Ateraction, a screen saver will est image acquired. OK Cancel
	Default Project Data Directory : Displays the default path with a browse button to allow you to choose a different path.			
	Stage cursor color: Click the eyedropper to select a new stage cursor color. Measurements default color: Click the eyedropper to select a new color for line and angle measurements.			
				ect a new color

Options Menu, All Selections Overview(Continued)				
Menu Selection	Description			
	Site/area of interest of default color for sites of	lefault color : Click the of interest and areas o	eyedropper to select a new f interest.	
	 Display 360° Scan Re When selected, dis When not selected 	otation: splays 360° scan rotat I, displays -180° to 180	ion.)° scan rotation.	
	Viewer Profile: Displays a list of viewing choices.			
	Viewer Profile:	Default Default Avizo Classic Maps		
	Choices include:			
	Default: See Default Profile for detailed inputs.			
	 Avizo: The inputs defined for the viewer are based on Avizo/Amira® software. See Avizo Profile for detailed inputs. 			
	 Classic Maps: The inputs defined for the viewer are based on the previous Maps version (v2.5). See Classic Maps Profile for detailed inputs. 			

	Options Menu, All Selections Overview(Continued)		
Menu Selection	Description		
	Viewer Fonts:		
	 Note: Changes to Viewer Font values affect the viewer scale bar and layer name annotations. These changes are applied to newly created annotations or measurements. Existing annotations and measurements displayed within the Maps project are not impacted by these font changes. 		
	The first field displays a list of available fonts for the annotation and measurement tool description. Select a font from the dropdown menu to designate is as the default font. Times New Roman ✓ Sylfaen Σψμβολ Tahoma Times New Roman		
	The second field displays a list of font sizes for the annotation and measurement tool description. Select a font size from the dropdown menu to designate it as the default font size. 14 ✓ 14 ✓ 15 16		
	 Bold: Click the B button to designate bold as the default font weight for the annotation and measurement tool description fonts. Italic: Click the <i>I</i> button to designate italic as the default typeface for the annotation and measurement tool description fonts. 		

Options Menu, All Selections Overview(Continued)			
Menu Selection	Description		
	Default Holder : Displays a list of possible sample holders. Select a holder to be used as a default for all new Maps projects.		
	Default Holder: none Histogram black tail: none 3mm Grid holder 8 Block Holder		
	14x30mm Round Quanta Moun		
	Histogram white tail: Trims a percentage of histogram pixels to eliminate noise.		
	Histogram black tail: Trims a percentage of histogram pixels to eliminate noise.		
	Screen Saver: Select the checkbox to display a screen saver when the application is busy for the specified number of minutes.		
	Improvement Program : Shows the client ID for tracking your anonymous usage data that the Maps application sends back to the Maps team. The "Allow usage logging" setting is activated by default. To opt out of the improvement program, select the slider and move it to the left to deactivate the setting (a dark grey slider is deactivated).		
Tile Set Templates	Displays the Tile Set Templates window. See Tile Set Templates for a full explanation of templates.		
	Tasks related to tile set templates include:		
	Default templates		
	Edit custom templates		
	Import and export tile set templates		

Options Menu, All Selections Overview(Continued)		
Menu Selection	Description	
Microscope Settings	These are system-specific.SEM/SDB Options MenuTEM Options Menu	
Snapshot Settings	These are system-specific.SEM/SDB Options MenuTEM Options Menu	
XPS Settings	These are system-specific. See TEM Options Menu.	
Cryo Settings	These are system-specific. See SEM/SDB Options Menu.	
Amira-Avizo2D Settings	See Amira-Avizo2D Connectivity.	

Help menu

This section describes options listed in the Help menu.


Help Menu Overview		
Menu Selection	Description	
Help (press F1)	Displays an interactive help system that contains release notes, installation information, and user guide topics. PDF versions of the content are accessible from within the help system.	
Generate Bug Report	Generates a bug report and opens the folder with the generated bug report.	
Open Bug Report Folder	Displays the folder where the generated bug reports are stored.	
Open Maps Min Help Desk	Opens the Maps Min Help Desk for online support (see Maps Min). You can request/renew a license or report bugs through the help desk.	
Open Application Logs (press Ctrl+L)	Displays the Log Viewer dialog.Image: Colspan="2">Image: Colspan="2">Image: Colspan="2">Image: Colspan="2">Image: Colspan="2">Image: Colspan="2">Image: Colspan="2">Image: Colspan="2">Image: Colspan="2"Image: Colspan="2"<	
About	Opens the About screen, which displays information about the software versions and a list of your installed license options with expiration dates. For descriptions of the license options, see Package options.	
License Terms	Displays the accepted end-user license agreement (EULA).	

Tomography menu

When Maps is installed with the Thermo Scientific[™] Tomography 5 Software application, the Tomography menu is available.

Tomography menu

Click **Tomography** to view the menu items, which include Import Presets and Import Batch Positions.

File	View	Tomography	Options	Help
		Import Pre	esets	
		Beta Import Bat	tch Positions	

Tomography Menu Overview		
Menu Selection	Description	
Import Presets	Imports Tomography microscope presets. The presets are stored as Maps templates to be used later for acquisitions and autofunctions.	
(Beta) Import Batch Positions	Imports batch positions from Tomography with relevant Exposure, Tracking, Focus, and Condition areas.	

Tomography menu with Maps Data Exchange Service

 Note: If you also have the Maps Data Exchange Service installed with the Tomography application, then the Tomography menu contains different options than if the service is not installed. For information on how to install the Maps Data Exchange Service, refer to the latest Maps Software Installation Guide.

Click **Tomography** to view the menu items, which include Import Presets, Synchronize With Tomography, and Send To Tomography.



Tomography Menu with Maps Data Exchange Service Overview		
Menu Selection	Description	
Import Presets	Imports Tomography microscope presets. The presets are stored as Maps templates to be used later for acquisitions and autofunctions.	
(Beta) Synchronize With Tomography	Toggles the automatic synchronization of data with Tomography between ON and OFF.	
	When synchronization is active, the following data are automatically exchanged between Maps and Tomography:	
	 Annotations are sent to Tomography, including areas of interest, sites of interest, and lamella sites. 	
	 Batch positions are imported from Tomography, including exposure, tracking, focus, and condition areas. 	
(Beta) Send To Tomography	Sends all annotations of given type to Tomography when live synchronization is turned OFF.	
	Select the annotation type from the submenu:	
	Send All Lamellae	
	Send All Areas of Interest	
	Send All Sites of Interest	

Toolbar

The common toolbar selections are described in the following table.

For system-specific selections, see:

- SEM/SDB Toolbar, Right Side
- Phenom Toolbar, Right Side

Topics in this section include:

Left Side, All Systems

Secondary Menu

Toolbar Descriptions

Left Side, All Systems

The left side toolbar appears as below on all systems.

Toolbar, Left Side, All Systems



Secondary Menu

If the window is resized too small to display all tools, some of them are stored in a secondary menu. Click the down arrow to access them.



Toolbar Descriptions

	Toolbar Descriptions
Tool	Description
Move & Resize Tool	Moves or resizes the tile grid with click and drag.

	Toolbar Descriptions(Continued)
ΤοοΙ	Description
Measurement Tool	Draws a line and displays the length as a callout.
Angle Measurement Tool	Draws a line and a third point; displays the angle as a callout.
Zoom to Selection	Zooms the field of view to the selected area.
Add Tile Set	Adds a tile set over a selected area in the viewer.
Add Circular Tile Set	Adds a circular tile set over a selected area in the viewer.
Add Freehand Tile Set	Adds a freehand drawn tile set over a selected area in the viewer.
Stamp Tile Set	Stamp a new tile set using the active template.

Toolbar Descriptions(Continued)		
Tool	Description	
Add Area of Interest	Displays the New Annotation dialog box for adding the drawn selected area of interest (AOI).	
	Ma New Annotation Name: Area of interest Focus: 5 mm Z: 100 µm AT: 0 OK Cancel The new AOI becomes a new Annotation layer in the Layer control.	
Save Image to File	Saves an image to file. See "Save Image to File" on page 69.	
Stitch Selection	Stitches a subset of tiles, as selected by the rectangle. See "Stitching" on page 99.	
Center View on Sample	Zooms out to the full sample view. By default, this view animates the pan and zoom function to orient you within the sample space. To deactivate the animation functionality, see the "Animate Pan & Zoom" under View menu.	
Center View on Selection (Ctrl+Shift+C)	Unavailable until a project is open. Pans and zooms the viewer so that the currently selected layer item is centered. By default, this view animates the pan and zoom function to orient you within the sample space. To deactivate the animation functionality, see the "Animate Pan & Zoom" under View menu.	
Go to Microscope FOV	Pans and zooms the view to match the microscope (electron beam) field of view (FOV). By default, this view animates the pan and zoom function to orient you within the sample space. To deactivate the animation functionality, see the "Animate Pan & Zoom" under View menu.	

	Toolbar Descriptions(Continued)	
Tool	Description	
Save Screenshot (F2)	Briefly displays a dialog box with the saved (automatically incremented) file name. Also available from File > Save Screenshot .	
Copy Screenshot to Clipboard	Copies the current image display to the system clipboard. You can then paste the image into other applications using paste commands.	
Global Alignment	Performs a 1-, 2-, or 3-point alignment to align all previously acquired data to the stage. It is useful if the sample was removed from the tool and you want to continue with it. See Coarse Alignment (1-, 2-, or 3-Point).	
Align to Holder (SEM/SDB only)	Performs an Automated Holder Alignment; requires a specific holder hardware. If no sample holder has been selected for the project, a dialog box appears to prompt you to select a holder. Ma Holder Selection × A sample holder must be set in the project before continuing. Please select a sample holder from the list below.	
	Holder: Single 22 mm coverslip holder Version: 1.0 Manufacturer: FEI Part number: 1101666 OK Cancel	

Layer/Holder Definition Control

This feature is only available on SEM/SDB systems (see Holder Tab).

Layer Tree Tabs

This section describes the tabs within the layer tree depending on the selected item such as a specific tile set, line, or angle. For example, the image below shows the Tile Set (20) and its associated tabs.

Topics include:

Image Tab

Tile Set Tab

Annotation Tab

Measurement Tab

Font Format Tab

Visualization Tab

Overview Viewer



Image Tab

The Image tab appears in the left panel when you select an image layer in the layer tree view. This tab displays properties of images, including imported images and stitched layers.



Control	Description
Name	Displays the name of the current image.

Control	Description	
Planes	Controls the image planes. Use the slider to change the current plane index and to show the plane in the main viewer.	
	You can also click the left and right arrows to navigate one plane at a time.	
	Note: The control is activated only for images with multiple planes.	
	 Note: If the image plane depth in the current image is available, then the Depth control is displayed instead of the Plane control (see Optional controls). 	
Current Plane	Displays the current plane index.	
Time Frames	Controls the image time frames. Use the slider to change the current frame index and to show the frame in the main viewer.	
	You can also click the left and right arrows to navigate one time frame at a time.	
	Note: The control is activated only for images with multiple frames.	
Current Time Frame	Displays the current frame index.	
Pixels	Displays the resolution of the current image.	
Pixel Size	Displays the physical size of a single pixel within the current image.	

Control	Description
Physical Size	Displays the physical size of the current image.
Show metadata	Opens a screen to display the metadata of a selected item in the layer tree (see Metadata Viewer). If there is not any metadata associated with the selected item, then this button is deactivated.

Optional controls

If the image plane depth in the current image is available, then the Depth control is displayed instead of the Plane control.

□ ≉ ⊘	
Name	DICOMDIR
Depth	
7.41 µm 🖪 🦲	3.67 mm
Plane	441 of 495 (3.27 mm)
Pixels	1900 px x 2048 px
Pixel Size	5.26 nm x 5.26 nm
Physical Size	10 μm x 10.78 μm
	Show metadata

Control	Description
Depth	Controls the plane depth of the current image. Use the slider to change the depth, to change the plane, and to show the plane in the main viewer. You can also click the left and right arrows to navigate one depth plane at a time.
Plane	Displays the current plane index and its depth.

Metadata Viewer

This topic describes the Metadata Viewer, which displays the information associated with the selected item from the layer tree such as images, tiles, and tile sets. For example, the metadata for a stitched image contains the original microscope information.



UI Item	Description
Title	Displays the name of the layer tree item associated with the metadata. The screen title is located in the upper-left corner of the viewer.
Metadata tabs	Displays the metadata of the element and is organized by the type of content. These tabs are located in the upper-left corner of the viewer.
Copy to clipboard	Copies the tab content to the clipboard, so that you can paste it to another text editor if needed. This button is located in the upper-right corner of the viewer.

Tile Set Tab

When you create a new tile set, the parameters for basic tile acquisition are populated with defaults.

III 🕂 🕂 🗄]	ĄŶĄ	۲		
Name		Tile Set	t (3)		
Acquisition Type		Electro	n		~
Tiles X, Y	a	10		10	
Tile HFW	G	1 mm			
Total HFW	G	9.1 mm	ı		
Resolution		2048 x	1768		~
Pixel Size		488.28	13 nm		
Dwell		300 ns			~
Frames		1			
Reduced Area			Select		
Image acquisition time: 00:00:04 Estimated acquisition time: 00:04:11 Elapsed acquisition time: 00:03:36 158 of 200 images acquired; 1.71 GB					
FROM MICROSCOPE			TO MICR	OSCOPE	

The selections are system specific:

- SEM/SDB Tile Set Tab
- TEM mode: TEM Tile Set Tab
- STEM mode: STEM Tile Set Tab

Annotation Tab

This Annotation tab contains the properties for the annotations within the main viewer. This tab appears in the left panel when you select either a *site of interest* or an *area of interest*.

Site of interest

Name Site of interest Notes		
Name Site of interest Notes		
Notes Color Ellipse Stage X -21.0649 mm Stage Y 11.6892 mm Source System SIMULATION Drive To	Name	Site of interest
Color Ellipse Stage X -21.0649 mm Stage Y 11.6892 mm Source System SIMULATION Drive To	Notes	
Ellipse Stage X -21.0649 mm Stage Y 11.6892 mm Source System SIMULATION Drive To	Color	C.
Stage X -21.0649 mm Stage Y 11.6892 mm Source System SIMULATION Drive To		Ellipse
Stage Y 11.6892 mm Source System SIMULATION Drive To	Stage X	-21.0649 mm
Source System SIMULATION Drive To	Stage Y	11.6892 mm
Drive To	Source System	SIMULATION
		Drive To

Property	Description
Name	Displays annotation name.
Notes	Displays annotation notes. Click in this field to add or edit these notes.
Color	Displays color of annotation. Click the color picker (that is, the eyedropper icon) to change the color.
Ellipse	Appears if the annotation is displayed with an ellipse form within the main viewer. Select this checkbox to change the value.
Stage X	Displays position of stage for X axis.
Stage Y	Displays position of stage for Y axis.

Property	Description
Source System	Displays name of source system.
Drive To	Click button to drive microscope field of view to annotation location.

Area of interest

If the annotation is an *area of interest*, then the opacity property is shown in addition to the properties described in the previous table.

Name	Area of interest
Notes	
Color	l de la companya de l
Fill On a site	
	Ellipse
Stage X	Ellipse
Stage X Stage Y	Ellipse -7.1668 mm 5.9121 mm
Stage X Stage Y	Ellipse -7.1668 mm 5.9121 mm
Stage X Stage Y Source System	Ellipse -7.1668 mm 5.9121 mm Scios 2
Stage X Stage Y Source System	Ellipse -7.1668 mm 5.9121 mm Scios 2 Drive To

Property	Description
Fill Opacity	Sets the level of opaqueness for the entire annotation. Click and move the slider to the right to increase the level of opaqueness, or click and move the slider to the left to decrease the level of opaqueness. The less opaque, the more transparent the annotation becomes. Transparent annotations are useful when overlaying a layer over another layer.

Measurement Tab

The Measurement tab appears in the left panel when you select a line or angle measurement layer. This tab displays properties of measurements.



Property	Description
Name	Displays measurement name.
Notes	Displays measurement notes. Click in this field to add or edit these notes.
Color	Displays color of measurement. Click the color picker (that is, the eyedropper icon) to change the color.
Fill Opacity	Sets the level of opaqueness for the entire measurement. Click and move the slider to the right to increase the level of opaqueness, or click and move the slider to the left to decrease the level of opaqueness. The less opaque, the more transparent the measurement becomes. Transparent measurements are useful when overlaying a layer over another layer.

Font Format Tab

Use the controls under the Font Format tab to set layer font properties as rendered within main viewer. This tab appears only when you select a site of interest or when you select a line or angle measurement layer.



Property	Description
Font Family	Displays a list of available fonts. Select a font from the dropdown menu to be displayed within the main viewer. Times New Roman Σψμβολ Tahoma Times New Roman
Font Size	Displays a list of font sizes. Select a font size from the dropdown menu to be displayed within the main viewer. 14 14 15 16
Bold	Click the B button to display the font weight as bold within the main viewer.
Italic	Click the <i>I</i> button to display the font style as italics within the main viewer.
Font Viewer	Provides a rendering of the selected font format properties. This viewer is located to the right of the other four properties.

Visualization Tab

The Visualization tab (formerly called "Results" tab) contains the correlation controls, which you can use to apply simple coloring to the tile set images when acquired. Some correlation controls appear unavailable if they are not applicable to the system type.

Correlation controls

III - AE II 😨	t Ø
Opacity	•
Additive Opacity	
Stitching Channel	Stem_HAADF 🗸 🗸
	ப்பகிலை histogram
int net	wt% at%
> ✓ Stem_HAADF	•
🗲 🔽 Pd (At)	
> Nb (At)	
🗲 🗹 Mn (At)	•

Correlation Controls Overview

Control	Description
Opacity	Sets the measure of opaqueness for the entire image. The less opaque, the more transparent it becomes. Transparent images are useful when overlaying on image over another.
Additive Opacity	Sets the control the opacity between fluorescence and regular channels. This is useful for the Analytics tile sets as they contain many non- overlapping channels. The Additive Opacity option is the default for EDS tile sets. Increasing the additive opacity makes fluorescence channels more visible.
Stitching Channel	Sets the channel used for stitching.
Show histogram	Shows/hides the histogram panel. See Histogram.

Correlation Controls Overview(Continued)				
Control	Description			
Quantification Modes	EDS channels acquired with Velox can have up to four different quantification modes. Click one of the activated buttons (shown in white) to select a quantification mode, which is used to visualize the elements. A deactivated button (shown in grey) indicates that the quantification mode was not calculated and is not available.			
	 Int net wt% at% Stem_HAADF The elements display data in each quantification mode as follows: int: the counts according to the raw integrated spectrum net: the counts according to the background corrected and fitted model wt%: the weight fractions at%: the atomic fractions 			



Histogram

Displays a histogram for the selected channel. Click a different channel name to select its histogram and associated color.

Selected Channel Histogram



Selected Channel Histogram Overview			
Interface Item	Description		
Selected channel name	Displays the selected layer item (a tile set or image) from the Layer control.		
Hide	Hides the histogram.		
Zoom	Expands the histogram area of images with less than 255 colors.		
Semi-log	When selected, changes the histogram intensity display to logarithmic.		
Invert	Inverts the intensity of the image, e.g., black to white.		
Auto	When selected, continuously adjusts the min and max for the histogram display.		
Current max position	Displays the current max slider position.		
Max slider	Any data above the max line is white. Gray levels display between the max and min sliders.		
Histogram data	Displays the histogram data for the selected channel.		
Min slider	Any data below the min line is black. Gray levels display between the max and min sliders.		
Current min position	Displays the current min slider position.		

Overview Viewer

The Overview viewer, located in the Overview tab, shows a lower-zoom view of your data, allowing you to see where the main Viewer field of view (FOV) is in relation to the whole project. See Viewer for more information about the main Viewer.



The yellow crosshair shows where the current FOV in the main Viewer is in relation to the entire data set. Double-click in the Overview viewer to move the FOV of the main Viewer.

There are two toolbar buttons:

Button	Description
5	Center view on sample. Click to snap the Overview viewer FOV over all the sample data in the project. This action moves the FOV to the center of the sample and adjusts the zoom, so that it is visible.
þ	Center view on selection. Click to snap the Overview viewer FOV to the currently selected data in the Layer tree. This action moves the FOV to the Layer tree and adjusts that zoom, so that it is visible.

There is one unique context menu selection:

Menu Selection	Description
Drive Stage Here	Moves and centers the stage to the selected tile (see the green cross in the center of the grid).

Layer Control

Use the controls to view existing layer items and create, delete, reorder, select, hide, and show layers.

Topics include:

Layer Icons

Change Layer Order

Progress of Running Layer Tasks

Annotation Context Menu

Layer Context Menu

Tile Set Context Menu

Image Context Menu

Nav-Cam Context Menu

Multiple Layer Context Menu



- When a check box is not selected, that layer item is not shown in the viewer.
- Double-click an item in the Layer control to center it in the viewer.

Layer Control Overview			
Control	Description		
Add Layer	Adds another Layer to the tree. When you have multiple layers, the selected layer will be green and the non-selected layers will be yellow.		
Add Tiles	You can only create a tile set if a microscope is connected.		
:#:	See Define a Tile Set for more information about adding tiles.		
Template	Displays the name of the active template for the connected microscope. In the example above, it is Default Template . See Tile Set Templates for more information about templates.		
Layer	Displays a list of existing tile sets, annotations (areas of interest), imported images (aligned and unaligned), sample holder, and fiducials for each layer.		
	See Layer Icons for icon descriptions.		

Layer Icons

Each created layer item has a unique icon.

Layer Control Icons

✓
🗹 🌐 EMTileSet7
🗹 🟳 Area of interest
🗹 📖 Line
🛃 📐 Angle
🗹 👜 ImportedImage
🗹 🗔 ImportedImage_aligned
👻 🔽 3mm Grid holder
🗹 🗔 Holder outline
Fiducials

Layer Icons			
lcon	Indicates the layer contains		
⊞	Complete tile set registration and acquisition.		
	Annotations (text callouts).		
	Annotations are directly created in the layer selected at the time of creation.		
	Line measurement.		
	Angle measurements.		
•	An imported image that is not yet aligned.		
	An imported image was aligned and stitched into this layer.		
::	A holder outline.		
\oplus	Fiducial information.		

Change Layer Order

Context Menu

Layers can be moved in the layer tree to change their position, depending on their Permissions.

Right-click on a layer in the layer tree (or click the ellipsis icon ... next to a layer) to access the context menu for moving the layer forward or backward.

Bring To Front
Bring Forward
Send Back
Send To Back

Layer Order Context Menu Options			
Menu Selection	Description		
Bring to Front	Moves the layer to the front or top-most layer.		
Bring Forward	Moves the layer forward one layer.		
Send Back	Moves the layer backward one layer.		
Send to Back	Moves the layer to the back or bottom-most layer.		

Move layers

Layers can be moved in the layer tree to change their position, depending on their Permissions. If allowed, you can click and move a layer up or down in the layer tree (or move multiple selected layers by pressing **Ctrl+left-click** before moving them). If a layer is locked, then it cannot be moved within the layer tree.

The mouse changes when dragging a layer to show what action is available:

- When dragging a layer, the cursor changes to a cursor with a rectangle below.
- When the dragged layer cannot be placed in a location, the mouse changes to a no-entry icon.

• When a layer is a valid drop location, the target group highlights with a rectangle :

✓ SnC_OptiPlan	
SnC_5kV_T2 (2) (stitched)	
₩ SnC_5kV_T2 (2)	
SnC_5kV_T2 (stitched)	
✓ Ⅲ SnC_5kV_T2	

• When dropping a layer between another layer, a divider appears:

👻 💟 SnC_OptiPlan	
🗹 🛄 SnC_5kV_T2 (2) (stitched)	
□	
SnC_5kV_T2 (stitched)	
✓ Ⅲ SnC_5kV_T2	×

Permissions

The table below shows if a parent layer can accept a child layer. When working with multiple layer selections, these permissions are applied to each layer selection.

Layer Order Permissions					
		Parent Layer			
		Layer Group	Annotation	Tile Layer	Image Layer
Child	Layer Group	Yes	No	No	No
Layer	Annotation	Yes	Yes	Yes	Yes
	Tile Layer	Yes	No	No	No
	Image Layer	Yes	Yes	Yes	Yes

Progress of Running Layer Tasks

- Acquisition progress is reported in the Job Control.
- Registration progress is reported by hovering the mouse over the tile set symbol.

-	🛃 Layer		
	т 🌐 🔽	ile Set (2)	· · · ·
V III Tile Set			
⊞		Running layer tasks:	
Name		Tile registration	

• When both Registration and Acquisition have completed, the green Tile Set icon appears.

🗾 🌐 Tile Set

Annotation Context Menu

Right-click on an annotation in the layer tree to access a context menu for centering and rotating the view, driving to, and deleting.

All of these menu commands, except **Drive To** and **Save Image to File...**, are common to most all other layer item context menus.



Annotation Context Menu Overview			
Menu Selection	Description		
Center View	Centers the view on the annotation.		
Center and Rotate View	Pans, zooms, and digitally rotates the viewer so that the selected annotation is centered and displayed at its natural point of rotation.		
Drive To	Drives the stage to the annotation. This menu selection is available only while Maps is running from the Microscope PC and thus connected to the microscope.		
Save Image to File	Saves an image from the region outlined by the annotation (see Save image to file under Viewer). This menu selection is only available for an <i>area of interest</i> annotation (see Area of interest under Annotation Tab).		
Delete	Deletes the annotation.		

Layer Context Menu

Right-click on any Layer selection to access a context menu.



- For descriptions of common selections to all systems, see Annotation Context Menu.
- For descriptions of Layer-specific selections, see the table below.

Layer Context Menu Overview		
Menu Selection	Description	
Copy Layer	Copies all tile sets in the layer to a new layer.	
Import Grid	Displays a standard Open window for browsing to a saved tile grid file.	
Import Images	Displays a standard Open window for browsing to a saved image file from another source, such as an optical microscope.	
	navigation guide for you to find areas of interest, like the Nav- Cam does.	
	See Import Images from Other Sources.	
Export To Project	Exports to a new or existing project.	
Alignment	Displays alignment choices. See Magnification Alignment.	
	Submenu selections include:	
	 Align: Aligns a single tile set or image. Similar to Global Alignment, but for a single item. See Coarse Alignment (1-, 2-, or 3-Point). 	
	 Align: Aligns a single tile set or image. Similar to Global Alignment, but for a single item. See Coarse Alignment (1-, 2-, or 3-Point). Fine Alignment: Use Fine Alignment to fine-tune the alignment of layers and tile sets visually. The Fine Alignment controls only display after clicking Next after a 1-, 2-, or 3-point alignment. See Fine Alignment for descriptions of the controls. 	

Tile Set Context Menu

Right-click on any tile set to access the context menu.

	Bring To Front		
	Bring Forward		
	Send Back		
	Send To Back		
	Show Grid Lines	Ctrl+G	— Unique selection for tile set
	Center View	Ctrl+Shift+C	
	Center and Rotate View		
1	Save and Set As Default Template		
(Save As Template		
	Copy Grid		
	Export Grid		
	Export To Project		
	Alignment		
	Drive To		
	Square Up		——— Unique selections for tile set
	Use For Magnification And Scan Rotation Alignment		
	Apply Settings To Microscope		
	Apply Microscope Settings To Grid		
	Check Magnification and Rotation Error		
	Stitch		
	Tile Registration		
	Split Channels		
	Create MIP Layer		
	Mark as done		
	Reset Acquisition		
\backslash	Open Z-Stack		
	Delete		

- For descriptions of common selections to all systems, see Annotation Context Menu.
- For descriptions of tile set-specific selections, see the table below.

Tile Set Context Menu Overview				
Menu Selection	Description			
Show Grid Lines (Ctrl+G)	Toggles to show/hide the tile grid lines onscreen.			
Save and Set As Default Template	After editing the defaults, saves the edits as the new default template. See Default templates for further instructions.			
Save As Template	Displays a submenu New Template , that allows saving the current tile layers template as a new template or overriding a previous template. See Create custom templates for further instructions.			
Copy Grid	Creates a grid above the one that is copied, and incrementally numbers the name. In this example, EMTileSet 5 is a copy of EMTileSet 4.			
Export Grid	Displays a standard Save As window for browsing to a folder in which to save the exported grid. Grids are saved in the *.gridx format.			
Export to Project	Copies the selected tile sets into a new or existing Maps project.			
Alignment	Displays a submenu of alignment choices. See Layer Context Menu for descriptions.			
Drive To	Drives to the selected tile set.			
Square Up	Rotates the stage to match the orientation of the tile set.			

Tile Set Context Menu Overview(Continued)		
Menu Selection	Description	
Use For Magnification And Scan Rotation Alignment	When selected, uses the offsets calculated in the acquired tile set to correct for misalignment between the scan axes and the stage axes by applying a small magnification and scan rotation. This will allow better acquisition and alignment of future tile sets. To activate this menu selection, all tiles must be successfully registered (that is, no tiles outlined in red). You can use a tile set of any size, but it is recommended that you use a tile set of at least (9) tiles.	
Apply Settings to Microscope	Applies all defined tile set settings to the microscope.	
Apply Microscope Settings to Grid	Applies all microscope settings to the tile set.	
Check Magnification and Rotation Error	Displays the calculated error of the microscope magnification and field-of- view rotation based on the tile registration results. To activate this menu selection, all tiles must be successfully registered (that is, no tiles outlined in red). You can use a tile set of any size, but it is recommended that you use a tile set of at least (9) tiles.	
Post processing	The following menu selections are related to post-processing functions. Stitch Tile Registration Split Channels Create MIP Layer Mark as done Reset Acquisition Open Z-Stack	
Stitch	Stitches the entire tile set.	
	Tile Set Context Menu Overview(Continued)	
-----------------------------------	---	
Menu Selection	Description	
Tile Registration	Launch registration on tile set. See	
Split Channels	Divides the individual channels of a tile set into separate tile sets.	
	i Note: This operation is permanent and cannot be undone.	
Create MIP Layer	Collapses the Z-Stack into a single image using the highest intensity pixels from the stack. This menu selection is enabled for tile layers with more than one plane.	
Mark as done	Sets the selected tile set that has started acquisition, but is not yet done. This option deactivates all the remaining non-acquired tiles and designates the acquisition job status as complete.	
Reset Acquisition	Clears all acquired image data in the selected tile set and unlocks the parameter fields in the Basic tab for editing.	
Open Z-Stack (SEM/SDB only)	Opens the Z-Stack Browser to view the focus stack of the tiles. See Z- Stack Browser Controls.	

Z-Stack Browser Controls

The Z-Stack Browser allows you to select and view individual image planes in the Z-Stack.

After acquiring a Z-Stack tile set, right-click on a single tile in the tile set and then click **Open Z-Stack** in the menu that appears.

Maps User Interface

Drive Stage Here		
Rotational Alignment	•	
Snapshot Here		
Snapshot Here w/HFW	•	
Add Tiles Here		
Add Site of interest		
Tile Set (4)	•	Open Z-Stack
Tile 3,3	•	Stitch
		Snap to stage position
Clear Digital Rotation		

The Z-Stack Browser appears on the right-side of the interface.

Z-Stack Browser Open with Controls on the Right



Z-Stack Browser Controls

	Z-Stack Browser Controls Overview
Interface Item	Description
Channel Selector	Selects the channel to be displayed in the multi-channel Z-Stack. The pulldown list contains the same choices as in the Microscope controls.
Histogram	Displays the same histogram as in the Microscope controls. See Histogram.
Close	Closes the Z-Stack Browser.
Plane Selector	Selects the plane from the Z-Stack to be displayed with the channel selection. Move the slider or use the mouse scroll wheel to step through the planes.
Current Plane Control:	Current Plane Select for Display Select for Stitching Reset

	Z-Stack Browser Controls Overview(Continued)
Interface Item	Description
Select for Display	Selects the plane to be displayed in the main viewer after you close the Z-Stack Browser.
	i Note: Only one plane can be displayed in the main viewer.
Select for Stitching	Selects the current plane to be used for the registration step for stitching. Choosing a plane with more image content is best.
Reset	Returns to the default gamma setting.
Go To	Sets the microscopes current focus to the value in the Z-Stack viewer control.
Gamma Control	Adjusts the gamma for the live image.
Tile Set View	Displays the selected tile (with a thick green outline) from the main viewer and its neighboring tiles.
	You can use the mouse wheel to zoom in and out.

-

Tile Registration

The tile registration feature allows you to re-run the registration on a previously acquired tile set.



4. Click Queue Job.

Ma Tile Registration - Define Parameters		
Re-running registration on Overview 84x tile layer will replace the current registration. Do you want to queue registration job?		
Select a channel:	ETD; SecondaryElectrons	~
Select a stitching profile:	Default	~
	Queue job	Cancel

5. A registration job processes the tile set.

Image Context Menu

Images are directly stored in the selected layer. Right-click on an image to access its context menu.



Maps User Interface



- For descriptions of common selections to all systems, see Annotation Context Menu.
- For descriptions of image-specific selections, see the table below.

Image Context Menu Overview		
Menu Selection	Description	
Export to Project	Copies the data to a new or existing project.	
Save Image to File	Saves the image to disk as a TIFF image.	

Nav-Cam Context Menu

The Nav-Cam image is stored in the selected layer. Right-click on a Nav-Cam entry to access a context menu.



[Bring To Front		
E	Bring Forward		
S	Send Back		
5	Send To Back		
(Center View	Ctrl+Shift+C	
C	Center and Rotate View		
E	Export To Project		
ļ	Alignment	•	
F	Persist Nav-Cam Alignment		—— Unique selection for Nav-Cam
[Delete		

This only applies to SEM/SDB systems (see SEM Nav-Cam Context Menu).

Multiple Layer Context Menu

This menu is only available when multiple layers are selected.

i Note: This menu is only available on SEM/SDB.

🔻 🗹 Layer			
🔽 🗔 Helios PFIB Nav	Cam		
🔽 🗔 Zn_K_map		Bring To Front	
🔽 🗔 Si_K_map		Bring Forward	
🔽 🗔 S_K_map		Send Back	
🔽 🗖 O_K_map		Send To Back	
🔽 🛄 Na_K_map			
🔽 🗖 Mg_K_map		Center View	Ctrl+Shift+C
V 🛄 K_K_map		Center and Rotate View	
V 🗖 Fe_K_map		Export To Project	
🔽 🗔 CI_K_map		Alignment	•
	•••	Marra Chanada	
		Greate MID Laure	
Name	Mg_K_map		
Plane	1	Delete	

Menu Selection	Description
Merge Channels	Merges the selected image layers together by combining all their channels together and creates one resulting image layer. This requires all selected layers to be image layers and have the same resolution, the same number of z planes, and the same time frames.

This only applies to SEM/SDB systems (see SEM Nav-Cam Context Menu).

Job Control

This topic describes the Job Control.

Summary toolbar

The summary toolbar is always visible in the user interface by default. Once a job is queued, the Run button becomes available in the summary toolbar.





- **Progress Bar**: This indicates the progress of an individual job (tile set acquisition, post processing, etc.).
- Remaining Time: This indicates the time required to complete all selected jobs.

Job Control panel

Use the Job Control panel to manage all of your queued jobs, including microscope acquisition jobs and data processing jobs. This panel contains an Acquisition tab for your microscope acquisition jobs and a Processing tab for your data processing jobs. Click the up arrow next to **Job Control** on the summary toolbar to open the Job Control panel.

• Note: To close the Job Control panel, click the user interface anywhere outside the control panel.

Microscope acquisition jobs

The Acquisition tab displays the current status and progress of the queued jobs related to microscope acquisition. Refer to the Acquisition Tab Overview table below for descriptions of each control within the Job Control panel.

Run queued microscope acquisition jobs

Click **Run** in the summary toolbar to begin job execution for the microscope acquisition jobs.

AQUISITION	PROCESSING	
		DESELECT ALL SELECT ALL
		Show completed jobs
Overview Tile	Set	00:09:58 🗹
Feature Detail	Tile Set	00:01:36 🗹
ICE detail		00:00:24 🗹
ESTIMATION DISK S	PACE	ESTIMATION TIME
1.07 GB / 2.43 TB		
POST ACQUISITI	ON	POST PROCESSING
Turn Off Bear	n When Done	Stitch Tile Sets When Done
Retract Detect	tors When Done	
Sleep When I	one	
Pump to HiVa	c When Done	
	Þ RUN	

Stop running microscope acquisition jobs

Click **Stop** in the summary toolbar at any time during job execution to stop running the microscope acquisition jobs.

Maps User Interface

AQUISITION	PROCESSING		
			DESELECT ALL SELECT ALL
			Show completed jobs
Overview Tile S	Set	49% Acquiring imag	es 00:01:16 🜌
Feature Detail	Tile Set		00:01:36 💆
ICE detail			00:00:24 🛃
ESTIMATION DISK S 649.06 MB / 2.45 TB	PACE		
POST ACQUISITI	ON	POST PROCES	SSING
Turn Off Beam	n When Done	💿 Stitch Tile	Sets When Done
Retract Detec	tors When Done		
Sleep When D	Done		
Pump to HiVa	c When Done		
	L STOP		0:03:16

Control	Description
Deselect All	Deselects all jobs for processing.
Select All	Selects all jobs for processing.
Show completed jobs	Displays completed jobs.

Acquisition Tab Overview(Continued)		
Control	Description	
Job list items	 Each job row lists: Job name Job progression Time estimation if job has not started. If estimation is not provided by job, ':' is displayed instead. Time remaining for job in progress. The value calculation is based on previous progress and appears after the job is started and the estimation is ready. Jobs are listed in the order they will be processed. Select the check box to the right of a job to add it to the processing queue. For non-running jobs, you can click and drag the job rows to reorder them. You can also use the Context Menu to reorder them. 	
Estimation disk space	Displays estimation disk space needed to complete all selected jobs and the available disk space. The field turns red if there is not enough space for the jobs.	
Estimation time	Displays total estimation time needed to complete all selected jobs. The field displays ':' if there is no estimation provided by selected jobs.	
Post acquisition	Post-acquisition actions are listed in this section. If selected, actions are performed after acquisition. For system-specific action, see: TEM Post-Acquisition Actions SEM/SDB Post-Acquisition Actions	
Post processing	Post-processing actions are listed in this section. If selected, actions are performed after processing.	
Stitch Tile Sets When Done	Automatically adds a stitching job to the end of the Processing queue for each completed acquisition job.	

Data processing jobs

The Processing tab displays the current status and progress of the queued jobs related to data processing such as stitching, registration, or maximum intensity projection creation. Since these jobs do not require the use of the microscope, they can run in parallel to the microscope acquisition jobs. The Processing tab has the same functionality as the Acquisition tab described in the Acquisition Tab Overview table, however, the Processing tab contains its own run and stop buttons.

Run data processing jobs

Click **RUN PROCESSING** in the summary toolbar to begin job execution for the data processing jobs.

ACQUISITION PROCESSING	
	DESELECT ALL SELECT ALL
	Show completed jobs
Stitch: Fracture Detail	00:00:43 🗙
Stitch: Sample Overview	00:03:54 🗙
Stitch: Interesting Feature	00:03:54 🗙
ESTIMATION DISK SPACE 0 bytes / 2.4 TB	ESTIMATION TIME 00:08:33
A JOB CONTROL RUN	

Stop running data processing jobs

Click **STOP PROCESSING** in the summary toolbar at any time during job execution to stop running the data processing jobs.

ACQUISITION PROCESSING	
	DESELECT ALL SELECT ALL
STOP PROCESSING	Show completed jobs
Stitch: Fracture Detail 23% Stitching tile [3,6]	00:00:36
Stitch: Sample Overview	00:03:54 🗙
Stitch: Interesting Feature	00:03:54 🗙
ESTIMATION DISK SPACE 0 bytes / 2.4 TB	

Processing Tab Overview		
Control	Description	
RUN PROCESSING	Starts execution of the data processing job. This button is available if any data processing jobs are queued, but the queue is not running.	
STOP PROCESSING	Stops the execution of the data processing jobs. This button is available only if the queue is running.	

Processing Tab Overview(Continued)		
Control	Description	
Spinner icon	The spinner icon is located next to the Run button in the summary toolbar. This icon appears and animates while the data processing queue is running to indicate that background work is active. You can click the spinner icon to open the Processing tab whenever this tab is closed and the spinner icon is present within the summary toolbar.	

Context Menu

For non-running jobs, you can right-click on any job row to open the context menu and change the processing order.

Bring To Front
Bring Forward
Send Back
Send To Back

- Bring To Front: Moves the selected job to the front of the queue. The job is processed as the first one.
- Bring Forward: Moves the selected job forward in the queue.
- Send Back: Moves the selected job backward in the queue.
- Send To Back: Moves the selected job to the back of the queue. The job is processed as the last one.

Job processing issues

When a job processes, you might encounter the following error and warning messages to indicate that there are processing issues.



- Error: This message indicates that the job execution is blocked, but it is not preventing other jobs from running.
- **Warning**: This message indicates that job execution needs attention, but it is not blocked.

You can also view the error and warning messages in the summary toolbar next to the job status.

^	JOB CONTROL		RUN	STEM mode not available	
^	JOB CONTROL	•	RUN	Column valves are closed.	

Messages are displayed according to the following strategy:

- Errors have priority over warnings.
- If more than one error (or warning) occurs during the same job, then the most recent job processing issue is displayed first.
- A job processing issue related to an unscheduled job is not displayed.

Viewer

This topic describes the Viewer in the Maps user interface.

Topics in this section include:

Tile Set Grid Color Key

Alt +Left-Click and Drag Selected Area

Save image to file

Viewer tile context menu

This example shows a 4 x 4 tile set grid.

- Double yellow lines indicate the overlap between tiles.
- The purple FOV (field of view) rectangle displays the HFW (horizontal field width).

i Note: The coloring and labels of the FOV rectangle may change depending on the context.

- The *scale bar* (not shown) in the bottom right corner shows the distance (for example, in microns) in relation to the current magnification of the displayed image.
- The **Delete** key works in Maps for any selected item. You will receive a Confirmation request.



Tile Set Grid Color Key

- Yellow = Unacquired tiles. Visually indicates the tile overlap.
- Grey = Tiles that have been disabled; they will not be acquired
- Lime Green = Acquired tiles that are awaiting registration
- No outline = Acquired tiles that have been successfully registered
- Red = Acquired; but these tiles failed registration and will require manual alignment (see Manual Alignment)
- **Bright Red** = Edges of groups of red tiles are highlighted with a brighter red, to show where the actual break in the connection is. Often, fixing a bright red connection (via Manual Alignment) will allow previously red tiles to become successful.

Alt +Left-Click and Drag Selected Area

Alt+left-click and drag a rectangle in the viewer to create a selected area. A menu appears with selections for that area.

Zoom to selection
Add tiles
Add area of interest
Save image to file
Stitch selection

Menu Selection	Description		
Zoom to selection	Zooms the field of view to the selected area.		
Add tiles	Places a new tile set in the set	lected area, using the Active Template.	
Add area of interest	Places a new Annotation laye	r in the Layer control.	
	Name	Area of interest	
	Notes		
	Color	Ċ\$	
	Fill Opacity		
		Ellipse	
	Stage X	-9.3604 mm	
	Stage Y	3.8232 mm	
	An area of interested is used t named, notes captured, color	o highlight regions of the sample. It can be selected, and it can also be an ellipse.	
Save image to file	Saves an image to file (see Save image to filebelow).		

Viewer Selected Area Overview

	Viewer Selected Area Overview(Continued)
Menu Selection	Description
Stitch selection	Stitches a subset of tiles, as defined by the rectangle area of interest (see Stitching below). This menu selection is deactivated and not available if a tile set is not selected.

Save image to file

When you use Alt+left-click and drag to create a new tile set, the menu selection **Save image to file** appears.



1. Click Save image to file. The Save Selection Options screen appears.



2. Either enter a desired resolution into the **Image Size** or **Pixel Size** field, or click the **Use full resolution** button to set the exported image width to the native pixel size.

Note: Full resolution exports for large selections of high-resolution data are not available for all image types.

- 3. Select the Add scale bar check box to place the viewer scale bar on the exported image.
- 4. Click the browse button (...) to display the browser window.
- 5. File type options include TIFF Image and Tiled TIFF Image. Select a file type, file name, and location and then click **OK**.
- There are limitations for the TIFF file type. If TIFF is desired, but the resolution of the selected area is too large for TIFF export, a warning appears. Click Shrink for Tiff and the exported image width in pixels is automatically reduced to the maximum resolution for TIFF export.



7. To save large resolution images, use the Tiled TIFF format.

i Note: Not all image processing software supports the Tiled TIFF format. We recommend that you use the TIFF format to ensure compatibility with external image processing packages.

8. The Save Selection Options window is still open. Click OK to close it.

Viewer tile context menu

Right-click on a tile within the tile grid to access a context menu. Alternatively, you can select the small menu button on the side of the selected tile to have the tile-specific menu appear.



The menu items outlined in purple below are system-specific; whereas, all other menu items are common to SEM/SDB and TEM systems. See the Common Viewer Tile Context Menu Overview table for descriptions of the common menu items, which includes common items in the main menu and associated submenus. For system-specific menu selections outlined in purple, click the following links.

- SEM/SDB Viewer Tile Context Menu
- TEM Viewer Tile Context Menu

Viewer Tile Context Menu for SEM/SDB

Drive Stage Here			
Rotational Alignment	·		
Snapshot Here			
Snapshot Here w/HFW	·		
Add Tiles Here			
Add Site of interest			
Tile Set (4) ►			
Tile 10,10 →		~	Acquire
Clear Digital Rotation			Open tile image folder
, , , , , , , , , , , , , , , , , , ,			Open source image
		Г	Preview Tile(s)
		L	Sub-tile >
			Open analytics data folder

Viewer Tile Context Menu for TEM



Common Viewer Tile Context Menu Overview		
Menu Selection	Description	
Drive Stage Here	Moves and centers the stage to the selected tile (see the green cross).	
Snapshot Here	Acquires a snapshot image centered on this stage location at the current horizontal field width.	
Add Tiles Here	Adds a new Tile Grid centered on the clicked location.	
Add Site of Interest	 Displays a New Annotation dialog box for creating the Site of Interest (SOI). An SOI is a specific stage location used to mark sample features. Z: Stage Z and T: Alpha Tilt can optionally be stored to be recalled later. The new SOI is created is the currently selected layer group. Note: This is different from the New Annotation dialog box for creating an Area of Interest, which marks an area of the stage. 	
Clear Digital Rotation	Resets the digital rotation to 0 degrees if rotation had previously been applied. This only changes the displayed rotation of the images and does not rotate the microscope stage.	
Tile Set submenu		
Tile Set	Displays the name of the selected tile set. Examples include: Tile Set, Tile Set (2), and Tile Set (3).	
Open Z- stack	Opens the Z-Stack Browser to view the focus stack of the tiles. See Z-Stack Browser Controls.	
Stitch	Initiates stitching for this tile set. See Stitching.	

Common Viewer Tile Context Menu Overview(Continued)				
Menu Selection	Description			
Snap to stage position	Moves the clicked position on the tile layer to the current stage position.			
Preview Corners	This item is only available in the Tile Set submenu before acquisition. After tuning Preview Corners, preview images appear in the corners of the tile set.			

Common Viewer Tile Context Menu Overview(Continued)			
Menu Selection	Description		
Preview Edges	This item is only available in the Tile Set submenu <i>before</i> acquisition. After running Preview Edges, preview images appear in center of each edge of the tile set.		
Tile X, Y	Displays the row (X) and column (Y) number for the selected tile.		
Tile submenu selections			
Tile	Displays the name of the selected tile. Examples include: Tile 1,1 and Tile 1,3.		
Open tile image folder	If activated, opens File Explorer to the directory that contains the tile image data.		
Open source image	If activated, opens the tile's source image file with the default image viewer (Windows Picture Viewer).		
Open analytics data folder	If activated, opens the tile's analytics data folder that contains the metadata files from Velox.		

Rotation Control in the Viewer

Grab the lollipop-looking control with the mouse to easily rotate tile layers (not for TEM) and annotation layers.



Micron Bar

The micron (scale) bar scales to the magnification.



Basic Operations

This section describes the basic process flow for Maps.

Topics include:

- Launch the application
- **Application Window**
- Create a Project
- Setting Options
- Define a Tile Set
- Tile Set Templates
- Acquire a Tile Set
- Import Images from Other Sources
- Stage Navigation
- Stitching

Launch the application

Click the desktop icon 🥮, or use the Windows Start menu to launch the Maps application.

When the End User License Agreement dialog opens, do the following:

- 1. Read the entire end user license agreement by using the vertical scroll bar as you read.
- 2. When finished reading the end user license agreement, click **I have read the EULA Terms** and Conditions.
- 3. If you are authorized to accept the license agreement terms and conditions, then click I confirm that I am authorized on behalf of the licensee of the software to and hereby accept the EULA Terms and Conditions.
- 4. If you do not want the Maps application to anonymously track usage and send the data back to the Maps team, then deselect the **Allow usage tracking** checkbox.
- If you agree with the end user license agreement terms and conditions, then click I Accept to proceed with the Maps application. If, however, you do not agree with the end user license agreement terms and conditions, then click I Decline to close the End User License Agreement dialog.



Application Window

The application window appears with the **Project History** window open.

Select a display mode from the icons below the **New** button:

• Thumb View: Displays thumbnail images of project data.

B Project History						
NEW IMPORT OPEN			REFRESH			
111						
Ð	VOLCANICSAND 7/23/2020 2:55:57 PM	PROJECT DEMO 1 7/23/2020 2:54:12 PM	PROJECT DEMO 2 ● = 7/23/2020 2:52:07 PM			
MULTI PROJECTS = 7/22/2020 10:00:32 AM						
			Quit Maps			

• List View: Displays a list of project data.

No Project History					
N	W IMPORT				REFRESH
88 (8				
		Project		Description	
	VolcanicSand				
	Project Demo 1				
•	Project Demo 2				
	Multi projects				
					Quit Maps

New Project

If this is your first project, then the list/thumb view will be empty and the **Open Project** button will be unavailable. Click **New** and proceed to **Create a Project**.

Import a Project

To import a project that was created on another system and does not appear in the project list, click **Import Project**. The project opens and its name is added to the project history list. See **Import a Project**.

Open a Project

To open an existing project, highlight a project name in the list and click **Open Project** and then proceed to **Open a Project**.

Yellow Exclamation Mark in Project List

This symbol appears if the project path is not found for that data, due to a disconnected network drive, USB drive, or other drive.

Refresh

Updates the status of any disconnected projects that have been reconnected properly.

Quit Maps/Cancel

- If you have not opened a project yet, then you can directly quit Maps by clicking Quit Maps.
- If a project is already opened, then this button is replaced by a **Cancel** button that allows you to close the **Project History** window.

Create a Project

Create a new project, import a project, or open an existing project using File menu selections or Project History buttons.

New Project

1. To create a new project, click File > New Project.



Alternatively, click **File > Open Project** and then click **New** within the **Project History** window.



2. When the **New Project** screen appears, enter the new into the **Project Name** field, enter a brief description into the **Project Description** field, select a holder from the Sample Holder dropdown menu, and then click **Create**.

Ma New Project			×
Project Name			
Project Description			
Project Path			
C:\Maps Projects			
Sample Holder:			
3mm Grid holder			~
កម្មមូមក្រ	Holder:	3mm Grid holder	
ā ā ā ā [Version:	1.1	
	Manufacturer:	FEI	
	Part number:	1046764	
		Create Canc	el

- 3. The default data directory is used unless you browse to a new location.
- 4. The new project name appears in the title of the Layer control panel and the Layer control shows a default Layer.
5. The purple rectangle represents the current field of view on the sample and the purple + is the current stage position.



6. The white outline shows the stage boundaries, which can be circular or rectangular shaped depending on the instrument.



Open a Project

To select an existing project, select File > Open Project .

File	View Options Help					
	New Project					
	Open Project					
	Import Project					
	Project Properties					
	Open Project Folder Ctrl+O					
	Import Grid					
	Import Images					
	Save Screenshot As					
	Save Screenshot F2					
	Quit					

The **Project History** window appears, showing all existing projects. See the list of options below to locate, open, and search for a project.

Basic Operations



 Note: The padlock located near the project title indicates that the current project cannot be opened because it is being used by the current instance or another Maps instance. The exclamation icon and a gray project title in italics indicates that the current project is not available (on a removable device).

- To open an existing project, double-click the thumbnail of the desired project and then click **OPEN**.
- To open the selected project in a new Maps Offline Viewer instance, double-click the thumbnail of the desired project and then click **OPEN IN OFFLINE VIEWER**.
- To search for an existing project, enter a project name or project description into the search bar and then click the **Search** icon.

• Click the down arrow menu of the selected project to view two submenus: **Open Project** Folder and Remove Project...



- Click **REFRESH** to refresh the status of the project list (for example, Project not found).
- Click the List View icon to switch to a list view of projects.

Ma Proj	ect History		×
N	EW IMPORT OPEN IN		REFRESH
G			
888	器 Search Q		
	Project	Description	
e	VolcanicSand		
	Earthworm	Earthworm testicles Array Tomo sample from Tilman Franke	
	CL mineral sample Maps 3.5	Mineral sample imaged with the Cathode Luminescence Detector (CLD) from James Ranney (Hillsboro NanoPort)	
	Welcome to Quattro Slim 3.5	A collection of various SEM samples from James Ranney (Hillsboro NanoPort)	
	Brain Tissue VS	VS standard mouse brain sample	
	Mouse Kidney	Mouse Kidney imaged in CorrSight from Todd Hanson	
	Gibeon Meteorite 062016	SEM imaging of a sample from the Gibeon Meteorite that fell during prehistoric times in Namibia. From Rick Passey.	
	BPAE - CorrSight	Bovine Pulmonary Artery Endothelial (BPAE) cells imaged in CorrSight from Todd Hanson	
	Welcome to Quattro Slim 3.5a	A collection of various SEM samples from James Ranney (Hillsboro NanoPort)	
	NP Blend	Nanoparticles including NiSi. LiNiMnO, CdS-ZnS, FeAI, all with some degree of segregation to make things more interesting. We usually use these to de	r
	Mouse Myoblasts - Correlative Workflow	Mouse Myoblast cells imaged in CorrSight and TEM, from Oliver Raschdorf	
	Spider	SEM z-stack of a common spider. From Edhouse.	
	Battery electrode	Correlative analysis on battery electrode	
	Zebrafish Embryo	Zebrafish Embryo Array Tomo dataset from Tilman Franke.	
	Tungsten Carbide Steel RIP 10-2016	Tungsten Carbine Steel sample imaged in SEM from Rick Passey (MSBU)	
	RamJetabradable	Ram Jet tip SEM data correlated with uCT from Bart Winiarski and Eric Goergen	
	Notre Dame	Red Trinitite from the original Trinity nuclear bomb test. Includes optical and SEM image data. From Rick Passey.	
	LLNP - Melt	Reflected optical and SEM images of a witness slip from Lawrence Livermore National Labs ignition testing via Rick Passey.	
	IN718_AM_Polishing_BW2	Propeller sample SEM data correlated with EBSD, uCT, reflective optical from Bart Winiarski and Eric Goergen	
	IC Repeating Features	Integrated circuit showing repetetive features, acquires and stitched by Maps.	
			Cancel

• Right-click a project name in the list view to access the submenu for opening a project folder and removing a project.



Import a Project

Importing a project is useful for passing projects between computers. Once imported, you can select it as an existing project.

1. To import (save) an existing project, select **File > Import Project**.



Alternatively, click File > Import Project and then click Import within the Project History window.



2. Browse to a Source directory. Name the project to be exported and then click Import.



3. The imported project opens and the file name is added to the project history list.

Ma Pro	ject History		×
~	IEW IMPORT OPEN OPE	N IN OFFLINE VIEWER REFRESI	н
888	器 Search Q		
	Project	Description	
	Project Demo 2000		۲
	VolcanicSand		
	Earthworm	Earthworm testicles Array Tomo sample from Tilman Franke	
	CL mineral sample Maps 3.5	Mineral sample imaged with the Cathode Luminescence Detector (CLD) from James Ranney (Hillsboro NanoPort)	
	Welcome to Quattro Slim 3.5	A collection of various SEM samples from James Ranney (Hillsboro NanoPort)	
	Brain Tissue VS	VS standard mouse brain sample	
	Mouse Kidney	Mouse Kidney imaged in CorrSight from Todd Hanson	
	Gibeon Meteorite 062016	SEM imaging of a sample from the Gibeon Meteorite that fell during prehistoric times in Namibia. From Rick Passey.	
	BPAE - CorrSight	Bovine Pulmonary Artery Endothelial (BPAE) cells imaged in CorrSight from Todd Hanson	
	Welcome to Quattro Slim 3.5a	A collection of various SEM samples from James Ranney (Hillsboro NanoPort)	
	NP Blend	Nanoparticles including NiSi. LiNiMnO, CdS-ZnS, FeAl, all with some degree of segregation to make things more interesting. We usually use these to der	
	Mouse Myoblasts - Correlative Workflow	Mouse Myoblast cells imaged in CorrSight and TEM, from Oliver Raschdorf	
	Spider	SEM 2-stack of a common spider. From Edhouse.	
	Battery electrode	Correlative analysis on battery electrode	
	Zebrafish Embryo	Zebrafish Embryo Array Tomo dataset from Tilman Franke.	
	Tungsten Carbide Steel RIP 10-2016	Tungsten Carbine Steel sample imaged in SEM from Rick Passey (MSBU)	
	RamJetabradable	Ram Jet tip SEM data correlated with uCT from Bart Winiarski and Eric Goergen	
	Notre Dame	Red Trinitite from the original Trinity nuclear bomb test. Includes optical and SEM image data. From Rick Passey.	
	LLNP - Melt	Reflected optical and SEM images of a witness slip from Lawrence Livermore National Labs ignition testing via Rick Passey.	
	IN718_AM_Polishing_BW2	Propeller sample SEM data correlated with EBSD, uCT, reflective optical from Bart Winiarski and Eric Goergen	
	IC Repeating Features	Integrated circuit showing repetetive features, acquires and stitched by Maps.	

Unsupported Project version of Converting From Maps 1.1 or 2.0 Software

Only projects of Maps version 2.5 or higher can be imported. The import process is only oneway. Older software projects must be converted. When a project is imported from Maps v1.1 and 2.0, a warning is displayed telling the user that the import is one-way and the project will no longer be compatible with v1.1 and 2.0.

If you atttempt to import older projects, following dialog will be displayed:



You are prompted to browse to a default location for storing project data.



We recommend that you do not use the C: drive, but you can use a supplemental data storage device such as a network drive, external USB drive, or additional internal hard drive.

Setting Options

Select new project defaults from the Options menu.

Application SettingsTileSet TemplatesDefault Autofunctions Templates (TEM Only)Microscope SettingsSnapshot Settings (SEM/SDB Only)SnapShot Settings (TEM Only)Animate Pan & ZoomNav-Cam Alignment (SEM/SDB Only)Running the Automated Holder Alignment (SEM/SDB Only)

Application Settings

Click **Options > Application Settings** to display a window for setting defaults. See **Options menu** for descriptions of the controls.

TileSet Templates

Click **Options > TileSet Templates** to display a dialog that shows a collection of templates and their respective settings. These settings are system specific:

- SEM/SDB Tile Set Tab
- TEM Tile Set Tab

Default Autofunctions Templates (TEM Only)

Click **Options > Default Autofunctions Templates** to display a dialog that shows a collection of templates and their respective settings. These settings are system and mode specific:

• TEM Default Autofunction Templates

Microscope Settings

Click **Options > Microscope Settings** to display the **Microscope Settings** window.

These are settings system-specific:

- SEM/SDB Options Menu
- TEM Options Menu

Snapshot Settings (SEM/SDB Only)

Click Options > Snapshot Settings to display the Microscope Settings window. See SEM/SDB Options Menu.

```
SnapShot Settings (TEM Only)
```

Click Options > Snapshot Settings to display the Snapshot Settings Window. See TEM Options Menu.

Animate Pan & Zoom

Click **Options > Animate Pan & Zoom** to allow a smoother transition of panning and zooming.

- Pan: To make a sweeping movement
- Zoom: To simulate movement rapidly away from or toward a subject using a zoom lens

i Note: This is a slower process and takes more memory.

Nav-Cam Alignment (SEM/SDB Only)

See SEM Nav-Cam Alignment

Running the Automated Holder Alignment (SEM/SDB Only)

See Automated Holder Alignment.

Define a Tile Set

This section describes how to define tile sets. Refer to Tile Set Templates to learn how to change default template settings.

Topics include:

Methods Define a Circular Tile Set Define a Freehand Tile Set Stamp a Tile Set Manually Adjust Tile Sets Tile Set Display with Grid Configure Tiling Parameters Methods

Define the tile set for your project using one of the three methods below:

• Click the Add Tiles button on the Layer control.



• Press Alt+click and drag a rectangle in the Viewer, and then click Add Tiles in the menu.

-
Zoom to selection
Add tiles
Add area of interest
Save image to file

• Click the Add Tile Set button in the tool bar, then click and drag an area in the viewer.



Maps creates a layer in the Layer control, depending on the type of microscope connected. You cannot create a tile set when not connected to a microscope.

The new name is added to the list of tile sets beneath Layer in the Layer control. The Tile Set tab displays default parameters, and a grid (green square) is centered on the stage location.

Define a Circular Tile Set

Click the Add Circular Tile Set icon on the tool bar, then click and drag an area in the viewer to create the tile set.



Example



Define a Freehand Tile Set

Click the Add Freehand Tile Set icon on the tool bar, then click and drag an area in the viewer to create the tile set.



Basic Operations

Example



Stamp a Tile Set

Go to the toolbar and click **Stamp Tile Set** to enter the stamp tile set mode. While in this mode, Maps creates a new tile set after each time you single-click in the Viewer. Maps creates tile sets using the active template, and it draws a tile set outline on the cursor location if it can determine the tile set size. To quit the stamp mode, either press **Esc**, right-click in the Viewer or select another tool from the toolbar.



Manually Adjust Tile Sets

Once a tile set is created, you can adjust the size of the tile set with the drag handles in the viewer.

When a tile set is selected, its outline turns green and drag handles appear as small circles in the corners and on the sides of the tile set border.



Click and drag within the tile set region to reposition it. This can be performed on multiple tile sets at once by pressing **Ctrl+left-clicking** each tile set to multiselect them.

Click and drag the handles to perform resizing actions on the tile set. While dragging, you will see the adjustment result in a light blue border as shown below.

Basic Operations

The screenshots below show resizing in process. The original border is green. The new border is light blue. The purple arrows show the direction that the border is dragged.



To rotate the tile set, click and drag on the small rotation control located on the side of the tile set. This can be performed on multiple tile sets at once by pressing **Ctrl+left-clicking** each tile set to multiselect them. This control is only visible for tile sets on systems that support rotation. See Rotation Control in the Viewer.

Tile Set Display with Grid

Two views are shown below. They look different in STEM Mode.

SEM/SDB View



TEM View



Configure Tiling Parameters

- Change the default *tiling* parameters to customize the grid for the data you will be acquiring.
- Modify the *imaging* parameters and resize the grid for the level of detail and size you need.

Tile Set Templates

Tile set templates are a set of pre-defined microscope settings designed to make your workflow more efficient. You can use the factory template, or create custom templates to match your use cases.

Template terms

Here is some terminology to help you understand tile set templates.

Template: A set of tile set settings that are applied to a newly-created tile set to help create defined tile sets more efficiently.

Custom template: A template created by a user.

Active template: The template that will be used when creating a tile set. To change the active template, click the **Template** list and then click the template you want to use. The example below shows a TEM with the factory template and one custom template.



You can see your more details about your templates under Settings > Tile Set Templates.

Default template: A template that will be set as the active template when the Maps software is launched. The default template is initially set to the factory template until changed. Any template can be set as the default template in the ! window.

Factory template: A pre-defined, immutable template that comes with the Maps Software. There is one factory template per microscope mode.

Templates on SEM

On a SEM tool, a tile set template will have the properties shown below saved and applied when used. This screen capture shows three templates: two factory templates and one custom template. The factory template is set as the default template.

SEM Tile Set Templates screen

Ma Tile Set Templates						×
Templates	Mode	Electron				~
	Factor	y Template	(Default) 🏦	Properties		
Holder Defaults	Factor	y Mineralogy Template		Name	Factory Templa	ate
	Custor	n		Tile Set Name	Tile Set	
				These settings a Grids. Items wi unchecked and microscope value	re used as defau th a check box they will instea at the time of cr	Its for new Tile may be left take on the eation.
				Beam Type	Electron	~
				Tile HFW		
				Columns	10	
				Rows	10	
				Overlap X	10%	
				Overlap Y	10%	
				Resolution		~
					16-bit image	
				Dwell		
				V Frames	1	
	Imp	ort Export all				Set As Default
						ОК

Templates on Phenom

On a Phenom tool, a tile set template will have the properties shown below saved and applied when used. This screen capture shows two templates: one factory template and one custom template. The factory template is set as the default template.

Phenom Tile Set Templates screen

🔤 Tile Set Templates									×
Templates	Mode	SEM							~
Holder Defaults	Factor	y Template	(Default)	Properties Name		Custom			
	Custor			Tile Set Name		Tile Set			
				Detector		BSD Full		~	
				HDR		OFF		~	
				Averaging		Live		~	
				Contrast	2	1.00			
				Brightness		0.00			
				High Voltage		5.3 kV		~	
				Intensity		Image		~	
				Vacuum		High		~	
				- Analytics					
				EDS Acquisition	n		2		
				Acquisition Pa	aram	eters			
				EDS Dwell			1 ms		
				Elements			C, Cu, Al, Ni, O, F, N		
				EDS Resolution	۱		Eighth (64x64)	~	
	Imp	oort Export all						Set As Defau	ult
								ОК	

Templates on TEM

On a TEM, only the settings shown in the **Tile Set Templates** screen are saved in **Factory Template (TEM)**.

TEM Tile Set Templates screen

Ma Tile Set Templates						2	×
Templates	Mode	ТЕМ					~
	Factor	y Template (TEM)	(Default) 前	Properties			
Holder Defaults	Base te	emplate		Name	Factor	ry Template (TEM)	
				Tile Set Name	Tile Se	et	
				These settings are TEM Tile Sets.	used as	s initial values for new	
				Tiles X, Y		3, 3	
				Overlap X, Y		20%, 20%	
				Resolution		Maximum	
				Exposure Time		1s	
				Focus Method		Fixed	
				Stage Settling Time	e	0 s	
			Image Correction				
	Imp	ort Export all				Set As Defau	lt
						ОК	

On a TEM, custom templates are also read-only after you set them up. You are prompted to define the settings you want when creating the template.

Ma Tile Set Templates							\times
Templates	Mode TE	М					~
	Factory Ter	mplate (TEM)	(Default)	Properties			
Holder Defaults	Base temp	late	Ē	Name	Base te	mplate	
				Tile Set Name	Tile Set	t	
				These settings are TEM Tile Sets.	used as	initial values for new	'
				Camera		BM-Ceta	
				Tiles X, Y		4, 5	
				Overlap X, Y		20%, 20%	
				Tile HFW		6.9264 µm	
				Total Area		23.5 µm x 29.1 µm	
				Resolution		512 x 512	
				Magnification		SA 5500 x	
				Pixel Size		13.8528 nm	
				Exposure Time		1 s	
				Intensity		0	
				Spot Size			
				Probe Mode		MicroProbe	
				Focus Method		Fixed	
	Import.	. Export all				Set As Def	ault
						OK	

Tasks associated with tile set templates include:

Default templates

Depending on what type of system you are using, you can either update settings in the Default Template or choose a custom template to be the new default.

Save new Default Template settings

This procedure works on all tool types. It overrides the Default Template settings.

- 1. Create a tile set with the settings you want in the new default tile set template. See Define a Tile Set.
- 2. Right-click on the tile set and then click **Save and Set As Default Template**.

Change Default Template

You can only directly edit the Default Template in SEM) but not in TEM. However, you can change the settings that Maps software uses to define default tile set templates.

Follow the steps from below.

Select a new default template

- 1. Create a tile set with the settings you want in the new default tile set template. See Define a Tile Set.
- 2. Open the **Tile Set Templates** window. On the **Options** menu, click **Tile Set Templates**.
- 3. Click the template you want to be the new default and then click Set As Default.

The (Default) label is displayed next to the new default template.

4. Click OK.

Create custom templates

Custom templates are collections of settings that you define and can reuse when you create tile sets.

- 1. Create a tile set with the settings you want in the custom tile set template. See Define a Tile Set.
- 2. Right-click the tile set. In the context menu, hover the mouse over **Save as Template** and then click **New Template**.

🔻 🔽 Layer							
🔽 🌐 Custom 2		**					
🔽 🌐 Custom 1				Bring To Front			
⊞ ₽ 5	Ø			Bring Forward Send Back Send To Back			
Name	Custom 2						
Tile Set Type	TEM		~	Show Grid Lines Center View	Ctrl+G Ctrl+Shift+C		
Camera	BM-Ceta			Center and Rotate View			
Tiles X, Y	3	3		Save and Set As Default Template			
Overlap X, Y	γ 20% 20%			Save As Template	•	New Template	
	800 рх	800 px		Copy Grid			
Tile HFW 6.9264 µm			Export Grid				
Total Area 18 μm x 18 μm			Export To Project				

- 3. On the New Tile Set Template window, type the Name of the new template.
- 4. Type the **Tile Set Name**. This becomes the default name of each new tile set created based on this template.
- 5. (SEM only) Update fields in the **Properties** section to the values you want in the custom template.
- 6. (SEM only) For the **Tile HFW**, **Resolution**, **Dwell**, and **Frames** fields, clear the check box if you always want Maps software to use the value from the microscope settings for that field.
- 7. Click OK. Maps software saves a new template based on the tile set settings.
- 8. To use the new template, click the **Template** list and then click the template you want to use.



Edit custom templates

After you Create custom templates, you can change or edit the settings to match your needs.

Edit a custom template (TEM)

You cannot directly edit custom templates on TEM microscopes, but you can override them.

- 1. Create a tile set with the settings you want in the custom tile set template. See Define a Tile Set.
- 2. Right-click the tile set. In the context menu, hover your mouse over **Save as Template** and then click the name of the template you want to update.

~			III TE	EMPLATE	Factory Template	(TEM) 🗸			
•	🔽 Laye	r							
	 ✓ 	⊞ Custo ⊞ Custo	r mo	Bring To) Front	¥:			
⊞	÷	1		Bring Fo Send Ba	orward ock				
Name				Send To	Back rid Lines		Ctr	rl+G	
Tile Set Type				Center \	/iew		Ctı	rl+Shift+C	
Tiles X, Y				Save and Set As Default Template					
Overla	Overlan X V			Save As Template				New Template	
				Copy G	id				Best template

Maps software updates the tile set template settings. All new templates created with that template use the new settings.

Edit a custom template (SEM)

- 1. Open the Tile Set Templates window. On the Options menu, click Tile Set Templates.
- 2. Click the template you want to edit.
- 3. Update fields in the **Properties** section to the values you want in the template.
 - If you clear the check box next to the **Tile HFW**, **Resolution**, **Dwell**, or **Frames** fields, then Maps software always uses the value from the microscope settings for that field.
 - Click **Reset** to reset the template to the default settings.
- 4. When you are done editing settings, click OK.

Delete custom templates

You cannot delete the template marked as (Default).

- 1. Open the Tile Set Templates window. On the Options menu, click Tile Set Templates.
- 2. Click the template that you want to delete.
- 3. Click the trash can icon.

Maps software deletes the template.

Import and export tile set templates

Maps software allows you to export tile set templates developed on one tool and import them onto another.

Import tile set templates

This procedure replaces all current templates with the imported set, including the default. You cannot recover the tile set templates replaced by the import process.

- 1. Obtain an XML file containing tile set templates and save it on your local machine. See Export tile set templates for instructions.
- 2. Open the Tile Set Templates window. On the Options menu, click Tile Set Templates.
- 3. Click Import.
- 4. Navigate to the XML file that contains your templates and then click **Open**.

Maps software replaces your current custom templates and the settings in the **Default Template** with the templates in the imported file.

Export tile set templates

You can export tile set templates from a Maps software installation on one tool to another. Once exported, you can import the templates on another tool using Import tile set templates.

- 1. Open the Tile Set Templates window. On the Options menu, click Tile Set Templates.
- 2. Click Export all.
- 3. Navigate to the place you want to save the XML file that contains your templates and then click **Save**.

Acquire a Tile Set

When ready to begin acquisition, click Start Execution located in the status bar next to No Job.

Tile images are registered for stitching as they are acquired and displayed in the grid. The tile sets are colored as they are registered to give immediate feedback on the stitching confidence. See Tile Set Grid Color Key.



Preconditions

Microscope preconditions must be met to proceed with tile set acquisition. For example, if the microscope is not enabled or an acquisition parameter is invalid, then a validation error appears in relation to the tile layer and you cannot start acquisition.

JOB QUEUE PROC	Essing 😑		
			DESELECT ALL SELECT ALL
			Show completed jobs
SEM Tile Set		🙁 Microscope is offline	
ESTIMATION DISK SPACE 0 bytes / 496.36 GB			ESTIMATION TIME
POST ACQUISITION		POST PROCESSING	
Turn Off Beam When D	one	Stitch Tile Sets When	n Done
Retract Detectors When	Done		
Sleep When Done			
Pump to HiVac When D	one		

Re-Adjust Acquired Tile Sets

When a tileset is acquired, you can adjust the size of the tile set with the drag handles in the viewer. See Manually Adjust Tile Sets for instructions.

When you resize an acquired tile set in such a way that some acquired tiles will be deleted (smaller in either dimension), you are asked to confirm the action:



- Click **Delete** to delete acquired tiles.
- Click Cancel to cancel the resize operation.

Import Images from Other Sources

You can import an external image for navigation purposes to find the area of interest.

1. Right-click on a Layer in the Layer control and then click **Import Images...** on the context menu that appears.



2. In the standard **Open** window that appears, browse to a saved image and then click **Open**. Most standard image formats are supported, including BMP, JPG, GIF, PNG, and TIFF. Various microscopy, scientific, and biological image data formats are also supported including LIF, MRC, RAW, and LEI as well as DICOM data with DICOMDIR, DIC, DCM, and other DICOM standard formats.

In the importing dialog that appears, you can cancel an image import or send the import to the background.



Sending the import to the background allows you to use Maps while the import is running. When the import is in the background, other image imports are unavailable.

You can track the background import with the status bar import control. Click the control to reopen the importing dialog.



The imported image is displayed to the right of the main viewer while the alignment workflow is in progress. See Manual Alignment.



If you imported an image with known stage position (like Nav-Cam images, SEM/SDB images), then you are finished. Otherwise continue to the next step.

3. Align the two images with a simple one-, two-, or three-point alignment procedure: Move the alignment points on the imported image to match the feature on the tile set image and then click **Next**. See Coarse Alignment (1-, 2-, or 3-Point) for detailed information.

After clicking **Finish**, the viewer is restored to full size and the newly imported image is displayed aligned to the stage. The imported image name is added to the Layers panel.

Stage Navigation

To move the stage to a location, do the following:

1. Double-click anywhere within the stage boundaries as shown by the white circle or square.



2. In the Overview Viewer, right-click anywhere within the stage boundaries as shown by the white circle or square and click **Drive Stage Here**.



Stitching

After a tile set with at least one tile has been acquired, the Stitch action is made available.

Initiate Stitching

Manual Alignment

Stitching Job Options

Stitching Profile Best Practices

Initiate Stitching

Stitch single tile set

Select the acquired tile set in the Layer control, and then do one of the following:

• Click **Stitch** in the tile set context menu.

The manual UI appears so that you can confirm the tile alignments.



• Or, left-click and drag a rectangle around the tile sets to be stitched and then click **Stitch Selection** in the stitch context menu.



Stitch selection stitches all tiles inside the rectangle for the selected tile set. Tiles that are partially inside the rectangle are included as well. See Alt +Left-Click and Drag Selected Area.

Stitch multiple tile sets

Select the acquired tile sets in the Layer control, and then click **Stitch** in tile set context menu. Note that the manual UI alignment of tiles is skipped, and the acquired tile sets will be stitched using the default stitching options (see **Stitching Job Options**).

Manual Alignment

Follow the guided instructions to confirm or edit the alignment of the tiles.



Stitching Manual Alignment Histogram

Displays the histogram for the Stitching Channel.

Stitching Manual Alignment Histogram



Interface Item	Description
Channel name	For SEM/SDB only: Displays the channel name, which is the detector name.
Apply	Applies changes to the histogram.
Auto	When selected, continuously adjusts the min and max for the histogram display.
Semi-log	When selected, changes the histogram intensity display to logarithmic.
Current Max position	Displays the current max slider position.
Histogram data	Displays the histogram data for the selected channel.
Current Min position	Displays the current min slider position.
Overlap Transparency

Overlap Transparency options are available for assigning a color for the **Center Tile** and its **Neighbor Tiles**.



Transparency is useful for seeing through the layers to line up the tiles in manual tile alignment.

Two Views

Manual stitching makes use of two views.

- The Tile Alignment View is the larger view:
 - This view displays the currently selected tile and it's top, bottom, left and right neighbors. The positions of the neighboring tiles are based on their individual registrations to the center tile. This is not the same as the final position when stitched, which will be a result of the optimization of all tile-to-tile registrations of the tile set.
 - When manual alignment is first initiated, the first tile that failed alignment is selected and displayed here.
 - The view can be panned and zoomed in the same way as the other views. The center tile is fixed, but you can click left-mouse+drag to move its surrounding tiles.
- The Project View is the smaller version of the standard viewer in the bottom right.
 - This is a miniaturized version of the standard Maps project view.
 - It can be panned, zoomed, and each tile can be selected with the mouse. When a tile is selected, it appears as the center tile in the Tile Alignment View, and it can then be manually aligned to it's neighbors.

Manual Alignment with Transparency Off



Manual Alignment with Transparency On



Stitching Alignment Status Color Key

Each tile outline is colored to indicate stitching alignment status.

- Green: The automatic tile alignment succeeded.
- **Orange**: One side of the tile alignment failed, but another succeeded. These tiles do not require manual alignment.
- **Red**: The tile failed to align with either neighbor, which means that there is not enough valid registration for a tile or one of its neighboring tiles to be properly positioned in the larger picture. Often multiple tiles can be red, but only one tile needs to be manually connected to the green tiles to achieve alignment for all. After realignment, the outline turns **blue**.

Realign Failed Tiles

Use the Next and Previous buttons to jump through the list of failed tiles.

Confirm and Stitch

When manual alignment is finished, click **Confirm and Stitch** to begin stitching. You will be prompted to confirm and click **Yes**.

Stitching Job Options

After you click **Yes** to the Confirmation prompt, the **Stitching Job Options** screen appears.

Some options are not available for all tile sets.

Ma	Stitching Job Options	×		
Fi	nal output size: 5734 x 3359 x 8-bit grayscale			
	Crop Image			
	Blend Tiles			
	Tile-to-Tile Contrast Normalization (Experimental)			
	Always Use Calculated Positions			
	Stitch to Project			
	Stitch to File			
Des	Destination file name:			
C:\	C:\Work\Maps data\scan\SampleTissue (stitched).tif			
	Specify output image width 1024 pixels			
	Add scale bar			
	Set as Default OK Cancel			

Stitching Job Options Overview

Interface Item	Description
Crop Image	Crops the outer edges of the image to make it rectangular. Otherwise, it will potentially have jagged edges where the tiles have been shifted.
Blend Tiles	Blends the overlapped regions of the tiles together to make it appear as one continuous image. In some rare cases, you may not want to blend the tiles.
Always Use Calculated Positions	Maps uses calculated relative positions between tiles even if they are outlined in red (that is, neighbor tiles in which Maps was not able to find a relative position with a high enough score).

	Stitching Job Options Overview(Continued)
Interface Item	Description
Tile-to-Tile Contrast Normalization (Experimental)	Removes differences in brightness and contrast between tiles in the stitched tile set. The example shows the <i>standard</i> output in the left image and the <i>normalized</i> stitched output in the right image.
Stitch to Project	Stitches the tile set into the project for display in the Viewer. The stitched output appears as a new layer in the Layer control. Image: Im
Stitch to File	Stitches the tile set into an image and exports the file to the specified folder location. The output file is formatted as either RAW or TIFF in the Destination File name field and then stored in the folder. The default folder location is the current project's data folder and the default file format is TIFF. When stitching to the TIFF image format, the image size limit is adjusted based on the following image bit depth: • 8-bit grayscale: 2 gigapixel, or about 46,340 pixels square • 16-bit grayscale: 1 gigapixel, or about 32,767 pixels square
Destination File Name	Specifies the path and file name for the stitched project.

Stitching Job Options Overview(Continued)		
Interface Item	Description	
Specify output image width	Determines the resolution of the stitched image. Enter the desired stitched output width in the pixels field. When not selected, stitches to full resolution.	
Add scale bar	When selected, a scale bar is burned into the final image. This option is not available for RAW image format.	
Set as Default	Current settings are stored and used as default values for all stitching jobs once you click Set as Default . You can edit the default stitching settings under Main menu > Options > Settings > Stitching .	
OK	Click OK to accept the default stitching job options and begin the stitching process.	
Cancel	Closes the window.	

Stitching settings

To edit the default settings for stitching jobs, navigate to the Stitching tab located under **Main menu > Options > Settings > Stitching**.

Ma Settings				×
Application	General			
Microscope Snapshot	Stitching Profile: Image Im	Default		~
XPS	 Tile-to-Tile Contrast Normalization (Experimental) Always Use Calculated Positions 			
Analytics	 Stitch to Project Stitch to File 			
Сгуо	For stitching to file only			
Stitching	Specify output image width 1024 pixels Add scale bar			
		Apply	ОК	Cancel

	Stitching Settings Overview
Settings Item	Description
Stitching Profile	Displays a list of stitching profiles. Default Default Natural Structures Repetitive Structures No Stitching
	 Choices include: Default: The settings in this profile are intended for stitching a broad range of samples. In general, it is recommended that you start with this profile. Natural Structures: This profile is intended for life science or geology samples. The main difference between this profile and the default profile is that this profile has more tolerance in its assessment of alignment validity. Use this profile if the default settings mark too many tiles as invalid while the alignment is okay. Repetitive Structures: This profile is intended for stitching of electronic devices, wafers, or repetitive structures. These settings force the algorithm to be more sensitive to edges and the algorithm uses a different method of assessing alignment validity (that is, since repetitive structures may mislead the image recognition algorithms). No Stitching: The profile will not attempt to align the images. This option prevents any kind of stitching. Use this option to save CPU resources when stitching is not desired, or if stitching will be done by another application.
Align stitched data to stage	For data stitched to project, the final image is aligned to the stage for more accurate measurements and navigation.
Crop stitched image	See Crop Image in the above table.

Stitching Settings Overview(Continued)		
Settings Item	Description	
Blend stitched tiles	See Blend Tiles in the above table.	
Tile-to-Tile Contrast Normalization (Experimental)	See Tile-to-Tile Contrast Normalization (Experimental) in the above table.	
Always Use Calculated Positions	See Always Use Calculated Positions in the above table.	
Stitch to Project	See Stitch to Project in the above table.	
Stitch to File	See Stitch to File in the above table.	
Specify output image width	See Specify output image width in the above table.	
Add scale bar	See Add scale bar in the above table.	

Troubleshoot: Contrast gradients in stitched images

Sometimes uneven illumination results in contrast gradients in stitched images. Individual tiles might appear even, but the stitched image shows contrast gradients (shadow squares). The image below is an example of this issue.



Solutions

- To remove the illumination gradients, use the Tile-to-Tile Contrast Normalization (Experimental) option.
- To resolve this issue on the TEM system, contact your Thermo Fisher Scientific application engineer and request the procedure.

Stitching Profile Best Practices

- Start with the Default profile and only change it if the default settings are not working well enough.
- Before starting a large job, to do a few small runs (e.g., 4 x 4 tiles) and visually inspect the result to find out what the best profile and acquisition settings are. Then use the same acquisition setting on the small runs as on the large run.
- If neither profile gives satisfactory results, try to improve the image quality via the acquisition settings; high quality image data has better chance to be successfully stitched. On SEM, consider using larger dwell times as higher image quality may reveal small features that facilitate image recognition
- If the image data is highly repetitive or very sparse, use a tile size such that there are enough distinct features present in each tile.

- When the success rate of the alignment is not sufficient, try to look at it from an image recognition point-of-view: the algorithm looks at the pixel data in the overlapping border regions; if there are not enough distinct features in that area, the alignment is very difficult. Also think about ambiguity. For example, if the overlapping area only contains a horizontal line, the vertical alignment is trivial but horizontally any position is a valid solution. The tile size, overlap, or the image resolution may need to be increased in order to get data that aligns well.
- On SEM/SDB, recommended overlap for systems with a mechanical stage (for example, Quanta FEG):
 - HFW > 50 μm overlap: 10%
 - HFW 20-50 μm overlap: 15%
 - HFW < 20 μm overlap: 20%

If stitching is still unsuccessful, consider increasing the overlap even more.

- On TEM, recommended overlap:
 - Low Magnification: 20%
 - Magnification 3000x-50000x: 10%-15%
 - Magnification > 50000: 15%-20%

Note: The amount of required overlap is strongly related to how well the system is aligned.

SEM/SDB

This section describes user interface elements and procedures that are specific to the SEM/SDB (Small DualBeam) systems.

Topics include:

SEM User Interface Elements SEM Analytics Cryo Maps Min Maps Min Maps XPS Rotational Alignment SEM and SDB Scan Rotation Stage Rotation Set sample tilt SEM Nav-Cam Alignment

SEM User Interface Elements

SEM Microscope Menu

The Microscope menu only appears for SEM/SDB systems.

Microscope	Options	Help	
Take Snapshot			
Import fr	om NavCa	m	
Get Imag	e from XT	UI	
Set Samp	ole Tilt		0.0°
Restore l	ast Sampl	e Tilt	

Maps Microscope Menu Overview Menu Selection Description Take Acquires a SEM image at the current stage position and HFW and places it Snapshot on the Maps viewer as a snapshot. The HFW can be changed from the SEM/SDB Tile Set Tab. Not available while under vacuum if the door mounted Nav-Cam is installed... Import from NavCam... Acquires and imports a Nav-Cam image, if the Nav-Cam is available. This top down image of the stage can be used for point and click navigation of the sample. See SEM Nav-Cam Alignment. Get Image Imports a SEM image or ColorSEM image (as a single color image), currently from XTUI... displayed in the xT UI without reacquiring the image and places it on the Maps Viewer as a snapshot. Set Sample Unlinks the stage and stores the current sample tilt to the current tilt of the Tilt microscope. When the sample tilt is set, there will be a check mark next to this option and the stage tilt that was set. See Set sample tilt for instructions. **i** Note: Setting sample tilt makes the linked focus unavailable.

	Maps Microscope Menu Overview(Continued)
Menu Selection	Description
Restore Last Sample Tilt	After you make the Set Sample Tilt unavailable, this menu option appears. Click Restore Last Sample Tilt to restore the last sample tilt value.

SEM/SDB Options Menu

- For descriptions of common selections to all systems, see Options menu.
- For descriptions of SEM/SDB-specific selections outlined in purple, see the table below. Note that microscope settings are system specific.



	Options Menu, SEM/SDB Overview
Menu Selection	Description
Microscope Settings	Min Settings Application Acquisition times Image: Display critical acquisition time warning Critical tile acquisition time [mm:ss]: Snapshot Optics Settings Cryo Stage Settings Cryo Stage Settings Beam wake up time [s]: Internal Magnification and Scan Rotation Alignment Magnification correction (Electron): 0% Clear Nav-Cam Clear Nav-Cam Alignment Nav-Cam Clear Nav-Cam Alignment
	 Application times: Display critical acquisition time warning: Select this check box to enable Maps to present the critical time warning as a dialog window that must be dismissed before proceeding. If this option is disabled, then the warning is displayed on the Layer Definition control, but it does not require acknowledgment. Critical tile acquisition time [mm:ss]: Maps displays a warning whenever the acquisition parameters selected for a tile set, result in a per-tile acquisition duration larger than this threshold.

Options Menu, SEM/SDB Overview(Continued)		
Menu Selection	Description	
Microscope	Optics Settings:	
Settings (continued)	 Use Degauss when changing focus: Select the check box to automatically Degauss when changing focus during queued acquisitions. Optics settle delay [ms]: Allows setting a delay (in milliseconds) before acquisition to allow the column action to settle. Without a delay, some 	
	image distortion can occur.	
	Stage Settings:	
	• Stage settle delay [ms]: Allows setting a delay (in milliseconds) before acquisition to allow the stage movement to settle. Without a delay, some image distortion can occur.	
	Electron Beam Settings:	
	• Beam wake up time [s]: Required wait time between commanding the beam to turn on and starting image acquisition. This period of time ensures that the beam has enough time to provide an acceptable signal. If this wait time is too short, then the first tile might have a weak signal and appear black.	
	Magnification and Scan Rotation Alignment:	
	 Magnification correction (Electron): Corrects the magnification factor for the current beam. If the correction is in use and has a value other than zero (that is, a positive or negative value), then the Clear button becomes available to clear the alignment value. 	
	• Scan rotation correction (Electron): Corrects the scan rotation alignment for the current beam. If the alignment is in use and has a value other than zero (that is, a positive or negative value), then the Clear button becomes available to clear the alignment value.	
	Nav-Cam Settings:	
	 Clear Nav-Cam Alignment: Clears the saved default Nav-Cam alignment that is applied to all acquired Nav-Cam images. See SEM Nav-Cam Alignment. 	
	• Door-mounted : Select this check box if the system's Nav-Cam is mounted to the exterior of the chamber door. Do not select the check box if there is a Nav-Cam inside the microscope's vacuum chamber.	

	Options Menu, SEM/SDB Overview(Continued)
Menu Selection	Description
Snapshot Settings	Displays the Snapshot Settings window and the scan settings that will be used to acquire preview images and snapshots.
(Not available for SEM	Ma Settings X Application Electron Beam
Offline)	Microscope Resolution: 1024 x 884 Dwell: 500 ns Frames: Ion Ream
	Analytics Interference Analytics 1024 x 884 Dwell: 500 ns Frames: 1
	Cryo
	Apply OK Cancel
Do Not Change Beam Mode	When this option is selected, it prevents the Maps application from switching the SEM into or out of UHR (immersion) mode. Setting this option to Disabled , allows a queued tile set to automatically activate UHR mode for image acquisition without manual user interaction in the xT UI. In most cases, leave this option set to its default value of Enabled .
	Warning: The Do Not Change Beam Mode feature exists to block the ability of Maps to perform an action with a risk of hardware damage. The SEM should never be put into UHR mode while a magnetic sample is loaded in the chamber, since this could damage equipment inside the microscope. Only make this feature unavailable if you are certain the sample is non-magnetic and you are operating in a use case that requires Maps to change lens modes without user intervention.

	Options Menu, SEM/SDB Overview(Continued)
Menu Selection	Description
Prefer Scan Rotation	If there is viewer rotation when creating a new tile set, the rotation will either be applied as stage rotation or scan rotation.
	 When selected: Uses scan rotation. When not selected: Uses stage rotation.

SEM/SDB Toolbar, Right Side

See Left Side, All Systems for common selections on all systems.

Tool Bar, Right Side, SEM/SDB



	Toolbar Descriptions, SEM/SDB
Tool	Description
Electron Beam (SEM)	Selects the Electron beam for imaging.
(Online version only)	
Ion Beam (FIB) (Online version only)	Selects the Ion beam for imaging.

Toolbar Descriptions, SEM/SDB(Continued)		
Tool	Description	
Import from Nav-Cam	Not available while under vacuum if the door mounted Nav- Cam is installed. Acquires and imports a Nav-Cam image, if the Nav-Cam is available. This top down image of the stage can be used for point and click navigation of the sample. See Nav-Cam Alignment (SEM/SDB Only).	
Take Snapshot	Acquires a SEM image at the current stage position and horizontal field width (HFW), and then places it in the Maps viewer as a snapshot image. The HFW can be changed from the Tile right-click menu selection: Snapshot Settings . See Options menu .	
Import Fluorescence Data	Opens dialog where fluorescence data from optical microscope can be selected and imported to the current Maps project. <i>Note: Requires Cryo features to be activated within the</i> <i>Settings.</i>	
Start Workflow	Opens workflow selection dialog, from where specific alignment workflow can be started.	

SEM/SDB Tile Set Tab

This topic describes options in the Tile Set tab for SEM systems.

Basic Tab



Tile Set Ta	ab, Basic Tab Overview
Control	Description
Name	Displays the name of the current tile set.
Acquisition Type	Specifies the beam to be used for imaging. Choices are: Electron or lon.
Tiles X, Y	Sets the number of tiles in the X and Y directions.

Tile Set Tab, Ba	sic Tab Overview(Continued)
Control	Description
Tile HFW	Specifies the horizontal field width (HFW) for each tile.
Total HFW	Shows the total HFW for all tiles.
Resolution	Specifies the image resolution. Choices are specific to each system type.
Pixel Size	Displays the physical size of a single pixel at the current acquisition settings. This is based on Tile HFW and Resolution.
Dwell	Specifies the amount of time the beam dwells on each pixel when images are acquired.
Frames	Sets the number of image frames.
Image acquisition time	Displays the estimated time required to acquire a single image in the tile set.
Estimated acquisition time	Displays the estimated remaining time required to acquire the whole tile set.
Elapsed acquisition time	Displays the actual elapsed acquisition time.
From Microscope	Applies the current microscope setting.
To Microscope	Applies these settings to the microscope.

Autofunction Tab

0

Note: The list of available autofunctions under the Autofunction tab may differ according to your Maps configuration and connected microscope.

× 🧕	TEMPL	ATE Factory Ter	nplate 🗸
👻 🗹 Layer			
🔽 🌐 Tile Set	AF		Si
⊞ ₹ [4]]	5 \$	Ø	
Name		Mod	e
Contrast Brightness		None	~
Focus		None	~
Lens Alignment		None	~
Focus After Lens Alignmer	nt	None	~
Stigmator Centering		Lower Edge	~ /
Stigmator		None	~
Focus After Stigmator		None	~
Final Contrast Brightness		None	~
Final Focus		None	~
		Cont	figuration

Autofunction	Description
Contrast Brightness	Optimizes the contrast and brightness of the image
Focus	Focuses the electron beam to produce a sharp image (see Focus Mode and Set interpolated focus)
Lens Alignment	Performs the lens alignment procedure
Focus After Lens Alignment	Performs another focus after lens alignment
Stigmator Centering	Performs the stigmator centering procedure
Stigmator	Compensates for any astigmatism
Focus After Stigmator	Performs another focus after stigmation

Autofunction	Description
Final Contrast Brightness	Performs one last contrast and brightness optimization
Final Focus	Performs on last focus optimization

Set autofunctions

- 1. Select the autofunctions to use during acquisition.
- 2. If needed, click **Configuration** to set up a selected autofunction and then click the **play** button to perform a test run of the selected autofunction.
- 3. In the **Mode** combo box, select the mode in which the autofunction should be performed. Available modes include the following:
- None The autofunction is not used.
- Every tile The autofunction is run *locally* at every tile.
- First tile The autofunction is run *once* at the first tile before the acquisition.
- Custom tile list... The autofunction is run over specified list of tiles. Select this option to create a new list.

Create a custom tile list

Click **Custom tile list** option and the selection workflow panel opens and displays instructions to guide you. In the viewer, select tiles where the autofunction should be performed and name the new list. You can use the **Select Nth tile** option in the right panel to quickly select tiles at the selected Nth internal. You can finish the selection with closing the workflow panel or with leaving the autofunction tab. The new list is inserted to the mode combo box and can be re-used by other autofunctions in this tile set. Use the **edit** button if you need to further edit any tile selection.



Advanced Autofunction Settings

Hover your mouse cursor icon over any field name for more information.

Ma Advanced Autofunction S	Settings (Tile Set AF)			×
Contrast Brightness	Focus	Mode	AutoFocus	~
Focus	Absolute Auto scan rotati	HFW	0 µm	
Stigmator	Dwel	Time	5 μs	
Focus After Stigmator	Line Integ Number of F	ration rames	1	
Final Contrast Brightness	Random Po	sition		
Final Focus	Relative	HFW	1.0	
	Reso Use 16-bit	lution mage	1024 x 884	~
	w) Step	1.000 µm	
	Copy settings from: Focus	After S	itigmator 🗸	Сору
		Res	et Reduced Area	Reset
				OK

Focus Mode

Select the Focus Mode to automatically focus before acquisition.

- AutoFocus: AutoFocus mode runs the auto focus routine at each tile before acquisition. This can result in better focus per tile, but adds considerably more time to the acquisition.
- LegacyAutoFocus: Runs the default autofocus algorithm provided by the microscope platform.

	Focus Mode	LegacyAutoFocus	~

Focus Tab

Select the Focus Strategy to focus the electron beam to produce a sharp image.

• None: The application does not touch the focus at all. You are free to change the focus using the xT UI during the run.

⊞ 📮 [AF]	5 \$ ⊘	
Focus Strategy	None	~
Focus Stack		
Top Focus Plane	Not set SET	
Bottom Focus Plane	Not set SET	
Number of Planes	2	
Total Stack Depth	Not set	
Plane Spacing	Not set	

• Fixed: Uses the same focus for every tile.

Focus Strategy	Fixed		~
Focus	10 mm	SET	
Use Focal Plane			

• **Interpolated**: Select three focus points on the sample. The focus plane is then defined by these points. Each tile's focus is set based on where it is located in this plane.



Set interpolated focus

- 1. Navigate to the first location.
- 2. Manually focus the image.
- 3. When image is in focus, click **Set**. The focus is retained for Point 1 and the **Set** checkbox is selected.
- 4. Repeat steps 1-3 for Point 2 (Set) and then for Point 3 (Set). Once all three Set checkboxes are selected, the tiles are ready to be acquired.

Note: These points are used to define a focus plane, so that all tiles use a focus as defined by this plane. Choosing points that are at opposite ends of the sample is important to get the best results.

Clear (unset) a focus point

To clear a focus point, simply toggle Set.

Focus Strategy	Interpolated		~
Point 1		SET	
Point 2		SET	
Point 3		SET	

Focus Stack

Focus stacking is a technique that allows you to acquire a stack of images on differing focal planes. The focus stack control has two states.

State 1

The first state is used to set up an explicit top and bottom working distance for the focus stack.

Next to the Top Focus Plane and the Bottom Focus Plane field, there is a Set button that grabs the current working distance from the XT platform position. You can change the number of planes to acquire in the Number of Planes text box. The Stack Depth is calculated from the setup values and the spacing between each image plane in the Plane Spacing field.

Focus Strategy	None		~
Focus Stack			
Top Focus Plane	10 mm	SET	
Bottom Focus Plane	9.1011 mm	SET	
Number of Planes	3		
Total Stack Depth	898.9008 µm		
Plane Spacing	449.4504 μm		

State 2

The second state of the focus stack control is the relative focus setup. This is the automatic state of control if you are using a Fixed, Interpolated, or Auto Focus autofunction. In this state, the focal planes are relative to focus values set from the fixed, interpolated, or auto focus.

In the second state, you can set the Number of Planes and the Total Stack Depth. After setting the planes and stack depth, you can see the relative top and bottom planes change. A small notification appears at the bottom of the control pane that notifies what the focus stack is based upon.

Focus Stack		
Top Focus Plane	+449.4504 μm	SET
Bottom Focus Plane	-449.4504 µm	SET
Number of Planes	3	
Total Stack Depth	898.9008 µm	
Plane Spacing 449.4504 µm		
The top and bottom focus planes are offsets from the Fixed working distance.		

Focus Tab Overview

Control	Description
Focus Stack	Perform the focus stack acquisition.
Top Focus Plane	The working distance for the top plane of the focus stack.
Bottom Focus Plane	The working distance for the bottom plane of the focus stack.
Number of Planes	The number of focus planes acquired in the focus stack.
Total Stack Depth	The physical depth of the total focus stack.
Plane Spacing	The physical spacing between each focus plane in the focus stack.

Advanced Tab

Use the controls on this tab to set advanced tile acquisition properties for each tile set.

⊞ 🕂 [AF]	-	٩ţ٩	Ø		
Line Integration	1				
Scan Interlacing		1	1		
Overlap X, Y		10%		10%	
		205 px		177 рх	
Stage X		440.58	74 µm		
Stage Y	. -	-881.1	749 µm		
Scan Rotation	L	0			
High Voltage				~	
Beam Current				~	
Beam Mode				~	
Drift Corrected					
16-bit Images					
Stage R		0			
Stage T					
Stage Z					
Order		Raster		~	
Maximum Run Time					
Display Acquisition Selection:					
Display 1		Di	splay 2		
Store Detector			Store Deteo	tor	
Display 3		Di	splay 4		
Store Detector			Store Detec	tor	
Stitching Display		Display	y 1	~	

Tile Set Tab, Advanced Tab Overview		
Control	Description	
Line Integration	Specifies the number of times each raster line is scanned and integrated to combat certain types of drift.	
Scan Interlacing	Specifies the number of times each frame is interlaced to combat certain types of drift.	
Overlap X, Y	Displays the percentage of overlap between tiles in the tile set, as well as the pixels.	
Stage X	Sets the X stage location of the center of the tile grid.	
Stage Y	Sets the Y stage location of the center of the tile grid.	
Scan Rotation	Allows manual adjustment of scan rotation for a tile set. Either enter a value in the edit box here or rotate the tile set in the viewer to update the scan rotation value.	
High Voltage	Selects value of the high voltage setting.	
Beam Current	Selects value of the total amount of electron or ion current striking the sample.	
Beam Mode	Selects the electron beam lens mode.	
Drift Corrected	If selected the correction of the image drift is used when the number of Frames is higher than 1.	
16-bit image	When selected, saves image as 16-bit (65K grayscale levels), else image is saved as 8-bit (256 grayscale levels).	
Stage R	Sets the stage rotation where the tiles are acquired. Existing tile sets rotate with the stage. New tile grids are created at the current rotation. Select the check box to enable the text field to change the default rotation.	
Stage T	Sets the stage tilt where the tiles are acquired. Select the check box to enable the text field to change the default tilt.	

Tile Set Tab, Advanced Tab Overview(Continued)			
Control	Description		
Stage Z	Sets the stage height (Z) to be stored and used for acquisition. This allows multiple tile sets to be acquired at different heights, that is sometimes necessary for using different detectors. If left unselected, the stage height is not changed.		
Order	 Determines the order in which the contained tiles are acquired. Options are: Raster: Left to right, then top to bottom. Serpentine: Alternating left to right pattern per row, running top to bottom. Spiral In: Iterating over tiles in a clockwise pattern starting with the outer most tiles moving inward. Spiral Out: Iterating over tiles in a clockwise pattern starting with the inner most tiles moving outward. 		
Maximum Run Time	Sets the maximum time spent acquiring data for this tile set. The tile set will run until this time limit has been reached, or if all the tiles have been acquired. The time limit entered is only used by this tile set and the countdown starts when this tile set starts collecting images.		
Display (1-4)	Select multiple displays to acquire images from multiple display channels.		
Store Detector	Selects the method for acquiring detector settings. You can set up different tile sets with different detectors and line them up for acquisition. When selected, a label displays the current setting.		
Stitching Display	Select the display used for stitching. Choose the display with the most detail of the sample for best results.		

SEM Tile Set Context Menu

Right-click on any tile set to access the context menu.

Alignment	Þ
Drive To	
Square Up	
Use For Magnification And Scan Rotation Alignment	
Apply Settings To Microscope	
Apply Microscope Settings To Grid	

- For descriptions of common selections to all systems, see Tile Set Context Menu.
- For descriptions of SEM/SDB-specific selections, see the two items outlined above as well as the table below.

Menu Selection	Description
Square Up	Rotates the stage to match the orientation of the tile set.
Use for Magnification And Scan Rotation Alignment	When selected, uses the offsets calculated in the acquired tile set to correct for misalignment between the scan axes and the stage axes by applying a small magnification and scan rotation. This will allow better acquisition and alignment of future tile sets.

SEM Nav-Cam Context Menu

The Nav-Cam image is stored in the selected layer. Right-click on a Nav-Cam entry to access a context menu.

Bring To Front	
Bring Forward	
Send Back	
Send To Back	
Center View	Ctrl+Shift+C
Center and Rotate View	
Export To Project	
Alignment	•
Persist Nav-Cam Alignment	
Delete	

- For descriptions of common selections to all systems, see Annotation Context Menu.
- For descriptions of SEM/SDB-specific selections, see the table below.

	Nav-Cam	Right-Click	Menu	Overview
--	---------	-------------	------	----------

Menu Selection	Description
Persist Nav-Cam Alignment	Keeps the current Nav-Cam alignment.

SEM/SDB Viewer Tile Context Menu

Right-click on a tile within the tile grid to access a context menu.

SEM/SDB

Drive Stage Here			
Rotational Alignment	•		
Snapshot Here			
Snapshot Here w/HFW	·		
Add Tiles Here			
Add Site of interest			
Tile Set (4)	•		
Tile 10,10 🔹 🕨		~	Acquire
Clear Digital Rotation			Open tile image folder
5			Open source image
		Γ	Preview Tile(s)
		L	Sub-tile 🕨
			Open analytics data folder

- For common selections, see SEM/SDB Viewer Tile Context Menu.
- For SEM/SDB-specific selections, see the four items outlined above as well as the table below.

	Viewer Tile Context Menu, SEM/SDB Overview
Menu Selection	Description
Rotational Alignment	Allows rotation of the stage to Place Point 1 , Place Point 2 , or Cancel . See Rotational Alignment.
View	er Tile Context Menu, SEM/SDB Overview(Continued)
-------------------------------------	---
Menu Selection	Description
Snapshot Here with HFW	Acquires a preview SEM image at the selected stage coordinate and current HFW. Submenu choices:
	2 μm 10 μm 50 μm 100 μm 200 μm 500 μm 1 mm This is the same functionality as using the Take Snapshot Image tool
	box button. See Tool Bar, Right Side, SEM/SDB.
Acquire/Queue for re-acquisition	 Displays whether or not the tile has already been acquired as defined below. You can toggle between the two tile acquisition states as needed; however, only the <i>selected</i> acquisition state appears in the menu. Acquire: This acquisition state displays if the tile <i>has not been acquired</i>. If you select the check box for this menu item, then Maps queues the selected tile to be acquired. Queue for re-acquisition: This acquisition state displays if the tile <i>has already been acquired</i>. If you select the check box for this menu item, then tile <i>has already been acquired</i>. If you select the check box for this menu item, then tile <i>has already been acquired</i>. If you select the check box for this menu item, then Maps queues the selected tile to be re-acquired and replaces the original image.
Preview Tile(s)	Acquires preview images for all selected tiles. Any tiles without pertinent image data can be disabled from image acquisition.

Viewe	er Tile Context Menu, SEM/SDB Overview(Continued)
Menu Selection	Description
Sub-tile	Displays a submenu for selecting the number of sub-tiles to be in a mini grid within an individual tile, using the selected number of images. 2 x 2 3 x 3 5 x 5 10 x 10 25 x 25 50 x 50 100 x 100 The sub-tile names are automatically added to the Layers control (TileSet: (3,4)).

SEM/SDB Post-Acquisition Actions

This section describes the post-acquisition actions for SEM/SBD microscopes.

POS	r acquisition
	Turn Off Beam When Done
	Retract Detectors When Done
\bigcirc	Sleep When Done
	Pump to HiVac When Done

Post-Acquisition Actions

Control	Description
Turn Off Beam When Done (Not available for SEM Offline)	Turns off the beam after all jobs are completed.

Post-Acquisition Actions(Continued)				
Control	Description			
Retract Detectors when Done (Not available for SEM Offline)	Retracts any insertable detectors after all jobs are completed			
Sleep When Done (Not available for SEM Offline)	Puts the system into Sleep mode after all jobs are completed.			
Pump to HiVac When Done	Returns the microscope to HiVac mode after completing a LowVac acquisition.			

SEM Analytics

With the Analytics plug-in application, the Maps software can acquire EDS data through a software application interface on a SEM/SDB microscope. EDS is an analytical technique that enables you to collect information about chemical composition and elemental analysis. This section describes how you can use Maps to collect this information in large area mapping scenarios.

Note: This plug-in is not to be confused with the Maps Min that can also acquire EDS data on SEM microscopes.

Preparation for SEM Analytics

Preconditions

A

To run Analytics, your system must include the following components:

- EDS detector
- SEM platform with ChemiSEM support and license
- Valid Maps license with XT Online and Analytics options

To verify that your Maps license includes the necessary options, on the Help menu, click About.

thermoscientific	Maps 3.	31
	Product	Version
	r logaci	
	Application	3.31.0.3619
	smaging Com Munic	7.3.0.2080
	Sem Plugin	3.31.0.3019
	VPS Plugin	3.31.0.3019
	Avizo20 Plunin	3 31 0 3619
	Avizo Bridge Plugin	3,31,0,3619
	ringe en age ringen	2.2.1.2.2.1.2
	Licensed Options	Expiration Date
	XT Online	Unlimited
	External Image Import	Unlimited
	 Stage Correlation 	Unlimited
	Color Correlation	Unlimited
	 Stitching 	Unlimited
	Array Tomography	Unlimited
		Or mining of

Indication that Analytics is available

If all of the preconditions are satisfied, you can observe the following in the status bar:

- Analytics button is visible
- The color indicator on the button is green



Analytics Workflow for SEM

Proceed with the workflow in the following order:

- 1. Define a Tile Set: Use the normal procedure to define a tile set, and then navigate to the tile set section and click the **Analytics** tab.
- 2. Configure Analytics for EDS Acquisition: Once you have defined the tile set, you must configure the parameters on the Analytics tab before you acquire the EDS tile set.
- 3. Acquire EDS Tile Set: When you acquire the EDS tile set, you can observe the status of the EDS acquisition and EDS count data during acquisition.
- 4. View Results of EDS Acquisition: Once an Analytics tile set has finished EDS acquisition, you can view the results in the layer tree, Visualization tab, and main Viewer.
- 5. Reprocess EDS Tile Set: You can reprocess the spectrum data that is saved within the tile set for a different selection of chemical elements and EDS maps.
- 6. Stitch EDS Tile Set: You can stitch EDS tile set in the same way as any other tile set.

Configure Analytics for EDS Acquisition

Note: If the **Analytics** tab is not visible, one or more Analytics preconditions is not satisfied. Refer to **Preparation for SEM Analytics** for more information.

Follow the steps below to configure the parameters on the Analytics tab.

1. Click the **Analytics tab**, and then select the **Perform EDS Acquisition** checkbox to activate the tile set for EDS acquisition.



- 2. Configure the EDS acquisition parameters.
- a. Observe the **Resolution** of the EDS maps. Note that EDS resolution is the same as the selected tile set resolution up to the maximum supported EDS resolution, which is 512 x 442 or 768 x 512 based on the aspect ratio of the tile set resolution.
- b. Enter the **Dwell** time and number of **Frames** to achieve satisfactory EDS counts during acquisition. Note that **Acquisition Time** and **Frame Time** are automatically computed.
- c. Locate and adjust the EDS Pixel Size parameter if necessary. Note that any changes to EDS Pixel Size parameter affect the HFW of the tile set. Adjust the Tile HFW parameter to match the specified EDS pixel size.

⊞ 📮 [AF]	₹	ھ لي	\oslash	
Perform EDS Acquisition				
Acquisition Parameters				
Resolution		512 x 442		
Acquisition Time	ê	3.62 s		
Dwell	G	500 ns		
Frames		2		
Frame Time		1.81 s		
EDS Pixel Size		1 µm		
Tile HFW		512 µm		
Analusia Davanatara				
		AL CH AN		
Count Mana		AI, CU, Ag		
Count Maps		✓ 089.29 MB		
Quant Maps		✓ 689.29 MB		
Quantification Type		Weight Percent	age	~
Segmentation Type		Object-based		~
Normalize Maps		2		
Enable EDS Reprocessing		🗹 45.94 MB		
	Repr	rocess		

3. Locate the **Elements** parameter under the Analysis Parameters section of the Analytics tab.

E - [AF]	≢ [ی لب	\oslash	
Perform EDS Acquisition		•		
Acquisition Parameters				
Resolution		512 x 442		
Acquisition Time	ê	3.62 s		
Dwell	6	500 ns		
Frames		2		
Frame Time		1.81 s		
EDS Pixel Size		1 µm		
Tile HFW	512 µm			
Analysis Parameters				
Elements		Al, Cu, Ag		
Count Maps		🗹 689.29 MB		
Quant Maps		🗹 689.29 MB		
Quantification Type		Weight Percenta	ge	~
Segmentation Type		Object-based		~
Normalize Maps				
Enable EDS Reprocessing		🗹 45.94 MB		
	Repr	ocess		

- 4. Click the ellipsis icon to open the Element Selector.
- a. Select the individual elements that you want to identify during EDS acquisition by selecting them with a mouse click.
- b. Optional: Associate a specific color with each selected element by right-clicking the element and selecting the desired color from color picker. The selected color is used for the EDS map generated for given element. If a color is not associated with the element, the color is automatically selected from the system palette to match the color that you typically see in the xT UI.



5. The acquired X-ray spectrum is processed and displayed as a color map for each selected element. There are two types of EDS maps that are currently supported: **Count Maps** and **Quant Maps**. To select the EDS maps associate with your specific type, do the following:

- a. For Count Maps, select one or more EDS maps to be generated after acquisition by selecting the corresponding checkbox. If you do not select any maps, then you must activate EDS reprocessing on the tile set so that processing can occur at a later point.
- b. For Quant Maps, specify the quantification parameters to be used for processing of the EDS spectrum:
 - Quantification Type: available values include Atomic Percentage and Weight Percentage.
 - Segmentation Type: available values include Object-based and Regular.
- c. For each selected type of map, observe the estimated disk size necessary for storing the maps after acquisition. If you anticipate that the estimated disk size will be exceded, consider deselecting some elements if they are not critical for the experiment.

6. Observe that the **Normalize Maps** checkbox is selected by default. This option allows you to normalize the intensity of EDS maps across the entire tile set. See **Reprocess EDS Tile Set** for more details.

7. Select the **Enable EDS Reprocessing** checkbox to include EDS reprocessing data in the tile set after acquisition. This allows you to reprocess the spectrum data later for a different selection of chemical elements and EDS maps. If the data is not saved, reprocessing will not be available for this tile set.

• Note: If you select this option, the disk size estimate is displayed next to the checkbox. Note that the exact size is not available before acquisition, so the estimate is refreshed during EDS acquisition based on the measured data.

Acquire EDS Tile Set

Follow the steps below to acquire the EDS tile set.

- 1. Follow the standard tile set acquisition procedure described in Acquire a Tile Set.
- 2. Observe status of the EDS acquisition in the following:
 - Progress bar and remaining time in the Acquisition job queue
 - Status notifications in the status bar
 - Remaining time in the status bar

ACQUISITION PROCES	SING			
			DESELECT ALL SELECT ALL	
			Show completed jobs	
Tile Set (2)	18%	Acquiring images	00:02:35 🜌	
ESTIMATED DISK SPACE 1.09 GB / 2.51 GB				
POST ACQUISITION		POST PROCESSING		
Turn Off Beam When Done		Stitch Tile Sets When D	Done	
Retract Detectors When Done				
Sleep When Done				
Pump to HiVac When Done				
V JOB CONTROL	P		2:35	

3. Click the **Analytics** button on the status bar. A dialog will then open for you to view data about incoming counts, tile counts, and tile set counts during the EDS acquisition.



View Results of EDS Acquisition

Once EDS tile set has finished acquisition and processing, you can view the results in the layer tree, Visualization tab, Analytics dialog, and main Viewer (see Layer Tree Tabs).

View results in the layer tree

In the layer tree, the tile set has a collection of layers within it that corresponds to each processed element and SEM image. The inner layers are represented as color channels, and are organized within Quant Maps and Count Maps groups.



- 1. Expand the group by clicking on the small arrow next to the group name. The layer group is expanded to show the list of corresponding element channels.
- 2. Select the checkbox on the left side of the layer to toggle visibility of the element channel or group of channels.
- 3. Click the colored circle on the right to change the color of the element channel.

View results in the Visualization tab

The Visualization tab of the tile set shows the list of color channels that corresponds to each processed element and SEM image. The channels are organized within Quant Maps and Count Maps groups.



- 1. Expand the group by clicking in the group header. The group is expanded to show the list of corresponding element channels.
- 2. Expand the individual element channels by clicking in the channel header. In the expanded control, you can modify the visualization properties of the channel.

- 3. Select the checkbox on the left side of the group or channel to toggle visibility of the element channel or group of channels.
- 4. Click the colored circle on the right to change the color of the element channel or group of channels.



View results in main Viewer

The overall result of the EDS acquisition is displayed in the main viewer by blending the color maps that correspond to each processed element and SEM image. You can control what maps are included in the blended result by toggling visibility of the element channels in the layer tree or the Visualization tab.



View quantification data and summary of counts in Analytics dialog

The following additional information is collected during EDS acquisition:

- Summary of incoming counts
- Quantification Data: This is included if you have selected the Quant Maps analysis on the Analytics tab.

You can observe this additional information by following the steps below:

- 1. Go to the status bar and click **Analytics** to open the Analytics dialog.
- 2. Select one of the tiles within the acquired EDS tile set in the main Viewer.
- 3. When the Analytics dialog opens, notice the name of the selected *tile set* and *tile* coordinates in the header of the Analytics dialog.

Tile Set Tile[2,3]				
EDS Coun	Q	uantification Data		
Counts Per Frame	5 089	Element	Intensity	At%
Tile Counts	10 178	Al		100 %
Tile Set Counts	43 060	💛 Cu		100 %
Average Counts Per Tile	21 530	Ag		3%
Detector St	atus			
Count Rate	17228 cps			
Detector Dead Time	11.18 %			_
Analytics 🌑 🗸	Scan rotation	0.0	🔨 🗹 Link	Stage rotation

4. View the quantification data measured on the selected *tile* in the Quantification Data panel.

- 5. The elements are ordered by maximum concentration on the *tile set*, and for each element you can view the following:
 - Element: This displays the name and color of the element.
 - Intensity gradient: The range of intensities within the Quant Maps on the selected *tile set*.
 - White marker on the intensity gradient: This marks the highest intensity that is associated with maximum element concentration on the selected *tile*.
 - Maximum concentration in At% or Wt%: This percentage is the maximum concentration for a given element measured on the selected *tile*. This percentage also indicates that there is at least one pixel on some part of the selected *tile* with that concentration level. Note that the unit of measurement (At% or Wt%) depends upon the selected Quantification Type on the Analytics tab.
- 6. View the summary of detected counts in the EDS Counts panel of the dialog.

Reprocess EDS Tile Set

The reprocessing functionality allows you to reprocess the spectrum data that are saved within the tile set for a different selection of chemical elements and EDS maps.

 Note: EDS reprocessing must be activated for the tile set in order for this functionality to be available. Refer to Configure Analytics for EDS Acquisition for more information.

You can include multiple changes, as described in the sections below, within the same reprocessing operation:

- Change the selection of elements, EDS maps, or normalization option.
- Observe that warning is displayed on the Analytics tab and the Reprocess button after each individual change: "EDS parameters have been changed. Reprocess the tile set to reflect the changes."
- Continue adjusting EDS parameters until you complete all intended changes.
- Click the Reprocess button to reprocess the tile set and have all changes applied to the tile set within the same EDS reprocessing job.

Change selection of elements

1. Locate the **Elements** parameter and click the ellipsis icon to open the Element Selector.

⊞ 📮 [AF] ∃	≝ 💷 👁 🗵
Perform EDS Acquisition	
Acquisition Parameters	
Resolution	512 x 442
Acquisition Time	
Dwell	Ω 1 μs
Frames	2
Frame Time	3.62 s
EDS Pixel Size	1.9531 µm
Tile HFW	1000 µm
Analysis Parameters	
Elements	Al, Cu, Ag
Count Maps	🗹 1.13 GB
Quant Maps	🗹 1.13 GB
Quantification Type	Atomic Percentage 🗸 🗸
Segmentation Type	Object-based V
Normalize Maps	
Enable EDS Reprocessing	S9.67 MB
	Reprocess

2. Observe the elements that are currently selected for EDS processing.

3. Choose new elements that you want to include in EDS reprocessing by selecting them with a left mouse click. Remove elements that you no longer want to include in EDS reprocessing by clicking the selected element with a left mouse click again.

4. Optional: Change the colors that are associated with each selected element by right-clicking on the element and choosing the desired color from color picker. You can also remove color that was previously associated with the element by clicking the Reset button in the color picker.

Change selection and parameters of EDS maps

- 1. Select or deselect **Count Maps** and **Quant Maps** checkboxes to include or exclude given type of EDS maps from EDS reprocessing.
- 2. For Quant Maps, you can also change **Quantification Type** and **Segmentation Type** that should be used for EDS reprocessing.

Analysis Parameters		
Elements	Al, Cu, Ag	
Count Maps	🖌 1.13 GB	
Quant Maps	🗹 1.13 GB	
Quantification Type	Atomic Percentage	~
Segmentation Type	Object-based	~
Normalize Maps		
Enable EDS Reprocessing	🗹 59.67 MB	
	Reprocess	

Normalize EDS maps

Select the **Normalize Maps** option to have the intensity of the EDS maps normalized across the whole tile set during EDS reprocessing.



The following occurs during the normalization process:

- The highest concentration of each element over the entire tile set is determined.
- The intesity maps are adjusted across all tiles to make the intensity ranges for given element relative to the global maximum concentration.
- Visually, this process ensures that the intensity values for a given element are consistent and comparable across different tiles.

Keep or remove EDS reprocessing data

Deselect the **Enable EDS Reprocessing** checkbox if you no longer want to include EDS reprocessing data in the tile set. This will save storage space necessary for the project, but you will no longer be able to reprocess the tile set. If you deselect the option, final reprocessing will be performed after which the EDS reprocessing data will be removed.

Reprocess the tile set

1. Reprocess the tile set by clicking **Reprocess** at the bottom of the Analytics tab.

Analysis Parameters			
Elements	Al, Cu, Ag		
Count Maps	🗾 1.13 GB		
Quant Maps	🗹 1.13 GB		
Quantification Type	Atomic Percentage	~	
Segmentation Type	Object-based	~	
Normalize Maps			
Enable EDS Reprocessing	🛃 59.67 MB		
Reprocess			
Warning: EDS parameters have been changed. Reprocess the tile set to reflect the changes			

2. Observe the status of the reprocessing job in the Processing job queue.

ACQUISITION PROCESSING			
STOP PROCESSI			Show completed jobs
Processing EDS for: Tile Set 💻	11%	Processing tile (2, 1)	00:03:20
ESTIMATED DISK SPACE			
0 bytes / 2.53 GB			
V JOB CONTROL NUN	\bigcirc		

Stitch EDS Tile Set

You can stitch an EDS tile set in the same way as any other tile set. Intensity maps for the selected elements are stitched during the process as well and available in the stitched result. SEM electron image is used as the default stitching channel, and it is used to align tiles based on the overlaps.

View structure of the stitched EDS layer

In the layer tree, the stitched EDS layer has the same structure as the original EDS tile set. The stitched EDS layer contains a collection of layers that correspond to each processed element and SEM image. The inner layers are represented as color channels, and these inner layers are organized within Quant Maps and Count Maps groups.



- 1. Expand the group by clicking on the small arrow next to the group name. The layer group is expanded to show the list of corresponding element channels.
- 2. Select the checkbox on the left side of the layer to toggle visibility of the element channel or group of channels.
- 3. Click the colored circle on the right to change the color of the element channel.

View quantification data in Analytics dialog

The stitched EDS layer also contains quantification data, if the following conditions are satisfied:

- The original tile set has Quant Maps analysis selected
- The original tile set is normalized

You can observe the quantification data by following the steps below.

- 1. Go to the status bar and click **Analytics** to open the Analytics dialog.
- 2. Select the stitched EDS layer in the layer tree, or select it in the main Viewer.
- 3. When the Analytics dialog opens, the name of your selected layer is displayed in the header and the normalized quantification data is displayed in the Quantification Data panel.

Tile Set (stitched))			
EDS Coun	ts	Q	uantification Data	
Counts Per Frame		Element	Intensity	At%
Tile Counts		🔵 AI	_	100 %
Tile Set Counts		💛 Cu		100 %
Average Counts Per Tile		Ag		100 %
Detector St	atus			
Count Rate	- cps			
Detector Dead Time	5.86 %			

The elements are ordered by maximum concentration in the Quantification Data panel, and for each element you can view the following:

- Element: This displays the name and color of the element.
- Intensity gradient: The range of intensities within the Quant Maps.
- White marker on the intensity gradient: This marks the highest intensity that is associated with the maximum element concentration in the stitched layer.
- Maximum concentration in At% or Wt%: This percentage is the maximum concentration for a given element measured on the stitched layer. This percentage also indicates that there is at least one pixel on some part of the sample with that concentration level. Note that the unit of measurement (At% or Wt%) depends upon the selected Quantification Type on the Analytics tab.

Cryo

The Maps Cryo workflow is a collection of controls to help create a list of sites that support the cryo lamella TEM prep workflow.

In the Cryo workflow, you can target a collection of cells on a frozen sample grid by adding a cryo lamella site at each cell of interest.

The Maps Cryo workflow consists of the following tasks:

- 1. (optional) Import and view fluorescence data.
- 2. Set sample tilt to make the cryo sample perpendicular to electron beam for imaging.

This procedure is required for the Cryo workflow because the cryo samples are on holders that are tilted. To image the sample correctly with the Maps software, the sample must be perpendicular to the electron beam at all times.

- 3. Add lamella sites.
- 4. Calculate the eucentric position of lamella sites.
- 5. Determine cryo milling position.
- 6. Perform GIS deposition. When you have a collection of lamella sites fully setup and ready for milling, you must then perform a GIS deposition on the sample to prepare the sample for milling.

GIS deposition is performed after setting up the sites because it is hard to see cells of interest with a layer of a deposited substance covering them.

After you complete the workflow, you can analyze the milled-in-place lamellae in a TEM.

Configure Cryo settings

The Cryo settings allow you to activate the Cryo features and then select from the options described below.



Turn on Cryo features

- 1. Navigate to the Cryo Settings tab. On the **Options** menu, click **Settings** and then click the **Cryo** tab.
- 2. Click **Enable Cryo** to activate the Cryo features. The Maps application alerts you if a system restart is required to proceed. When the Maps software is in Cryo mode, then all the Cryo controls are available.

Choose a directory to watch

This option can speed up attaching files to selected lamella sites within Maps software. You can set this to the directory where xT saves images. When files are added to that directory and a lamella site is selected in the Maps software, you are prompted to attach the files to the selected lamella site timeline. See Cryo lamella timeline for more about the timeline.

- 1. Navigate to the Cryo Settings tab. On the **Options** menu, click **Settings** and then click the **Cryo** tab.
- 2. Select **Watch Directory** and then browse to a folder to choose a directory for Maps software to watch for new files.

Choose a lamella site color

- 1. Navigate to the Cryo Settings tab. On the **Options** menu, click **Settings** and then click the **Cryo** tab.
- 2. Click the eye dropper to choose the initial color used for a lamella site in the viewer when it is created.

Default lamella site color:

Choose a default import directory for fluorescence data

The Leica Microsystems' Application Suite X (LAS X) software can export fluorescence data specifically for importation into the Maps software. The fluorescence data is a collection of images and a special XML file with "{CLEM_Positions_TFS}" in its name within the same directory. If you use this feature, then you can choose a default directory for Maps software to look for Leica fluorescence data import files.

- 1. Navigate to the Cryo Settings tab. On the **Options** menu, click **Settings** and then click the **Cryo** tab.
- 2. In the **Default Fluorescence Import Directory** field, browse to and select the directory where your fluorescence data is stored. This directory is used as the starting directory when you **Import and view fluorescence data**.

Choose a default import file type for fluorescence data

The open file dialog defaults to a file filter for the selected fluorescence data type. Select the data type that will be imported most often.

Options include:

- **iFLM**: Integrated fluorescence light microscope, available on some Thermo Fisher Scientific systems.
- Leica: Fluorescence data from the Leica Microsystems' LAS X software.

Import and view fluorescence data

You can import fluorescence data and then view it within the Optical Data Browser.

Import procedure

When the Maps Cryo features are turned on, a command for importing fluorescence data is added to the Maps toolbar.



Prerequisites:

- Requires Cryo features to be activated within the Settings.
- Have fluorescence data available.
- 1. Click the Import Fluorescence Data icon within the Maps toolbar.



- 2. Select an image to import, and then Click Open.
- 3. The Maps software imports the fluorescence data.

View fluorescence data

When you select multilayer fluorescence data, hover your mouse cursor over the slider and move your mouse scroll wheel up/down to browse through the image layers displayed in the Optical Data Browser. Alternatively, you can use the Page Up and Page Down keys on your keyboard to browse through the image layers. When the slider has focus, use the Arrow keys on the keyboard to change the value by 1 and use the Page Up and Page Down keys to change the value by 10%.

Maps displays the actual position of the image planes in microns and adds tick marks in the browser's main viewer to display individual image planes for more accurate navigation. If Z coordinates are available within the imported file, then you can view the minimum and maximum depth values at each end of the Depth control slider. If Z coordinates are *not available* in the imported file, then a plane number value appears for the Plane control.



Import and view Arctis data

You can import and view Arctis data in the Maps software application on TEM systems. The Arctis data include fluorescence images and information about milled lamellae.

Note: This Cryo functionalilty is specific to TEM systems and is not available on SEM systems.

Import procedure

When the Maps Cryo features are turned on, a command for importing Arctis data is added to the Maps toolbar.



Prerequisites:

- Requires Cryo features to be activated within the Settings.
- Have Arctis data available.
- 1. Click the Import Arctis Data icon within the Maps toolbar.



- 2. Select Arctis data to import, and then Click **Open**.
- 3. The Maps software imports the Arctis data.

View individual lamella sites

You can view individual lamella sites by expanding the corresponding lamella group in the layer tree. You will find electron images, optical images, and a lamella annotation within the group. When selecting the lamella annotation, you can see information about the milling angle, lamella thickness, and rating.



Add lamella sites

There are two ways to add lamella sites in Maps software. One way is to find a spot on images in the Maps software using the viewer. The other is to drive to the spot within xT and create the site based on the current microscope field of view position.

Prerequisites:

• Requires Cryo features to be activated within the Settings.

Add a lamella site in Maps software

This is the preferred/recommended method when you already have data in Maps and can find the desired site in the viewer.

1. In the Maps viewer, navigate to the location where you want to make a lamella site.

2. In the Maps viewer, right-click where you want to create a new lamella site and then click Add Lamella Site Here.



 \oslash Name Lamella ð Mapping Position Drive To Update Ø **Eucentric Position** Calculate Update **Milling Position** Store Angle Update Milling Angle Not Set Success Failed

Maps software creates a new lamella site at that location and adds it to the active layer.

Add a lamella site using the current stage position.

Use this method when you do not have enough data in the Maps software to see the desired site from within the viewer. This method uses xT to find and navigate to the site of interest.

1. In xT, find and drive to the site of interest and center it in the field of view (FOV).

2. In the Maps viewer, right-click anywhere and then click Add Lamella Site On FOV Position.

Drive Stage Here	
Rotational Alignment	•
Snapshot Here	
Snapshot Here w/HFW	•
Add Lamella Site Here	
Add Lamella Site On FOV Position	
Add Tiles Here	
Add Site of interest	
No tile set selected	•
No tile selected	•

Maps software creates a new lamella site in the center of the current microscope field of view and adds it to the active layer.



Calculate the eucentric position

Maps software can help you determine the eucentric position of a lamella site for the Cryo workflow. *Eucentric position* refers to a position where the sample can be tilted, and the area of interest does not move out of the field of view or go out of focus.

Prerequisites:

- Requires Cryo features to be activated within the Settings.
- At least one lamella site has been created.
- Stage tilt set for current sample. See Set sample tilt for instructions.

This procedure can be performed only for lamella sites in the Cryo workflow. If you have calculated the eucentric position and you want to change it, then you can refine it using the same procedure.

To calculate or refine the eucentric position

- 1. Click the lamella site in the Layer tree to select it.
- 2. In the Lamella Site pane, next to Eucentric Position, click Calculate.



If the eucentric position has already been calculated for this site, then click **Refine Eucentric Position**.



3. Maps software prompts you to center the feature under the electron beam in xT. When finished, return to Maps software and click OK.

- 4. To calculate or refine the eucentric position of the site, perform a series of stage tilts, and recenter the feature at each tilt position.
 - a. On the dropdown list, click the tilt increment you want to use to tilt the stage. You can change the tilt step at any point while you are calculating. Starting with smaller tilt steps might make this task easier.



- b. Use the arrow buttons on either side of the dropdown list to tilt the stage up or down by the selected tilt step.
- c. Once the stage is done tilting, recenter your feature in xT.
- d. Click Calculate Eucentric Position.

Repeat the previous step as needed to increase the accuracy of the eucentric position. If the eucentric position is calculated correctly, then the feature stays centered in xT view window as you change the stage tilt.

5. Once you see the feature stays in the center of the display in xT between tilts of at least +/-10 degrees, you can close the **Eucentric Position** dialog.

The eucentric position is now calculated and stored within the lamella site in Maps software.

Name	Lamella		i.
Mapping Position	Drive To	Update	Ø
Eucentric Position	Drive To	Update	\oslash
	Refine Eucentric Position		
Milling Position	Store Angle	Update	()
Milling Angle Not Set			
Success		Failed	

Determine cryo milling position

The *milling position* is the stage tilt and position to be used when milling the lamella. You must determine this position before you can Perform GIS deposition.

Prerequisites:

- Requires Cryo features to be activated within the Settings.
- At least one lamella site has been created and the eucentric position has been calculated.

The yellow warning symbol indicates that the milling position needs attention.

SEM/SDB

Name	Lamella		d d
Mapping Position	Drive To	Update	\oslash
Eucentric Position	Drive To	Update	\oslash
	Refine Eucentric Position		
Milling Position	Store Angle	Update	()
Milling Angle Not Set			
Success		Failed	

To determine the milling position:

- 1. Select the desired lamella site.
- 2. Click Drive To for Eucentric Position.
- 3. In xT, change the stage tilt to the angle at which you want to mill a lamella.
- 4. In Maps software, click **Store Angle** to finish setting up the milling position. Maps software stores the angle. The lamella site is now fully setup.
5. The green check mark symbol confirms that you have successfully determined the milling position.

Name	Lamella		i de la constante da la consta
Mapping Position	Drive To	Update	Ø
Eucentric Position	Drive To	Update	\bigotimes
	Refine Eucen	tric Position	
Milling Position	Drive To	Update	\oslash
Milling Angle	48.0°		
Success		Failed	

Perform GIS deposition

While creating sites, the sample tilt is set and the stage is unlinked. xT requires the stage to be linked when performing deposition. To make sure you do not accidentally perform any unlinked moves with the Maps software during deposition, you must lock the Maps user interface until you are done with deposition in xT.

Prerequisites:

- Requires Cryo features to be activated within the Settings.
- At least one lamella site must be fully setup and ready for milling.

1. On the Microscope menu click **GIS Deposition**.

Note: This command only locks out the Maps user interface. It does not actually perform any deposition.



- 2. Using the xT software, perform a GIS deposition.
- 3. When deposition is finished, return to the Maps software and click **OK** to resume using the Maps software.

After deposition, the sample is ready to perform lamella milling in a TEM.

Cryo menu items

When Maps software is in Cryo mode, new commands are available on the Microscope menu.





Menu item	Description
Set Sample Tilt	Unlinks the stage and stores the current sample tilt to the current tilt of the microscope. When the sample tilt is set, there is a check mark next to this option and the stage tilt that was set. See Set sample tilt for instructions.
Restore Last Sample Tilt	Restores the previous sample tilt that was set. This command is only visible when the sample tilt is not set.
GIS Deposition	Sets up the Maps software for GIS deposition. See Perform GIS deposition for instructions.

Cryo lamella site toolbar

The cryo lamella site toolbar provides acces to controls and actions relevant to the Cryo workflow. It is displayed only when the Cryo features are turned on as described in Configure Cryo settings.

	thermo	oscientífic		_ O X
+ w k	Q			* 🕷 🌂 🛍 o 💡 💈
Lamella sites	🗸 🔺 Lamella	~ ▶	≯ +⊭ [≯] +⊭ ⊕	

Toolbar item	Description
Lamella sites 🗸	Dropdown menu containing layer types that can be used in this toolbar.
▲ Lamella ✓ ▶	List of fully set up lamella sites. Changing the lamella selection changes the selected site in the layer tree. Use the arrow buttons on either side of the list to select the next or previous lamella site in the collection.
и к л к	Drives to the selected lamella site eucentric position. See Cryo lamella sites for more information about lamella sites.

Toolbar item	Description
[™] + ^ĸ	Drive to the selected lamella site mapping position. See Cryo lamella sites for more information about lamella sites.
•	Starts the process to calculate the eucentric position. See Calculate the eucentric position for instructions.

Cryo lamella timeline

The timeline is a collection of notes, files, and actions that have happened or been attached to a selected lamella site. By default, the timeline automatically tracks property changes as text notes. This control can help you identify what has happened during the setup of the selected lamella site.



Add a note to the timeline

You can add simple text notes to lamella sites. The text appears on the timeline.

- 1. Click the site to which you want to add a simple text note.
- 2. Click Add Note.



3. Type a text note in the dialog and then click OK.

Add a file to the timeline

You can attach files to lamella sites. The file is copied into the Maps project, so it is still available if you move the Maps project to another directory. In Cryo Settings, you can Choose a directory to watch.

- 1. Click the site to which you want to add a file.
- 2. Click Add File.



3. Browse to the file you want to attach, click it, and then click **Open**.

Cryo lamella sites

A *cryo lamella site* is a special annotation layer defined on a SEM tool to mark a specific location on a cryo sample where you want to mill a lamella. This layer contains special functionality and positioning. Each lamella site is used to mark cells of interest and to determine the angle at which to mill the lamella.



Lamella sites have the following properties in Maps software.

Lamella site status

The status symbol indicates the setup state of the lamella. This status symbol is displayed in the layer tree.

User interface item	Description
🗹 🔎 Lamella	The grey status symbol means the lamella site is in the initial state. It has been created, but nothing else has been done to it.
🖌 🄁 Lamella	The orange status symbol means the eucentric position has been calculated and stored for the lamella site.
🗹 🔂 Lamella	The green status symbol means the milling position and angle have been stored. The site is ready to be used for deposition and milling.

Lamella site properties

The following properties are displayed when you click on a lamella site.

Property	Description
Name Lamella	Name of the lamella site.
it.	Color of the lamella site. Click this to change the color of the lamella site shown in the viewer.

Lamella site positions

Positions listed in the Lamella Site property pane. A position has a yellow exclamation mark

when it is not set up and green check once it is set up. You must set up a lamella site before you can Perform GIS deposition.

Depending on the lamella site status, there are a different options per position. Notice the change in the buttons and the status indicators. A lamella site is considered fully setup when all

three positions have a green check end and them.

Initial Eucentric position calculated Setup complete Ø \bigotimes \oslash Lamella ð Name Lamella ð Lamella 1 Mapping Positio Mapping Position Drive To Update Drive To Updat ntric Positio Drive To Update ntric Position Drive To Updat Refine Eucentric Po Refine Euce Milling Angle Not Set Milling Position Store Angle Update Milling Position Drive To Milling Angle Failed Not Set Milling Angle 48.0°

Mapping Position

The *mapping position* is the XYZ position where the lamella was initially added. See Add lamella sites for instructions.

- Drive To Centers the Maps viewer on that position.
- Update Changes the mapping position to the current microscope field of view location.



Initial setup

Eucentric Position

The *eucentric position* is the location where the stage can be tilted through a range of angles and the cell of interest the lamella site is marking does not move in X, Y, or Z in the field of view. It is not initially set up and must be calculated manually. See Calculate the eucentric position for instructions.

- Calculate Click if the position has not been calculated.
- Drive To Centers the Maps viewer on that position.
- Update Changes the mapping position to the current microscope field of view location.
- Refine Eucentric Position Click to re-calculate the position.

Eucentric position calculated

Name	Lamella		ð
Mapping Position	Drive To	Update	Ø
Eucentric Position	Drive To	Ø	
	Refine Eucent	ric Position	
Milling Position	Store Angle	Update	()
Milling Angle I	Not Set		
Success		Failed	

Milling Position

The *milling position* is the stage tilt and position to use when milling the lamella. See Determine cryo milling position for instructions.

- Drive To Centers the Maps viewer on that position.
- Update Changes the milling position to the current microscope field of view location.

Setup complete

Name	Lamella		đ
Mapping Position	Drive To	Update	\bigotimes
Eucentric Position	Drive To Update		\bigotimes
	Refine Eucen	tric Position	
Milling Position	Drive To	Update	\oslash
Milling Angle	48.0°		
C		E-ilad	

Maps Min

Maps Min includes a workflow spanning multiple software applications to collect and analyze EDS and mineralogy data from geological samples. Software includes Thermo Scientific Maps, Thermo Scientific[™] Nanomin (called "Nanomin" within this user guide), and Bruker ESPRIT. Hardware includes the Support PC and the Microscope PC.



Get started

To get started in Maps Min, see Configure Mineralogy Acquisition.

Online support

Go to the Help menu and click **Open Help Desk** to access online support from within the Maps application. Alternatively, click https://thermofisher-

asg.atlassian.net/servicedesk/customer/portal/16 to gain access outside of Maps. Click Report Maps Min Bug for mineralogy issues.

Configure Mineralogy Acquisition

Before you proceed with the Maps Min workflow, you must first configure the Maps software for mineralogy acquisition on the Microscope PC with the Nanomin software and Bruker ESPRIT software on the Support PC.

Prerequisites:

- Set up the Support PC according to the installation instructions.
- There must be a shared folder on the Support PC that the Microscope PC can access.

Note: This task needs to be performed only once.

Procedure

On Support PC

• Note: Only one user can be logged in at any given time on the Support PC. If you need to switch users, then first log off from your user account before another user logs into the Support PC.

- 1. Confirm that the required services are running.
 - a. Thermo Scientific Nanomin Service

(E)	Windows Task Manager					
File	e Options View Help					
Ap	pplications Processes Services Perfo	ormance	Networking U	sers		
	Name	PID	Description	Status	Group	
	MSDTC MSiSCSI msiserver	2656	Distributed Microsoft i Windows I	Running Stopped Stopped	N/A netsvcs N/A	
	NanominService.Host	3520	Thermo Sci	Running	N/A	
	napagent		Network A	Stopped	NetworkSe	
	Netlogon		Netlogon	Stopped		

b. Thermo Scientific Bruker Service x64 and Thermo Scientific Bruker Service x86 (two blue icons appear in the task tray).



2. Perform the Energy Calibration procedure to ensure that the Bruker EDS detector is properly calibrated.

On Microscope PC

- 1. Launch the Maps application.
- 2. When prompted for a default Maps project directory, select a location that is local on the Microscope PC.
- 3. When the project history window appears, click New to create a new project.
- 4. Go to the Menu bar, and click **Options > Settings > Mineralogy**.
- 5. Go to the Bruker Detector Type dropdown menu, and then select the detector that is installed on your microscope system.

- 6. Go to the **Mineralogy section >> Data Export >> Final Data Directory** field, and click the ellipsis [...] to search for the shared folder location on the Support PC. Select this folder location to designate it as the location where the final mineralogy data will be saved to and processed from during the mineralogy workflow.
 - Note: At this time, you may also specify a Staging directory location. Go to the Mineralogy section >> Data Export >> Staging Directory field and enter a nonnetwork location on the Microscope PC. The staging directory location must NOT be a network location since Maps will store files in this location during acquisition to improve reliability of the transfer to the Final Data Directory.

Ma Settings		×
Application	- Image Settings	
XPS		
Mineralogy	 Please make sure that the Bruker Service has been installed and is running on the Bruker PC. You can find the installer on the install media Maps came on. 	
Velox	Bruker Detector Type: XFlash 6 Series 🗸	
Сгуо	Bruker Alignment Aspect Ratio 8:7 🗸	
Stitching	XRay Batch Size: 384	
Internal	Data Export Staging Directory: 2	
	D:\MineralogyStagingArea	
	Final Data Directory: 🕜 😃	
	Apply OK Canc	el

7. Click OK to save the settings.

8. Wait for a solid, green status icon to appear in the status bar. This status icon indicates that Maps is connected and ready for Bruker acquisitions.



9. Continue to Mineralogy Workflow now that the configuration is complete.

Energy Calibration

Correct calibration of the energy-channel (or gain) is a required for reliable qualitative and quantitative analysis results.

Prerequisites:

- The Bruker ESPRIT application must be installed.
- Dead time value is bellow 10%. If it is too high, decrease the spot size on the microscope.

Note: It is important to check the calibration on a weekly basis and adjust it if necessary to compensate for spectrometer drift.

Note: Calibration must always be performed after switching on the signal processing unit. This action is not required when using the standby mode, but it is highly recommended.

Procedure

Refer to the ESPRIT v2.2 user interface shown below for an overview of the procedure before you begin. The user interface slightly differs depending on the version of your ESPRIT software package.

1. Click EDS, and then click the Spectrometer tab to select the spectrometer workspace.

Note: Since system factor calibration is accessible through the calibration buttons in every workspace that includes spectra acquisition, do not confuse *energy calibration* with *system factor calibration*.

- 2. Move the calibration sample into the analysis position and set the beam current (spot size) to produce an intermediate count rate. Any element in the mid-energy range (that is, elements Mn-Zn) in a given sample can be used, provided that it is contained in a reasonable concentration. It is recommended that you use single element standards, such as pure copper, to avoid peak misidentification, increase speed, and improve accuracy.
- 3. Under the Settings to calibrate panel:
 - a. Pulse throughput: Select the check boxes of all the pulse throughput settings. The list of available speeds in this option is dependent on the license.
 - b. Energy range: Select **20kV** for the energy range to be calibrated for the spectrometer.
 - c. Detector segments: Select all detectors that are physically present on the system.
- 4. Under the periodic table, click the element of interest and then click K- α . It is recommended to use K- α lines whenever possible.
- 5. Under the Calibration panel, click **Precise** which is sufficient in most use cases, and then click **Start**.
- 6. When prompted to save the new calibration data, click **Yes**. Note that the storage of calibration data is not user specific, so changes in calibration affect all users of the spectrometer.

ESPRIT v2.2 software package



Mineralogy Workflow

The Maps Min workflow explains how to use of Maps in tandem with Nanomin to acquire EDS data from a sample. It also explains how to generate a Microsoft[®] Office Excel[®] report with the acquired data and how to publish the data to the web-based Maps Min Reporting application. You can use the Maps Min Reporting application to filter, combine, and view the data online as desired. You can also export reports as Excel files similar to the Nanomin application.

Note: The Maps Min workflow is only available on SEM systems and requires an active Mineralogy license. See Package options for information about licenses.

Prerequisites:

- The Maps software is configured for mineralogy acquisition on the Microscope PC (see Configure Mineralogy Acquisition).
- The Nanomin software is configured on the Support PC (see Configure Mineralogy Acquisition).
- The Maps Min Reporting application is installed on the Support PC.
- You have loaded your sample and aligned your holder in Maps.

Note: You can repeat this workflow for each mineralogy sample if needed.

Set up mineralogy tile set

- 1. Complete the procedure in Set Up Mineralogy Tile Set.
- 2. Repeat this task for as many tile sets as needed.
- 3. (Optional) Use the Mineralogy Measurement Template to quickly create multiple tile sets at once for repeated work.

Acquire mineralogy data

- 1. Start the Job Queue.
- 2. Wait for acquisition and processing to finish.

Open mineralogy data

- 1. Complete the procedure in Open Mineralogy Data.
- 2. Diagnose the data.

Export mineralogy report

- 1. Complete the procedure in Export Mineralogy Report.
- 2. Analyze data in report.

Publish mineralogy data

- 1. Complete the procedure in Publish Mineralogy Data.
- 2. Analyze data in Maps Min Reporting.

Set Up Mineralogy Tile Set

This procedure explains how to set up a tile set in Maps, which performs a mineralogy EDS acquisition with the Bruker ESPRIT software and detector.

Prerequisites:

- Maps must be configured for mineralogy acquisition (see Configure Mineralogy Acquisition).
- xT software must be installed and running on the Microscope PC.

Note: If you have Mineralogy tile set templates that you created with an earlier version of Maps 3.28, then you will need to recreate them.

Note: Repeat this procedure for every data sample of interest to capture EDS information. Refer to Mineralogy Acquisition Properties for more information about each acquisition parameter.

Procedure

- 1. Create a tile set with one of the toolbar controls (see Toolbar).
- 2. Select the new tile set.
- 3. Click the Mineralogy Acquisition tab, and then select the **EDS Acquisition** check box to designate the tile set for mineralogy acquisition.

⊞	.	[AF]	.,	1	¢۴	Ø	
Perform	n EDS Acq	uisition			 		

4. Click the ... to select the desired elements to acquire EDS data.



5. In the **Select Elements** window, click on the elements you want to select, and then click **OK**. If you do not want to view elements during acquisition, then only click **OK** and continue to the next step.

Ma Select Elements									×								
Select the elements you want to identify during acquisition.																	
н																	He
u	Be												С		0	F	
Na	Mg											AI	Si	Р	S	α	Ar
К	Ca	Sc	Ti	V	Cr	Mn	Fe	Co	Ni	Cu	Zn	Ga	Ge	As	Se	Br	Kr
Rb	Sr	Y	Zr	Nb	Mo	Тс	Ru	Rh	Pd	Ag	Cd	In	Sn	Sb	Te	I	Xe
Cs	Ba		Hf	Ta	w	Re	Os	lr	Pt	Au	Hg	Π	Pb	Bi	Po	At	
		La	Ce	Pr	Nd	Pm	Sm	Eu	Gd	ТЬ	Dy	Но	Er	Tm	Yb	Lu	
		Ac	Th	Pa	U	Np	Pu	Am	Cm		Cf	Es	Fm	Md		lr	
													ОК			Cance	ł

- 6. Complete the Define Background BSE procedure.
- 7. Click the **New** button next to the Segmentation Settings dropdown menu to open the Segmentation Wizard. Use this wizard to set up the segmentation settings for acquisition and post processing. The wizard also helps you to define the background BSE.



8. Optional: In the Recipe dropdown menu, select which recipe you want to process the tile set with post acquisition in Nanomin.



9. Optional: To stop acquisition after a specific number of acquired particles, select the **Particle Count Termination** check box and enter a positive value in the dedicated field.



10. Optional: To acquire only a sub-set of particles based on the size metric:

a. Click **Particle Size Metric**, and then select one of the available values from the dropdown menu.



b. Enter the minimum and/or maximum particle size value in the dedicated fields. If needed, you can deactivate each field by deselecting the associated check box.

Particle Size Metric	Long Axis 🗸 🗸
Minimum Particle Size 🛛 🧹	1 µm
Maximum Particle Size 🛛 🧹	100 µm

11. Optional: To perform Bright Phase Search acquisition, select the **Use BSE Search Ranges** check box and follow the **Optional:** Bright Phase Search Acquisition Workflow procedure.

Use	BSE Search Ranges	2	
1	Note: When you activa sends to Nanomin for p	te Use BSE Search Ranges, the uncalibrated in	ne BSE image that Maps nage.

- 12. Select an Acquisition mode:
 - Grid: This mode performs a simple serial scan to collect X-rays.

Acquisition Mode	Grid	\mathbf{v}
EDS Dwell Time	8 ms	
XRay Spacing	10 µm	

After you select **Grid** mode from the dropdown menu, enter the **X-Ray Spacing** value to determine the distance between each location to collect X-rays. This is essentially the pixel size for the resulting EDS image.

• Centroid (beta): This mode performs some image processing on the BSE electron image to detect grains and attempts to perform a single X-ray scan on each grain to improve acquisition performance.

Acquisition Mode	Centroid (beta)	~
Process Edge Particles		
Grain Size	3 µm	

After you select **Centroid (beta)** mode from the dropdown menu, enter a **Grain Size** value to determine the grains of interest. Maps ignores any grains that are smaller than the Grain Size value; therefore, increasing the Grain Size value improves acquisition time, but may result in missing measurements on grains. The Process Edge Particles check box is deselected by default since there are known issues with edge particles while in Centroid mode. If you keep the default, then the particles that fall on acquisition tile edges are ignored during Nanomin processing.

Caution: The Centroid mode is still in beta and may have varying results. It is recommended that you only use a resolution of 512x442 for tile sets.

13. Enter the **EDS Dwell** time. This is the time spent collecting X-rays per EDS pixel. The default is 8 ms. Longer dwell times result in more acquired X-rays and higher quality results.



- 14. Optional: Fill out the Tags fields based on the sample associated with the tile set. These tags are used for sample tracking and generating reports. For more information on what the tags represent, see Mineralogy Acquisition Properties.
- 15. Optional: Repeat this procedure to create additional tile sets if needed.
- 16. Return to the Mineralogy Workflow to continue.

Optional: Bright Phase Search Acquisition Workflow

The Bright Phase Search Acquisition workflow obtains mining data for particles containing bright phase, which results in faster acquisition and results within the web-based Maps Min Reporting application.

Workflow overview

The Bright Phase workflow utilizes user-defined standards to calibrate the incoming BSE images. The workflow then compares the BSE pixel values against another set of userdefined mineral search ranges to acquire EDS on particles of interest at a different resolution and dwell time. The goal of this workflow is to minimize data size and reduce acquisition time when only a subset of minerals is of interest.



Prerequisites:

- The Maps software is configured for mineralogy acquisition on the Microscope PC (see Configure Mineralogy Acquisition).
- The Nanomin software is configured on the Support PC (see Configure Mineralogy Acquisition).
- You have loaded your sample and aligned your holder in Maps.
- The xT software is installed and running on the Microscope PC.

Procedure

a

1. Set up standard sample locations with the Mineralogy Standards Workflow. You are required to have standard sample locations set up to continue with this procedure.

Note: To ensure that the data are accurate, these standard sample locations must be from a sample holder that is currently loaded.

2. Begin the Set Up Mineralogy Tile Set procedure. Before you select an acquisition mode, select the Use BSE Search Ranges check box to opt for Bright Phase Search acquisition.

3. Click the Mineralogy Acquisition tab and set the acquisition settings for **Target** and **Non-Target**. The Target settings are acquisition scan settings for particles that contain pixels within defined BSE search ranges, and the Non-Target settings are acquisition scan settings for all remaining particles.

Note: Skipping Non-Target particles is not currently supported. As a workaround, you can lower the EDS Dwell Time to 1 ms to speed up acquisition of Non-Target particles.

4. Click Add Search Range to add a new BSE search range.



5. When the **Add New Search Range** window appears. Click one of the target phases in the dropdown menu (or type a target phase in the Search box), and then click **Add**.

👪 Add New Search Range	×
Select a target phase to create a BSE search range for	
Q	
Abelsonite	•
Abenaküte-(Ce)	
Abhurite	
Abswurmbachite	
Acanthite	
Acetamide	
Achavalite	
Actinolite	
Adamite	
Adelite	
Aeoirine-auoite	
Selected: Add Cance	

6. Your selected target phase will be added under **Search Ranges**. It will be populated with a search range based on the phase's average atomic weight. You can change the upper and lower threshold values of the search range by adjusting the left and right sliders bordering the histogram. Alternatively, you can click the up and down arrows of the two text boxes to change the upper and lower threshold values. The upper and lower threshold values range from zero to one hundred percent (that is, 0%-100%).



7. (Optional): Repeat this procedure to create an additional search range if needed.

8. Return to Set Up Mineralogy Tile Set to select an acquisition mode and continue setting up the mineralogy tile set.

Optional: validate procedure

Once tile set acquisition is complete, validate that the particles were successfully acquired with the Target settings.

1. Select the stitched result layer, and click the Visualization tab.



2. Deselect all channels except the calibrated channel. As shown in the example below, the calibrated channel displays "(Calibrated)"at the end of the channel name.

□ ‡ Ø	
Opacity	•
Additive Opacity	•
	L Show histogram
•	
>	•
> 🗆 si	•
> 🗖 Ag	•
> 🗷 BSD; None (Calibrated)	ê •
> BSD; None	•

3. Expand the Calibrated channel, and then click Add Color.



4. Review the BSE search range setup in the Bright Phase procedure above, and convert the upper and lower threshold values into 16-bit values (that is, pixels). For example, the 84%-94% range for gold would be converted to 55,049 and 61,603 pixels, respectively.

5. In the solid color pane, move the upper and lower threshold values to match the converted values that you calculated in the previous step. These updated threshold values define your BSE search range.



6. View the acquired image to locate the red pixels, which represent your BSE search range. This means that if a particle contains red pixels, then the particle was acquired with the Target scan settings.



Note: If none of the particles in the acquired image contain red pixels, then it is recommended that you acquire another tile set with a different BSE search range values.

Mineralogy Standards Workflow

The Mineralogy Standards workflow explains how to set up standard sample positions used within the Optional: Bright Phase Search Acquisition Workflow.

1. Click the Start Workflow icon located on the right side of the Toolbar.



2. In the Select Workflow window, click Mineralogy Standards Setup, and then click Begin.

Ma Select Workflow	×				
Please select a workflow					
Alignment Wizard					
Global Alignment Wizard					
Global Focal Reference Plane					
Debug Layout Test Workflow					
Mineralogy Standards Setup					
Array Tomography					
	•				
	^				
Begin Car	ncel				

3. The **Workflow: Mineralogy Standards Setup** window appears on the right side of the main screen.

Wor	kflow: Mineral	ogy Standards Setup	×
1 Using XT, driv controls to d position. Repe the workflow	ve to a standard on esignate which sta eat this for 2 more s to save the standard	the sample holder. Use the l andard it is and to save its standards then click the check ds positions.	below stage box in
Standard 1			~
Drive To	Location	Save Current Stage Locatio	on
Standard 2			~
Drive To	Location	Save Current Stage Locatio	on
Standard 3			~
Drive To	Location	Save Current Stage Locatio	n
2 Ensure all 3 s the check ma workflow.	standard locations and standard locations and states the states of the s	are saved and valid before cli ottom right to complete the	icking setup
)	
	Setup Standa	rd Locations	

- 4. Use the xT interface to drive the stage to a standard location on the sample holder.
- 5. Select the standard location from the **Standard 1** dropdown menu.
- 6. Click Save Current Stage Position.
- 7. Repeat steps 4-6 to set up Standard 2 and Standard 3.
- 8. Click the check mark icon located in the lower right corner of the workflow window to save the standard positions.

Segmentation Wizard

This procedure explains how to use the Segmentation Wizard, which you use to set the segmentation settings. Maps uses these settings during acquisition and classification to find particles and grains.

Prerequisites:

- Maps must be configured for mineralogy acquisition (see Configure Mineralogy Acquisition).
- xT software must be installed and running on the Microscope PC.

Note: Repeat this procedure for every data sample of interest to capture EDS information. Refer to Mineralogy Acquisition Properties for more information about each acquisition parameter.

Procedure

- 1. Create a tile set with one of the toolbar controls (see Toolbar).
- 2. Select the new tile set.
- 3. Click the Mineralogy Acquisition tab, and then select the **EDS Acquisition** check box to designate the tile set for mineralogy acquisition.



4. Click the **New** button next to the Segmentation Settings dropdown menu to start the Segmentation Wizard. The Segmentation Wizard automatically starts the workflow, which includes defining the BSE background threshold, setting up the particle segmentation, setting up the grain segmentation, and saving your new segmentation settings.

Segmentation Settings	Default	✓ New
-----------------------	---------	-------

- 5. Define the BSE background threshold. This step determines the areas of the image that will not have any X-rays collected. Once a desired region is excluded from X-ray collection, move to the next step in the workflow.
 - a. Click Drive To Tile Set to move the stage to the center of the current tile set.
 - b. Click **Acquire Snapshot** to acquire an image. When prompted to apply tile set settings to the microscope, click **Yes**. This step is required to ensure that the image captured mimics the image that will be acquired by the tile set.





c. Once Maps acquires the image, the image appears in the main panel of the workflow and a threshold histogram appears in the right panel.

d. Adjust the slider control for the threshold histogram to designate the areas of the image to be excluded from the collected X-rays. Continue adjusting the slider control until all the designated areas are highlighted in red.



e. Once a desired region is excluded from X-ray collection, move to the next step in the workflow.
- 6. Set up the particle segmentation. In this step, you to test the particle segmentation with the BSE image taken in the previous step to ensure that the segmentation settings perform as desired.
 - a. Click the Strength slider control to determine how aggressive you want the particle splitting. When you increase the strength, the number of found particles increases.
 - b. Click **Run Particle Splitter** to test the particle segmentation. Found particles appear in color, so you can identify them. Click **Show/Hide Particles** to toggle the visibility of the particles on and off so that you can compare them with the BSE image.



c. When you are satisfied with your particle results, move to the next step in the workflow.

- 7. Set up the grain segmentation. In this step, you test the grain segmentation results from within the found particles of the particle segmentation. Note that this step is only available when the acquisition mode is set to **Centroid (beta)**.
 - a. Adjust the Grain size slider control to determine how granular you want the segmentation. Each grain appears in color, so you can identify them. Click **Show/Hide Particles** to toggle the visibility of the grains on and off so that you can compare them with the BSE image.



b. When satisfied with the grain results, move to the next step in the workflow.



8. Click **Save** to save your new segmentation settings, so that you can reuse them with other tile sets. This step completes the segmentation workflow.

Mineralogy Measurement Template

The Maps Min software enables you to quickly set up multiple mineralogy tile sets at once with a customized CSV template file specific to your use case. Although a customized CSV template file is optional, it is recommended that you create one and import it into Maps.

A CSV file is a comma-separated values text file, which allows you to save data in a tabular format. Each line of data in the CSV file is a record, and each record consists of one or more fields separated by commas. In this instance, Microsoft Office Excel automatically opens the provided CVS file template and organizes the data. Each row in the Excel spreadsheet represents one mineralogy tile set record, and the columns contain the tile set properties associated with each mineralogy tile set record. Note that each mineralogy tile set record contains 2-10 tile set properties.

Create custom CSV template file

- 1. Navigate to the Support Files folder in the Mineralogy Installer.
- 2. Locate the file entitled *MineralogyMeasurementTemplate.csv* and copy/paste it to the desktop.
- 3. Double-click the desktop file copy to open the file as a standard Excel spreadsheet.
- 4. To create the first tile set record, refer to the table below and do the following in the first row of the spreadsheet.
 - a. Enter a value into the Hole Position column.
 - b. Enter a value into the Template column.
 - c. (Optional): Enter values into the other columns as desired.
- 5. To create more tile set records, repeat steps 4a-4c for each additional row.
- 6. Click File, select Save As, rename the file, and then click Save.
- 7. Continue to Import custom CSV template file.

Column Header	Data Type	Required Field	Notes
Hole Position	Integer	Yes	This field represents which hole in the holder that you designate for the tile set creation (for example, 1 = hole 1, 2 = hole 2, 3 = hole 3, etc.).
Mineralogy Project Name	String	No	Use this field to designate a parent directory name for the acquisition results. This can help organize your sample data on disk.
Replicate	String		
Sample Code	String	No	Use this field to name the created tile set.
Sample Description	String	No	
Sample Name	String	No	
Sample Weight	Double	No	
Size Fraction	String	No	
Size Fraction Weight %	Double	No	
Survey Name	String	No	
Survey Description	String	No	
Template	String	Yes	This field represents the name of the <i>tile</i> <i>set template</i> within Maps. The tile set template must exist prior to importing your custom CVS template file into Maps.

Import custom CSV template file

After you have created your custom CSV template file, you can then import it into the Maps.

Procedure

- 1. Confirm that a project is open and that Maps has a Bruker connection.
 - a. View the Maps status bar located at the bottom of the main screen. A green status icon next to the Bruker label indicates that Maps is connected to the Bruker label, and a green status icon next to the Nanomin label indicates that maps is connected to the Nanomin Service (refer to the Main Screen for more information on the microscope connection states).



- b. If the green status icon is not present, then go to On Microscope PC and configure the Maps software for mineralogy acquisition on the Microscope PC with Nanomin.
- 2. Click File, and then select Import Mineralogy Measurement Template.



- 3. When Maps prompts you, select your custom CSV template file from the desktop to import it into Maps. Maps uses this template to create a tile set for each mineralogy tile set record in the file.
- 4. Return to the Mineralogy Workflow to continue.

Open Mineralogy Data

This procedure explains how to open and view acquired mineralogy data in Nanomin.

Prerequisites:

- Nanomin must be installed and licensed.
- You must have access to a Maps project that has acquired mineralogy data.

Note: This task can be performed anytime and repeated as needed.

Procedure

- 1. Launch the Nanomin application.
- 2. Click **File** from the toolbar, and then click **Open Maps Tile Set** from the dropdown menu to open the directory dialog.



3. Locate the Maps project file that contains the mineralogy tile set to open. If there is more than one mineralogy tile set within the selected Maps project, then Nanomin prompts you to select a mineralogy tile set to open.

- 4. Double-click the desired mineralogy tile set to open it.
 - If Nanomin has not yet classified the mineralogy tile set, then Nanomin prompts you to select a recipe to process the tile set. Select a recipe to continue.
 - If Nanomin has classified the mineralogy tile set, then classification automatically runs.
- 5. The mineralogy tile set opens and appears in the viewer. You can then view the mineral and elemental layers for the tile set.
- 6. Continue to Export Mineralogy Report.

Export Mineralogy Report

Nanomin can generate a mineralogy report containing a list of calculated results from one or more Maps Min tile sets. Follow the steps below to export a mineralogy report that can be read in Microsoft Office Excel.

Prerequisites:

- The Nanomin application must be installed.
- You must have a Maps Min tile set acquired and classified by Nanomin.

Procedure

- 1. Launch the Nanomin application.
- 2. Click **Data** from the toolbar, and then click **Export Reports to Excel** from the dropdown menu to open the directory dialog.



3. Locate the Maps project that contains the tile sets to be included in your report and then double-click it.

Include	☑ Thumbnail	Measurement Details	Metadata	SolidSolution	
		Tile Set (3) Mineralogy/old/2019-10-19 Garth Particle Scan\Garth Particle Scan_3.11 - Copy (Recipe: All Minerals Classified: 10/18/2019 6:07:23 AM Measured: 10/18/2019 4:05:43 AM	Survey: Sample: Size Fraction: Replicate:		
-		Tile Set (2) Mineralogy/old/full centroid 5um sample Recipe: Everything Classified: 10/25/2019 12:18:58 PM Measured: 10/24/2019 1:48:17 AM	Survey: Sample: Size Fraction: Replicate:		
Z		Tile Set (2) Mineralogy/old\new centrold Recipe: Everything Classified: 10/21/2019 6:19:46 PM Measured: 10/22/2019 2:18:53 AM	Survey: Sample: Size Fraction: Replicate:		

4. Click the check boxes of the tile sets to designate their data for export.

- 5. Click Export.
- 6. When prompted, enter a file name, select a file location, and then click **Save**.
- 7. The report automatically opens after the file is saved. See Mineralogy Report Data for more information.

Publish Mineralogy Data

Nanomin can publish one or more Maps Min tile sets data to use with the web-based Maps Min Reporting application. Follow the steps below to publish data to the Maps Min Reporting server.

Prerequisites:

- The Nanomin application must be installed.
- You must have at least a Maps Min tile set acquired and classified by Nanomin.

Procedure

- 1. Launch the Nanomin application.
- 2. Click **Data** from the toolbar, and then click **Publish Multiple Measurements** from the dropdown menu to open the directory dialog.



- 3. When prompted, enter the URL address of the Maps Min Reporting server and then click **Next**.
- 4. When the Reporting Server Selection screens opens, enter the reporting PC name.

Reporting Server Selection
Please enter the name of the PC hosting the reporting software package.
Reporting PC Name reporting-spc
Connection validated 💉
Next Cancel

5. Locate the Maps project that contains the tile sets that you want to publish to the reporting server, and then double-click the project.

6. Select the checkboxes in the Publish column to designate the desired tile set data for publishing, and then click **Publish**.

		Publish			
Directory D:\	Mineralogy projects\Project Test				
Preview	Dataset	Metadata	Acquisition Date	Last Modified Date	Publish
	Test 201 Recipe: Copper Sulfides Measurement Mode: Grid	Survey: Survey A Sample: Sample A Size Fraction: +25 Replicate: Replicate A Description: Tile Set	6/13/2023 4:49 PM	6/14/2023 10:46 AM	
	Tile Set Recipe: Bauxite Measurement Mode: Grid	Survey: Survey A Sample: Sample A Size Fraction: +25 Replicate: Replicate B Description: Tile Set	7/27/2023 10:44 AM	8/9/2023 1:44 PM	
				Publish	Cancel

7. Verify and edit the selected dataset's information, and then click **Publish**.

				Publish				
Dataset	Project	Survey	Sample	Sample Weight	Size Fraction	Size Fraction Weight %	Replicate	Status
Tile Set	Test 1 ~	Survey A 🗸	Sample A 🗸 🗸	100	+25 ~	87	Replicate A 🗸 🗸	
Tile Set	Test 1 v	Survey A 🗸	Sample A 🗸 🗸	100	+25 ~	87	Replicate B 🗸 🗸	
							Publish	Cancel

8. The data are automatically published to the Maps Min Reporting server.

SEM/SDB

References

The Maps Min workflow includes supporting information as listed below.

Topics include:

Define Background BSE Mineralogy Acquisition Properties Mineralogy Report Data

Define Background BSE

The Background BSE is a percentage that represents the darkest percent of pixels in the BSE image. Properly defining the Background BSE ensures that X-rays in any of these dark pixel locations are excluded from the acquisition, which optimizes EDS acquisition time.

Prerequisites:

- Maps Min must be configured (see Configure Mineralogy Acquisition).
- A Maps tile set must be created and selected (see Set Up Mineralogy Tile Set).
- The Mineralogy Acquisition tab must be selected in the tile set's property pane and the EDS Acquisition check box must be selected.

Procedure

1. Click the gear icon next to the Background BSE parameter.

Background BSE	9.41%	۲
----------------	-------	---

- 2. Follow the beginning of the Segmentation Wizard procedure to define the background BSE.
- 3. Once you define the background BSE and move to the next step in the workflow, Maps closes the workflow and updates the Background BSE in the setup tab.

Mineralogy Acquisition Properties

There is a collection of acquisition properties for you to complete within Mineralogy Acquisition >> Mining Tag for each mineralogy tile set. Refer to the three tables below for property names and descriptions.

Property	Description
Acquisition Mode	 Determines the image processing mode used for collecting X-rays. Choices include: Grid: This mode performs a simple serial scan to collect X-rays. When this mode is selected, you must provide an X-Ray Spacing value. This value is the distance between X-ray scan locations. Centroid (beta): This mode performs some image processing on the BSE electron image to detect grains and attempts to perform a single X-ray scan on each grain to improve acquisition performance. When selecting this mode, you must provide a Grain Size value to determine the minimum grain size when performing the image processing. Note: Centroid acquisition may have some known issues
Background BSE	Percentage that represents the darkest pixels in the BSE image. Maps software does not acquire X-rays in any of these dark pixel locations. The reason for this functionality is to optimize acquisition time by not measuring resin or dark materials that are not of interest. See Define Background BSE for details.
EDS Acquisition	Determines if the tile set performs an EDS acquisition. Select the check box to activate.

 Table 1 - Acquisition Parameters

Property	Description
EDS Dwell	Time spent collecting X-rays per EDS pixel. The default is 8 ms. Longer dwell times result in more X-rays being acquired and higher quality results.
Elements	The selection of elements displayed in Maps. Each selected element is its own channel in the tile set result.

Table 2 - Mining Tags

Property	Description
Mineralogy Project Name	This is a label that represents the name of the project containing a collection of mineralogy surveys. The mineralogy project name typically includes the customer name, date, and/or business case.
Survey Name	This is a collection of related sample measurements, which normally represents a set of samples taken from a plant or ore body at a particular time.
Sample Name	This represents a single instance of sample material being taken from a particular location. For example, the concentration stream of a flotation cell. This entity represents the raw and unsized material. This may optionally be further subdivided into multiple size fraction ranges if you have mechanically separated this material for analysis.
Sample Weight (kg)	This represents the weight of a single instance of sample material being taken from a particular location. The default value is 100 kg.
Size Fraction (µm)	This is a label that represents the size range of particles for this sample, assuming that the original sample has been sorted into size ranges. For samples that are not sorted, leave this field blank (null in the database) indicating that the sample is not sized.

Property	Description
Size Fraction Weight %	If the original sample is split into size fractions, this field indicates the proportion of total sample mass that exists within the current size range. This is a percentage, because the sum of all size fractions equals 100%. If they do not equal 100%, then you can conclude that there are samples missing or a data entry mistake has occurred.
Replicate	This represents the facility for making duplicate physical epoxy pucks, or thin sections for analysis, that share an identical size fraction, survey, and sample. This is typically done when searching for rare minerals to increase the reliability of statistics derived from them. For example, since platinum minerals are rare (parts per billion) you might only see 1-5 grains per physical block. Companies typically create 10,20,30 replicates from the original source material to increase the number of grains they find, so that they have decent aggregate data such as grain size histograms.

Table 3 - Post Processing

Property	Description
Recipe	Name of the Nanomin processing recipe classified with the tile set after acquisition. If Nanomin has not yet classified the mineralogy tile set, then Nanomin prompts you to select a recipe to process the tile set. If Nanomin has classified the mineralogy tile set, then classification automatically runs.

Mineralogy Report Data

The mineralogy report contains a collection of calculated results for the measured sample. The table below contains descriptions of the information that is stored in each tab of the Microsoft Office Excel report. Refer to this table when analyzing the data in your report.

Tab	Description
Deportment	This shows an inversion calculation of the elemental assay and modal mineralogy weight data. For a specified element (for example, Cu) in the Elemental Assay report, it shows what minerals are contributing to the weight of that selected mineral. It can be used in tailings analysis to determine whether a valuable element is worth recovering, or whether the mineral contributing to that signal is too difficult to process.
Deportment Pivot Table	This shows the elemental deportment in a pivot table where you can easily select an element of interest (and one or more samples) to see a report of that element alone.
Elemental Assay	Shows a bulk computed assay derived from the minerals identified in the Modal Mineralogy Weight report (including for region-of-interest). Minerals are converted into their equivalent chemistry from their theoretical textbook chemistry. However, if a specific mineral has a large number of pixels in the image, the chemistry is not taken from the textbook definition but instead from an EDS elemental quantification of the spectrum obtained for that mineral. Data here can be compared to other bulk analytical techniques, such as XRF, LIBS, wet chemistry, etc.
Grain Size Distribution	This records the grain size distribution for individual minerals found in the samples exported. They are recorded in a histogram.
Grain Size Distribution Data	This records the grain size distribution for all mineral grains across all measurements. You should use the filters at the top of the spreadsheet to select a mineral of interest.

Tab	Description
Liberation Histogram	This shows a liberation report for all minerals and samples found. Liberation is computed as the total area of a mineral within each particle, and the report shows a summary view for all particles in the sample.
Locking Details	This is a detailed breakdown of the locking summary report. For each target mineral, it shows the percentage of the other minerals found in those particles as their weight percentage in binary or ternary classes.
Locking Summary	This is a refinement of the liberation report in which the percentage of minerals that are liberated or not liberated are shown. For particles that have two minerals, they are shown in the binary column. Particles with three minerals are shown in the ternary column. Particles with four or more minerals are shown in the complex column.
Modal Mineralogy Area	Shows the bulk mineralogy by area for the entire sample for one or more measurements. It also shows the mineralogy for all regions-of-interest defined in the exported measurements. Data here can be compared to other AM systems from other vendors.
Modal Mineralogy Weight	Based on the Modal Mineralogy Area, but shows the bulk results by weight instead of by area. Thus, heavy minerals (for example, pyrite or gold) have a higher effect than light minerals (for example, quartz or carbon). Data here can be compared to other bulk analytical techniques such as XRD, FTIR, RAMAN, or other AM systems.
Particle Size Distribution 1	This records a non-cumulative particle size distribution as a histogram across all measurements.
Particle Size Distribution Data	This records the cumulative particle size distribution for all particles across measurements.

Maps XPS

Maps XPS allows you to view XPS data and prepare sites, lines, and areas of interest for post processing within the Avantage software. To get started, follow the procedure below:

Get started

- 1. Go to the **Options** menu.
- 2. Click Settings, and then click the XPS tab.

Ma Settings		×
Application	Enable XPS	
XPS	Select this option to enable these Avantage XPS data features:	
Mineralogy	- Toolbar button for XPS data import - Toolbar button to create line of interest for line scan	
Velox	- Export to Avantage	
Сгуо	Use Theme Colors for Charts	
Stitching		
Internal		
	Apply OK	Cancel

- 3. Click **Enable XPS** to activate the XPS features. The Maps application alerts you if a system restart is required to proceed.
- 4. Click OK.
- 5. Proceed to the XPS Workflow.

XPS Workflow

In the Maps XPS workflow, you can view imported XPS data and create annotations for post processing with the Avantage software. During post processing, the Avantage software uses the exported sites of interest for point measurements, converts the lines of interest to line scans, and converts the areas of interest to snap maps.

Prerequisites:

- You must enable XPS features within the Maps application settings (see Maps XPS).
- You must have a file containing XPS data that was prepared by the Avantage software (version 6.1 or newer).
- You must have an External Image Import license (see Package options).

Workflow procedure

- 1. Import XPS Data that are prepared in Avantage software.
- 2. Initialize XPS Holder (optional).
- 3. View XPS Layer Types within the Maps application.
- 4. Create and acquire tile sets (or snapshots).
- 5. Create Annotations based on acquired data from the microscope.
- 6. Export XPS to Avantage new annotations to an XML file.
- 7. Import the XML file to the Avantage software for post processing.

Import XPS Data

When the Maps XPS features are enabled, the Maps application adds an icon for importing XPS data to the SEM/SDB toolbar. If the XPS features are not enabled, then the icon does not display within the SEM/SDB toolbar.

- 1. Go to the right side of the SEM/SDB toolbar.
- 2. Click the Import XPS Data icon.



- 3. Select a file to import (file type *.tfs.xml).
- 4. Click **Open** to start the data import.
- 5. Proceed to View XPS Layer Types.

Initialize XPS Holder

If the imported XPS data contain information about the XPS holder, then the XPS holder initialization workflow starts automatically at the end of XPS data import. If desired, you can abort the initialization workflow anytime after it starts (see Abort holder initialization).

1. Maps locates the imported images that contain the fiducials and then populates them in the right panel.



- 2. Maps processes the fiducials and places reference points for each one. If Maps processes at least three fiducials, then it prompts you to check the fiducial positions.
 - If the status of each fiducial within the right panel indicates success (that is, indicated by a green status icon), then check for false positives by confirming that the blue crosshair locations in the fiducial and small reference screens match. If the crosshair locations match for each successful fiducial, then click Yes to initialize the holder and go to step 5 to initiate the automated holder alignment. If the crosshair locations do *not* match for each successful fiducial, then click No and continue the remaining steps to perform a manual placement for the false positives.

placement		
Wo	rkflow: Holder initializatio	n X
Find locations of t located to initializ	the holder fiducials. At least 3 fiduc te the holder.	ials must be
Bottom right Reference:		
Upper left		
Bottom left		-6-
💽 Upper right		
Bottom right		
	Find fiducials	
	Initialize holder	
	Abort	
	•	
	Fiducial locations	~

• If the status of each fiducial within the right panel indicates failure (that is, indicated by a red status icon), then click **No** and continue the remaining steps to perform a manual placement for the failed fiducials.

Wo	rkflow: Holder initialization	×
Find locations of I located to initializ	the holder fiducials. At least 3 fiducials must l re the holder.	he
Bottom right Reference:		
Upper left		
Bottom left		
Upper right		
Bottom right		
	Find fiducials	
	Initialize holder	
	Abort	
	•	
	Fiducial locations	~

- 3. Go through failed fiducials in the right panel and manually set the correct fiducial positions by double-clicking within the fiducial screen to move the blue crosshair so that it is in the same location as shown in the small reference screen. The small reference screen is located to the left of the fiducial screen. If needed, you can use the mouse wheel to drag and zoom to precisely set the fiducial position.
 - Note: Sometimes the Maps algorithm produces false-positives and places fiducial positions incorrectly (that is, the blue crosshair locations in the fiducial and reference screens do not match). If this occurs, you can manually change the fiducial positions defined by the automated workflow.

SEM/SDB



4. After you properly set at least three fiducial positions, click **Initialize holder** at the bottom of the panel.



5. After holder initialization completes, click Yes to start the automated holder alignment.



Alternatively, you can later start the holder alignment from the main toolbar by clicking the **Align to Holder** icon. Refer to the **Automated Holder** Alignment section for more information.



Abort holder initialization

1. Click Abort at the bottom of the right panel.



2. When the application prompts you, click Yes to confirm.



3. If you abort the holder initialization and would like to start the workflow again, navigate to the main menu and click **XPS** >> **Initialize Holder**. The Initialize Holder submenu is only visible if you have XPS data with an XPS holder that has not yet been initialized. Once you initialize the holder, this submenu disappears.

Ma	Maps 3.1	8				
File	View	Microscope	Options	Help	XPS	
		li li	nitialize Hold	er		

View XPS Layer Types

After importing the XPS data into the Maps application, the Maps project contains the following new layer types for you to view.

- Measurement point
- Line scan
- Snap map



Measurement point

Maps displays the measurement point as an elliptical area. This layer type contains multiple spectra, and you can select which spectrum to display in the left panel.



At the bottom of left panel, there are two options:

• Open Peak Table: Select this option to open the peak table (if present) in a separate window.



• Open VGP File: Select this option to open the VGP file associated with the measurement point in the Avantage software. You must have the Avantage software installed to use this feature.



Line scan

Maps displays the line scan as thin rectangular area. This layer type contains multiple spectra, and you can select which spectrum to display in the left panel.



SEM/SDB

Additionally, you can select a profile as shown below to display.



Snap map tab

The Snap Map tab contains the default properties.

Name Snap Map Pixels 100 px x 100 px Pixel Size 30 µm x 30 µm Physical Size 3 mm x 3 mm Show metadata	□ ⋧ Ø	
Pixels 100 px x 100 px Pixel Size 30 µm x 30 µm Physical Size 3 mm x 3 mm Show metadata Technique Snap Map	Name	Snap Map
Pixel Size 30 µm x 30 µm Physical Size 3 mm x 3 mm Show metadata Technique Snap Map	Pixels	100 px x 100 px
Physical Size 3 mm x 3 mm Show metadata Technique Snap Map	Pixel Size	30 µm x 30 µm
Show metadata Technique Snap Map	Physical Size	3 mm x 3 mm
Technique Snap Map		
	Technique Snap Map	
Open VGP File	Ор	en VGP File

Property	Description
Name	Displays Snap Map name.
Pixels	Size of the image in pixels.
Pixel Size	Physical size of one pixel.
Physical Size	Physical size of the whole image.
Show metadata	This property is deactivated and not available since Snap Map does not include metadata.
Technique	Technique used to create the data in XPS. For example, Snap Map is the technique in this case.

Visualization tab

Maps imports the Snap Map as a multichannel image layer.



Create Annotations

With the XPS data imported into the Maps application, you can create new annotations for your lines of interest, areas of interest, and sites of interest.

Toolbar



Create line of interest annotation

- 1. Go to the SEM/SDB toolbar. The Line of Interest icon only appears in the toolbar when the Maps XPS features are enabled.
- 2. The Line of Interest icon is located in the middle of the toolbar (see image above). Click the Line of Interest icon.
- 3. Click and drag your mouse cursor in the Maps Viewer to draw a line.
- 4. Enter a name for the annotation in the dialog that opens, and then click OK.
- 5. The Line of interest tab opens in left panel. Update the Line of Interest properties under this tab.
- 6. (Optional) Add notes if needed.
- 7. Click the color picker to designate the annotation color.
- 8. Set the fill opacity (see details in the table below).

/ / Ø	
Name	Line
Notes	This is line of interest
Color	Ċ.
Fill Opacity	•

Property	Description
Name	Displays annotation name.
Notes	Displays annotation notes. Click in this field to add or edit these notes.
Color	Displays color of annotation. Click the color picker (that is, the eyedropper icon) to change the color.

Property	Description
Fill Opacity	Sets the level of opaqueness for the entire annotation. Click and move the slider to the right to increase the level of opaqueness, or click and move the slider to the left to decrease the level of opaqueness. The less opaque, the more transparent the annotation becomes. Transparent annotations are useful when overlaying a layer over another layer.

Create area of interest annotation

- 1. Go to the SEM/SDB toolbar. The Area of Interest icon is present in the toolbar even if the Maps XPS features are not enabled.
- 2. The Area of Interest icon is located in the middle of the toolbar (see image above). Click the Area of Interest icon.
- 3. Click and drag your mouse cursor within the Maps Viewer to draw a rectangular area.
- 4. Update the properties as defined and described in the Annotation Tab.

Create site of interest annotation

- 1. Right-click in the Viewer to access the Viewer Right-Click Menu.
- 2. Select Add Site of interest.
- 3. Update the properties as defined and described in the Viewer.

Export XPS to Avantage

After you have created create new annotations for your sites of interest, areas of interest, and lines of interest, you can export them to an XML file and then later import the XML file into the Avantage software for post processing.

Export procedure

1. Go to the right side of the SEM/SDB toolbar.

Note: Alternatively, you can start the export with the Main screen method and then skip to step 5.
2. Click the Workflow icon.



3. Select Export to Avantage from the menu.



4. Click Begin.

5. A single-step workflow panel opens on the right side of the export screen. This panel contains all annotations that you have created in the project.

Workflow: Export to Avantage				
	Site of interest	Point	~	
	Site of interest (2)	Point	~	
	Line	Line	~	
	Site of interest (3)	Point	~	
	Site of interest (4)	Point	~	
	Line (2)	Line	~	
	Area of interest	Area	~	
		Area		
		Line		

Here you can do the following:

- All of your annotations are preselected, but you can deselect any of them if desired.
- You can double-click an annotation to center and zoom Maps on it.
- You can also select how these annotations will be exported.

Note: Lines are exported *as lines*, sites are exported *as points*, and areas are exported *as either areas or lines*. If you export an area as a line, then the area is collapsed into a central line from within the area.

6. When the File Save dialog opens, select a file system location, enter the name of the exported file, and then click **Save**.

Main screen method

- 1. Go to the File menu.
- 2. Select Export to Avantage from the menu.



Rotational Alignment

This topic gives instructions for using the Rotational Alignment feature.

Caution: Take care when using this feature with detectors inserted. Also, verify the clearance of installed stage hardware manually using the xT UI before performing the rotation in Maps.

- 1. Right-click where point 1 is to be placed.
- 2. On the context menu, click Rotational Alignment and then Place Point 1.



- 3. Right-click where point 2 is to be placed.
- 4. On the context menu, click Rotational Alignment and then Place Point 2.



- 5. Select the rotation that you want to use.
 - Stage Rotation: Uses stage rotation to align the image.
 - Scan Rotation: Uses scan rotation to adjust the image.

The microscope rotates to make the line between Points 1 and 2 horizontal and zooms the viewer to the line.



SEM and SDB Scan Rotation

This function rotates the scan and aligns the image. It has no effect on the stage movements and is solely a scan coil function, but is used to orient the image relative to mechanical rotation and detector direction.

			40 mm	179.1*
Scan rotation 179.1	• ^ 🗸	Link	Stage rotation 30.0	•

Scan Rotation Descriptions, SEM and SDB				
Control	Description			
Scan rotation	Displays and controls the current scan rotation in the microscope. Click the up arrow to assign the viewer rotation value to the scan rotation text box.			
Link	Select the Link checkbox to link the scan rotation with the viewer rotation. When linked, changes to the scan rotation angle automatically rotate the viewer and vice versa. This is indicated by the <i>green</i> rotational dial in the main viewer. When unlinked, the rotational dial is white.			
Stage rotation	See Stage Rotation.			

Stage Rotation

Type a value into the text box and press Enter to rotate the stage in the microscope.



Set sample tilt

Setting the sample tilt allows Maps software to perform stage moves with the stage unlinked. Having this tilt set lets Maps know how to drive the stage in x and y while maintaining the same height of the sample. This is vital when performing imaging within the Maps software.

Before you perform this procedure, confirm that the sample is perpendicular to the electron beam. This preliminary step adjusts the Maps stage to keep the sample at the same distance from the beam; thus, allowing navigation and tiling at stage tilt.



To set the sample tilt:

Before the sample tilt is set, Maps software displays the current stage tilt in degrees next to the menu command.

- 1. On the Microscope menu, click Set Sample Tilt.
- 2. Maps software asks if you want to set the current sample tilt to the current stage tilt and it will unlink Z from the working distance. Click **Yes**.

The Sample Tilt is now set. A check mark is displayed next to the menu command along with the stage tilt that was set.

Task Result

With the sample tilt set, you can now image the tilted sample with Maps software.



SEM Nav-Cam Alignment

For systems that have the Nav-Cam, Maps is able to acquire Nav-Cam images and align them. Maps is also able to save the Nav-Cam Alignment to be automatically applied to all future Nav-Cam images acquired with Maps.

Align the SEM Nav-Cam

- 1. Click the Import from Nav-Cam tool bar icon .
- 2. When the Nav-Cam image appears in Maps, right-click on its name in the Layer control.
- 3. On the context menu, click Align.

👻 🔽 Layer		·	12
Image: Scios 2 NavCarr Image: Scios 2 NavCarr <	Bring To Front Bring Forward Send Back Send To Back	6	
	Center View Center and Rotate View	Ctrl+Shift+C	100
	Export To Project		
	Alignment	•	Align
	Persist Nav-Cam Alignment		Fine Alignment
	Delete		Clear Alignment
			COLUMN SHE

- 4. Follow the alignment wizard to perform the coarse point alignment (1-, 2- or 3-Point) as described in Coarse Alignment (1-, 2-, or 3-Point).
- 5. Follow the alignment wizard to perform fine alignment adjustments as described in Fine Alignment.
- 6. Test the Nav-Cam alignment by navigating to features on the sample and acquiring preview images.

If the Nav-Cam image is not sufficiently aligned, then perform the alignment procedure again.

- 7. When the Nav-Cam image is sufficiently aligned, right-click on its name in the Layer control.
- 8. On the context menu, click Persist Nav-Cam Alignment.



All subsequent Nav-Cam acquisitions using Maps will use this alignment.

Clear the Nav-Cam alignment

This clears the alignment for an individual image.

- 1. Right-click the Nav-Cam image in the Layer control.
- 2. On the context menu, click Alignment and then Clear Alignment.

Alignment	Align
Persist Nav-Cam Alignment	Fine Alignment
Delete	Clear Alignment

Clear the saved default Nav-Cam alignment

This clears the alignment for all acquired Nav-Cam images.

- 1. On the **Options** menu, select **Settings -> Microscope**.
- 2. Click Clear Nav-Cam Alignment.

Acquisition times						
Display critical acquisition	time warning	Critical tile acquisition time [mm:ss]:	15:00			
Optics Settings						
Use Degauss when changing	ng Focus					
Optics settle delay [ms]:	1000					
Stage Settings						
Stage settle delay [ms]:	1000	Stage alignment (Electron): 0°	Clear			
Nav-Cam						
Clear Nav-Cam Alignment Door-mounted						
- From Microscope / To Microscope						
Use position alignment for	"To SEM"					

Phenom

This section describes user interface elements and procedures that are specific to the Phenom desktop SEM systems.

Topics include:

Phenom Connectivity

Phenom Settings

Phenom Operational States

Phenom User Interface Elements

Phenom Connectivity

When Maps starts for the first time, or when you have configured Maps to ask for a connection (see Phenom Settings), then the connection dialog opens as shown below.

Ma Connect to Pheno	m			×
	Device	Phenom-L (12.148.118.229)		∨ ∂
	is Phenom at the si	tart		
			Connect	Continue offline

Connect microscope

- 1. Click the **Device** dropdown menu and select a Phenom microscope from the list of microscopes. If you do not see the desired microscope in the list, then do the following:
 - a. Click Refresh and then select the desired microscope from the refreshed list.
 - b. If the refreshed list does not contain the desired microscope, then enter the IP address of the desired microscope in the **Device** field.
- 2. *Optional*: If you added a new Phenom microscope to the list of microscopes (see Phenom Settings), then select the **Remember this Phenom** checkbox. Maps will then add the new Phenom microscope to the list of known microscopes, so that it is a menu option in the future.
- 3. *Optional*: If you want Maps to automatically connect to your selected microscope upon startup, then select the **Automatically reconnect with this Phenom at the start** checkbox.
- 4. Select one of the following options to complete this procedure:
 - Click **Connect** to initiate the connection between Maps and the selected Phenom microscope.
 - Click **Continue offline** if you do not want to connect to a microscope. Maps will continue to perform in the offline mode.

Connection states

The connection state icon is located in the lower-right corner of the Maps main screen. There are four color-coded connection states as described below.

States	Description
	A green icon indicates that Maps is connected to the microscope with full use and control.
	A blue icon indicates that Maps is connected to the microscope, but the connection is read-only since the microscope is locked by some other process (es). With this status, you can still use Maps, but you cannot use or control the microscope.
0	A yellow icon indicates that Maps has lost the connection to the microscope. Note that the lost connection might be temporary since Maps automatically attempts to reestablish the connection while in this status. With this status, you can still use Maps, but you cannot use or control the microscope.
	A red icon indicates that Maps is not connected to the microscope. With this status, you can only use Maps in the offline mode.

View microscope details and disconnect

If you want to view the microscope details and/or disconnect the microscope from Maps, perform the procedure described below.

Ma Phenom instrument				×
	State	Connected		
	ID	XL-Sim-6.1		
	Address	12.148.118	.229	
	Туре	PhenomXL		
	Software version	6.1.1.rel.43	7984a.25336	
			Disconnect	Close

- 1. From the Main screen, click **Phenom->Phenom instrument** to open the Phenom instrument dialog. This dialog displays the connection state and microscope details.
- 2. Click **Disconnect** to disconnect the Phenom microscope. When disconnected, Maps switches to the offline mode. *Note that Maps cannot reconnect or connect to another microscope when it is in the offline mode.*

Phenom Settings

The Phenom settings allow you to define the Maps connection startup behavior and manage the known Phenom microscopes.

Ma Settings					×
Application	Maps startup				
XPS	Ask for connection	XL-Sim-6.1.1-rel (12.148.1)	18.229)	~	
Phenom	Start offline				
Сгуо	- Known Phenom microscopes -				
Stitching	Instrument ID	IP	User		
	MVE049379-10002-L	12.183.245.133	maps		<u>ش</u>
Internal	XL-Sim-6.1.1-rel	12.148.118.229	PhenomSimulationAcces	s	世
			Apply	ОК	Cancel

Configure settings

- 1. Navigate to the **Options** menu, click **Settings**, and then click the **Phenom** tab.
- 2. Click one of the following Maps startup options:
 - Ask for connection: The connection dialog appears when Maps starts (see Phenom Connectivity).
 - **Connect to microscope**: Maps automatically connects to the selected Phenom microscope when Maps starts.
 - Start offline: Maps starts in the offline mode without a Phenom microscope connection.
- 3. Under Known Phenom microscopes, review the list of stored Phenom microscopes. This list displays the instrument ID, IP address, and user name associated with the microscope.
 - To add a new microscope to the list, refer to the connection procedure under Connect microscope.
 - To remove a microscope from the list, click the **trash can** icon next to the microscope's user name. When a warning dialog prompts you to confirm the removal, click **OK** to save the change and close the warning dialog.
- 4. Click **Apply** to save your settings, or click **OK** to save your settings and close the Settings screen.

Phenom Operational States

This topic describes the operational states of the connected Phenom microscope. Maps can use and control a Phenom microscope only when the following criteria is met:

- The microscope is operational.
- The microscope is in the operational mode, which is the SEM mode.

If this criteria is not met, then Maps will prompt you to take corrective action through the Phenom instrument dialog.

Corrective actions

With each of these operational states you have the two options. You can either bring the microscope to the operational mode, or disconnect the microscope and continue with Maps offline.

Note: Note that Maps cannot reconnect to the microscope when it is in the offline mode.

Operational state: The sample is unloaded and the microscope door is open

- Click Close the door and move to SEM to bring the microscope to the operational mode.
- Click **Continue offline** to disconnect and switch Maps to the offline mode.

Ma	Phenom instrument					
		Sample is u	nloaded			
			Continue offline	Close the door and move to SEM		

Operational state: The microscope is on standby

- Click Activate and move to SEM to bring the microscope to the operational mode.
- Click **Continue offline** to disconnect and switch Maps to the offline mode.

Ma	Phenom instrument						
		Instrument is o	n standby				
			Continue offline	Activate and move to SEM			

Operational state: The sample is under NavCam

- Click **Move to SEM** to bring the microscope to the operational mode.
- Click **Continue offline** to disconnect and switch Maps to the offline mode.

Ma	Phenom instrume	nt		
		Sample is under NavCam		
			Continue offline	Move to SEM

Phenom User Interface Elements

Phenom Toolbar, Right Side

See Left Side, All Systems for common selections on all systems.

Toolbar, right side, Phenom



Toolbar Descriptions, Phenom		
ΤοοΙ	Description	
Import Nav- Cam	Imports a Nav-Cam image or acquires a new one if none is available, and imports the image to project. This top- down image of the stage can be used for point and click navigation of the sample. See Nav-Cam Alignment (SEM/SDB Only).	

Toolbar Descriptions, Phenom(Continued)		
ΤοοΙ	Description	
Take Snapshot	Acquires an image at the current stage position and horizontal field width (HFW), and then places the image in the Maps viewer as a snapshot image. The image is acquired using the current parameters set in the microscope.	

Phenom Tile Set Tab

This topic describes options in the Tile Set tab for the Phenom microscope.

Basic tab



Tile Set Tab, Basic Tab Overview		
Control	Description	
Name	Displays and sets the name of the current tile set.	
Tiles X, Y	Sets the number of tiles in the X and Y directions.	
Tile HFW	Specifies the horizontal field width (HFW) for each tile. Values between 10 µm and 1 mm are always available. Other values can be used too, but their usage is limited to certain High Voltage and Working Distance settings. Confirm that your desired value is available for the acquisition. Limited values are indicated with the following warning icon:	
Total HFW	Shows the total HFW of the whole tile set.	
Resolution	Specifies the image resolution. Select the resolution from the list or enter a custom resolution.	
Pixel Size	Displays the physical size of a single pixel at the current acquisition settings. This is based on Tile HFW and Resolution.	
Averaging	 Sets averaging, or the number of image frames. Select a value from the list of preset options, or enter a custom value (number 1-255). The preset options are: Live (1) Medium (4) High (16) Best (32) 	
Contrast	Sets contrast. The value is a number between zero and one hundred (that is, 0.0-100.0).	
Brightness	Sets brightness. The value is a number between zero and one (that is, 0.0-1.0).	

Tile Set Tab, Basic Tab Overview (Continued)		
Control	Description	
Image acquisition time	Displays the estimated time required to acquire a single image in the tile set.	
Estimated acquisition time	Displays the estimated remaining time required to acquire the whole tile set.	
Elapsed acquisition time	Displays the actual elapsed acquisition time.	
From Microscope	Applies the current microscope settings to the tile set.	
To Microscope	Applies the tile set settings to the microscope. In some instances, this operation might require additional time to process as shown in the dialog below:	
	Ma Apply settings	
	Applying tile set settings to microscope, please wait	
	Setting acquisition parameters	

Focus tab

Select the Focus Strategy to focus the electron beam to produce a sharp image.

• None: The application does not touch the focus at all.



• Fixed: Uses the same focus for every tile.

III 📮 [AF]		Ø
Focus Strategy	Fixed	~
Focus	5.5 mm	SET

To set the fixed focus:

- 1. In the microscope, navigate to the sample location and manually focus the image.
- 2. Click Set to retain the working distance in the tile set.

• Interpolated: Select three focus points on the sample. The focus plane is then defined by these points. Each tile's focus is set based on where it is located in this plane.



To set up the interpolated focus:

- 1. Navigate to the first location.
- 2. Manually focus the image.
- 3. When the image is in focus, click **Set**. The focus is retained for Point 1.
- 4. Repeat steps 1-3 for Point 2 and then for Point 3. Once all three points are set, the tiles are ready to be acquired.
- 5. Choosing points that are at opposite ends of the sample is important to get the best results.
- 6. Position of the points is visible in the viewer when the tile set is selected:



Autofunction tab

NAME	MODE
Auto Focus	First tile 🗸 🗸
Auto Contrast Brightess	LeftEdge 🗸 🗸
Final Auto Focus	None 🗸

Autofunction	Description
Auto Focus	Focuses the electron beam to produce a sharp image
Auto Contrast Brightness	Optimizes the contrast and brightness of the image
Final Auto Focus	Performs on the last focus optimization

Set autofunctions

- 1. Select the autofunctions to use during acquisition.
- 2. In the **Mode** combo box, select the mode in which the autofunction should be performed. Available modes include the following:
- None: The autofunction is not used.
- Every tile: The autofunction is run *locally* at every tile.
- First tile: The autofunction is run once at the first tile before the acquisition.
- Custom tile list...: The autofunction is run over specified list of tiles. Select this option to create a new list.

Create a custom tile list

Click **Custom tile list** option and the selection workflow panel opens and displays instructions to guide you. In the viewer, select tiles where the autofunction should be performed and name the new list. You can use the **Select Nth tile** option in the right panel to quickly select tiles at the selected Nth internal. You can finish the selection with closing the workflow panel or with leaving the autofunction tab. The new list is inserted to the mode combo box and can be re-used by other autofunctions in this tile set. Use the **edit** button if you need to further edit any tile selection.



Advanced tab

Use the controls on this tab to set advanced tile acquisition properties for each tile set.

▦	.	[AF]		لمب	ĄŶ	\oslash	
Over	ар Х, Ү			10%		10%	
				96 px	96 px 60 px		
Stage	X			-932.4	-932.439 µm		
Stage	Υ			4.411	2 mm		
Scan	Rotation			0 °	0°		
Detector		BSD Full			~		
HDR		OFF			~		
High Voltage		5.3 kV			~		
Intensity		Imag	Image		~		
Vacuum		Medium (10pa) 🛛 🗸 🗸		~			
Acquisition Order		Raster		~			
							_
	FROM	MICROSCO	DPE		TO MIC	ROSCOPE	

Tile Set Tab, Advanced Tab Overview

Control	Description
Overlap X, Y	Displays the percentage of overlap between tiles in the tile set, as well as the pixels.
Stage X	Sets the X stage location of the center of the tile grid.
Stage Y	Sets the Y stage location of the center of the tile grid.
Scan Rotation	Allows manual adjustment of scan rotation for a tile set. Either enter a value in the edit box here, or rotate the tile set in the viewer to update the scan rotation value.

	Tile Set Tab, Advanced Tab Overview (Continued)
Control	Description
Detector	Selects the detector for the acquisition. Available detectors may be different for different microscopes.
	Note: The SED detector requires High vacuum. If the SED detector is selected, then the vacuum control is set to High and deactivated.
	Activation of the SED detector might need some time, and the Maps application must wait for the activation while the tile set settings are applied to the microscope. During this waiting period, an estimate of remaining time is displayed in the progress dialog.
	Ma Apply settings
	Applying tile set settings to microscope, please wait
	Enabling SED (remaining time: 1 min : 19 sec)
	Cancel
	Note: You can abort the SED activation by clicking Cancel . When aborted, the SED detector is not set in the microscope.
HDR	Specifies whether HDR (High Dynamic Range) should be applied to images.
	Note: If HDR is applied, it also sets the image format to 16-bit.
High Voltage	Sets the high voltage setting. Select a value from the list of presets, or enter a custom value.
	The preset options are:
	• 5 kV
	• 15 kV
	• 20 kV

	Tile Set Tab, Advanced Tab Overview (Continued)
Control	Description
Intensity	 Sets beam intensity. Select a value from the list of preset values, or enter a custom value. The preset options are: Low (2.1) Image (3.3) Point (4.3) Map (5.1)
Vacuum	 Specifies the required vacuum pressure during acquisition. Options are: High: 0.1 Pa or 1.0 Pa (according to the specification of the connected Phenom microscope) Medium (10 Pa) Low (60 Pa) If the tile set settings are applied to a microscope, then the microscope might need some time to reach the required pressure. The Maps application waits until the pressure level stabilizes. During this waiting period, current pressure and an estimate of remaining time are displayed in the progress dialog.
	Ma Apply settings Applying tile set settings to microscope, please wait Setting High vacuum (pressure: 5.4 Pa, remaining time: 2 min : 27 sec) Cancel Note: If you need to abort the waiting period for the pressure stablization, click Cancel.

	Tile Set Tab, Advanced Tab Overview (Continued)
Control	Description
Acquisition Order	 Determines the order in which the contained tiles are acquired. Options are: Raster: Left to right, then top to bottom. Serpentine: Alternating left to right pattern per row, running top to bottom. Spiral In: Iterating over tiles in a clockwise pattern starting with the outer most tiles moving inward. Spiral Out: Iterating over tiles in a clockwise pattern starting with the
	inner most tiles moving outward.
From Microscope	Applies the current microscope setting to the tile set.
To Microscope	Applies the tile set settings to the microscope. In some instances, this operation might require additional time to process as shown in the dialog below:
	Ma Apply settings
	Applying tile set settings to microscope, please wait
	Setting acquisition parameters

EDS tab

⊞ 🕂 [AF] 🗄							
EDS Acquisition							
Acquisition Parameters							
EDS Dwell	1 ms						
Elements	Al, C, O						
EDS Resolution	Eighth (57x57) 🗸						

	EDS Tab Overview
Control	Description
EDS Acquisition (checkbox)	When selected, activates the acquisition of X-ray spectrums that are processed into separate element channels.
EDS Dwell	Sets the minimum time to acquire X-rays per tile. Values between 1 millisecond and 10 seconds are allowed.

	EDS Tab Overview (Continued)
Control	Description
Elements	Defines the elements that will be processed from the acquired spectrum. A color channel will be added for each element that you select: Image: ROI Image: ROI Image: ROI Image:
	additional elements. Note that any elements that you deselect will have their channels deactivated, but they will remain in the layer tree unless you manually delete them. See Element Selector for more information.
EDS Resolution	Sets the resolution of the EDS map, this is directly proportional to the tile set resolution

Phenom

Element Selector

To select the elements that you want to identify during acquisition, do the following:

1. Go to the **Elements** field and click its ellipsis icon ... to open the Element Selector as shown below.

M	Se	lect Ele	ments															×					
					5	elect th	e elem	ents vo	u want	to ider	ntifv du	ring ac	auisitio	n.									
							Right	click e	lement	to cha	nge its	color.											
1	H																		0		8	8	
	Li	Be											В								R	60	•
	Na	Mg									_			ζ.		H					G	242	•
	к	Са	Sc	Ti	v	Cr	Mn	Fe	Co	Ni	Cu	Zn	Ga					_			в	233	Ç
	Rb	Sr	Y	Zr	Nb	Mo	Тс	Ru	Rh	Pd	Ag	Cd	In	Ľ					Ж	Cance	H	Rese	ŧ
	Cs	Ва		Hf	Та	w	Re	Os	Ir	Pt	Au	Hg	П	Pb	Bi	Ро	At	Rn					
			La	Ce	Pr	Nd	Pm	Sm	Eu	Gd	ть	Dy	Но	Er	Tm	Yb	Lu						
			Ac	Th	Pa	U	Np	Pu	Am	Cm	Bk	Cf	Es	Fm	Md	No	Lr						
															ок	T	Can	cel					

- 2. Right-click each element that you want to identify during EDS acquisition. This action selects the individual elements.
- 3. Optional: To associate a specific color with each selected element, right-click the element and select the desired color from the color picker. The selected color is used for the EDS map generated for given element. If a color is not associated with the element, then the color is automatically selected from the system palette to match the color that you typically see in the xT UI.
- 4. Click OK.

Phenom Scan Rotation

This function rotates the scan and aligns the image. It has no effect on the stage movements and is solely a scan function, but is used to orient the image relative to the detector direction.



	Scan Rotation Descriptions, Phenom
Control	Description
Scan rotation	Displays and controls the current scan rotation in the microscope. Click the up arrow to assign the viewer rotation value to the scan rotation text box.
Link	Select the Link checkbox to link the scan rotation with the viewer rotation. When linked, changes to the scan rotation angle automatically rotate the viewer and vice versa. This is indicated by the <i>green</i> rotational dial in the main viewer. When unlinked, the rotational dial is white.

Phenom Post-Acquisition Actions

This section describes the post-acquisition actions for Phenom microscopes.



The table below contains the post-acquisition actions. You can activate any combination of these actions or all of them depending on your use case. If you activate all of them, then the Maps application performs them in the following order: Pump to HiVac When Done, Move to NacCam When Done, Go to Standby When Done.

	Post-Acquisition Actions
Control	Description
Pump to HiVac When Done	Returns the microscope to HiVac mode after all jobs are completed. Select this action This can be useful if you used LowVac or MedVac mode for the acquisition.
Move to NavCam When Done	Moves the sample to the NavCam after all jobs are completed.
Go to Standby When Done	Puts the microscope in the standby operational state after all jobs are completed.
TEM

This section describes user interface elements and procedures that are specific to the TEM, EFTEM, and STEM modes on TEM systems.

Topics include:

TEM User Interface Elements Define a Tile Set for TEM TEM Tiling Corrections Tiling Corrections Quality Magnification Alignment Use Interpolated Focus Method STEM Acquisition with Multiple Detectors STEM Intermediate Images Automated Image Filtering TEM Analytics

TEM User Interface Elements

TEM Options Menu

- For descriptions of common selections to all systems, see Options menu.
- For descriptions of TEM-specific selections outlined in purple, see the table below. Note that microscope settings are system specific.



		Options Menu Over	view	
Menu Selection	Description	1		
Microscope Settings	Ma Settings Application Microscope Acquisition Data Exchange XPS Analytics Amira-Avizo2D Cryo Stitching Autofunction Ignore fa autofunct set acquis desired, y Notificatio Autofunct specifies error dialo	Autofunctions Ignore failed autofunctions Autofunction failed dialog timeout Stage Use Backlash correction for Stage Stage Axis Angle Correction * Doptics STEM Scan Orientation Correction * Beam Blanker Unblank Setting Time Beam Blanker Unblank Setting Time iled autofunctions: Wh ion errors during tile set sition automatically con you can view failed autofunction automatically con you can view failed autofunction panel. They are lister tion failed dialog timeof the timeout period for a og. Tile set acquisition of modify the default value	30 s 0.00 0.00 100 ms Apply en you select this ch t acquisition are igno tinues without an err function errors within d as pipeline warning out: When you deseler r Settings >> Microso utomatically closing continues after the er	X OK Cancel A A A A A A A A A A A A A

	Options Menu Overview(Continued)
Menu Selection	Description
	Stage settings:
	• Use Backlash correction for Stage: When this check box is selected (default), stage moves are more accurate because the moves are performed in a way to reduce backlash and wind-up effects. Usually there is no need to turn this correction off, and this switch is mostly used for diagnostics. One might consider temporarily turning off the correction to speed up low-magnification overviews (at the cost of accuracy) as the corrected move takes a bit longer.
	• Stage Axis Angle Correction: This is a small correction angle that allows for compensating for mechanical imperfections, if present, in the orthogonality between the X and Y axis of the stage. The default value is 0. Only adjust this value if there is a recurring, significant misalignment between tiles in TEM and STEM modes.
	 Optics settings: STEM Scan Orientation Correction: This is a small correction angle for compensating imperfections in the calibration of the STEM scan rotation. The default value is 0. Only adjust this value if there is a significant systematic misalignment between tiles in STEM mode. Beam Blanker Unblank Setting Time: After blanking or unblanking the beam, it takes a short time for the optics to settle. The default values are sufficient for most systems, but on Tecnai systems without a fast beam blanker, a value of 1-3 seconds is recommended. An insufficient setting time manifests itself only in STEM mode, resulting in tile images where the top part starts out black and unstable.

	Optio	ons Menu Over	view(Contin	ued)
Menu Selection	Descriptior	1		
Acquisition	STEM Mode			
Settings (STEM Mode)	Displays the Detector, Dwell Time, and Resolution Settings to be used for acquisitions of snapshot images when in STEM mode. This screen also displays the Default Autofunctions Templates. These dropdown menus enable you to select a template for autofunctions with default imaging parameters that differ from the imaging parameters of tile images. See the TEM Default Autofunction Templates for details.			
	The default settings for Eucentric Template and Focus Template contain default autofunction parameters for new tile sets.			
	Ma Settings			×
	Application	EFTEM STEM	ТЕМ	
	Microscope	Snapshots Detector: HAAD	F 🗸	
	Acquisition	Dwell Time: 10 µs		
	Data Exchange	Resolution: 512 x	512 🗸	
	XPS	Default Autofunctions Templa Eucentric Height Template:	tes	None (STEM)
	Analytics	Focus/OptiSTEM Template:		None (STEM) V
	Amira-Avizo2D	 Eucentric Height Default Settin Final Stage Angle ° 	ngs 15.0	
	Сгуо	Max. Z-Height Deviation	250 nm	
	Stitching	Focus Default Settings Coarse Step	3 um	Include Coarse Search
		Fine Step	600 nm	
		Number Of Steps Dwell Time	10 2 μs	
				Apply OK Cancel

	Optic	ons Menu Over	view(Contin	ued)
Menu Selection	Description	1		
Acquisition Settings (TEM Mode)	TEM Mode Selects the E Snapshots an such as binni Templates. T autofunctions parameters of for details. The default s	Exposure Time to re always acquire ing 1. This screen These dropdown is with default ima of tile images. See settings for Eucer	use for the ac ed with the ma n also display menus enable oging paramet e the TEM De ntric Template	equisition of snapshot images. Eximum camera resolution, as the Default Autofunctions be you to select a template for ers that differ from the imaging fault Autofunction Templates and Focus Template contain
	default autofu	EFTEM STEM Snapshots Camera: Current Exposure Time: 1 s Resolution: Maximu	TEM TEM	× sets.
	XPS Analytics Amira-Avizo2D	Default Autofunctions Templat Eucentric Height Template: Focus Template: Eucentric Height Default Settin Final Stage Angle °	ngs 15.0	Search Map None (TEM) Use Image Filtering
	Cryo Stitching	Max. Z-Height Deviation Focus Default Settings Desired Defocus Iterate to Focus Method	250 nm 0 μm -5 μm Objective Lens	Use Auto Stigmate Use Three Image Method
				Apply OK Cancel

	Optic	ons Menu Over	view(Contin	ued)
Menu Selection	Description	1		
Acquisition Settings (EFTEM Mode)	EFTEM Mod Selects the E Snapshots an such as binn Templates. T autofunctions parameters of for details. The default s default autofu	e Exposure Time to re always acquire ing 1. This screet These dropdown s with default ima of tile images. Se settings for Eucer unction paramete	use for the a ed with the ma n also display menus enable aging parame e the TEM De ntric Template ers for new tile	cquisition of snapshot images. aximum camera resolution, vs the Default Autofunctions e you to select a template for ters that differ from the imaging efault Autofunction Templates e and Focus Template contain e sets.
	Max Settings Application Microscope Acquisition Data Exchange XPS Analytics Amira-Avizo2D Cryo Stitching Stitching	EFTEM STEM Snapshots Curren Camera: Curren Exposure Time: 1 s Resolution: Maxim Default Autofunctions Template: Focus Template: Focus Template: Maxim Eucentric Height Default Setting Max. Z-Height Deviation Focus Default Settings Desired Defocus Iterate to Focus Method	TEM t um tes 15.0 250 nm 0 μm -5 μm Objective Lens	None (EFTEM) None (EFTEM) None (EFTEM) Use Image Filtering Use Auto Stigmate Use Three Image Method See Use Image Filtering

	Options Menu Overview(Continued)
Menu Selection	Description
Data Exchange Settings	Missettings X Application GRPC server Port: 500 GRPC Server version 3.0.0.69 is running. Start Acquisition Start Data Exchange GRPC client Server response : XPS Analytics Test Connection Stitching Stitching
	 GRPC server: Port: Shows the port number that the clients use to communicate with the server. This is used for diagnostics purposes, in case there is a conflict with another GRPC-based server. Server status: Shows the current version and status of the server. Possible states are Not installed, Running, Stopped. Start: Starts the server if it is not running. The server must be installed. Stop: Stops the server.

Options Menu Overview(Continued)		
Menu Selection	Description	
	 GRPC client: Server response: Shows the response from the server when the connection test is in progress. Shows the result of the test when the test is complete. 	
	• Test Connection : Tests the connection to the server while the server is running. This connection test is used for diagnostics purposes. It performs actions that Maps client does when connecting to the server and shows the result of each of the actions.	
Tile Set Templates	On TEM, you cannot edit custom tile set templates. See Tile Set Templates for full template instructions.	
Alignment Corrections	Displays the corrections that currently apply for metrology analysis.	
Tiling Corrections	Shows dialog with the list of tiling corrections that are currently available. See TEM Tiling Corrections topic for more details.	
Automatic Histogram Stretching	Automatically equalizes the minimum and maximum of associated tile set histograms during acquisition.	

TEM Toolbar, Right Side

See Left Side, All Systems for common selections to SEM/SDB systems.

Toolbar, right side, TEM, EFTEM, STEM modes



٦	oolbar TEM, EFTEM, STEM mode Descriptions
Tool	Description
Run Eucentric Height autofunction at current position	Runs Eucentric Height autofunction at current stage position. Default autofunction template for Eucentric Height autofunction will be used to set acquisition parameters on the microscope. Click Stop to cancel operation.
Take Snapshot	Acquires an image at the current stage position and magnification and places it on the Maps viewer as a snapshot image. See TEM Options Menu.
Magnification Alignment	Aligns tile sets from one magnification to another to correct small shifts. See Magnification Alignment.
Import Fluorescence Data	Opens dialog where fluorescence data from optical microscope can be selected and imported to the current Maps project. Note: Requires Cryo features to be activated within the Settings.
Import Arctis Data	Opens a dialog for you to select and import Arctis data into the current Maps project. See Import and view Arctis data. Note: Requires Cryo features to be activated within the Settings.

Toolbar TEM, EFTEM, STEM mode Descriptions(Continued)		
Tool	Description	
Toggle Automatic Synchronization	Toggles automatic synchronization of data with Tomography ON and OFF. Note: Requires Thermo Scientific Maps Data Exchange Service to be installed and running.	
Start Workflow	Opens workflow selection dialog, from where specific alignment workflow can be started.	

TEM Tile Set Tab

Basic group, TEM and EFTEM modes

🌐 🕂 🏝	** 📀		
Name	Tile Set (2)		
Tile Set Type	тем		
Camera	BM-Ceta 🗸 🗸		
Tiles X, Y	3 3		
Overlap X, Y	20% 20%		
	800 рх 800 рх		
Tile HFW	10.0251 µm		
Total Area	26.1 µm x 26.1 µm		
Resolution	4096 x 4096 🗸 🗸		
Magnification	SA 3800 x 🗸 🗸		
Pixel Size	2.5063 nm		
Exposure Time	1s		
0 of 0 tiles acquired; 19.48 KB			
FROM MICROSCOPE	TO MICROSCOPE		

Tile Set Tab, TEM and EFTEM Modes, Basic Group Overview		
Control	Description	
Name	Displays the name of the current tile set.	
Tile Set Type	Indicates that the acquisition mode for this tile set is TEM, EFTEM, or STEM.	
Camera	Selects the camera to use for the acquisition of tile images. Unsupported cameras are also listed and indicated in the camera dropdown list. These unsupported cameras may work with Maps, but compatibility is not guaranteed.	
Tiles X, Y	Sets the number of tiles in the X and Y directions.	
Overlap X, Y	Displays the percentage of overlap between tiles in the tile set as well as the pixels.	
Tile HFW	For TEM, this is not editable. It is determined by the magnification, camera, and other optical settings.	
Total Area	Displays the total HFW for all of the tiles.	
Grid Squares Mask	Selects the detected grid layer to apply as a mask for the current layer.	
Resolution	Specifies the image resolution. The choices are specific to each system type.	
Magnification	Selects the magnification to use for the acquisition of the tile images	
Pixel Size	Displays the physical size of a single pixel at the current acquisition settings. This is based on Tile HFW and Resolution.	
Exposure Time	Specifies the exposure time on the camera for each tile acquisition.	
Number of Tiles Acquired	Displays the number of tiles and the total disk space required to acquire the current tile set. A message appears if there is not enough space on disk for the acquisition.	

Tile Set Tab, TEM and EFTEM Modes, Basic Group Overview(Continued)		
Control	Description	
From Microscope	Reads the current optical settings from the microscope and stores them in the acquisition settings of the selected tile set.	
To Microscope	Sets the optical parameters of the acquisition settings of the selected tile set to the microscope.	

- **i** Note: From Microscope/To Microscope controls:
 - Camera settings are not read from or written to the microscope; camera settings only apply during acquisition.
 - Not all the optical acquisition settings can be edited directly. Some optical acquisition settings require clicking **From Microscope** to update the values.

Focus group, TEM and EFTEM modes

⊞	Ŧ	[AF]	-	ţţţ	Ø		
Focus Method Fixed V					~		
Nom	Nominal Defocus 0 µm						
Objective Lens 89.4448 %							
Stage Z						0 µm	Get
	FROM		DE		TOM	ICROSCORE	
	PROM M	nenosee			IO M	ICROSCOPE	

Tile Set Tab, TEM and EFTEM Modes, Focus Group Overview				
Control	Description			
Focus Method	 Specifies whether the tile set is acquired with at all. The options are: None: The focus on the microscope is left Fixed: The tile set is acquired at a fixed for tile set with the From Microscope button Defocus for contrast formation. Interpolated: You select three focus point plane is defined by these points. Each tile is located in this plane. See Use Interpolated the interpolated focus method procedure. 	et as-is. t as-is. tocus value to be stored in the plus an optional relative ats on the sample and a focus of ocus is set based on where it ated Focus Method to perform transference focus 3 oscope		
Nominal Defocus	 When Focus Method is set to Fixed, specifier to apply when acquiring the tiles. Note: Sets the focus to the microscope defocus is only applied during acquisiti top of the Focus setting that was stored 	e, not the defocus. The ion as a relative value on d with the tile set.		

Tile Set Tab, TEM and EFTEM Modes, Focus Group Overview (Continued)				
Control	Description			
Objective Lens	When the Focus Method is set to Fixed , this is the objective lens strength setting to apply when acquiring the tile set. To update this value, click From Microscope .			
Stage Z	When Focus Method is set to Fixed, specifies the Z coordinate at which the tile set is acquired.Click the check box to retrieve the current value from the instrument.			
From Microscope	Reads the current optical settings from the microscope and stores them in the acquisition settings of the selected tile set.			
To Microscope	Sets the optical parameters of the acquisition settings of the selected tile set to the microscope.			

Autofunction group, TEM and EFTEM modes



Use the controls in this group to define the Eucentric Height and Focus autofunction parameters for each TEM tile set. The Eucentric Height autofunction parameters are initially set at the creation of the tile set and filled with the values of the selected template, or with the current values on the instrument if the default template is used. After creation of the tile set, click **From Microscope** to update the values to the current state on the microscope. Some settings can be entered directly. For each autofunction, for example, you can specify which tiles of the tile set to run the autofunction. The list of values is specific for each autofunction.

Tile	Set Tab, TEM and EFTEM Modes, Autofunction Group Overview
Control	Description
Name	 Available autofunctions are listed in this column. Eucentric Height: The Eucentric Height autofunction will adjust the sample height so that the sample can be tilted and the area of interest remains in the field of view. For the current tile set, you can select a custom autofunction template that is different from the default autofunction template.
	 Focus: The Focus autofunction will adjust the focus for a clear, sharp image. For the current tile set, you can select a custom autofunction template that is different from the default autofunction template. For both autofunctions, the options are:
	 The default autofunction template is the template that you defined under the Tile Set Templates settings.
	 The other options include a dynamic list of templates that you define as the user.

Control	Description					
Mode	Mode selection of the autofunction.					
	Note: For the Eucentric Height autofunction, only the <i>None</i> and <i>First Tile</i> modes are available.					
	Choices from the dropdown menu include the following: • None: Autofunction does not run on the tile set					
	 First Tile: Autofunction runs only on the first tile within the tile set. 					
	• Every Tile: Autofunction runs on every tile within the tile set.					
 Custom Tile List: This allows you to select a custom collection or where autofunction runs. Select your collection of tiles by pressin Ctrl+right-click, or by using the Select Nth tile option in the right 						
	With Management Image: Margine Management Image: Margine Management Image: Margine Margine Management Image: Margine Ma					
	Ci Tim A DOS CONTROL MANY					

Tile Set Tab, TEM and EFTEM Modes, Autofunction Group Overview(Continued)

Configure autofunctions

You can configure parameters for each autofunction by selecting **Autofunction**, and then clicking **Configuration** or the gear icon.

Ma Advanced Autofunction Sett	tings (Tíle Set)		×
Eucentric Height	Final Stage Angle •	15.0	
Focus	Maximum Z-Height Deviation	250 nm	
	🚥 Use Image Filtering		
			Reset
		Ok	Cancel

Tile Set	Tile Set Tab, TEM and EFTEM Modes, Autofunction Configuration					
Autofunction	Control	Description				
Eucentric Height	Final Stage Angle°	Specifies the final stage angle. The default value is set to 15.0 degrees and accepts values from 1.0 to 90.0 degrees.				
	Maximum Z- Height Deviation	Specifies the maximum accepted Z-height difference measured between negative and positive final stage angles. The default value is set to 250 nm and it accepts values from 250 nm to 2 μ m. Autofunction processes more quickly if you use a high value for this control, but the result will be less accurate.				
	(Beta) Use Image Filtering	When activated, automatic image filtering is applied to cross-correlation algorithm.				

Autofunction	Control	Description
Focus	Desired Defocus	Desired result of Focus Autofunction - expressed as a difference against optimal focus.
	Iterate to	Specific focus level to which Focus Autofunction iterates - expressed as a difference against optimal focus.
	Focus Method	Method used to adjust current focus. Possible values are Objective Lens and Stage Z Height .
	Use Auto Stigmate	When enabled, performs astigmatism correction.
	Use Three Image Method	When enabled, uses three image method to determine cross-correlation and peaks, as the function computes focus adjustment.
	(Beta) Use Image Filtering	When activated, automatic image filtering is applied to cross-correlation algorithm.

Autofunction error dialog

If you experience an autofunction error (and the **Ignore autofunction errors** option is not selected under Settings >> Microscope), then the following error dialog appears. This error dialog closes after the specified time (defined in the **Autofunction failed dialog timeout** option under Settings >> Microscope), and the tile set acquisition job continues.



343 | Original Instructions

Choices include:

- Yes {28s}: Click to ignore the error dialog and continue with the image acquisition. The value in brackets indicates the total number of seconds required to perform the image acquisition.
- **Ignore**: Click to prevent these error dialogs from appearing in the future. If you opt to ignore future error dialogs, you can still view failed autofunction errors within the TEM Notification Panel. They are listed as pipeline warning messages.
- No: Click to stop the acquisition job.

Advanced group, TEM mode

The advanced parameters are initially set at the creation of the tile set and filled with the settings of the selected template (or filled with the current settings on the system if the default template is used). After creation of the tile set, click **From Microscope** to update the settings to the current state on the instrument. For some settings, you can enter values directly into the user interface.

III 🕂 [AF] 🗉	¢۴	Ø				
BM-Ceta Noise Reduction						
Frames Summed	4					
- Optics						
Illuminated Area		35.508	9 µm			
Spot Size		6	~			
Probe Mode		Micro	Probe 🗸			
C2 Aperture		None	~			
 Correction 						
Use image corrections						
Use Tiling Correction						
Stage						
Settling Time		0 s				
FROM MICROSCOPE		TO MICRO	DSCOPE			

Tile	Set	Tab.	TEM	Mode.	Advanced	Group	Overview
		,		,			

Control	Description
Noise Reduction	Specifies the Noise Reduction setting associated with the BM-Ceta camera. See the <i>Ceta 16M Application Instructions</i> for details.
Frames Summed	Specifies the Frames Summed setting associated with the BM-Ceta camera. See the <i>Ceta 16M Application Instructions</i> for details.

Tile Set Tab, TEM Mode, Advanced Group Overview (Contin- ued)			
Control	Description		
Illuminated Area	Specifies the beam intensity to use during the tile set acquisition.		
Spot Size	Specifies the spot size to use during the tile set acquisition.		
Probe Mode	Specifies the probe mode, such as the NanoProbe or MicroProbe, to use during the tile set acquisition.		
C2 Aperture	Specifies the C2 aperture size to use during the tile set acquisition. Select a value to activate, or select None to deactivate.		
Use image corrections	When this setting is activated, a small image correction is applied to the acquired tile images resulting in improved tile set accuracy. Using this setting slightly affects the raw data recorded by Maps. If this result is undesired, turn off the image correction. This results is less accurate navigation.		
	• Note: When the correction is activated, the tiles are slightly smaller. Also, with the correction activated you can notice a small difference between the purple field-of- view rectangle and outline of the tiles.		

Tile Set Tab, TEM Mode, Advanced Group Overview (Contin- ued)			
Control	Description		
Use Tiling Correction	When this setting is activated, tile alignment is corrected during acquisition according to a matching tiling correction. If matching correction is not available, then by default Maps performs an uncorrected acquisition. See TEM Tiling Corrections.		
Settling Time	Specifies the duration of time in which acquisition is paused before each tile, which allows the stage to settle after moving to the target position.		
From Microscope	Reads the current optical settings from the microscope and stores them in the acquisition settings of the selected tile set.		
To Microscope	Sets the optical parameters of the acquisition settings of the selected tile set to the microscope.		

Advanced group, EFTEM mode

The advanced parameters are initially set at the creation of the tile set and filled with the settings of the selected template (or filled with the current settings on the system if the default template is used). After creation of the tile set, click **From Microscope** to update the settings to the current state on the instrument. For some settings, you can enter values directly into the user interface.

III 🕂 (AF) 🗐	<u>ا</u> به)		
EF-Falcon				
Camera Mode	Linear		~	
- Ontics				
Intensity		0.400		
Spot Size		б	~	
Probe Mode		MicroProbe	~	
C2 Aperture		None	~	
Insert Slit		Yes	~	
Slit Width		10.0 eV		
Correction				
Use image corrections				
Stage				
Settling Time		0 s		
FROM MICROSCOPE	тс) MICROSCOPE		

Tile Set Tab, EFTEM Mode, Advanced Group Overview

Control	Description
Camera Mode	Specifies the Camera Mode setting associated with Falcon camera.
Intensity	Specifies the beam intensity to use during the tile set acquisition.
Spot Size	Specifies the spot size to use during the tile set acquisition.
Probe Mode	Specifies the probe mode, such as the NanoProbe or MicroProbe, to use during the tile set acquisition.

Tile Set Tab, EFTEM Mode, Advanced Group Overview (Continued)					
Control	Description				
C2 Aperture	Specifies the C2 aperture size to use during the tile set acquisition. Select a value to activate, or select None to deactivate.				
Insert Slit	When this setting is set to Yes , energy-selecting slit is inserted before tile set acquisition. When the value is set to No , energy-selecting slit is retracted before tile set acquisition.				
Slit Width	Specifies the width of the energy-filtering slit.				
 Note: This setting is applied only when the energy-filtering slit is set to be used for the acquisition (Insert Slit settings is set to Yes). 					
Use image corrections	When this setting is activated, a small image correction is applied to the acquired tile images resulting in improved tile set accuracy. Using this setting slightly affects the raw data recorded by Maps. If this result is undesired, turn off the image correction. This results is less accurate navigation.				
	• Note: When the correction is activated, the tiles are slightly smaller. Also, with the correction activated you can notice a small difference between the purple field-of-view rectangle and outline of the tiles.				
Use Tiling Correction	When this setting is activated, tile alignment is corrected during acquisition according to a matching tiling correction. If matching correction is not available, then by default Maps performs an uncorrected acquisition. See TEM Tiling Corrections.				
Settling Time	Duration for which acquisition is paused before each tile, to let the stage settle after moving to the target position.				

Tile Set Tab, EFTEM Mode, Advanced Group Overview (Continued)			
Control	Description		
From Microscope	Reads the current optical settings from the microscope and stores them in the acquisition settings of the selected tile set.		
To Microscope	Sets the optical parameters of the acquisition settings of the selected tile set to the microscope.		

STEM Tile Set Tab

Basic group, STEM mode



The Set Tad, SIEM Mode, Basic Group Overview		
Control	Description	
Name	Displays the name of the current tile set.	
Tile Set Type	Displays the microscope mode for the tile set, which is either STEM, TEM, or EFTEM (see TEM Tile Set Tab for TEM or EFTEM modes).	
Tiles X, Y	Sets the number of tiles in the X and Y directions.	
Overlap X, Y	Displays the percentage of overlap between tiles in the tile set as well as the pixels.	
Tile HFW	For STEM, this is not editable. It is determined by the magnification, camera, and other optical settings.	
Total Area	Displays the total HFW for all of the tiles.	
Grid Squares Mask	Selects the detected grid layer to apply as a mask for the current layer.	
Mode	Selects the probe mode of the beam: Microprobe, Nanoprobe, or LM. Note that to use LM STEM, it is strongly recommended to first enable LM STEM via the TEM User Interface.	
Scan Field Of View	Defines the size of the scan area and magnification for the acquisition of each tile.	
Resolution	Specifies the image resolution. The choices are specific to each system type.	
Pixel Size	Displays the physical size of a single pixel at the current acquisition settings. This is based on Tile HFW and Resolution.	
Dwell Time	Specifies the amount of time the beam dwells on each pixel when images are acquired.	
Scan Rotation	Defines the orientation of the scanning, which determines the orientation of the image. The tile set orientation is adjusted accordingly. Typical use of this property is to line up the acquired images to a feature on the sample.	

ΤE	Μ	

Tile	Tile Set Tab, STEM Mode, Basic Group Overview (Continued)			
Control	Description			
Selected Detectors	Specifies the detectors to use for the acquisition of tile images. You are required to select at least one detector. Choices depend on which detectors are available on the system.			
Default Detector	Specifies the default detector for the acquisition. This detector is used for situations in which only one detector is required for acquisition such as within stitching and autofunctions. Choices only include the detector(s) that you selected for acquisition under Selected Detectors.			
Frame Time	The estimated time for each tile acquisition.			
Number of Tiles Acquired	Displays the number of tiles and the total disk space required to acquire the current tile set. A message appears if there is not enough space on disk for the acquisition.			
From Microscope	Reads the current optical settings from the microscope and stores them in the acquisition settings of the selected tile set.			
To Microscope	Sets the optical parameters of the acquisition settings of the selected tile set to the microscope.			

i Note: From Microscope/To Microscope controls:

- Camera settings are not read from or written to the microscope; camera settings only apply during acquisition.
- Not all the optical acquisition settings can be edited directly. Some optical acquisition settings require clicking **From Microscope** to update the values.

Focus group, STEM mode

				* * * * *			
▦	.	[AF]	-	ĄŶ	Ø		
Focus	s Method					Fixed	~
Inten	sity Lens					0.5560 %	
Obje	ctive Lens					79.5483 %	
Stage	żΖ						Get
	FROM N	AICROSCO	OPE		то м	ICROSCOPE	

	Tile Set Tab, STEM Mode, Focus Group Overview					
Control	Description					
Focus Method	Specifies whether to acquire the tile set with specific focus settings or none at all. There the following options:					
	None: The focus on the microscope is left as-is.					
	 Fixed: The tile set is acquired at a fixed focus value to be stored in the tile set with the From Microscope button. 					
	• Interpolated: You select three focus points on the sample to define a focus plane. Each tile focus is set based on where it is located in this plane. See Use Interpolated Focus Method to perform the interpolated focus method procedure.					
	Focus Method Interpolated 🗸					
	Intensity Lens 0.5560 %					
	Objective Lens 79.5483 %					
	Stage Z Get					
	Interpolated Focus					
	Set Focus 1 Set Focus 2 Set Focus 3					

Tile Set Tab, STEM Mode, Focus Group Overview (Continued)			
Control	Description		
Intensity Lens	When the Focus Method is set to Fixed , specifies the beam intensity percentage value at which the tile set is acquired. This lens focus setting is set during tile set acquisition. To update this value, click From Microscope .		
Objective Lens	When the Focus Method is set to Fixed , specifies the objective lens percentage value at which the tile set is acquired. This objective lens setting is set during tile set acquisition. To update this value, click From Microscope .		
Stage Z	When Focused Method is set to Fixed , specifies the Z coordinate at which the tile set is acquired.		
From Microscope	Reads the current optical settings from the microscope and stores them in the acquisition settings of the selected tile set.		
To Microscope	Sets the optical parameters of the acquisition settings of the selected tile set to the microscope.		

Maps 3.32.1 User Guide 355 | Original Instructions

Advanced group, STEM mode

Use the controls in this group to set advanced tile acquisition properties for each STEM tile set. The advanced parameters are initially set at the creation of the tile set and filled with the settings of the selected template or with the current settings on the instrument if the default template is used. After creation of the tile set, click **From Microscope** to update the settings to the current state on the microscope. Some settings can be entered directly.

⊞	.	[AF]	t1 ∳≬∮	Ø		
- Opti	Optics					
Spot	Size			8	~	
Came	era Lengt	h		14 m	102	
	5					
- Dete	ectors					
			Gain		Offset	
BF			50%		00%	
DF2			34.3%	5	6%	
DF4			53%	9	1%	
HAA	DF		25%	5	0%	
- Corre	ection —					
Use image corrections						
Stage	e					
Settlir	ng Time			0 :	5	
	FROM N	AICROSCO	PE	то міс	CROSCOPE	

	Tile Set Tab, STEM Mode, Advanced Group Overview
Control	Description
Spot Size	Specifies the spot size to use during the tile set acquisition.
Camera Length	Specifies the STEM camera length, which is typically used to tune the contrast of the images.

Tile Set Tab, STEM Mode, Advanced Group Overview (Continued)		
Control	Description	
Detectors	 Specifies the individual detector settings for the detectors available on the system. You can only change settings of the detectors selected for STEM acquisition. Choices for each detector include: Gain: Specifies the contrast setting of the detector, which represents the dynamic range of the detected signal (that is, it defines the amplification of the detector signal). Offset: Specifies the brightness setting of the detector, which represents the null level of the detected signal (that is, it defines the offset of the detector signal). 	
Use image corrections	When this setting is enabled, a small image correction is applied to the acquired tile images resulting in improved tile set accuracy. Using this setting slightly affects the raw data recorded by Maps. If this result is undesired, turn off the image correction (that is, at cost of slightly less accurate navigation).	
	• Note: When the correction is enabled, the tiles are slightly smaller. Also, with the correction enabled you can notice a small difference between the purple field-of-view rectangle and outline of the tiles.	
Settling Time	Specifies the amount of time it takes for the stage to settle.	
From Microscope	Reads the current optical settings from the microscope and stores them in the acquisition settings of the selected tile set.	
To Microscope	Sets the optical parameters of the acquisition settings of the selected tile set to the microscope.	

Autofunction group, STEM mode

Use the controls in this group to define the Eucentric Height and Focus autofunction parameters for each STEM tile set. The Eucentric Height autofunction parameters are initially set at the creation of the tile set and filled with the values of the selected template, or with the current values on the instrument if the default template is used. After creation of the tile set, click **From Microscope** to update the values to the current state on the microscope. Some settings can be entered directly. For each autofunction, for example, you can specify which tiles of the tile set to run the autofunction. The list of values is specific for each autofunction.

Name			Mode
Eucentric Height		None	
Default	~	None	*
Focus		Neee	
STEM 1um	~	None	~
Beea OptiSTEM			
STEM 1um	~	None	~
			Configuration

	Tile Set Tab, STEM Mode, Autofunction
Control	Description
Name	 Available autofunctions are listed in this column. Eucentric Height: The Eucentric Height autofunction will adjust the sample height so that the sample can be tilted and the area of interest remains in the field of view. For the current tile set, you can select a custom autofunction template that is different from the default autofunction template. Focus: The Focus autofunction will adjust the focus for a clear, sharp image. For the current tile set, you can select a custom autofunction template that is different from the default autofunction template autofunction template. (Beta) OptiSTEM: The OptiSTEM autofunction adjusts both the focus and
	condenser astigmatism (A1/C1). For the current tile set, you have the option to select a custom autofunction template that is different from the default autofunction template. This feature is only available on 2.15 and 3.15 TEM Server versions.
	 For all autofunctions, the options are: The default autofunction template is the template that you defined under the Tile Set Templates settings.
	 The other options include a dynamic list of templates that you define as the user.

	Tile Set Tab, STEM Mode, Autofunction (Continued)	
Control	Description	
Mode	Mode selection of the autofunction.	
	Note: For the Eucentric Height autofunction, only the <i>None</i> and <i>First Tile</i> modes are available.	
	Choices from the dropdown menu include the following:None: Autofunction does not run on the tile set.	
	• First Tile: Autofunction runs only on the first tile within the tile set.	
	• Every Tile: Autofunction runs on every tile within the tile set.	
	 Custom Tile List: This allows you to select a custom collection of tiles where autofunction runs. Select your collection of tiles by pressing Ctrl+right-click. 	
	Image: Control of Cont	

Configure autofunctions

You can configure parameters for each autofunction by selecting Autofunction, and then clicking Configuration or the gear icon.

Ma Advanced Autofunction Settings (OptiSTEM 1um)			×
Eucentric Height	Final Stage Angle *	15.0	
Focus	Maximum Z-Height Deviation	250 nm	
Bea OptiSTEM			
			Reset
		Ok	Cancel

Autofunction	Control	Description
Eucentric Height	Final Stage Angle°	Specifies the final stage angle. The default value is set to 15.0 degrees and accepts values from 1.0 to 90.0 degrees.
	Maximum Z-Height Deviation	Specifies the maximum accepted Z-height difference measured between negative and positive final stage angles. The default value is set to 250 nm and it accepts values from 250 nm to 2 μ m. Autofunction processes more quickly if you use a high value for this control, but the result will be less accurate.
Focus	Include Coarse Search	Allows using the coarse search method.
	Coarse Step	Specifies the coarse step value.
	Fine Step	Specifies the fine step value.
Autofunction	Control	Description
--------------------	-----------------------------	--
Autorunction	Control	Description
	Number of Steps	Specifies the number of steps for focus autofunction.
	Use Tree Image Method	Specifies if the tree image method is used.
(Beta) OptiSTEM	Algorithm	This read-only field specifies the OptiSTEM algorithm. Currently, only the A1/C1 algorithm is supported. This feature is only available on 2.15 and 3.15 TEM Server versions.

Autofunction error dialog

If you experience an autofunction error and the Ignore autofunction errors option is not selected under Settings >> Application, then you can view the following error dialog. The error dialog automatically closes after the time specified in the Autofunction failed dialog timeout option under Settings >> Microscope, and the tile set acquisition job continues.



Choices include:

- Yes {28s}: Click to ignore the error dialog and continue with the image acquisition. The value in brackets indicates the total number of seconds required to perform the image acquisition. This option is the default.
- Ignore: Click to prevent these error dialogs from appearing in the future. If you opt to ignore future error dialogs, you can still view failed autofunction errors within the TEM Notification Panel. They are listed as pipeline warning messages.
- No: Click to stop the acquisition job.

TEM Notification Panel

You can view additional information about the progress of the tile set acquisition jobs (and potentially any other job) in the Notification Panel, which is located under Job Control >> Notifications.

The Notification Panel contains these three sections.

- Filters: This section is located at the top of the screen and contains three filter buttons: Notifications, Warnings, and Errors. Toggle each button to show/hide the notification, warning, and error messages associated with each selected pipeline. If needed, click Clear All to delete the list of all pipelines and their associated messages.
- **Pipelines**: This section is located in the middle of the screen and contains a list of pipelines. A pipeline typically corresponds to a single job, and each pipeline indicates when it was started. If the pipeline is closed/ended, then is also indicates the time it finished and job duration.
- **Pipeline messages**: Click a single row to select a pipeline. Once selected, you can view the list of associated messages for the pipeline at the bottom of the screen.

JOB QUEUE PROCESSING ONOT		PROCESSING ON	NOTIFICATIONS	
i 5	9 NOTIFICAT	IONS 0 WARNINGS	B 3 ERRORS CLEAR ALL	
Na	me	Started Finished	Duration	
Tile S	et 2021/	01/04 08:30:44 2021/01/04 08:31:33	3 00:48	
Tile S	et 2021/	01/04 08:31:45 2021/01/04 08:32:58	8 01:12	
Tile S	et 2021/	01/04 08:34:53 2021/01/04 08:36:35	5 01:41	
	Name	Details	Date	
0	Tile Set	Preparing microscope	2021/01/04 08:34:53	1
0	Tile Set	Normalizing lenses	2021/01/04 08:34:53	
0	Tile Set	Performing initial auto functions	2021/01/04 08:35:07	L
0	Tile Set	Performing Focus autofunction	2021/01/04 08:35:07	
8	Tile Set	Focus failed. Preconditions not me	et. 2021/01/04 08:35:07	
0	Tile Set	Microscope set up	2021/01/04 08:35:36	
0	Tile Set	Preparing for tile [1,1]	2021/01/04 08:35:36	
0	Tile Set	Acquiring tile [1,1]	2021/01/04 08:35:36	
0	Tile Set	Saving image(s) of tile [1,1]	2021/01/04 08:35:42	
	Tile Set	Prenaring for tile (2.1)	2021/01/04 0R-25-42	
~ J	OB CONTROL	► RUN		

TEM Default Autofunction Templates

See the TEM Options Menu to open Settings >> Acquisition. This screen enables you to assign a template to each autofunction and mode (TEM, EFTEM, or STEM), which then applies to all projects. You can select any autofunction template from this screen except for the default factory template. If a template is not selected for an autofunction and mode (displayed as **None** in the dropdown menus), then it is not possible to run an autofunction.

TEM			

Ma Settings							×
Application	EFTEM	STEM	TEM				
Microscope	Snapshots Camera:	Current					
Acquisition	Exposure Time:	1 s					
Data Exchange	Resolution:	Maximun	1				
·····	Default Autofunction	s Templates					
XPS	Eucentric Height Temp	plate:		TEM 7k			~
Analytics	Focus Template:			TEM 70	k		~
Analytics	Furnantzia Uniaht Dafe						
Amira-Avizo2D	Eucentric Height Dera	ault Setting	15.0		🔽 💷 lise ima	ne Filtering	
	Man 7 Height Deviation	1	250		ose inte	gernering	
Сгуо	Max. 2-Height Deviation	on	250 nm				
	- Focus Default Setting	js					
Stitching	Desired Defocus		0 µm		Use Auto Stig	jmate	
	Iterate to		-5 µm		Use Three Im	age Method	
	Focus Method		Objective Lens	~	📃 📴 Use Ima	ge Filtering	
					Apply	ОК	Cancel

Template parameters

There are two parts to the autofunction parameters.

- Common imaging parameters (such as magnification, resolution, and pixel size): These parameters are defined by the autofunction template and used during the tile set autofunction run.
- Autofunction-specific parameters (such as final stage angle, use auto stigmate, and iterate to): These parameters are initially determined by the tile set template, but can be changed for each tile set instance. These settings are available under the configuration button under the Autofunction tab.

Follow these guidelines to achieve the best results from the autofunctions during your workflows:

- Either manually create tile set templates, or import templates from the Tomography application to be used as default autofunction templates.
- For manual template creation, create a new tile set and set parameters to be used for a specific autofunction run.
- Parameters suitable for autofunction templates can vary based on the specimen or tool type. For the Eucentric Height autofunction, it is recommended that you use a low SA magnification range; whereas, for the Focus autofunction, it is recommended that you use a higher SA magnification range.
- Consider selecting (Beta) Use Image Filtering if you want to improve the reliability of the default cross-correlation algorithm (depends on your sample type). See Automated Image Filtering for more information.
- If you right-click your mouse cursor directly on the tile set in a layer tree, then it is possible to save the tile set as a template. This template can then be aligned to a specific autofunction by selecting it from the Focus Template dropdown menus listed under Settings >> Acquisition.

Focus autofunction example

- 1. Create a new tile set (manually or import it).
- 2. Set the magnification to 120 kx and other parameters suitable for the Focus autofunction.
- 3. Right-click your mouse cursor directly on the tile set in layer tree and select **Save as Template**, select, **New Template**, and then enter **Focus Template** as the name.
- 4. Go to the main screen, click the **Options** menu, and then select **Default Autofunctions Templates**.

- 5. From the Focus Template dropdown menus, select your new focus template and close the dialog.
- 6. Create a new tile set and activate Focus Autofunction under that Autofunction tab.
- 7. Set additional parameters (such as desire defocus for, focus method, etc.) for the Focus autofunction through the **Configuration** button under the Autofunction tab.

TEM Viewer Tile Context Menu

 Drive Stage Here				
Take Snapshot				
Add Tiles Here				
Add Site of interest				
Tile Set	٠			
Tile 1,1	•	\	Acquire	
 Clear Digital Rotation			Open tile image folder	
Rotate Viewer to FOV			Open source image	
Rotate Viewer to Tile Set			Open analytics data folder	
			Send to Amira-Avizo2D	

Right-click on a tile within the tile grid to access a context menu.

- For common selections, see the Common Viewer Tile Context Menu Overview.
- For TEM-specific selections, see the table below. and Define a Tile Set for TEM.

	Viewer Tile Context Menu, TEM Overview
Menu Selection	Description
Take Snapshot	Takes a single snapshot at the current microscope location.

Vie	Viewer Tile Context Menu, TEM Overview (Continued)					
Menu Selection	Description					
Acquire/Queue for re-acquisition	Displays whether or not the tile has already been acquired as defined below. You can toggle between the two tile acquisition states as needed; however, only the <i>selected</i> acquisition state appears in the menu.					
	• Acquire: This acquisition state displays if the tile has not been acquired. If you select the check box for this menu item, then Maps queues the selected tile to be acquired.					
	• Queue for re-acquisition: This acquisition state displays if the tile has already been acquired. If you select the check box for this menu item, then Maps queues the selected tile to be re-acquired and replaces the original image.					
	For TEM, turning off a tile acquisition state will only be remembered while the tile is inside the stage boundary.					
	See Define a Tile Set for TEM for more information.					
Send to Amira- Avizo2D	If activated, sends the tile image to Amira-Avizo2D.					
Rotate Viewer to FOV	Digitally rotates the viewer to line up the display to the orientation of the current field-of-view (as indicated by the purple rectangle).					
Rotate Viewer to Tile Set	Digitally rotates the viewer so the current tile set is aligned with the screen. See Viewer Tile Context Menu for TEM.					

TEM Post-Acquisition Actions

This section describes the post-acquisition actions for TEM microscopes.

POST ACQUISITION	POST PROCESSING
Close Column Valves when Done	Stitch Tile Sets When Done

	Post-Acquisition Actions					
Control	Description					
Close Column Valves When Done	Closes the column valves after all jobs are completed. Select the check box to close the column valves after all jobs are completed.					
Switch off Emission when Done	Applies only to systems with thermionic guns, such as a Talos LaB6. Select the check box to turn off the gun emission for any queued acquisition to preserve the filament. In order to turn the gun emission on again after it is switched off, you must use the TEM UI, which is outside of Maps.					

Define a Tile Set for TEM

This section is a supplement to the general operations on tiles and tile sets. Tiles for TEM cannot be acquired when they are situated outside the stage boundary. That is why defining, resizing and moving a tile set on TEM is different from those actions on other systems. It also affects enabling and disabling **Acquire** property for a tile.

Methods for defining a tile set

A tile set can be defined the same way as described in Define a Tile Set. When creating a tile set, Maps software creates a layer in the Layer control. Tiles in that layer that are created inside the stage boundary have **Acquire** available by default. Tiles in that layer that are created outside the stage boundary cannot be acquired (**Acquire** command is unavailable).

Manually adjust tile sets

Once a tile set is created, you can adjust the size of the tile set with the drag handles in the viewer as described in Define a Tile Set.

Acquire menu command

	Drive Stage Here				
	Take Snapshot				
	Add Tiles Here				
	Add Site of interest				
	Tile Set	•			
	Tile 1,1	•	N	Acquire	
	Clear Digital Rotation			Open tile image folder	
Г	Rotate Viewer to FOV			Open source image	
L	Rotate Viewer to Tile Set			Open analytics data folder	
				Send to Amira-Avizo2D	

The **Acquire** menu command is described in **TEM** Viewer Tile Context Menu. When a tile Acquire property is turned off, and the tile is moved inside the stage boundary, then the **Acquire** command remains unavailable. When the tile Acquire property is turned off, and the tile is moved outside the stage boundary, and then back inside the stage boundary, then the **Acquire** command is made available.

Automatically deactivate a tile after acquisition

When a tile set has been acquired, the tile set cannot be moved again. Also, the **Acquire** menu option is not available when the tile has been acquired.

TEM Tiling Corrections

On TEM systems, you will observe imperfections in tile alignment during tile set acquisition. The imperfections vary across different magnifications and result in visible artifacts along the edges of individual tiles. The sources of these errors include magnification error, rotation error, and imprecision of stage movement. You can use the Tiling Correction beta feature to compensate for these errors and achieve better tile alignment during acquisition.

• Note: This functionality is intended to compensate only for relatively small errors that remain on an aligned and calibrated system. It is recommended that you always ensure that your system is properly calibrated and aligned before using Tiling Corrections.

Stitching verses tiling correction

There are two tiling correction mechanisms available in Maps for TEM systems: stitching and tiling correction.

Stitching creates a new image layer from a given tile set, in which individual tiles are moved so that the overlapping portions of neighboring tiles match each other. While the relative shift between neighboring tiles is computed and stored in tile registration data already during acquisition, stitching is a post-processing action and can be performed only after the acquisition is finished.

Tiling correction, on the other hand, takes advantage of the fact that the tile alignment error is static for a given set of optical settings. This means that the alignment error can be compensated by correcting tile rotation and magnification *during* tile set acquisition. The correction is computed from the same tile registration data as stitching, but instead of using the correction only on a single layer it is saved for subsequent acquisitions of all tile sets that match a given set of optical settings.

Create tiling correction based on existing tile set

1. Click the acquired tile set that has misaligned tiles. For example, the tiles in the following 3x3 tile set are misaligned since the cross-grating pattern is broken along the edges of neighboring tiles.



- 2. Right-click on the selected tile set.
- 3. In the context menu, click Beta Save As Tiling Correction.



4. When the Create Tiling Correction screen appears, review the correction parameters and quality score generated by Maps.

Ma Create Tiling Correction					
%Hfw	3.10				
∆Rotation	3.34°				
Quality		High			
(OK	Cancel			

5. Click **OK** to create the tiling correction.

Note: If the quality score is Medium or Low, move your mouse over the quality indicator to view the tool tip. The tool tip contains a short description of why the quality is lower than High and it provides suggestions for quality improvement. See Tiling correction quality for more information.

View available tiling corrections

When a tiling correction is created, Maps stores a record of the correction and associates it with optical settings that match the original tile set. Click **Options > Tiling Corrections** to view the records in the Tiling Corrections screen.

The tiling corrections are stored in a table per instrument mode within the Tiling Corrections screen, so that subsequent acquisitions benefit from the tiling corrections. Be aware that the tiling corrections are not part of the project; instead, they are part of the Maps global settings.

Ma Tiling Co	rrections						×
EFTEM	TEM						
						D	elete All
Mode	Magnification	Probe Mode	Camera	%Hfw	ARotation	Quality	
LM	210x	MicroProbe	BM-Ceta	1.81	0.22°	High	×
LM	700x	MicroProbe	BM-Ceta	1.92	-0.18°	High	×
нм	11000x	MicroProbe	BM-Ceta	3.10	3.34°	High	×
HM	150000x	NanoProbe	BM-Ceta	2.75	4.56°	Medium	×
Export	Import					ок	ancel

Correction Record Overview

Interface Item	Description
Mode	Indicates the magnification mode (values include: LM, HM).
Magnification	Specifies the magnification value.
Probe Mode	Indicates the probe mode (values include: MicroProbe and NanoProbe).
Camera	Displays the Camera name.

	Correction Record Overview(Continued)				
Interface Item	Description				
%Hfw	Specifies the horizontal field width correction that Maps applies during acquisition.				
ΔRotation	Specifies the rotation correction that Maps applies during acquisition.				
Quality	Indicates the quality score (values include: High, Medium, and Low). Maps uses a cross-correlation algorithm to determine the quality score. See Tiling correction quality.				

There are additional controls in the Tiling Corrections screen that you can use to import, export, or delete corrections.

	Tiling Corrections Table Overview				
Control	Description				
Export	Exports the corrections for TEM and EFTEM instrument modes to a file. Select the instrument tab (EFTEM or TEM), and then click Export .				
Import	Imports the corrections for TEM and EFTEM instrument modes from a file. Select the instrument tab (EFTEM or TEM), and then click Import .				
Delete All	Removes all corrections for the selected instrument mode (EFTEM or TEM).				
OK	Saves the changes made to the EFTEM and TEM tables, and then closes the window.				
Cancel	Cancels the changes made to the EFTEM and TEM tables, and then closes the window.				

Use tiling correction during acquisition

To use tiling correction, you will need to create a new tile set with parameters that match one of the saved tiling corrections and then acquire the tile set. For example, consider the correction that is highlighted below (parameters: HM, 11000x, MicroProbe, BM-Ceta, 3.10%, 3.34°, High).

Mode	Magnification	Probe Mode	Camera	%Hfw	A Rotation	Quality	
LM	210x	MicroProbe	BM-Ceta	1.81	0.22°	High	×
LM	700x	MicroProbe	BM-Ceta	1.92	-0.18°	High	×
НМ	11000x	MicroProbe	BM-Ceta	3.10	3.34°	High	×
HM	150000x	NanoProbe	BM-Ceta	2.75	4.56°	Medium	×

Create a TEM tile set

- 1. Click the TEM Tile Set tab, and then do the following:
 - a. Select **BM-Ceta** from the Camera dropdown menu.
 - b. Select **SA 11000x** from the Magnification dropdown menu.

⊞ ₽	[AF]	-	Ŷ\$	۲		
Name			Corre	cted Tile S	et	
Tile Set Type			TEM			
Camera			BM-C	ieta		~
Tiles X, Y			3		3	
Overlap X, Y			20%		20%	
			100 px		100 px	
Tile HFW			5.4871	μm		
Total Area			14.3 µ	m x 14.3 µr	m	
Resolution			512 x	512		~
Magnification			SA 11000 x			~
Pixel Size			10.974	13 nm		
Exposure Time			1s			
	0 of 9	images	acquired;	18.24 ME	B	
FROM	MICROSCO	PE		TO MI	CROSCOPE	

- 2. Click the Advanced tab.
 - a. Select MicroProbe from the Probe Mode dropdown menu.
 - b. Notice that the Use Tiling Correction checkbox is automatically selected.

III 🕂 🚝 🔳	₩ 📀
BM-Ceta Noise Reduction Frames Summed	4
Optics Intensity Spot Size Probe Mode C2 Aperture	0.590 5 ~ MicroProbe ~
Correction Image corrections Image corrections Image correction Image correction Stage Settling Time	0.
FROM MICROSCOPE	0 s TO MICROSCOPE

3. Acquire the tile set. If needed, refer to Acquire a Tile Set.

4. View the results in the main Viewer. Notice that the tiles are better aligned as a result of the tiling correction applied during acquisition.



Tiling correction quality

Tiling correction is computed from cross-correlation of overlapping portions of neighboring tiles. The quality and precision of the computation depends on several factors, including:

- Image intensity and focus
- Size of tile overlap
- Distinct features in the overlap portion of the tiles
- Stitching profile that matches sample type

- Note: As a general rule, it is recommended to use tile sets with the following parameters as a basis for tiling corrections:
 - Dimensions at least 3x3, with at least 20% tile overlap
 - All tile overlaps are successfully registered (no red outlines in the tile set)
 - Image has sufficient intensity and is in focus

You can observe the quality score of the tiling correction in the following screens:

- The Create Tiling Correction screen (when you are creating a new tiling correction)
- The Tiling Corrections screen (when you review available corrections)

See Tiling Corrections Quality for additional instructions on how to resolve issues related to quality.

Tiling Corrections Quality

This topic describes how you can resolve issues that negatively impact the quality and precision of tiling corrections.

i Note: For a general description of tiling corrections and the quality of tiling corrections, see TEM Tiling Corrections.

The following issues cause the quality score of a tiling correction to be lower than desired:

- Failed tile registrations
- Insufficient size of tile overlap
- Insufficient size of the tile set
- Large error margin in a portion of the tile set
- Large overall size of the correction

Failed tile registrations

You can recognize failed tile registrations by one or more red outlines visible in the acquired tile set. The failed registrations are caused by inability to compute cross-correlation on overlapping portions of neighboring tiles. This is usually due to poor imaging conditions, or when the tile set is being partially placed on a portion of the sample with no visible details.

For example, the tile set below has several failed tile registrations because the affected tiles are placed on a grid bar. To resolve this issue, reposition the tile set so that all tiles contain a sufficient level of detail and then acquire the tile set again.



It is recommended to use a tile overlap of at least 20% for tile sets that are used as a basis for tiling correction. The overlapping regions of neighboring tiles should ideally contain distinct features that the cross-correlation algorithm can use to compute necessary correction. A larger size of tile overlap gives the algorithm larger areas to work with, which is desirable.

For example, the following tile set has only 5% tile overlap, causing the quality indicator to report a low quality score. To resolve the issue, reset the tile set, increase the size of tile overlap to at least 20%, and then acquire the tile set again.



Insufficient tile set size

It is recommended to use a tile set of at least 3x3 size as a basis for tiling correction. This size gives the cross-correlation algorithm a sufficient number of overlapping regions to compute overall correction from since they are equally distributed across both columns and rows.

For example, the following tile set has only 3 tiles in a single row, thus causing the quality indicator to report a low level of quality. To resolve the issue, reset the tile set, increase its size to at least 3x3, and then acquire the tile set again.



Large error margin

Error margin is the difference between the average correction computed from the whole tile set and a local correction that is computed from a specific tile overlap. When the corrections computed for different tile overlaps differ by a large extent from each other, the quality indicator reports a large error margin.

Increased error margin can be caused by one of the following issues:

- Different imaging conditions in one portion of the tile set compared to the rest of the tile set.
- Recurring pattern in the specimen that causes false peak in the cross-correlation, so the correction targets the pattern rather than a distinct feature in the specimen.
- Lower signal-to-noise ratio such as low-dose conditions in which the signal counts are so low that recurring patterns and noise in the dark, and gain reference images, are interpreted as features of the specimen.
- Imprecision of mechanical stage movement at high magnifications.

For example, the following tile set contains a recurring grid pattern. The default stitching profile is not able to detect false peaks in cross-correlation, thus causing a large error margin in a portion of the tile set.



To resolve the issue in this specific use case, do the following:

1. Right-click the tile set to open the context menu and then select **Tile Registration** to open the Tile Registration - Define Parameters screen. See - topic for details.

2. Select **Repetitive Structures** from the stitching profile dropdown menu.

Ma Tile Registration - Define Parameters							
Re-running registration on Atlas 210x tile layer will replace the current registration. Do you want to queue registration job?							
Select a channel:	Default	~					
Select a stitching profile:	Repetitive Structures	~					
	Queue job Cano	el					

3. Click **Queue job** to rerun the tile registration with the selected stitching profile.

4. Create a tiling correction for the tile set based on the updated stitching profile (see Create tiling correction based on existing tile set).

Ma Create Tiling Correction					
%Hfw	3.63				
ΔRotation	0.49°				
Quality		High			
	ОК	Cancel			

5. View the improved quality for the tile set with the recurring grid pattern. As shown below, the tile alignment is properly aligned with the high-quality tiling correction compared to the original low-quality tiling correction.



Large correction size

Tiling corrections are intended to compensate only for relatively small errors that remain on a well aligned and calibrated system. If the computed size of the correction is too large, then the quality indicator will display a lower quality score. To resolve the issue, ask the system owner to verify and rerun the magnification calibrations on the microscope.

Magnification Alignment

On TEM systems, you will observe imperfections in the tile set alignment from one magnification to another. These imperfections result in small shifts when changing the magnifications. You must always properly calibrate and align the system, but small shifts may still remain since changes in the optics can lead to additional shift. To compensate for the small shifts that may remain, you can use the drag-and-drop functionality of Maps to align tile sets.

• Note: This functionality is available for both TEM and STEM mode, but the remainder of this section explains how it works for TEM tile sets. For STEM mode the functionality is similar, with the small difference that the magnification is a derived value (inferred from the Field of View).

Magnification Alignment versus Manual Alignment

There are two alignment mechanisms available in Maps: Magnification Alignment and Manual Alignment.

Magnification Alignment moves the data in the Maps canvas and *stores this alignment for subsequent acquisitions* at the same magnification.

Manual Alignment only aligns the data within the Maps canvas. Any subsequent acquisition will not have access to information about the alignment.

Align a Tile Set and Its Magnification

1. Click the acquired tile set that is slightly misaligned.

For example, the higher magnification tile set within the green border is misaligned with respect to the lower magnification.



2. Right-click on the selected tile set.

- 3. Do one of the following:
 - In the context menu, click Alignment and then click Magnification Alignment.

	Bring To Front				
	Bring Forward				
	Send Back				
i	Send To Back				
~	Show Grid Lines	Ctrl+G			
	Center View	Ctrl+Shift+C			
	Center and Rotate View				
	Save and Set As Default Template				
	Save As Template	•			
	Copy Grid				
	Export Grid				
	Export To Project				
	Alignment	•		Align	
	Alignment Drive To			Align Fine Alignment	
	Alignment Drive To Stitch	,	<u>s</u>	Align Fine Alignment Magnification Alignment	
	Alignment Drive To Stitch Tile Registration	,		Align Fine Alignment Magnification Alignment Clear Alignment	
	Alignment Drive To Stitch Tile Registration Split Channels	,		Align Fine Alignment Magnification Alignment Clear Alignment	
	Alignment Drive To Stitch Tile Registration Split Channels Create MIP Layer	,		Align Fine Alignment Magnification Alignment Clear Alignment	
	Alignment Drive To Stitch Tile Registration Split Channels Create MIP Layer Mark as done			Align Fine Alignment Magnification Alignment Clear Alignment	
	Alignment Drive To Stitch Tile Registration Split Channels Create MIP Layer Mark as done Reset Acquisition	,	22	Align Fine Alignment Magnification Alignment Clear Alignment	
	Alignment Drive To Stitch Tile Registration Split Channels Create MIP Layer Mark as done Reset Acquisition Open Z-Stack	,	22	Align Fine Alignment Magnification Alignment Clear Alignment	
	AlignmentDrive ToStitchTile RegistrationSplit ChannelsCreate MIP LayerMark as doneReset AcquisitionOpen Z-StackRecipe Processing			Align Fine Alignment Magnification Alignment Clear Alignment	

• On the right side of the tool bar, click Magnification Alignment

On the right side of the tool bar, click Workflow. See Workflow.
 The Workflow Magnification Alignment panel appears.



4. Reposition the tile set using the directional buttons in the Workflow Magnification Alignment panel.



You can reposition the tile set by clicking and holding the left-mouse button while hovering over the tile set.



5. Click the check mark within the alignment workflow panel to complete the current alignment, or click the X to cancel it.

The results of the alignment are twofold:

- The tile set with the existing data is realigned within the Maps canvas.
- An *alignment correction* record is stored for the magnification associated with the tile set. All subsequent acquisitions are automatically aligned based on this record. See Alignment Corrections Table.

TEM

Alignment Corrections Table

When a tile set is aligned, Maps stores a record of the magnification associated with the tile set. Click **Options > Alignment Corrections** to view the record in the **Alignment Corrections** table.



The **Alignment Corrections** are stored in a table per instrument mode, so that subsequent sessions benefit from the alignments. Note that the corrections are not part of the project; they are part of the Maps global settings.

Ma A	lignment C	Correction	s				\times
STE	м	TEM					
						Dele	te All
	Mode	;	Magnificatio	ΔΧ		ΔΥ	
N	lanoprobe	10	000x	-5 µm	-5 µm		×
E	xport	Imp	ort		ОК	Car	ncel

	Alignment Corrections Table Overview
Interface Item	Description
Import	Imports all the alignment corrections for TEM and STEM.
Export	Exports the alignment corrections for TEM and STEM.
Delete All	Removes all alignment corrections for the currently selected mode (STEM or TEM).
Cancel	Cancels the changes made to the tables and then closes the window.
OK	Saves the changes made to the tables and then closes the window.

Propagation of Alignment Corrections

The magnification alignments automatically propagate to higher magnifications that do not have an entry in the Alignment Correction table. However, individual magnifications can be aligned at any time, overruling the propagated alignment. With this approach, only a minimum number of alignments is required and it allows for refinement when needed. An additional benefit of this approach is that it naturally follows the typical Maps workflow from low-to-high magnification, allowing you to realign during the workflow when needed.

- In TEM, alignment corrections for low-magnification mode (LM) do not automatically propagate to the high-magnification modes (Mi/Sa/Mh), because on TEM systems these modes are mostly uncorrelated.
- In STEM, alignment corrections only propogate within the same mode (LM/NanoProbe/Microprobe).

The concept of the propagation is demonstrated in the example described below.

Example: Assuming you have an Alignment Corrections table as shown below and the system supports the magnifications as listed in Example of Propagation.

Example of Propagation

(Mode	Magnification	Correction	
ions	ſ	LM	50x	none	
ficat		LM	100x	none	<u> </u>
L.	-11	LM	500x	(-3 um, 4 um) 🛛 📍	ĺ –
N N		LM	1000x	(-3 um, 4 um)	- 2
ġ	Ц	LM	1800x	(-3 um, 4 um) 🗸	
	r	Mi	2000x	none	ר <u>ר</u> ן
S		Mi	2500x	none	-3
catio		SA	3000x	none	
gnifi		SA	5000x	(1 um, -2 um)	10
N	-	SA	12000x	(1 um, -2 um) 🗸	ΙŪ
눹		SA	20000x	(0.7 um, 0.2 um) 📍	1 I
-		SA	50000x	(0.7 um, 0.2 um)	
		Mh	100000x	(0.7 um, 0.2 um)	0
	L	Mh	200000x	(0.7 um, 0.2 um) 🗸	J
			-	-	•

In the example, the propagation works as follows. Numbers used below correspond to the annotation in the table above.

- 1. The first magnifications have no alignment correction.
- The LM 500x magnification has an alignment correction, which propagates to the next magnifications LM 1000x and LM 1800x, but not to the high magnification modes (3, 4 and 5).
- 3. The first high magnifications do not have an entry in the alignment corrections table, hence they have no correction.
- 4. The SA 5000x is the first high magnification with a correction, which propagates only to the next magnification SA 12000, because SA 20000x has its own entry in the alignment correction tables (5), which overrules the propagated correction of SA 5000x.
- 5. The correction of SA 20000x propagates to all the next magnifications.

Working with Magnification Alignments

- You can refine the alignment at any point in time, even when job execution is paused during tile set acquisition.
- A magnification alignment automatically applies to higher magnifications, so not all magnifications need realignment. See Propagation of Alignment Corrections.

- You can align multiple tile sets simultaneously by selecting a number of tile sets before starting the alignment.
- The alignment corrections are stored on disk and remain active for subsequent sessions.
- View which alignment corrections are present in the Alignment Corrections Table.
- If alignment corrections apply, you should see the field-of-view rectangle shift when changing the magnification on the system as Maps tracks the current field-of-view. Tracking includes the alignment corrections.
- Magnification Alignment can be used to compensate for shifts between STEM and TEM mode. This allows you to use a LM TEM tile set as reference for navigation while in NanoProbe STEM.

Alignment Tips

- The magnification alignments are for Maps only. They determine how Maps controls the stage and how acquired images are positioned in the Maps canvas without affecting the optics of the microscope. This mechanism can safely coexist with other applications that use different means for realignment.
 - The FOV visualization automatically incorporates the image beam shift set by other applications, such as with Thermo Scientific Tomography. This allows Maps to be used side-by-side with other applications, since it ensures that the FOV is the same in both applications. See Alignment with Other Applications.
- Magnification Alignment is related to Align Layer (see Layer Context Menu) and Global Alignment functions (see Toolbar), but serves a different purpose.
 - Global Alignment realigns existing data to the coordinate system of Maps. This functionality is typically used for imported data.
 - Magnification Alignment targets misalignments on the instrument itself.
- Magnification Alignment is a mechanism to compensate for imperfections in the optical alignment of the microscope. It is global in nature and applies to all positions. Small position-dependent shifts might remain due to other factors.

• Performing a Magnification Alignment for a magnification changes the current alignment state. That means that tile sets that have already been acquired before the adjustment were using a different Magnification Alignment. If you would re-acquire such tile sets, the images may be shifted due to the new alignment. This situation is detected and he system displays the following warning:



Alignment with Other Applications

- The FOV visualization automatically incorporates image beam shift set by other applications, such as Thermo Scientific Tomography. This allows Maps to be used side-by-side with other applications, since it ensures that the FOV is the same in both applications.
- Image beam shift is a mechanism that compensates for imperfections in the optical alignment of the microscope. The image beam shift changes the current alignment state. This means that tile sets that were already acquired before changing the image beam shift used a different magnification alignment. If you reacquire these tile sets, the images might shift due to the new alignment. The new alignment is not detected automatically, and no warning is displayed.

Note: With some applications, such as Thermo Scientific Tomography, the image beam can shift regularly.
Focus Method

Use Interpolated Focus Method

Focus M	lethod	Interpolated	~	
Intensity	y Lens	0.5560 %		
Objectiv	e Lens	79.5483 %		
Stage Z			Get	
- Interp	polated Focus			
	Set Focus 1	Set Focus 2	Set Focus 3	

This topic explains the procedure for the interpolated focus method.

- 1. Navigate to the first location.
- 2. Manually focus the image.
- 3. When the image is in focus, click **Set Focus 1**.

The current focus will be remembered for that point and the button is now selected.

4. Repeat for the other two **Set Focus** buttons.

Once all three are checked, tiles are ready to be acquired.

Note: These points will be used to define a focus plane. All tiles will use a focus as defined by this plane. For best results, choose points at opposite ends of the sample.

To unset a focus point, click Delete Focus for that point.

STEM Acquisition with Multiple Detectors

This topic describes how to set up tile set STEM acquisition with multiple STEM detectors.

Create STEM acquisition with multiple detectors

Do the following to set up STEM acquisition with multiple detectors, do the following.

1. Define a new STEM tile set (see Define a Tile Set for TEM).

2. Click the Tile Set tab, and then select the checkbox of more than one detector within the **Selected Detectors** section. Detector choices depend on which detectors are available on the system.



3. Select a default detector from the **Default Detector** dropdown menu. The default detector is used for operations that require only one detector (for example, autofunctions and stitching). The acquired images from the default detector automatically display in the main Viewer.

4. Click the Advanced tab and locate the Detectors section. Edit the values of the **Gain** and **Offset** fields as desired (see Advanced group, STEM mode). You can only edit the parameters of the detectors that you previously selected for acquisition. Unselected detectors are deactivated.

🌐 🕂 🕂] ≋ Ø)				
Optics						
Spot Size		5 🗸				
Camera Length	2	205 mm				
Detectors						
	Gain	Offset				
BF-S	31.25%	67.2%				
DF-S	31.25%	67.2%				
HAADF	24.22%	43.71%				
Correction						
Use image corrections						
Stage						
Settling Time		0 s				
FROM MICROSCOPE	то	MICROSCOPE				

Result of multiple detector acquisition

The result of STEM acquisition with multiple detectors is a multichannel tile set in which each channel corresponds to each of the selected detectors. The channel with the default detector is always displays as the first channel and its acquired image is automatically displayed in the main Viewer.



To view images from other detectors, deselect the tile set channels above the selected tile set channel in the layer tree. In the example below, you can view the BF-S image because this tile set channel is selected and the HAADF tile set channel above it is deselected.



STEM Intermediate Images

This topic describes the intermediate images during STEM AutoFocus and the STEM Eucentric autofunction. Intermediate images are visual observations that display in the right-side panel of the Main Screen, which appears when autofunction starts and then disappears when autofunction stops.

STEM Autofocus

You can observe the intermediate images acquired during the AutoFocus process to determine if the parameters used for AutoFocus are acceptable. The same feature can be found also in STEM Tomography software.

The AutoFocus image contains the defocus value for the image. A graph of the contrast quality appears directly below the intermediate image screen, and the contrast quality increases and then decreases within the range of different defocus values. When the graph reaches the maximum value for quality along the Y axis, then the image is focused.



Quality is defined as a unitless number that is based on the measured contrast of the image.



STEM Eucentric Autofunction

You can observe intermediate images captured at different angles, typically between -15° and 15°, during the Eucentric autofunction measurement. You can also change the angle value by configuring the Final Stage Angle of the Eucentric autofunction (see Configure autofunctions).

A cross correlation image displays under the positive and negative images. The cross correlation image contains a crosshair cursor at the center of the image, as well as a circle marker at the position of the highest cross correlation peak.



Automated Image Filtering

This topic describes automated image filtering, which you can use to improve image shift calculation associated with TEM and EFTEM autofunctions.

Image shift calculation

TEM and EFTEM autofunctions are often implemented by calculating the shift between two images. For example, the Eucentric Height autofunction determines the eucentric height by minimizing shift between positive and negative tilt images. The image shift calculation typically uses the cross-correlation of the Fourier transform for the images. Depending on the specimen and the required imaging conditions, there are multiple factors that can decrease the reliability and accuracy of the image shift calculation.

- Recurring patterns in the specimen can cause false peaks in the cross-correlation at various shifts. The peaks are related to the pattern rather than a distinct feature in the specimen.
- The signal-to-noise ratio of the images influences the quality of the cross-correlation. Especially in low-dose conditions, the signal counts can be so low, that recurring patterns and noise in the dark and gain reference images can be interpreted as features of the specimen. Because the reference images are identical in all acquired images, the crosscorrelation image may show a peak at zero shift.
- A Fourier transform is based on the assumption that the image can be extended periodically in all directions. In practice, however, the image is finite and has edges. This results in artifacts in the Fourier transform, such as horizontal and vertical lines (for example, when the image is dark on one edge and bright on the opposite edge). These artifacts can have a negative impact on the accuracy of the image shift calculation.

To overcome these factors, you can use automated image filtering to do the following:

- Improve the sharpness and relative height of the peak in the cross-correlation by enhancing features of a specific size range.
- Reduce artifacts introduced by the edges of the image.
- Decrease or eliminate the peak at zero shift.

Automated image filtering

With automated image filtering, the Maps software application selects the following set of filters and tunes them automatically by using recommended default values.

• Band Pass filter: Enhances features of size range between 1/200 of FOV width and 1/20 of FOV width.

For example, when running automated eucentric height at 7500X magnification with FOV width of 4 μm , enhances features between 20 nm and 200 nm in size.

- Hanning Window filter: Reduces artifacts introduced by the edges of the image.
- Suppress Zero Peak filter: Decreases or eliminates the peak at zero shift.

Configure automated image filtering

You can configure automated image filtering as the default setting *for all new tile sets* through the application settings. If desired, you can also override the default setting for all new tiles sets and configure the automated image filtering *for a specific tile set* through the tile set tab.

To configure the default setting in TEM mode (or EFTEM mode) for all new tile sets.

Settings				×
Application	EFTEM S	ятем тем		
Microscope Acquisition	Snapshots Camera: Exposure Time:	Current 1s		
Data Exchange	Resolution:	Maximum		
XPS	Default Autofunctions Eucentric Height Templ Focus Template:	i Templates late:	None (TEM) None (TEM)	ž
Marjus	Eucentric Height Defa	ult Settings		
Amira-Avizo2D	Final Stage Angle*	15.0	🖉 🔤 Use Image Filtering	
Слую	Max. 2-Height Deviatio	n 250 nm		
Stitching	Focus Default Settings Desired Defocus Iterate to	0μm -5μm	Use Auto Stigmate Use Three Image Method	
	Focus Method	Objective Lens	V 🔲 🔤 Use Image Filtering	
			Apply OK	Cancel

- Click the TEM Options menu >> Settings... menu >> TEM tab (or EFTEM tab) >> Acquisition tab.
- 2. Under the Eucentric Height Default Settings section, click the **Use Image Filtering** check box.
- 3. Under the Focus Default Settings section, click the (Beta) Use Image Filtering check box.
- 4. Click OK.

To override the default setting and configure the automated image filtering *for a specific tile set* through the tile set tab (TEM and EFTEM modes).



- 1. Under the tile set tab, click the Autofunction tab.
- 2. Click **Configuration...** to open the Advanced Autofunction Settings (Tile Set) screen.
- 3. Click the Eucentric Height tab within the Advanced Settings (Tile Set) screen.
- 4. Click the (Beta) Use Image Filtering check box.
- 5. Click the Focus tab within the Advanced Autofunction Settings (Tile Set) screen.
- 6. Click the (Beta) Use Image Filtering check box.
- 7. Click OK.

TEM Analytics

With the Analytics plug-in, the Maps software can acquire EDS data through a software application interface with Thermo Scientific Velox[™].

Note: This plug-in is not be confused with Maps Min or Generic Detector Interface (GDI).

Preparation for TEM Analytics

Preconditions

To run Analytics, your Maps installation must include the following components:

- At least one software package from a provider for EDS acquisition
- An Analytics plug-in with the correct license

To verify that your license includes the Analytics plug-in option, on the Help menu, click About.

thermoscientific	Maps 3.	20
	Product XPS Plugin Anlytics Plugin Avizo Bridge Plugin IQM Plugin Offline Plugin Cryo'temPrep Plugin	Version 1.000 3.200.2111 3.200.2111 3.200.2111 3.200.2111 3.200.2111
	Licensed Options Color Correlation Project Editing External Image Import Stage Correlation Stage Correlation TEM Online	Expiration Date Unlimited Unlimited Unlimited Unlimited Unlimited Unlimited
	Josennay TEM Online Analysics	Unimited Unimited

Analytics Workflow for TEM

The Analytics workflow explains how to set up and activate the Velox software on your TEM system before you define a tile set. It also explains how to configure Analytics, acquire a tile set, and view the results of the acquisition.

Proceed with the workflow in the following order:

- 1. Set Up Velox Software: Before you acquire EDS images in Maps with Velox, you must set up the Velox software.
- 2. Activate Velox Software: To perform the initial configuration for Analytics, you must activate Velox in Maps.
- 3. Define a Tile Set: Use the normal procedure to define a tile set, and then navigate to the tile set section and click the **Analytics** tab.
- 4. Configure Analytics for Acquisition: Configuration depends upon which EDS acquisition software you have, either Velox or Bruker. Follow the instructions for your specific EDS acquisition software.

- 5. Acquire a Tile Set: Use the normal procedure to acquire an Analytics tile set.
- 6. View Results of Acquisition: Once an Analytics tile set has finished acquisition, you can view the results in the layer tree and the Results pane.

Set Up Velox Software

Before you acquire EDS images in Maps with Velox, you must set up the Velox software.

1. Launch the Velox software application to the Acquisition and Processing screens.



2. In the Acquisition screen, under the Optics pane, click **STEM**.

Ve Acquisi	tion - Velox										
File Edit	t View Optics	DPC H	lelp								
₽ ₽ ₽		r1	μ.	🕹 TEM	HAADF	DF4	¥		-	+	512 x 512 🔻
₹ <u></u>	*	ij	· (STEM	DF2	BF		O		+	100 ns 🔻 🐃
Layout 🔻	Annotations 🔻			Optics	Det	ectors		S	TEM Ima	ging 🔻	

3. Within the Detectors pane, click HAADF.

Ve Acquisi	ition - Velox										_ []	×
File Edit	t View Optics	DPC F	Help									
	× .	r1	<u> </u>	🕹 TEM 🄇	O HAADF	DF4	*		=	- +	512 x 512 👻	
€∰	×	ij		STEM	DF2	BF		_►	<u>ہ</u>	- +	100 ns 🔻	»
Layout 🔻	Annotations 🔻			Optics	Det	tectors			S	TEM Imaging	17	

4. Within the STEM Imaging pane, click the Scanning button to pause.

Ve Acquisition - Velox										
File Edit	t View Optics	DPC I	Help							
E D.₽.	×.	r1	<u> </u>	🕹 TEM	O HAADF	😳 DF4	* ≡ :	_	- +	512 x 512 🔻
₹∰		ij		E STEM	DF2	O BF		<u> </u>	- +	100 ns 🔻 »
Layout 🔻	Annotations •			Optics	Dete	ectors		STE	M Imaging	ı •

5. Within the STEM Imaging pane, change the resolution to match the tile set resolution in the Maps application.

Ve Acquisition - Velox												
File Edit	t View Optics	DPC H	lelp									
		r1	<u> </u>	📥 TEM	O HAADF	DF4	*	=		-	+	512 x 512 👻
₹ <u></u>		ij	(TT)	STEM	DF2	BF		►	0		+	100 ns 🔻 »
Layout 🔻	Annotations •			Optics	Dete	ectors			ST	TEM Ima	iging 🔻	

6. Within the SI pane expander, select Auto stop.



7. In the Processing screen, click to the elements of interest under the Periodic Table (Spectrum Imaging) pane.



8. Go to Activate Velox Software.

Activate Velox Software

To perform the initial configuration for Analytics, you must activate Velox in Maps.

- 1. Under the **Options** menu, click **Settings**.
- 2. Click the Velox tab, and then select the Use Velox check box.
- 3. Click OK.

Ma Settings		×
Application	EDS Detector Provider	
Microscope	Image Settings	
Acquisition	Default Additive Opacity:	
Data Exchange		
XPS		
Amira-Avizo2D		
Velox		
Сгуо		
Stitching		
	Apply OK	Cancel

4. Restart the Maps software application.

5. After restarting Maps, a green status icon appears next to the Velox label on the status bar to confirm success.



Configure Analytics for Acquisition

Most Analytics details are set up in the software packages from the EDS acquisition providers. If you are using Velox software for EDS acquisition, then follow the setup instructions below. If you are using Bruker software for EDS acquisition, then follow the instructions under Set up Bruker Analytics.

Set up Velox Analytics

1. Select the EDS Acquisition check box.



2. If the Velox software did not provide the required fields, then Maps shows a warning icon and activates the Resolution dropdown menu. Select the same value here that you selected when you set up the Velox software.



3. Select 1-4 quantification modes as desired.



 Note: The tile set resolution must match the resolution of the image returned from the Velox application, so verify they are the same *before* acquisition begins. For example, resolution width: 512, and resolution height: 512.

The tile set is now set up and ready for acquisition. If you change parameters on the tile set (specifically anything related to resolution), then you might need to run the setup again. See Configure Analytics for Acquisition.

Set up Bruker Analytics

Follow the procedure in the Configure Mineralogy Acquisition section.

View Results of Acquisition

Once an Analytics tile set has finished acquisition, you can view the results in the layer tree and the Results pane.

Layer tree

In the layer tree, the tile set has a collection of layers within it that corresponds to each element and STEM image.



- Click the colored circle on the right to change the color of the element layer.
- Select the check box on the left side of the layer to toggle visibility.

Results tab

The Results tab of the tile set shows the list of channels for the tile set, which corresponds to each of element and STEM image. You can expand each any of these channel controls by clicking in the channel header.



In expanded control, you can modify the channel and view the metadata collection of the given element and STEM image. An example of an expanded channel follows.

Opacity Mdditive Opacity				•	•
Stitching Channel			Stem_D	F-S	~
			sha Sha	ow histogra	ım
int	net	wt	%	at%	
–					
🗸 🎽 Stem_DF-S					
Opacity					•
Gamma	- i		·		Reset
Saturation	Additive				Add Color
Ť					
ImageWidth	512 px				
ImageHeight	512 px				
ImageName	Image 1				
Detector	DF-S				
Optics					
PixelSizeX	0.02329308	28861061			
PixelSizeY	0.02329308	28861061			
Dwell	2000 ns				
Frames					
StemMagnification	8600				

Sample Holders

This section describes the sample holders that are available for Maps.

When a holder is selected and applied to the project:

- The outline displays in green on the viewer. The holder outline is red until it is aligned.
- The name, drawing, version, and so on are displayed on the Holder tab.
- A new Layer is added to the Layer Control. This is where you choose to show or hide the Holder outline and Fiducials. The Fiducials only show after image acquisition. The default is to not show them. Select the check box to view your fiducial images.

Refer to your application specialist for procedural information for each sample holder type.

Holder Tab

When a holder is selected and applied to the project:

- The outline displays in red on the viewer if automated holder alignment is supported. If automated alignment is not supported, then the outline displays in green.
- The name, drawing, version, and so on are displayed on the Holder tab.
- A new Layer is added to the Layer Control. This is where you choose to show or hide the Holder outline and Fiducials. The Fiducials only show after image acquisition. The default is to not show them. Enable the check box to view your fiducial images.

• The Left , Middle, and Right quick access buttons for the sample area appear adjacent to **Right Slide**.

✓ ✓ Single 22 mm coverslip holder	
🗾 ⊡ Holder outline	
Holder:	Single 22 mm cov
Right Slide Version:	1.0
Manufacturer:	FEI
Part number:	1101666
Opacity	
Sample areas: Template Factory	y Template 🗸 🗸
Right Slide	;;;;;;;;;;;;;;;;;;;;;;;;;;;;;;;;;;;;;;
Saved holder alignment	Save Clear
Saved acquisition parameters	Reset

- Left: Centers the view on the sample area
- Middle: Drives the stage to the center of the sample area
- Right: Creates a new tile set that covers the sample area

Holders Supported for Automated Holder Alignment

When selected, the holder outline displays in red on the viewer, if automated holder alignment is supported. See Automated Holder Alignment.

Topics include:

3 mm Grid Holder

Single 22 mm Coverslip Holder

Double 22 mm Coverslip Holder

3 mm Grid Holder

This is a holder for 3 mm diameter grids that can be loaded inside a transmission electron microscope (TEM).

3 mm Grid Holderwith Upper and Lower Fiducials Marked



Single 22 mm Coverslip Holder

This is a single 22 mm coverslip holder.



Double 22 mm Coverslip Holder

This carrier holds up to two samples.



Holders Supported for Graphical Overlay

Maps can draw a graphical overlay to assist with navigation. The graphical overlay can be aligned manually to other image layers, but the automated holder alignment is not supported on these layers.

Note: When selected, the holder outline displays in green on the viewer, indicating that automated holder alignment is not supported.

Topics include:

Cryo Corrsight Shuttle Holder Cryo AquilosSEM Shuttle Holder Slide 75 mm x 25 mm Holder 12xthin-section Quanta Mount Holder 14 x 30 mm Round Quanta Mount Holder

Cryo Corrsight Shuttle Holder

This holder can be connected to a cooling system for imaging samples at cryogenic temperatures.



Cryo AquilosSEM Shuttle Holder

This holder is used for Cryo applications. It can be connected to a cooling system for imaging and sample preparation at cryogenic temperatures.



Slide 75 mm x 25 mm Holder

Holder project 🕂 🛶 📐 Q | 🎬 🏶 🐵 🖻 🖻 🕄 🖂 💥 🗗 🔍 🖻 ت ش 🤇 Setting TEMPLATE Factory Template 🔽 Layer 🗾 🗔 Holder outline 🖸 🧭 Holder: Slide 75mm x 25n Version: Manufacture Part number Opacity • Template Factory Template ~ Saved holder align Sa ed acquisition p 20

This holder represents a slide with dimensions 75 mm x 25 mm.

12xthin-section Quanta Mount Holder

This mount holder is designed to handle up to 12 thin mineralogy samples. The mount holder includes a Faraday cup; meets gold, quartz, and copper standards; and has a magnesium calcium position.



14 x 30 mm Round Quanta Mount Holder

This mount holder is designed to handle up to 14 samples of 30-mm resin embedded mineralogy samples. The mount holder includes a Faraday cup; meets gold, quartz, and copper standards; and has a magnesium calcium position.



Automated Holder Alignment

This section describes the procedure for automated holder alignment (AHA) setup, calibration, and operation for a particular supported holder on SEM/SDB systems. It is a guideline for using Maps and the automated holder alignment, and can also be used for SEM/SDB-only as well as multi-system correlative Maps projects.

i Note: Automated Holder Alignment is not supported for Maps on TEM systems.

Topics include:

Holders Supported for Automated Holder Alignment

Automated Holder Alignment User Interface

Execute Holder Alignment

SEM/SDB Holder Calibration

Correlating Projects

Problems with Alignment

Holders Supported for Automated Holder Alignment

When selected, the holder outline displays in red on the viewer. They appear in green if automated holder alignment is *not* supported. See Sample Holders.

- 3 mm Grid Holder
- Double 22 mm Coverslip Holder

All supported holders must be manually calibrated on a given system before Automated Holder Alignment is possible. This is a one-time procedure. After manual calibration is complete, little to no manual action will be required.

Automated Holder Alignment User Interface

Click the **Align to Holder** icon in the tool box and select the appropriate holder from the **Sample Holder** dropdown menu to cause the Automated Holder Alignment user interface to appear next to the viewer.

Automated Holder Alignment UI



Automated Holder Alignment UI Overview

Interface Item	Description
Buttons at Bottom (Before Starting Alignment):	Run Alignment Retry Apply Close
Run Alignment	Moves the stage to the position where it expects to find the fiducial, auto- focuses, and attempts to match the fiducial.

Automated Holder Alignment UI Overview(Continued)	
Interface Item	Description
Apply	Applies the alignment to the project if the alignment is not immediately applied when the automated procedure completes. This can be used if the automated alignment completes but manual adjustment is needed.
Close	Closes the alignment UI and returns to the main viewer before the automated procedure has begun.
Buttons at Bottom (During Alignment):	Continue Retry Skip Abort
Continue	On SEM systems, the automated holder alignment requires manual optimization of the image (contrast, brightness, focus, and so on). After the stage is automatically moved to the location of the first fiducial and the prompt for image optimization has appeared, use Continue to proceed with fiducial matching.
Retry	Autofocuses and attempts to match the fiducial. If the match is successful, the automated alignment procedure will continue to the next fiducial.
Skip	Skips the current fiducial matching procedure and proceeds to the next fiducial. This can be useful if Maps encounters difficulty matching, however, the best practice is to optimize the image for successful automatic matching or to manually select the fiducial position.
Abort	Cancels the alignment procedure currently in progress and returns to the main viewer.
	Automated Holder Alignment UI Overview(Continued)
-------------------------------	---
Interface Item	Description
Fiducial Window	Vpper left Reference: Initial: N/A Acquire
Green/Red Ball	Indicates if the fiducial was found: Green = Found; Red = Not Found.
Fiducial Name	Displays the name of the fiducial. In this example, Upper Left.
Reference	Displays the reference image.
Initial	Displays the result of the previous alignment procedure for a project that has already been automatically aligned.
Current Fiducial Window	Displays the fiducial image acquired during the current automated holder alignment procedure.
Acquire	Acquires an image of the specified fiducial. Maps will move to the expected location for the specified fiducial and attempt to match.

	Automated Holder Alignment UI Overview(Continued)
Interface Item	Description
Fiducial Markers for	The fiducial marker is placed on the fiducial image as a point of reference. Maps will align these reference points during alignment.
Matching	• Light Blue markers are used in the reference images to display the reference point that the automated procedure expects to find. Light blue markers are also automatically placed on the fiducial image when a match is successful.
	 Light green markers denote the reference points from the previous alignment procedure and appear in the initial fiducial images.

Execute Holder Alignment

When using a sample holder that has already been calibrated on your microscope, the following procedure describes how to execute the alignment sequence.

1. Click Align to Holder icon.



2. Select the same holder that was just calibrated, in this case **Double 22 mm coverslip holder**, and then click **OK**.

Ma Holder Selection			\times
A sample holder n Please select a san	nust be set in the project nple holder from the list b	before continuing. pelow.	
Sample holder: Double 22 mm coverslip hold	er		~
Left SlideRight	Holder:	Double 22 mm coverslip holder	
	J Version:	1.0	
	Manufacturer:	FEI	
	Part number:	1043363	
		OK Cancel	

3. Click Run Alignment. The Alignment proceeds automatically.

We	orkflow: Holder Alignment		×
Find locations of t located for the ali	the holder fiduciols. At least 2 fiducia gnment.	ls must be	
Upper left Reference:	N/A Acquire		
Upper left		N/A	
Bottom left		N/A	
Upper right		N/A	
	Run Alignment Retry Apply Close		
	Fiducial locations		~

- If the automated alignment is able to match all fiducials successfully, the holder calibration is complete.
- If the alignment failed to recognize any of the fiducials, then perform the calibration procedure (see SEM/SDB Holder Calibration

SEM/SDB Holder Calibration

This topic explains the procedure for calibrating a SEM/SDB holder.

Mount the Coverslip Holder and Adaptor to the Stage

Follow this procedure to mount the Double 22 mm coverslip holder and adaptor to the stage.

Top View of the Holder



The holder only fits in the adaptor in one position.

1. Withdraw the clamp and place the holder with the cut off side facing the clamp in the holder. The holding device (encircled in red) fits the hole (encircled in red) in the adaptor.

Bottom View of the Holder



6 inch Stage Adaptor for Double 22 mm Coverslip Holder



2. If the holder is positioned, then release the clamp.

Holder Positioned on Adaptor



3. Place the holder and adaptor on the stage. This can only be done in one orientation. The adaptor has three screws (encircled in red) and two pins (encircled in green) that fit nicely into the stage.

Adaptor Base with Screw and Pin Positions



Stage Screw and Pin Placement Indicators



4. Tighten the screws so the stage cannot move.

- 5. Set the Z coordinate to 0 using the xT UI to make sure the holder will not touch the pole piece when inserting.
- 6. Close the chamber.
- 7. Press Pump.

Note: Refer to your application specialist for procedural information for each sample holder type. See Sample Holders for descriptions of the other types of holders.

SEM/SDB Preconditions

Before performing calibration on a Maps supported holder using an SEM/SDB system, verify that the following preconditions are satisfied:

- Home stage
- SEM mode: lowest magnification
- ETD
- BSE
- 10 kV
- 100 pA (Use higher beam current of about 1 nA if you have difficulty matching automatically)
- Eucentric
- Link stage to WD
- Optimize contrast such that the fiducial holes are black and the background is white

Find the Fiducials and Store their Coordinates

This procedure is only necessary when starting Maps for the first time on a particular SEM/SDB, using a particular holder.

Using the xT UI, find the upper right (UR), upper left (UL), and bottom left (BL) fiducials and store the coordinates on the Stage tab.



The fiducial orientation must match the reference image.

The fiducial next to the chamfer/cut off (encircled in red) is called upper right.

Upper Right Fiducial



The upper right fiducial should be present at coordinates:

- X: 9.5 mm
- Y: -35.0 mm

The upper left (UL) fiducial should be present at coordinates:

- X: 9.5 mm
- Y: 35.0 mm

The bottom left (BL) fiducial should be present at coordinates:

- X: -9.5 mm
- Y: 35.0 mm

i Note: Depending on the system type, X and Y coordinates might be reversed.

- If you are entirely unable to find the fiducials on the holder, perform a large overview tile for help in navigation.
- If the system is equipped with a NavCam, use the NavCam navigation image to assist with finding the fiducials.

Calibrate the Holder

1. Load the appropriate sample holder. In this example, it is Double 22 mm coverslip holder.

Note: Refer to your application specialist for procedural information for each sample holder type. See Sample Holders or descriptions of the other types of holders.

- 2. Create a new project or open a project that uses the loaded holder type. See Create a Project for more information about creating projects.
- 3. Click the Align to Holder tool box button.



4. Select the appropriate holder from the dropdown and then click OK.

Ma Holder Sele	ction			×
	A sample holder me Please select a sam	ust be set in the project ple holder from the list b	before continuing. below.	
Sample holder:	Double 22 mm coverslip holde	er		~
	_Left SlideRight \$	Holder:	Double 22 mm coverslip holder	
γ		Version:	1.0	
	։ լ լլ	Manufacturer:	FEI	
+		Part number:	1043363	
			0//	
			OK Cance	



5. In the fiducial alignment controls that appear, click **Run Alignment**.

The stage drives to the position where it expects to find the upper left fiducial. Maps displays a prompt to adjust microscope settings for optimal fiducial image.

6. Click Close.



- 7. Go to the xT UI and find the stored position for upper left fiducial. Select the upper left coordinates and then click **GoTo**. The stage navigates to the stored coordinates of the upper left fiducial.
- 8. Optimize the image (set correct focus and contrast/ brightness) of the upper left fiducial using the xT UI. The optimal image will have completely black fiducial holes against a completely white background.

9. When the image has been optimized, return to the Maps UI and click Continue.

i Note: If you are having trouble matching, make sure the:

- Contrast / brightness matches reference
- Fiducial orientation matches reference
- HFW is 1.5 mm
- Sample is clean with clearly distinguishable features and no dust / contamination



During calibration, Maps does not know the orientation of the holder. This can cause false positive matches, as shown in the diagram below. This is normal and these false positives will not occur after the holder has been successfully calibrated.



10. The Microscope Settings box appears again if Maps still does not recognize the acquired fiducial. Click **Close**.



11. Optimize the fiducial image again to match the reference image and then click Retry.





12. If Maps is not able to automatically match the fiducial, double-click to place the blue cross hair on the same fiducial hole that contains the red crosshair in the Reference image.

The red button turns green if Maps recognizes the fiducial.

🛑 Upper left 🔵 Upper left

If Maps does not automatically navigate to the second fiducial, then click Continue.

- 13. Maps automatically attempts to match the second fiducial. If the match is not successful, then repeat the calibration for the second fiducial. Once the second fiducial is recognized, Maps will automatically navigate to the final fiducial (for holders with more than two fiducials).
- 14. Since the automated alignment required manual adjustment, the Calibration prompt appears. Click **Yes** to save the holder calibration.



• Note: If the calibration prompt does not appear, Maps has low confidence in the calibrated fiducial positions and will not allow you to save the potentially bad calibration values. This can occur if there is a large unexpected rotation between fiducials. Perform the calibration a second time, and make sure fiducials are well centered and are oriented properly to match the reference.

Verify Alignment

- 1. On the File menu, click Project Properties.
- 2. In the Sample holder list, click none.

none	~
none	
3mm Grid holder	
8 Block Holder	
14x30mm Round Quanta Mount	
Corrsight Shuttle	
Aquilos Shuttle	
Leica Shuttle	
Single 22 mm coverslip holder	
12xthin-section Quanta Mount	
Slide 75mm x 25mm	
Double 22 mm coverslip holder	

3. Click **OK** to clear the holder from the project. Click **OK** again at the prompt.

Ma Project properties				×
Project Name:				
Demo project				
Project Path:				
C:\Maps Projects\Demo project				
Disk Space Usage:				
281.25 MB (294,914,519 bytes)				
Total Number of Images:				
25 tiles, 2 imported images				
Project Description:				
Sample Holder:				
none				~
	Holder:			
	Version:			
	Manufacturer:			
	Part number:			
		Ok	Apply	Cancel

4. Proceed to Execute Holder Alignment.

Correlating Projects

Prerequisites for running a multi-tool correlative experiment with Maps using Automated Holder Alignment:

- Sample is mounted on a supported sample holder
- Holder to be used has been calibrated on all systems

i Note: Manual correlation can be performed on non-supported sample holders using the Maps Global Alignment procedure.

To perform a correlative project:

- 1. Create a new project in Maps on the first system.
- 2. Run the AHA (Automated Holder Alignment).
- 3. Acquire the desired sample data.
- 4. Close Maps.
- 5. Make a copy of the entire project directory and move it to the second system's PC.
- 6. Install the supported holder into the second system.
- 7. Open Maps on the second system and Import the project.
- 8. Run the AHA.
- 9. Acquire the desired sample data.
- 10. Use alignment to adjust the position of sample data, as desired.
- 11. Export the results by stitching or saving images to file.

Problems with Alignment

Misalignment between data in a Maps project can be due to one or more of the following contributing factors:

- Stage positioning error
- Stage drift
- Modified scanning parameters
- Modified column parameters
- Microscope modality change
- Sample moved relative to holder
- Sample changed physically during sample prep
- Misalignment between imaging modes/microscope objectives
- Improper holder calibration or false positives during Automated Holder Alignment

Manual Alignment

This section describes manual alignment workflows.

Alignment is used in Maps to adjust project data to the current stage position. For example, this is useful after a sample has been removed and re-installed into a system in a different orientation.

Alignment can be performed on imported images, tile sets, stitched tile set data, layers, and entire projects (Global Alignment). The procedure is identical when aligning each of these different objects.

It is worth noting that objects are never aligned to each other in Maps, but instead are aligned to the current stage position. If two objects do not match up, align the first to the stage, then align the second to the stage. To maintain alignment between objects, place all the aligned objects into a separate layer. Then, you can perform alignment on the Layer and all objects in the layer will be moved in unison.

For Offline alignment where there is no stage present, simply choose one image or tile set as a standard and align all other data to this standard. When the project is reopened on a system, a single Global Alignment will correctly adjust all the project data.

There are four levels of alignment:

- Align: Aligns a single tile set or image. Selected from the Tile Set Context Menu.
- Align Layer: Aligns an entire layer, including all tile sets and images at once. Selected from the Layer Context Menu.
- **Global Alignment**: Performs an alignment to align all previously acquired data to the stage. Selected from the Toolbar.
- **Magnification Alignment**: Performs an alignment to align TEM-acquired layers based on magnification. Selected from the Layer Context Menu.

Topics include:

Coarse Alignment (1-, 2-, or 3-Point)

Fine Alignment

Coarse Alignment (1-, 2-, or 3-Point)

Use 1-, 2-, or 3-point alignment to match up a previously acquired image with the current stage position.

- 1 point: Adjusts the data using translation only.
- 2 point: Adjusts the data using translation, rotation, and scale.
- 3 point: Adjusts the data using translation, rotation, scale, and orthogonality.

Workflow

The workflow described below relates to aligning a single tile set. The same procedure is used for all alignments in Maps.

Coarse Alignment from Tile Set

- 1. Acquire a tile set as described in Acquire a Tile Set.
- 2. Right-click on the tile set in the Layer control and click **Alignment > Align**.

	Bring Forward				
	Send Back				
	Send To Back				
~	Show Grid Lines	Ctrl+G			
	Center View	Ctrl+Shift+C			
	Center and Rotate View				
	Save and Set As Default Template				
	Save As Template				
	Copy Grid				
	Export Grid				
	Evenet To Designt				
	Export to Project				
	Alignment	•		Align	
	Alignment Drive To	×		Align Fine Alignment	
	Alignment Drive To Stitch	•	5	Align Fine Alignment Magnification Alignment	
	Alignment Drive To Stitch Tile Registration		Ē	Align Fine Alignment Magnification Alignment Clear Alignment	
	Alignment Drive To Stitch Tile Registration Split Channels		ē	Align Fine Alignment Magnification Alignment Clear Alignment	
	Alignment Drive To Stitch Tile Registration Split Channels Create MIP Layer		E	Align Fine Alignment Magnification Alignment Clear Alignment	
	Alignment Drive To Stitch Tile Registration Split Channels Create MIP Layer Mark as done	•	E	Align Fine Alignment Magnification Alignment Clear Alignment	
	Alignment Drive To Stitch Tile Registration Split Channels Create MIP Layer Mark as done Reset Acquisition		Ð	Align Fine Alignment Magnification Alignment Clear Alignment	
	Alignment Drive To Stitch Tile Registration Split Channels Create MIP Layer Mark as done Reset Acquisition Open Z-Stack	•	đ	Align Fine Alignment Magnification Alignment Clear Alignment	
	Alignment Drive To Stitch Tile Registration Split Channels Create MIP Layer Mark as done Reset Acquisition Open Z-Stack Recipe Processing		đ	Align Fine Alignment Magnification Alignment Clear Alignment	

Coarse Alignment from Workflow

1. Click the workflow button, and then click **Alignment Wizard** and **Begin**. The workflow panel opens and lists layers for alignment.



2. Select an acquired tile set to align.

Workflow: Alignment Wizard	×
EMTileSet 3	
EMTileSet 2	
EMTileSet	
orrsight	
40x_SDgb_01	
LA Tile Set 2	
40x_TLC_01	
LA Tile Set	
M - STEM	
STEM tileset 2	
✓STEM tileset	
TEM tileset 2	
TEM tileset	
••••	
Select Layers	>

3. Click the right arrow button at the bottom of the workflow panel to go to the next step.

4. Select a Tile Set to which you want to align to (such as the Tile Set as acquired in step 1).

		Workflow: Alignment Wizard	×
Sele	ct which laye	rs to which needs to be aligned by checking them.	
	All	Clear	
	- Sam	nle Array	
	- Dam	Tile Set	
		Tile Set 3	
		Tile Set_2	
		Tile Set_1	
	-	Sample Array_Frames	
		Sample Array_Frames_Frame 3	
		Sample Array_Frames_Frame 2	
		Sample Array_Frames_Frame 1	
	- SEM		
		EMTileSet 3	
		EMTileSet 2	
		EMTileSet	
	Corre	sight	
		40x_SDgb_01	
		LA Tile Set 2	
	<u>~</u>	40x_TLC_01	
		STEM tileset 2	
		STEW theset 2	
<		Select Layers to align to	>

- Click the right arrow button at the bottom of the workflow panel to go to the next step.
 The Tile Set that was selected in step 1 appears in the workflow panel.
 The Tile Set that was selected in step 2 appears in the Viewer.
- 6. Follow the instructions in the workflow panel.



7. Follow the guided instructions in the workflow panel to find the same distinct feature in both windows and place Point 1, 2, or 3 in each window. You might have to zoom in close on a feature and grab a new preview image with more resolution to make the best match.



8. (Optional) Using the **Mirror Stage Axis** or **Mirror Image** controls below the viewer in the workflow panel, you can mirror the image horizontally or vertically. Use **Swap Stage Axes** to swap the X and Y stage positions. It is common to use combinations of these buttons when aligning optical images. If you select the checkbox for the **Link Rotation** option, the left and right image viewers will rotate together. Deselect the checkbox for the **Link Rotation** option to rotate them independently.



- 9. Select 1, 2, or 3 point alignment.
- 10. Click the right arrow button at the bottom of the workflow panel to enter Fine Alignment.

Fine Alignment

Use Fine Alignment to fine-tune the alignment of layers and tile sets visually.

Note: When you start a second alignment on a layer, the Maps application remembers and reuses the alignment points from the previous alignment.

Fine alignment user interface

The Alignment Wizard appears when you do either of the following:

- Click the right arrow button at the bottom of the workflow panel after a 1-, 2-, or 3-point alignment.
- Right-click the Tile Set in the Layer control and then click Alignment > Fine Alignment.



	Fine Alignment Overview			
Interface Item	Description			
Viewer Window	The work area within the Alignment Wizard. You can drag alignment points into the viewer during the fine alignment (see Fine alignment mouse controls) process. This window is empty when it opens.			
Back Arrow	Returns to the selection of alignment points in the coarse alignment workflow. From there, click Cancel if you want to undo the coarse alignment.			
Finish Check Mark	Closes the Fine Alignment workflow panel.			

Fine alignment mouse controls

During fine alignment, you can:

- Drag an alignment point into the viewer by left-clicking the alignment point with your mouse and holding the mouse button down as you drag the alignment point into the viewer.
- To pan within the viewer, left-click your mouse and then hold down the mouse button as move your mouse up, down, left, or right.
- Rotate the viewer by clicking the viewer and holding down the **Shift** button.
- Add a new alignment point, or reposition an existing alignment point, by positioning your mouse within the viewer and right-clicking your mouse to open the viewer's context menu. Click the menu option associated with your desired task.

Correlation

This section describes the correlation controls you can use to apply simple coloring to the tile set images when they are acquired. This can be used to highlight some features for demonstration or evaluation purposes.

Topics include:

Visualization Tab

Segmentation

Visualization Tab

The Visualization tab (formerly called "Results" tab) contains the correlation controls, which you can use to apply simple coloring to the tile set images when acquired. Some correlation controls appear unavailable if they are not applicable to the system type.

Correlation controls

III - (AE) II 😨	
Opacity	•
Additive Opacity	—— — —
Stitching Channel	Stem_HAADF 🗸
	<u> </u>
int net	wt% at%
🔪 🗹 Stem_HAADF	•
🗲 🔽 Pd (At)	
> Nb (At)	
📏 🔽 Mn (At)	•

Correlation Controls Overview

Control	Description
Opacity	Sets the measure of opaqueness for the entire image. The less opaque, the more transparent it becomes. Transparent images are useful when overlaying on image over another.

Correlation Controls Overview(Continued)	
Description	
Sets the control the opacity between fluorescence and regular channels. This is useful for the Analytics tile sets as they contain many non- overlapping channels. The Additive Opacity option is the default for EDS tile sets. Increasing the additive opacity makes fluorescence channels more visible.	
Sets the channel used for stitching.	
Shows/hides the histogram panel. See Histogram.	
EDS channels acquired with Velox can have up to four different quantification modes. Click one of the activated buttons (shown in white) to select a quantification mode, which is used to visualize the elements. A deactivated button (shown in grey) indicates that the quantification mode was not calculated and is not available.	
 Int net wt% at% Stem_HAADF The elements display data in each quantification mode as follows: int: the counts according to the raw integrated spectrum net: the counts according to the background corrected and fitted model wt%: the weight fractions 	



Histogram

Displays a histogram for the selected channel. Click a different channel name to select its histogram and associated color.

Selected Channel Histogram


Selected Channel Histogram Overview		
Interface Item	Description	
Selected channel name	Displays the selected layer item (a tile set or image) from the Layer control.	
Hide	Hides the histogram.	
Zoom	Expands the histogram area of images with less than 255 colors.	
Semi-log	When selected, changes the histogram intensity display to logarithmic.	
Invert	Inverts the intensity of the image, e.g., black to white.	
Auto	When selected, continuously adjusts the min and max for the histogram display.	
Current max position	Displays the current max slider position.	
Max slider	Any data above the max line is white. Gray levels display between the max and min sliders.	
Histogram data	Displays the histogram data for the selected channel.	
Min slider	Any data below the min line is black. Gray levels display between the max and min sliders.	
Current min position	Displays the current min slider position.	

Segmentation

Image segmentation is the process of partitioning a digital image into sets of pixels, called segments. The goal of segmentation is to simplify and/or change the representation of an image into something that is more meaningful and easier to analyze.

Add Color

In Maps, images and tile sets can be segmented such that each segment is assigned a color. Segments are defined using the image histogram and a min/max slider that determines the regions of the image to be included in the segment. Additionally, Maps automatically calculates the number of pixels within the segment as a percentage of the total number of pixels and displays the result.

Segmentation, One Color





Segmentation, Two Colors

1. Click Add Color to add an additional image segment.

~ 🛛	π	•
Opacity 🖉		•
Gamma		Reset
Satura	ation Fluorescence	Add Color

2. Choose a color from the **Color Selector** and select the **Solid** check box. When not selected, the color will be a gradient. See Segmentation, One Color with Gradient.



Correlation

- 3. Adjust the **Max** and **Min** sliders to change the range of the histogram data to be assigned to the segment.
- 4. Click Add Color again to add another image segment and get segmentation with two colors.
- 5. Adjust the **Max** and **Min** sliders to select a different range of the histogram data to be assigned to the second image segment.

Segmentation, One Color with Gradient

When the **Solid** check box is *not* selected, the histogram fill color is a gradient from light to dark.

Segmentation, One Color with Gradient

Segmentation Controls Overview



Interface Item	Description
Remove	Removes the segment from the image and returns to normal display.
Color Picker	Displays the color picker for selecting a color, then colors the area between the Min and Max sliders with the selected color.
Solid	When selected, the data between min and max on the Segmentation histogram will be one color instead of a gradient.
%	Calculates the number of pixels between min and max as a percentage of the entire image.

Segmentation Controls Overview

Amira-Avizo2D Connectivity

This section describes how to set up the connectivity between Maps and Amira-Avizo2D, so that you can do the following:

- Create a recipe in Amira-Avizo2D from an image sent by Maps
- Define post-processing options in Amira-Avizo2D

Prerequisites

Maps with Amira-Avizo2D is designed to run on a system with resources equal to or greater than the following:

- 8 GB of physical RAM
- Maps must run in offline mode, or with a SEM or TEM microscope
- Analyzer is installed by Amira-Avizo2D product installer (minimum version 2021.1).

Configuration options

There are four configuration options for connecting Amira-Avizo2D to Maps: None, Local, Direct, or NAT. Select the option based on your network setup, and then refer to the corresponding section below for detailed instructions.

- 1. None: Use the None option if connectivity with Amira-Avizo2D is not required.
- 2. Local: Use the Local option if Amira-Avizo2D is installed on the same PC as Maps. This will typically be for Maps Offline installations on a workstation.



- 3. Direct:
 - Use the Direct option if Amira-Avizo2D is installed on a different PC on the same network. This may be applicable if Maps is on the Microscope PC and Amira-Avizo2D is installed on the Support PC, or if Maps Offline is installed on a different PC other than Amira-Avizo2D.



- Also, use the Direct option for microscopes with a G5 network if a managed switch is used instead of a Support PC.
- 4. **NAT**: Use the NAT option if Amira-Avizo2D is installed on a different PC on the customer network and the Microscope PC is only connected to the Support PC on the private microscope network.



Local setup

- 1. From the Maps PC, navigate to **Maps Options>>Settings... menu item** to open the Amira-Avizo2D settings.
- 2. Click the Amira-Avizo2D tab.
- 3. Select Local from the Connection Type field.
- 4. Configure the shared folder path as indicated in the table below.

Amira-Avizo2D Local Settings	
Interface Element	Description
Maps - Amira- Avizo2D folder	Enter the path to the shared folder for the image and recipe data. See Configure shared folder for details.
Connection time out	Determines the duration time required for Maps to connect with Amira- Avizo2D before Maps cancels the connection. The duration time is measured in seconds.

5. From the same PC, launch the Analyzer application (see Start Analyzer application).

Direct setup

- 1. From the Microscope PC, navigate to **Maps Options>>Settings... menu item** to open the Amira-Avizo2D settings.
- 2. Click the Amira-Avizo2D tab.
- 3. Select **Direct** from the Connection Type field.

4. Configure the shared folder path as indicated in the table below.

Amira-Avizo2D Direct Settings		
Interface Element	Description	
Shared folder path from	Enter the path to the shared folder for the Maps software PC. See Configure shared folder for details.	
	Note: If the shared folder is on the Maps PC and the application encounters an I/O error while processing a recipe, consider moving the shared folder to the Amira-Avizo2D PC. See Define Shared Folder for instructions.	
Shared folder path from Amira- Avizo2D PC	Enter the path to the shared folder for the Amira-Avizo2D PC. See Configure shared folder for details.	
Hostname of Amira- Avizo2D PC	Enter the hostname of the Amira-Avizo2D PC. See Locate hostname for details.	
Connection time out	Determines the duration time required for Maps to connect with Amira- Avizo2D before Maps cancels the connection. The duration time is measured in seconds.	

5. From the Amira-Avizo2D PC, launch the Analyzer application (see Start Analyzer application).

NAT setup

It is required that you first set up the Support PC by following steps indicated in Maps settings to connect Maps with Amira-Avizo2D.

- 1. From the Microscope PC, navigate to **Maps Options>>Settings... menu item** to open the Amira-Avizo2D settings.
- 2. Click the Amira-Avizo2D tab.
- 3. Select **NAT** from the Connection Type field.

4. Configure the shared folder path as indicated in the table below.

Amira-Avizo2D NAT Settings		
Interface Element	Description	
Shared folder path from Maps PC	Enter the path to the Support PC shared folder for the Maps software PC. See Configure shared folder for details.	
Shared folder path from Amira- Avizo2D PC	Enter the path to the Support PC shared folder for the Amira-Avizo2D PC. See Configure shared folder for details.	
Hostname of Support PC	Enter the hostname of the Support PC. See Locate hostname for details.	
Hostname of Amira-Avizo2D PC	Enter the hostname of Amira-Avizo2D PC. See Locate hostname for details.	
Connection time out	Determines the time duration required for Maps to connect with Amira-Avizo2D before Maps cancels the connection. The time duration is measured in seconds.	

- 5. Click Setup Support PC.
- 6. Click **OK**, and then follow steps indicated to set up the Support PC.



7. From the Amira-Avizo2D PC, launch the Analyzer application (see Start Analyzer application).

Configure shared folder

The shared folder is used as an image and recipe exchange location between Maps and Amira-Avizo2D (see **Define Shared Folder** for details). Configure the shared folder the way it can be accessed by both Amira-Avizo2D and Maps applications. The shared folder must be mapped as drive according to the following connection options:

- Local: On the same computer as Maps if both Amira-Avizo2D are hosted on the same computer.
- **Direct**: Either on the Amira-Avizo2D computer and drive mapped to the Maps computer, or on the Maps computer and drive mapped to the Amira-Avizo2D computer.
- NAT: On the Support PC, drive mapped on both the Microscope PC and the Amira-Avizo2D computer.

Test connections

Test the Maps and Amira-Avizo2D connectivity after configuration of the hardware and shared folder to verify the connection. Go to the Connection with the Amira-Avizo2D field, and then click **Test Connection** to test the connection between Maps and Amira-Avizo2D.

When testing the connection to determine whether or not the images can be analyzed with Amira-Avizo2D, there are five possible results.

• **Connection**: If the Maps and Amira-Avizo2D connection and shared folder are correct, then the following message appears. Click **OK**.



The message indicates settings supported by an Amira-Avizo2D instance connected to Maps (see Recipe processing specification).

• No Connection: If the Amira-Avizo2D host is incorrect while in Direct or NAT mode, or the Analyzer application is not launched in Local or Direct mode, then the following message appears.



Solution: Click **OK**, and then check that the hostname is entered correctly and that Analyzer is running. If not, see Configuration options and Start Analyzer application.

 No Connection: If the Analyzer application is not launched in NAT mode, then the following message appears.



Solution: Click **OK**, and then check that Analyzer is running. If not, then see **Start Analyzer** application.

• **No Connection**: If the selected Amira-Avizo2D version does not correspond to the running application, then the following message appears.



Solution: Click **OK**, and then check that the Amira-Avizo2D version matches with the running application. If not, then select the matching version.

• **Connection without shared folder**: If Amira-Avizo2D connects and the shared folder is not set up correctly, then the following message appears.



Solution: Click **OK**, and then check that the entries for the shared folder point to the same location from both PCs. If not, see **Configure shared folder**.

Start Analyzer application

It is required that you run the Analyzer application on the Amira-Avizo2D PC to use Amira-Avizo2D capabilities with Maps. The Analyzer application is included with the Amira-Avizo2D software package.

Click the **Analyzer for MAPS** desktop shortcut, which launches the Analyzer application with the graphical user interface (GUI) used for recipe processing.

Process images with recipe

Once the connection to Amira-Avizo2D is successful and you have launched the Analyzer application, then the system is ready to process images with Amira-Avizo2D. To begin this process, you must open, create, or import a Maps project from the shared folder. You can use either a SEM/SDB or TEM tile set; SEM/SDB or TEM acquired tile set; or stitched image. Note that the recipe processing specification differs depending on the software version that is defined in the Amira-Avizo2D settings (see Recipe processing specification).

SEM/SDB or TEM Tile Set

- 1. On a newly created tile set, open the Amira-Avizo2D section.
- 2. Configure recipe processing options according to your use case:

Interface Element	Description	
Recipe file path	Select the .hxisp file used for recipe processing. This file must be located in the shared folder.	
Measure group	Select the measure group that will be processed on Amira-Avizo2D from all measure groups available in the A2DAnalyzer application. See the Amira-Avizo2D documentation for more information on Measures.	
	• Select the Refresh button to reload measure groups available in Amira- Avizo2D. Refresh is useful if measure groups were added or removed on A2DAnalyzer in the meantime.	
	• By default, a None entry is selected and it will not produce any .csv measure output in the layer directory.	
	Note: When measure groups cannot retrieved, the default measure group list contains a Custom measure group. Select the Custom measure group only when it is available in the A2DAnalyzer application, otherwise recipe processing fails.	

3. Finalize your tile set configuration and run the Job Queue. While the tile set is acquired, the resulting layer tiles are updated once they are processed by Amira-Avizo2D.



•	There is no recipe tile result for 'Z:\Shared Kronos' MD4\Tile_001-001-000000_0-000.s0001_e00.tif'. Something went wrong during tile recipe processi	\A2d\LayersData\Layer\Tile Set ing in Amira-Avizo2D application.
Don	ot show this message again	ОК

SEM/SDB or TEM Acquired Tile Set

- 1. Right-click on the tile set acquired on the layer tree, and then select **Recipe Processing**.
- 2. Configure recipe processing options according to your use case:

SE	M or TEM Acquired Tile Set Recipe Processing Overview	
Interface Element	Description	
Recipe file path	Select the .hxisp file used for recipe processing. This file must be located in the shared folder.	
Measure group	Select the measure group that will be processed on Amira-Avizo2D from all measure groups available in the A2DAnalyzer application. See the Amira-Avizo2D documentation for more information on Measures.	
	• Select the Refresh button to reload measure groups available in Amira- Avizo2D. Refresh is useful if measure groups were added or removed on A2DAnalyzer in the meantime.	
	• By default, a None entry is selected and it will not produce any .csv measure output in the layer directory.	
	Note: When measure groups cannot be retrieved, the default measure group list contains a Custom measure group. Select the Custom measure group only when it is available in the A2DAnalyzer application, otherwise recipe processing fails.	

- 3. Click **Queue Job** to create a new job in the job control.
- 4. Run the Job Queue to start processing.
- 5. When the job is finished, a tile set is created containing the processing results from Analyzer.

Stitched image

- 1. Right-click the stitched image on the layer tree, and then select **Recipe Processing**.
- 2. Configure recipe processing options according to your use case:

	Stitched Image Recipe Processing Overview
Interface Element	Description
Recipe file path	Select the .hxisp file used for recipe processing. This file must be located in the shared folder.
Measure group	 Select the measure group that will be processed on Amira-Avizo2D from all measure groups available in the A2DAnalyzer application. See the Amira-Avizo2D documentation for more information on Measures. Select the Refresh button to reload measure groups available in Amira-Avizo2D. Refresh is useful if measure groups were added or removed on A2DAnalyzer in the meantime. By default, a None entry is selected and it will not produce any .csv measure output in the layer directory.
	 Note: When measure groups cannot be retrieved, the default measure group list contains a Custom measure group. Select the Custom measure group only when it is available in the A2DAnalyzer application, otherwise recipe processing fails.

- 3. Click Queue Job to create a new job in the job control.
- 4. Run the Job Queue to start processing.
- 5. When the job is finished, an image layer is created containing the processing results from Analyzer.

Recipe processing specification

The recipe processing specification differs depending on the connected A2DAnalyzer instance (that is, the version and licensing). To confirm A2DAnalyzer supported settings, test the connection between Maps and Amira-Avizo2D.

• Test connection to A2DAnalyzer 2021.1 licensed without PRO extension



• Test connection to A2DAnalyzer 2021.1 licensed with PRO extension



Recipe processing settings overview		
Settings	Description	
Maximum resolution	Indicates the maximum resolution in pixels supported by the connected A2DAnalyzer instance.	
Input type	Indicates the recipe input types supported by the connected A2DAnalyzer instance.	
	Choices include:	
	 Single: Recipe supports only a single input image and generates a single output image. As a result, applying a recipe on a multichannel layer produces a multichannel layer with the same number of channels. 	
	• <i>Multiple</i> : Recipe supports multiple input images and generates a single output image. As a result, applying a recipe on a multichannel layer produces a single channel layer.	
Combined CSV file	Determines if a measure as a recipe parameter produces a single .csv file for all processed images named 'CombinedMeasure.csv' in the result layer folder. If not activated, Maps produces a .csv file for each processed image in the result layer folder.	
Default unit	Default unit supported by the connected A2DAnalyzer instance, which is used to exchange images.	

Define Shared Folder

A shared folder is necessary to allow data exchanges. This folder can be located in any place, as long as it is accessible to both software applications. It can be located on one of the two computers.

- 1. To share a folder, determine the folder you want to share.
- 2. Right-click on the folder and click Show Properties.
- 3. Click **Share** in the **Properties** dialog box.

src Properties
General Sharing Security Previous Versions Customize
Network File and Folder Sharing
Not Shared
Network Path: Not Shared
Share
Advanced Sharing
Set custom permissions, create multiple shares, and set other advanced sharing options.
Advanced Sharing
OK Cancel Apply

4. Select Everyone for Name, click Read/Write for Permission Level, and then click Share.

3 File Sharing	
Choose people on your network to share with	
Type a name and then click Add, or click the arrow to find som	neone.
.,,-	
	▼ Add
Name	Permission Level
Administrators	Owner
La Chuquillanqui, Sami (Sami.Chuquillanqui@fei.com)	Read/Write 💌
A Everyone	Read/Write 🔻
I'm having trouble sharing	
the name of the state states	

5. Click Done.

🚱 <u>8</u> File Sharing	
Your folder is shared.	
You can <u>e-mail</u> someone links to the	ese shared items, or <u>copy</u> and paste the links into another program.
Individual Items src \BOD1151\src	^
B Constituted Into these China	
Show me all the network shares on t	his computer.
	Done

6. Right-click on the shared folder and then click **Map Network Drive** to define the shared folder as a network drive to simplify access to this folder.

	08
G + Network + bod1151 +	■ 4 ₂ Search bod1151
Organize Search active directory Network and Sharing Center View remote printers	\$* CI (
Cropbox (Personnelle) Secont Places Maps, Data Share	Share Open
Documents Git Music	Open in new window Ajouter à la liste de lecture de VLC Lire avec VLC
E Pictures Videos	KD#3 • Git Init Here Git Gui
Compare systems (0)	Git Bach What is locking this folder?
도 및 en (, Sec 803) (2) 로 vol 5 (, Vac 8030) (8) 로 schward 5 (, Vac 8043) (12) 로 public (, Vac 8043) (12) 도 public (, Vac 8043) (12)	Share with Get Clone Get Clone Get Clone
♀ holdingsres (\\hbscmitg) (\\\) ♀ MapcShare (\\bodhasBL) (\k) ♀ holdingsres (\\bodhasB2 storage) (\\)	Scan for threats Shared Felder Synchronization
S≥ public (1/bod043) (2:) ♥ Natwork	Restore previous versions Map network drive
sec (\\bod1153) Offline availability: Not availabile Share Offline status: Online	Copy Create shortcut Properties

7. Select the network drive and click Finish.

9	😪 Map Ne	etwork Drive
	What ne Specify the	twork folder would you like to map? drive letter for the connection and the folder that you want to connect to:
	Drive:	T: •
	Folder:	\\bod1151\src
		Example: \\server\share
		Reconnect at logon
		Connect using different credentials
		Connect to a Web site that you can use to store your documents and pictures.
		Finish Cancel

Locate hostname

A hostname is a label assigned to a device (that is, a "host") on a network. It distinguishes one device from another on a specific network, or over the Internet.

Do the following to locate the hostname for your computer.

- 1. Click the Windows Start icon
 on the left end of the taskbar to open the Start menu.
- 2. Click the Settings cogwheel icon it to open the Settings menu.
- 3. Click the System icon to open the system settings screen.



- 4. Navigate to the left-side panel and click About.
- 5. Scroll down to the Device specifications section and locate the computer model (for example, Precision 7540).
- 6. View the Device name, which is also the hostname for your computer.

Device specifications

Precision 7540 Device name COMPUTER-12345

Tomography

This section describes how to use Tomography with the Maps software.

Topics include:

Import Tomography Presets Import Tomography Batch Positions View and Visit Tomography Batch Positions Send Annotations to Tomography (Beta)

Import Tomography Presets

This topic describes the Import Presets option under the Tomography menu.

 Note: For this menu option to work, you must have both Maps and the Tomography 5 Software installed. Also, you must have used Tomography at least once, and be aware that the import will only work for the instrument modes that have been used for Tomography (EFTEM, STEM, and TEM).

The following settings, which are available in both Tomography and in Maps, are stored as templates when you use the import option.

- Atlas
- Exposure
- Template
- Overview
- Search
- Template

Import presets

- 1. Navigate to the TEM Options menu and click Import Tomography presets....
- 2. If not all the instrument modes are available, then a message appears to indicate which instrument modes are missing. Click **Yes** to continue.



3. If a template already exists for Tomography presets, then the following warning appears to indicate that it will be overwritten by the import. Click **Yes** to continue.



4. The following message appears in the import is successful. Click Yes to continue.



View stored templates

Each successful import is stored as a template, and you can inspect or alter them from within the Tile Set Templates screen. Note that the imported presets end with (from Tomo) as shown below.

Node TEM				`
Factory Template (TEM) (Default)	Properties			
Exposure (from Tomo)	Name Atlas (from Tomo)			
Search / Template (from Tomo)	Tile Set Name Tile Se		t	
Overview / Positioning (from Tomo) Atlas (from Tomo)	These settings are used as initial values for new TEM Tile Sets.			
	Camera			
	Tiles X, Y			
	Overlap X, Y			
	Tile HFW			
	Total Area			
	Resolution		4096 x 4096	
	Magnification			
	Pixel Size		195.3125 nm	
	Exposure Time			
	Illuminated Area			
	Spot Size			
	Probe Mode			
	Focus Method			
	Defocus			
	Stage Settling Time	e		
	Image Correction			

Import Tomography Batch Positions

This topic describes how batch positions are imported from Tomography to Maps. How you import batch positions depends on whether or not the Maps Data Exchange Service is installed on the system.

Note: For the import operation to work, you must have both Maps and the Tomography 5 Software installed. Also, you must have imported the Exposure preset from the Tomography 5 Software. If you want to use automatic synchronization instead of the file-based exchange to import the positions, then the Maps Data Exchange Service must also be installed.

Import and update batch positions

- 1. Navigate to the **Tomography menu**, and then click **Import Batch Positions...** to open the system file dialog.
- 2. Click the file you want to import, and then click **Open**.
- 3. Maps imports the Tomography batch positions after successfully importing a new top layer. You can view the imported positions in the Viewer as shown below.



- 4. Hover your mouse cursor over an imported position and right-click to open the context menu.
- 5. Select **Update Tomography batch positions** to update the current list of batch positions. Positions are reloaded from the batch positions file last used. If the file is no longer available, the system file dialog opens so that you can select a different file to import. All existing positions that are not included in the new file will be deleted.

Synchronize batch positions with Maps Data Exchange Service (beta)

- 1. Confirm the following preconditions are met.
 - a. Maps is running
 - b. Tomography is running and a session has been created
 - c. Tomography is connected to the Maps Data Exchange Service
 - d. One or more batch positions have been created
- To activate the automatic synchronization of data with Tomography, either click Synchronize With Tomography under the Tomography menu, or click the Toggle Automatic Synchronization icon in the TEM Toolbar, Right Side. Once activated, the arrows in the toolbar icon turn green.



3. Maps creates a new layer that represents the active Tomography session. It also imports batch positions from the active session. You can view the imported positions by expanding the layer, as shown below.



4. While the automatic synchronization is turned ON, any batch position changes in the active Tomography session are automatically propagated to Maps (end-user interaction is not required).

View and Visit Tomography Batch Positions

This topic describes how you can view and visit Tomography batch positions after they are imported from the Tomography 5 Software application.

View batch positions

You can view the batch positions in the Viewer from two different modes, zoomed-out or zoomed-in.

Zoomed-out mode

When you view the batch positions from the Zoomed-Out mode, you can see an overview of all the lamellae locations (indicated in green).



Zoomed-in mode

When you view the batch positions from the Zoomed-In mode, you can see the locations of the tracking area (indicated in green), exposure area (indicated in yellow), and focus area (indicated in purple) for each individual lamella.



Visit batch positions

You can move the stage to the batch position or zoomed-in location of a lamella's areas for tracking, focus, and exposure. To move the stage, click **Drive To** under Position properties as shown below.



Send Annotations to Tomography (Beta)

This topic describes how annotations are sent from Maps to Tomography.

Note: For the operation to work, you must have both Maps and the Tomography 5 Software installed. Also, Maps Data Exchange Service must be installed.

Preconditions

- 1. Before sending the annotations, confirm the following preconditions are met.
 - a. Tomography is running, a session has been created, and the session is connected to the Data Exchange Service
 - b. Maps is running, a project has been created, and the project is connected to the Data Exchange Service
- 2. Create one or more annotations in Maps. The annotations may include any of the following:
 - a. Sites of interest
 - b. Areas of interest
 - c. Lamella sites

Synchronize all annotations automatically

- 1. To activate the automatic synchronization of data with Tomography, either click **Synchronize With Tomography** under the **Tomography menu**, or click the **Toggle Automatic Synchronization** icon in the TEM Toolbar, Right Side.
- 2. The arrows in the toolbar icon turn green once the automatic synchronization is activated. Maps sends all annotations to Tomography and then displays them in the Tomography viewer.



3. While the automatic synchronization is turned ON, any batch position changes in the active Tomography session are automatically propagated to Maps (end-user interaction is not required).

Send all annotations of given type manually

- 1. Click Send To Tomography under the Tomography menu.
- 2. Click an annotation type from the submenu to send all annotations of the selected type from the current Maps project to Tomography.



3. To update the annotations of your selected type at a later time, repeat the previous steps.

Send selected annotations manually

- 1. Select annotations in the layer tree of the current project.
- 2. Hover your mouse cursor over an imported position and right-click to open the context menu.

✓ 🔮 📰 TEMPLATE	Factory Template (TEM)	
👻 🗹 Annotations		
Site of interest		Pring To Front
Area of interest		Bring Forward
🔽 🔎 Lamella site		Send Back
Tomography Session		Send To Back
		Center View Ctrl+Shift+C
Name	Site of interest	Center and Rotate View
		Drive To
Notes		Export To Project
Color		Alignment
	Ellipse	Delete
Source System	Talos Beta	Send To Tomography

- 3. Click **Send to Tomography** to send selected annotations from the current Maps project to Tomography. Any annotations that have been previously sent, are replaced by the current annotations.
- 4. To update the annotations at a later time, repeat the previous steps.
Array Tomography

This section describes how to create an Array Tomography (AT) job using Maps software. Maps AT functionality requires an active AT license.

Topics include:

About Array Tomography

Array Tomography Workflow

Extended Controls

About Array Tomography

AT is an imaging technique in which 3D image data is reconstructed from serial recordings on serial sections that are spread out on a large sample carrier.

Samples

Samples for AT are produced by serial sectioning of a plastic-embedded sample in a microtome. Sections are deposited on ITO-coated glass substrates (that is, 22 x 22mm cover glasses) or on wafers.

Depositing samples is done manually or by a tape collection device (that is, ATUMtome). The resulting AT samples differ fundamentally.

Classic and Tape Ribbons



Sample Sections and Tile Set Arrays

The main purpose of the AT feature is to easily align high-resolution imaging on areas of interest on all sample sections. The AT feature set consists of a concept called a *Tile Set Array*. The Tile Set Array sets up local coordinate systems around the individual sample sections to create a collection of tile sets across *all* sample sections.

Terminology

- Section: Defined as a single slice of the overall sample placed on the stage. Alternatively used to refer to the outline of a given sample slice that is used in Maps.
- Tile Set Array: A collection of tile sets that span all or some of the sample sections. There will be one tile set per section initially and they are all placed in the same relative location within their section. Tile set arrays are used as the source images for creating a 3D data set. Most tile set settings are shared across the whole array. All the tile sets created within this array are linked. If the position, rotation, or imaging parameters of one tile set changes, then the rest of the tile sets will be updated.

Array Tomography Workflow

The AT workflow contains two branches to set up and capture images for an AT reconstruction:

- A branch for mapping out all sections on your sample manually.
- A branch for running an algorithm to auto detect the sections.

The following diagram shows the steps in the worklfow for both branches.



Start the Array Tomography Workflow

- Step 1: Define and Place Sections
- Step 2: Define Section Order
- Step 3: Preview Sections (Optional)
- Step 4: Refine Positions
- Step 5: Create Tile Set Arrays
- Step 6: Acquire Tile Set Arrays

Start the Array Tomography Workflow

An Array Tomography (AT) job is a workflow used within the Maps workflow panel. The workflow panel is a control that contains step-by-step processes that help make complicated jobs easier. The workflows available depend on your product license.

1. Click the **Start Workflow** icon located on the right side of the Toolbar.



2. In the Select Workflow window, click Array Tomography, and then click Begin.



3. The **Workflow: Array Tomography** window appears on the right side of the main screen. To navigate through the steps within the workflow, use the forward and back arrows located at the bottom of the window. To exit the workflow, click **X** at the top-right corner of the panel.

	Workflow: Array Tomography
1	Define section template
	Zoom into any of the sections so you can see the whole section. Rotate the viewer to square up the section with shift+drag. Then click the Define Section Template button and drag an outline around the whole section.
	Define Section Template
	Section Template Image
	No template image defined yet.
2	Run Section Finder
	Select Layers and Tile Sets to search, then click the Find Sections button. The Section Finder will automatically find sections based on the Template Image. Optionally click the cog wheel button to change the Matching Threshold.
	No search images available
	Please acquire or import some images to search with
	Find Sections 0 ⁰
3	Define sections manually (optional)
	Sections can be defined manually if the section finder was not used or if they were missed. To add section outlines, click the Place Sections button or hold down 's' and click in the viewer. To remove section outlines, (Ctrl-) select them and hit the Delete key. Be careful to not delete images!
	Place Section Outline (s)
	Move View To Next Predicted Location
	Section Definition

Step 1: Define and Place Sections

The first step of the workflow is to define a section template to pattern match across the entire sample for section placement.

Workflow procedure

- 1. Define the section template.
 - a. Use the mouse scroll wheel to zoom into one section of the sample.
 - b. Press **Shift+left mouse button** to rotate the viewer until the section aligns horizontally and vertically within the viewer.
 - c. Click **Define Section Template**, and then draw a rectangle around the sample slice in the viewer. Maps displays this image selection in the **Section Template Image** box within the workflow panel. Array Tomography then uses this image as the search template within the section finder.



- 2. Use the section finder to automatically search for regions of overview images that match the **Section Template Image** for fast section placement. *Note that if you have 10 sections or less in your workflow, then it is faster to perform this step manually.*
 - a. In the Searchable Layers pane, select the desired overview image layers for the Array Tomography search.

Select images to search in: All Cle		Clear		
•	▶ NavCam			
•	v (Overview Tile Sets		
🔽 Tile Set				
🗹 Tile Set (2)				
		Find Sections		° 🖻

- b. *Optional*: You can click the double gear icon to open a dialog to set a matching threshold for the section finder. This icon is located below the layers pane.
- c. *Optional*: You can click the trash can icon to either clear all sections if (a) you would like to start over, or (b) the section finder did not function as expected with your defined parameters. This icon is located below the layers pane.

d. Click **Find Sections** to start the automated section finder. When the automated section finder finishes searching, the viewer displays the outlined sections.



- e. The section finder might place an extra section outline where a sample section does not exist (this is called a "false positive"). If this occurs, then delete the false positive as follows:
 - a. Click the section outline in the viewer.
 - b. Press the **Delete** key to remove the extra section outline.

- 3. Add any remaining sections manually.
 - a. Confirm that the section requiring an outline is correctly aligned within the viewer. The software application typically aligns the section within the viewer.
 - b. Click Place Section Outline (s) to generate a preview outline of the missing section or press the letter s on your keyboard as a shortcut. Next, select the Move View To Next Predicted Location check box to auto-orient the view based on the two previously placed section locations and rotation.

3	Define sections manually (optional)
	Sections can be defined manually if the section finder was not used or if they were missed. To add section outlines, click the Place Sections button or hold down 's' and click in the viewer. To remove section outlines, (Ctrl-) select them
	and hit the Delete key. Be careful to not delete images!
ſ	and hit the Delete key. Be careful to not delete images!
ſ	and hit the Delete key. Be careful to not delete images! Place Section Outline (s)

- c. Click in the viewer to place the outline.
- 4. Once you place all the sections, click the forward arrow at the bottom of the workflow panel to continue to Step 2: Define Section Order.

Step 2: Define Section Order

The second step of the workflow defines the order of the sections.

Ordering tool

An ordering tool allows you to draw a line through a collection of sections within the viewer to determine the order. Sections closest to the start of the line are first in the order; whereas, sections near the end of the line are at the end of the sequence.

There are two available ordering tools for you to use:

- Freeform Tool: Use this ordering tool to draw a custom line in the viewer like a pencil tool used in graphical editors.
- Line Tool: Use this ordering tool to draw a line between two points by dragging a line in the viewer.

Workflow procedure

The viewer displays the newly ordered sections once you define their order. To define the order of the sections, do the following workflow procedure:



- 1. Click the user interface button of your chosen ordering tool to activate it.
 - Freeform Tool: Click the Freeform Tool (f) button.
 - Line Tool: Click the Line Tool (s) button.
- 2. Starting with the first section in the sequence, use the ordering tool to draw a line in the viewer tool to indicate section order. Note that the *unordered sections* are highlighted in brown and the *ordered sections* are highlighted in green.



Note: Use the keyboard shortcut to quickly switch ordering tools anytime during this step of the workflow. Press the letter **f** on your keyboard to activate the Freeform Tool; press the letter **s** on your keyboard to activate the Line Tool.

- 3. Repeat the previous step until all sections appear in the ordered list. If you want to delete the current order and start over, then click **Reset**. The number of unordered sections is displayed at the bottom of the panel for your quick reference.
- 4. After you have ordered all sections, click the forward arrow at the bottom of the workflow panel to continue to Step 3: Preview Sections (Optional).

Step 3: Preview Sections (Optional)

The third step of the workflow is optional, but this step is highly recommended since it will increase the accuracy of the section placement on your sample for more aligned final data capture. Proceed with this workflow step to acquire high-resolution images for each section, so that you can then perform section position refinement.

Note: To skip this step and Step 4: Refine Positions in the workflow, DO NOT acquire any additional section previews. Be sure to delete all acquired section previews and existing section preview tile sets. Once you have deleted these items, click the forward arrow at the bottom of the workflow panel to continue to Step 5: Create Tile Set Arrays.

Workflow procedure

- 1. Determine a section in the viewer to acquire and use for your section preview tile sets.
- 2. *Optional*: Select the desired tile set to use for previews in the viewer, and then click **Center View** to zoom in on and match the rotation of the tile set.

1	Define First Preview Image		
	Select a section in the viewer to use	when setting up section previews.	
	Center View	Create Tileset	

3. Click **Create Tileset** to define the section preview template, which is a tile set that will be used for all other section preview tile sets (including their properties).

- 4. Edit the section preview template to capture a quality image of the section. Note that you can edit the section preview template like other tile sets.
 - a. Set up the recommended parameters, so that the section preview template has good image contrast and recording speed:
 - Use a single tile with a 100-1000 nm pixel size
 - Set the dwell time to 2 µs
 - Use a minimum total resolution of 2K for the tile set
 - b. Click **Acquire** to acquire, or click **Reset** to reset the acquisition results of the section preview template.



5. Click **Acquire Previews** to create and acquire all remaining section previews based on the section preview template.



6. *Optional*: Click **Delete Previews** to delete all section previews except the template section preview.

 Note: If you encounter workflow issues while acquiring section previews, repeat Steps 1-5 to recreate the template section preview and delete all section previews. Array Tomography

7. Once you have acquired all section previews, click the forward arrow at the bottom of the workflow panel to continue to Step 4: Refine Positions.

Step 4: Refine Positions

The fourth step in the workflow will help increase the accuracy of the positions of the section outlines based on the preview images acquired during Step 3: Preview Sections (Optional) of the workflow.

Workflow procedure

1. Click View First Section, and then click Define Refinement Region.

1	Define a region in the first section to refine section positions with.		
	View First Section Define Refinement Region		
	Refinement Template Image		

 In the viewer, drag the mouse to create a yellow border box to define the refinement region. The Maps software application will then use this region to begin refining the section positions. The refinement region image cutout is displayed in the Section {section #} Result box.



- 3. Click Run Refinement to begin the refinement process.
 - Refinement progress is displayed in the workflow panel. The example below shows the refined result for a selected section. To stop the refinement process, click **Stop**.



• During the fine refinement sequence, Array Tomography places green rectangles in the viewer where the refinement process found an appropriate match. The refinement region image displayed and the offset results both update based on the incoming results. Refinement results appear based on the order of the sections.



- After the refinement process completes, you can browse sections using the slider, input a number in the input box, or press the play button to navigate through all sections.
- Adjusting the **Target-Result** slider will adjust the relative transparencies of the current result and the previous result.

4. If there are no failures during the refinement run, then the workflow skips this step. If there are failures, then a dialog box appears after the refinement completes. The dialog box displays the names of the sections that require your attention.



Failed sections are highlighted red in the viewer.

- a. To update or fix failed sections, navigate within the viewer to the failed section cutout and select it.
- b. Click **Adjust Result**. The cutout color should change to blue, indicating that a manual adjustment is in progress.
- c. Move and rotate the cutout in the viewer to correct its placement.
 - Move the slider control below the Section Result image to the Target side. This will overlay the current result on top of the previous result with varying transparency.
 - To help target a center spot, match the location of the green crosshair displayed within the blue box of the viewer to the red crosshair displayed in the Section Result image.

- d. There are several choices once the manual adjustment has been made.
 - Save
 - ^o Use the Save option to save the current manual adjustment for the section.
 - The section cutout will update to green in the viewer, thus indicating that it has been refined.
 - All other sections have their refinement status retained.
 - Save and Continue
 - Use the Save and Continue option to save the current placement of the section cutout. Array Tomography starts running the refinement again from this section onwards, thus redoing any previous refinement for sections following the saved section cutout.
 - The section cutout will update to green in the viewer, thus indicating that it has been refined.
 - Cancel
 - Use the Cancel option to cancel any current adjustments to the section.
 - The section cutout will update to its color in the viewer to the color it had prior to starting the adjustment.
 - Trash can
 - Use the Trash Can option for sections that were damaged during sample prep and will result in inadequate final data.
 - ° The Trash Can option marks the section to be deleted.
 - The section cutout will be updated to a light grey color in the viewer.
 - If the refinement is re-run, this section will be ignored from the run.
 - When applying the refinement results, this section will be deleted from the project along with its associated preview tile set.
- e. Repeat the previous steps for any remaining failed sections.
- 5. After all sections have passed refinement, check how well the refinement placed itself by using the slider bar above the Section Result image box. Do the following to check all refined sections:
 - a. Move the slider from left to right to show the final results in the Section Result image region.
 - b. Use the Play button to the right of the slider bar to automatically animate the sequence.

6. Click **Apply Refinement Results** to apply the refinement offsets to the section outline placement. Or, click **Reset All Results** to start from the beginning of this step.

3	Apply results to update all existing section positions accordingly.	
	Apply Refinement Results	Reset All Results
0	Note: Applying the refinen	ent results is a one-time action is not reversible and it
	will automatically progress you to the next step in the workflow.	

7. When you are satisfied with the section outlines placement, click the forward arrow at the bottom of the workflow panel to continue to Step 5: Create Tile Set Arrays.

Step 5: Create Tile Set Arrays

In the fifth step of the workflow, you create and edit the tile set arrays.

Workflow procedure

1. Click Create Tile Set Array.



This step creates a new tile set array. Each tile set is placed in the center of a section.

i Note: There is an exception to the tile set linking functionality in the array for one of the imaging parameters. The Focus Strategy for all tile sets within the array are shared. However, if the Focus Strategy is set to Fixed, then the actual Fixed Value is not shared between each tile set. This allows you to manually focus and store the found focus position for each tile set for samples on which autofocus does not work.

2. Once the tile set array is created, you can adjust the properties of the tile set array with the edit controls provided. To learn more about these edit controls, see Extended Controls.

2 Use the controls below to edi	t the selected tile set array.
These controls can also be found in array layer is selected and can be u	n the left panel of Maps when a tile set used outside of the workflow.
Section Span Slider	Total Available Sections 5
< 1 >	● < 4 >
Current Edit Mode	6 6 6
View Controls	Section 0001
	Center View On Section 0001

3. *Optional*: After final acquisition, select the **Stitch during acquisition** checkbox and a folder to output each stitched tile set to a designated folder. When activated, this functionality outputs a series of images to your specified folder after acquisition. These images are all fitted to be the exact same size and are aligned, so they can be reconstructed into a 3D volume within another software application. If data has already been acquired, then the **Finalize Stitch Results** button becomes activated and you can click it to manually export the images.

Stitch during acquisition Stitch All Output Directory:		
D:\Data Sets\Projects\Array Tomography Sample 1		
Finalize Stitch Results		

4. When you have created and edited all tile set arrays as desired, click the forward arrow at the bottom of the workflow panel to continue to Step 6: Acquire Tile Set Arrays.

Step 6: Acquire Tile Set Arrays

In the sixth step of the workflow, you acquire the tile set arrays.

Workflow procedure

1 Choose an order of acquisi grouped by section.	tion for acquiring the tileset arrays,
Û Į	
1. Array Tile Set 001	
	lead area the Area waited was the
2 Verify that all tile sets are p "Update Queue" button to sch	aced correctly. Once verified, use the redule all tileset arrays for acquisition.
Upd	ate Queue
3 In order to begin acquisition queue.	n, press the RUN button near the job

- 1. Double-check that all desired tile sets have been created.
- 2. Order the tile sets.
 - a. Click a tile set to select it.
 - b. Click the arrow buttons to move it either up or down.
- 3. Click Update Queue to add all tile sets to the acquisition queue, grouped by their section.
- 4. Click the **RUN** Button on the bottom status bar to begin final acquisition.



Extended Controls

Create and manipulate AT data sets with tools described in this topic.

Most of the work for AT occurs within the Start the Array Tomography Workflow. These extended controls allow you to create and manipulate AT data sets outside of the workflow.

Layer property pane

Some layers types that are associated with an AT job have extended controls in the layer property pane in the Array Tomography section. The options change depending on what you select in the layer tree.

Tile set array

Click the tile set array group layer. Controls appear in the **Array Tomography** pane that allow you to edit the tile set array.

>	
Name	Array Tile Set 001
Opacity	
Section Span Slider	Total Available Sections 5
	< 2 >
Current Edit Mode	6 6 6
View Controls	Section 0001
	Center View On Section 0001

- Name The name of the selected tile set array.
- **Opacity** Controls the level of transparency for the tile set array. Click the slider knob and move it to the right for low transparency, or move it to the left for high transparency. By default, the slider knob is on the extreme right so that the tile set array is completely visible.

- Section Span Slider Controls which sections the tile set array spans across. The left value is the section index that the tile set array starts on, the right value is the section index where the tile set array ends.
- **Current Edit Mode** Changes which tile sets within the selected array change when one tile set from the array is dragged or rotated in the viewer.



• View Controls - Center the viewer on the displayed section or move the view from one section to another while keeping the same relative field of view. This can help you inspect the results of a tile set array.

Section outlines

The layer group that contains all the section outlines. It includes **View Controls** described above and displays the number of **Total Available Sections**.

۵ 🗲	
Name Opacity	Section Outlines
Total Available Sections	4
View Controls	Section 0001
	Center View On Section 0001

Section

The Section layer includes **View Controls** described above and shows the **Section Index** and **Section Rotation** of the selected section.

Section Index	2
Section Rotation	303.63
View Controls	Section 0002
	Center View On Section 0002

Layer tree context menu

Tile set array



The tile set array layer group in the layer tree has some extended options in its right-click context menu.

- Copy Tile Set Array Creates a complete copy of the whole tile set array.
- Reset Acquisitions Resets the acquisitions of any tile sets within the tile set array.
- Enqueue Tile Set Array Queues up the whole tile set array for acquisition.
- Dequeue Tile Set Array Removes the whole tile set array from acquisition.
- Stitch All... Prompts you to start a Stitch All operation. This has the following outcomes:
 - All stitched images are saved to one output directory.
 - All images are saved with the same dimensions. This makes it easier to create image stacks.
- **Delete** Deletes the whole tile set array. When you select this option, a message dialog appears to summarize all the items designated for deletion. Verify the summary, and then click **Delete** to confirm.

Create a new tile set array

- 1. To create a new tile set array outside of the workflow:
- 2. Press Alt and drag to select an area that is completely contained within a section in the viewer.



3. Right-click and then click Create Tile Set Array here.

This creates a new tile set that spans all selected sections.

Known issues

The following known issues have been identified in the Maps software application.

Maps 3.32.1 known issues

Maps 3.30.2 known issues

Maps 3.29.1 known issues

Maps 3.28.1 known issues

Maps 3.27.1 known issues

Maps 3.24 known issues

Maps 3.22 known issues

Maps 3.20 known issues

Maps 3.18 known issues

Maps 3.14 known issues

Maps 3.13 known issues

General known issues

Amira-Avizo2D known issues

SEM known issues

TEM known issues

Mineralogy known issues

Maps 3.32.1 known issues

Channels for removed elements are not always properly deleted

When reprocessing of EDS channels starts before everything has been drawn to the layer from previous reprocessing or acquisition, the element maps might not be properly finished.

- To prevent this issue, wait until everything is drawn to the element channels.
- If the issue occurs, reprocess the layer again. It is recommended that you first remove count and quant maps, and then initiate the reprocessing again from the very beginning.

Maps 3.30.2 known issues

Mineralogy: Maps freezes when simultaneously updating Mineral Reference Editor and running a mineralogy acquisition with a recipe (also affects all previous Maps versions)

If you update Mineral Reference Editor (MRE) while running an acquisition with a configured MRE recipe, then the Maps user interface freezes during the X-ray acquisition steps. To prevent this issue, perform all necessary Mineral Reference Editor updates prior to initiating a Maps acquisition. If you encounter the user interface freeze, click **Stop** to pause the acquisition and then click **Run** to resume the acquisition.

Maps 3.29.1 known issues

TEM: Data exchange between Maps and Tomography does not work

The data exchange does not work with the version 3.0.0.72 of the Maps Data Exchange Service, which is included in the Maps 3.29.1 installation package (*ThermoScientificMapsDataExchangeService3.0.0.72.exe*).

Workaround: Use the *ThermoScientificMapsDataExchangeService3.0.0.70.exe* file, which is included in the Maps 3.27.1 installation package.

Mineralogy: Loading time for Mineral Association report

The Mineral Association report in Reporting may take some time to load due to the complexity of the analysis involved. However, to provide a better user experience, a loading icon is displayed to indicate that the report is loading. Additionally, a warning icon is displayed next to the report selection to alert you about the loading time.

Mineralogy: Reports data cannot be displayed after installation

When upgrading from Maps Min Reporting 3.28 to 3.29, you might encounter a problem in which the data are not displayed on the Reporting web page for a few minutes after the installation process successfully completes. To address this issue, it is recommended that you uninstall the Maps Min Reporting application and then reinstall it as a temporary workaround.

Mineralogy: Nanomin cannot open exported report in Excel (affects Maps 3.29.1 and any other Maps version on a Windows 11 system)

Nanomin might crash when attempting to open Excel with newly exported reports in Excel format on Windows 11. This issue occurs specifically when Excel displays the "Your privacy matters" dialog with Microsoft 365 (Office), or when the Excel license is not activated. To resolve this issue on Windows 11 systems, follow the steps listed below on the next page.

- 1. Ensure that your Excel license is activated and up to date.
- 2. Follow these steps to deactivate any Microsoft 365 (Office) privacy matters dialog that is triggered by Excel:
 - a. Open any Office 365 application, such as Word or Excel.
 - b. Click the "File" tab in the upper left corner.
 - c. Select **Options** from the drop-down menu.
 - d. In the Options window, navigate to the "Trust Center" tab.
 - e. Click Trust Center Settings...
 - f. In the Trust Center window, select Privacy Options.
 - g. Click Privacy Settings...
 - h. Uncheck the box that says, "Turn on optional connected experiences".
 - i. Sign out of the Microsoft application that you have open.
 - j. Open the Microsoft 365 (Office) application on your computer and sign out.
 - k. Restart all Office applications.
 - I. Sign back into your Microsoft 365 (Office) account.
 - m. Open a Microsoft application and log in. The privacy matters dialog should no longer appear.

Maps 3.28.1 known issues

TEM: Maps Data Exchange Service installer is missing in 3.28.1 installation package

The *ThermoScientificMapsDataExchangeService3.0.0.70.exe* file is not present in the Maps 3.28.1 installation package.

Workaround: Use the *ThermoScientificMapsDataExchangeService3.0.0.70.exe* file, which is included in the Maps 3.27.1 installation package.

Mineralogy: Centroid mode slows down Maps

When running the Centroid mode only or when using the BSE search ranges option within the Maps Min workflow, the Maps application could hang if you use high-resolution tiles, or if you use very fine particles or grains. The particles split into very fine particles if you use a high-strength setting during particle segmentation, and the grains split into very fine grains if you use a small grain size setting during grain segmentation.

The following can help you to prevent this issue:

- Use 512x442 for BSE resolution
- Use a lower strength setting for particle segmentation
- Use a larger grain size setting for grain segmentation

Maps 3.27.1 known issues

Inability to recover data if rare error occurs during splitting of tile set channels

If a rare error occurs during the splitting of tile set channels, data loss can occur in Maps and the data would not be recoverable. Also, if a Phenom EDS or Mineralogy tile set is split through the Split Channels feature, then the backing Mineralogy or Phenom data will be lost. This specific data loss would result in not being able to reprocess a Phenom EDS tile set and the data loss could also break Mineralogy tile sets.

Mineralogy: Distinct Particle Size Distribution report results between Nanomin and Reporting

Nanomin and Reporting use two different methods to calculate the Particle Size Distribution report results. Nanomin uses a calculated pixel area that encompasses all pixels within the particle, including those within porous regions; whereas, Reporting, uses a calculated pixel area equivalent to the sum of all grain pixel areas of the particle.

Phenom: Compatibility limitation (affects Maps 3.26.1 and Maps 3.27.1)

Maps requires the Phenom server version 6.6 or later.

Phenom: Remote connections not supported for EDS acquisition

Maps must be running on a computer directly connected to the Phenom instrument when performing EDS acquisition. Remote connections are not supported due to slow performance.

Maps 3.24 known issues

TEM: Lamella site imported from SDB might show incorrect milling angle

The milling angle of a lamella on SDB systems is computed based on several parameters, including the sample pre-tilt angle (which depends on the sample shuttle type), the stage tilt angle, and the angle of the ion beam. The sample pre-tilt angle is currently assumed to always be 35 degrees on TEM systems. If the shuttle type with a different pre-tilt angle was used on the SDB system, then the milling angle that is shown on the TEM system will not be correct.

This known issue will be resolved in an upcoming release of Maps. A workaround is not available.

Maps 3.22 known issues

TEM: Failure to acquire EDS tile set (affects Maps 3.19-3.22)

Maps encounters an error during acquisition of EDS tile set. The error description is the following: "Error encountered during job execution: Error trying to acquire an image: Index (zero based) must be greater than or equal to zero and less than the size of the argument list." The error is caused by a timing issue and thus appears to show up randomly.

Workaround: Restart Maps and Velox, and then perform the acquisition again to resolve the error. If the error persists, decrease the amount of data that is exchanged between Velox and Maps during acquisition (for example, by lowering the resolution of the acquired images and EDS maps).

Maps 3.20 known issues

Mineralogy: Segmentation Wizard hang

When running the Segmentation Wizard, the Maps application will hang if you use very fine particles or grains. The particles split into very fine particles if you use a high strength setting during particle segmentation, and the grains split into very fine grains if you use a small grain size setting during grain segmentation.

The following can help you to prevent this issue from happening:

- Use a lower BSE resolution
- Use a lower strength setting for particle segmentation
- Use a larger grain size setting for grain segmentation

Maps 3.18 known issues

SEM: Quattro acquisition failures

Maps may encounter an error during the first acquisition of a tile set or snapshot. The error description is "Error encountered during job execution: Job execution failed: Image grab action failed. The following error occurred: No description available. The error occurred was: No description available:" To resolve this error, close Maps and restart the xT Microscope Server.

Mineralogy: Data migration performance

Performing data migrations from Maps Min 3.14 or 3.1 can be slow. Terminating the process can corrupt the underlying database. Allow the migration process to complete to preserve data integrity.

Mineralogy: Mineral classification requires free space on C: drive

If the system C: drive is low on space, then Nanomin will report that the C: drive is out of space during classification even if the data is stored on a different drive. Ensure the C: drive has at least 10 GB of free space.

Maps 3.14 known issues

TEM: 4D-STEM acquisitions fail

If the camera is selected for STEM mode when acquiring EDS data in the Maps application with Velox, the acquisition fails and displays the following error message: "Error trying to acquire an image: Camera cannot be enabled in STEM mode during remote acquisition." When you receive this error message, deselect the camera and then restart Maps to continue acquisition.

TEM: Templates and camera replacement

If the camera is replaced with a different camera, then the tile set templates may no longer be valid and thus require manual updating.

Maps 3.13 known issues

SEM: Failure to acquire square resolutions

When acquiring square resolutions on Windows 7 microscopes (for example, 8096 x 8096), the acquisition may fail with an unhandled exception. If this error is encountered, either use a non-square resolution, or install the platform update for Microsoft Windows 7 SP1 KB2670838.

General known issues

Error message: "A generic error occurred in GDI+."

After installing the application and starting it for the first time, this message might appear.

Workaround: Restart the PC to resolve the issue.

Cannot select a tile set in the viewer

Workaround: Hide other items in the layer tree that are in front of the desired tile set. Alternatively, select the tile set in the layer control.

Error message: "There was a problem preparing the tiles for stitching. Please restart the application as soon as possible."

Workaround: Restart the application. If this does not resolve the problem, then restart the PC.

Stitching errors occur when stitching a tile set

See the "Stitching Best Practices" section of the Maps user guide.

Workaround: Optimize your tile set parameters and image quality, or try a different stitching profile.

Maps appears to hang while recalculating tile positions

For very large (> 100 GB) high-resolution tile sets, recalculation of tile positions can be time consuming.

Workaround: Allow a few minutes for this procedure to complete.

Maps does not open on the correct monitor

Workaround: Move the Maps window to the desired screen and click the X button located in the top right corner to close Maps. The next time Maps is opened, it opens to the correct monitor.

Newly-created tile set does not appear in the viewer

Workaround: Ensure the tile set and associated layer are enabled in the layer control. Also, the **Hide Annotations** feature (press Ctrl+H) can be used to make all Maps viewer graphics enabled or not enabled.

Projects from Maps v1.0 cannot be imported

Maps v1.0 projects are not supported.

Workaround: You can import Maps v2 and v1.1 projects.

Quit is not enabled in the Maps File menu

Workaround: If the automated holder alignment is running, Abort and Close the holder alignment UI. It should then be possible to quit Maps.

Amira-Avizo2D known issues

Sending two images in a row without performing the image calibration results in unexpected behavior from Analyzer. It is recommended that you always perform the image calibration after you send an image to Amira-Avizo2D.
SEM known issues

AutoFocus only runs in Quad 1

Even when acquiring multiple Quads or when Quad 1 acquisition is disabled, AutoFocus only runs in Quad 1. You cannot define a tile set that will AutoFocus on a different quad. Workaround: If you select AutoFocus or InitialAutoFocus in the Focus Setting property of a tile set, then the detector that you use for Quad 1 imaging must be the one that you intend to use for the AutoFocus operation.

Cannot use Line Integration with 8192 x 8192 or 4096 x 4096 image resolution

An error occurs when you use try to use Line Integration to capture a tile set with 8192x8192 or 4096x4096 resolution on the Helios PFIB microscope.

Workaround: If you need to use line integration, then select a different image resolution.

i Note: This issue only occurs on the Helios PFIB product.

SEM: Focus stack can result in blurry images

On Quanta and Prisma E systems, acquiring a focus stack can result in a blurry stripe along the top of the images.

There is no workaround. Avoid focus stack acquisition on these systems.

Line Integration and Scan Interlacing are ignored when acquiring multiple detector channels

When you use multiple quads on the **Advanced** tab of a SEM tile set, Maps ignores any values in the **Line Integration** or **Scan Interlacing** settings.

Workaround: If you need to use Line Integration or Scan Interlacing, then you can only specify a single detector channel in your tile set.

Unable to queue jobs that involve detector insertion/retraction

Maps does not insert or retract detectors.

Workaround: Manually insert retractable detectors.

Some tiles are not in focus when running tile sets with AutoFocus enabled

Workaround: Try modifying the AF parameters using the microscope control software. Alternatively, you can reacquire the tiles with bad focus using a different focus strategy.

Rotation is not supported when using Set Sample Plane

Workaround: Do not change stage R while using this feature.

Fast multi-detector tile acquisitions error message: "This detector combination is not supported for multiple detector acquisitions"

Fast acquisitions could be, for example, 1024 resolution with 50 ns dwell.

Workarounds:

- Use single-detector acquisition.
- Use a longer dwell time.
- Use a larger resolution.

SEM tile sets from earlier versions

Due to architectural changes, SEM tile sets from versions prior to Maps 3.6 can no longer be manipulated or acquired. This data can still be imported, visualized, and stitched. To acquire new SEM data into these projects, create new tile sets.

When acquiring a STEM tile set that is using auto focus and the auto focus routine fails on a tile, no image for that tile is acquired. This can result in some tile sets having blank tiles.

SEM: Lens Alignment auto-function is not available on Quanta FEG

This auto-function does not support the Quanta FEG.

SEM: PrismaE xT version 16.1 intermittently fails when using square resolutions

Square resolutions (such as 2048 x 2048 and 40960 x 40960) fail intermittently with an error: "Cannot move stage in X axis. It is either unavailable or locked." Do not use square resolutions on PrismaE with xT version 16.1. This issue is resolved in xT 16.2.

TEM known issues

After you select Ceta or Falcon I/II camera in PEOUI, it is not selected in Maps

When both Ceta and Falcon I/II cameras are not inserted, the Microscope Field of View (FOV) might show an unsupported camera.

Workaround: In PEOUI, insert the Ceta or Falcon I/II camera.

For large magnifications, the Microscope FOV might not be located at the exact position of the last acquired tile

At high magnifications, there can be a small differences between the tile positions and position of the Microscope FOV rectangle when the tile is being acquired. This small offset originates from small mechanical errors in the stage positioning.

Workaround: The FOV rectangle is the most accurate representation since it tracks the stage measuring system.

Snapshot image differs from a tile image.

When a snapshot image is taken at the same location and with equal settings that a tile was acquired with, the snapshot image differs from the tile image because the histogram of a snapshot is based on a single image; whereas, for a tile, the histogram is based on the entire tile set.

Unclear with which spot size a tile set is acquired when spot size is clipped

For example, when you select a spot size of 1 in Maps and acquire the tile set, the microscope clips the spot size to 5 for X-ray safety. In this case, the tile set is acquired with spot size 5. However, if you look at the tile set parameters it seems that it was acquired with a spot size of 1.

Workaround: Set the tile set to be acquired with a spot size that is not going to be clipped.

Tiles are misaligned

This can happen for a number of reasons. Maps uses the instrument alignments and calibrations to determine tile size and orientation. If these are off, then the tile set will not be properly aligned. Regular system alignments and calibrations are recommended as described in the system documentation.

Workarounds (contact your application support engineer for details):

- If the instrument is not at eucentric height and/or far out of focus, then adjust Z to set the stage at eucentric height and/or focus.
- If there are inaccurate lens alignments and magnification calibrations, redo the alignments and calibrations for the applicable magnification.

If the stage is non-orthogonal: Some TEM stages have a slight angle between the X and Y stage axis. This can be checked by performing pure X and Y moves using PEOUI and TIA. Then measure the angle. Maps has a provision to take this into account, but this requires adjustment of one of the configuration files.

When a tile set with invalid values is acquired, it is unclear with which settings it is acquired

When one or more parameters of a tile set are out of range, this is indicated by a red box. It is still possible to acquire such a tile set. The values used during acquisition are the last valid values that you entered.

Workaround: Enter valid values, so that a red box is not shown before acquisition.

When C2 or C3 lens is switched off, a tile set can be acquired with an unrelated intensity value

On a system with three condenser lenses, you can switch off either C2 or C3. Maps does not detect which lens is switched off. You can create a tile set when C2 (or C3) is switched off and then acquire it when C3 (or C2) is switched off, which results in an unrelated intensity value.

Workaround: Complete tile set creation and acquisition before switching to another lens.

Maps might prevent or fail acquisition in other applications

Maps does not check if other applications use a detector or camera. The detectors and cameras are retracted, except the one that the tile set is going to be acquired with. As soon as Maps starts acquisition, it detects that another application is using the detector or camera (and does not start acquisition). This behavior might prevent acquisition from being done in the other applications.

Workaround: Do not start acquisition of a tile set when another application is using a detector or camera.

Intensity or illuminated area is not set to the microscope

When you switch off C2 or C3 on a system with three condenser lenses, the intensity is not set to the Microscope for a tile set that was created when you switched on both C2 and C3.

Workaround: Switch on both C2 and C3 and then click **To Microscope**. When you switch on C2 and C3 on a system with three condenser lenses, the illuminated area is not set to the Microscope for a tile set that was created when you switched off C2 or C3.

Workaround: Switch off C2 or C3 and then click **To Microscope**.

Errors occur when TEM server, TIA and Digital Micrograph are not running

Before Maps can be started, the TEM server, TIA, and Digital Micrograph (when applicable) must be running. When Digital Micrograph, TIA, or the TEM server crashes or is stopped, Maps might crash, show error messages, or behave incorrectly. Examples include the following:

- TEM plug-in generated an exception during initialization.
- Object reference not set to an instance of an object.
- FOV size incorrect.
- Maps crash during acquisition as soon as TIA crashes or is stopped.

Workarounds:

- When using the Camera ocx in PEOUI, the camera combo box is empty, TIA has crashed, or is not running. Restart TIA and ensure that the camera combo box lists your cameras, and then restart Maps.
- When the TEM server has crashed or is not running, restart the TEM server, start TIA, and then restart Maps.
- For camera usage that depends on Digital Micrograph, the startup order is:
 - 1. TEM server
 - 2. TIA
 - 3. Digital Micrograph

When TEM server is restarted, restart TIA and then Digital Micrograph. When TIA is restarted, restart Digital Micrograph.

• After changing the high-tension value or executing camera calibrations, restart Maps.

Maps becomes unresponsive

This can happen when the stage is not working properly.

Workaround: In PEOUI, verify that the stage is working properly and restart Maps.

Maps becomes unresponsive when deleting a large amount of large tile sets

When you delete a large amount of large tile sets, Maps becomes unresponsive during this operation. It can take extra time to delete large tile sets. Do not end the program, or you will interrupt the operation and need to start over.

When Maps is running offline, sometimes Intensity is displayed instead of Illuminated Area

When Maps is running offline, Maps does not know whether the TEM tile set was acquired on a system with two or three condenser lenses. Maps incorrectly shows Intensity for a TEM tile set that is acquired on a system with three condenser lenses.

The disk space usage estimate may be incorrect

The total disk space required is only an estimate and might vary. Please allow extra space for large data sets.

Images that include part of a grid bar are slightly darker

In a tile set, the images that contain a part of a grid bar are slightly darker due to a difference in image contrast.

Workaround: Avoid images with grid bars.

New tile sets are deselected in the Job Queue after a Magnification

Alignment

When Magnification Alignment is performed on a tile, empty tile set jobs in the **Job Queue** might become deselected causing them to not get acquired.

Workaround: Go to the Job Queue and select the check boxes of the applicable tile sets.

Microscope FOV is not updated when the Stage Axis Angle Correction is updated

When the Stage Axis Angle Correction is updated in the **Settings** menu, the microscope Field of View (FOV) is not updated.

Workaround: Change the magnification to update the microscope FOV for magnification, or move the stage to update the microscope FOV for location.

Microscope FOV is not updated when the STEM Scan Orientation Correction is updated

When the STEM Scan Orientation Correction is updated in the **Settings** menu, the microscope FOV is not updated.

Workaround: Change the STEM FOV or magnification.

Maps software hangs when acquiring on a microscope with a Gatan camera or DigitalMicrograph

On a microscope with a Gatan camera or Gatan DigitalMicrograph running, Maps software can sometimes freeze during acquisition.

Workaround: Close Maps software and restart Maps.

In EFTEM mode, images can be acquired only with maximum camera resolution

In EFTEM mode, images can be acquired only with maximum camera resolution, such as binning 1.

In EFTEM mode, stitched images are incorrect

When tile sets acquired in EFTEM mode are stitched, the size of the stitched output is incorrect and tiles are incorrectly joined.

Mineralogy known issues

Mineralogy: The reporting page is not reachable due to WSL

The reporting page may fail to work properly if the Windows Subsystem for Linux (WSL) is not started.

This issue can occur in situations where the WSL service is not running or encounters an error during the startup process.

To resolve this issue, please follow the steps outlined below:

- 1. Open a command window as an administrator.
- 2. Type the command: tasklist /svc /fi "imagename eq svchost.exe" | findstr LxssManager and note the returned PID.
- 3. Run the Task Manager as an administrator.
- 4. Go to the "Details" tab in Task Manager.
- 5. Search for the svchost.exe process containing the previously noted PID.
- 6. Right-click svchost.exe and select "End Process Tree."
- 7. Return to the command window opened as an administrator and type **'wsl -shutdown'**, followed by **'wsl'**.
- 8. Ensure that a command line is returned without any errors.

By following these steps, you can resolve the issue related to the reporting page not working due to WSL not being started. This workaround will help restart the necessary services and ensure the proper functioning of the reporting page.

Mineralogy: Bruker Esprit error on long running acquisition

During the process of running a long acquisition with the Bruker Esprit tool, there is a known issue where an error message may appear stating 'An expected error occured on server. Consider to save your current and restart Esprit'. In some cases, this error message can cause the acquisition process to stop unexpectedly on the Support PC. A workaround is currently not available.

Release Notes

Be sure to read the release notes for the latest software version of the Maps software to learn about new features and changes in functionality.

Maps 3.31.1 Release Notes

Maps 3.30.2 Release Notes

Maps 3.30.1 Release Notes

Maps 3.29.1 Release Notes

Maps 3.28.1 Release Notes

Maps 3.26.1 Release Notes

Maps 3.25.1 Release Notes

Maps 3.31.1 Release Notes

This document specifies changes made to the Thermo Scientific[™] Maps 3.31.1 software application.

New Features

SEM Analytics

Maps can now acquire EDS data through a software application interface on a SEM/SDB microscope. This functionality is available on systems with ChemiSEM support, including Apreo ChemiSEM, Axia ChemiSEM, and Quattro ESEM.

See SEM Analytics for more information.

TEM: Support for TEM 3.22 and TEM 2.22

Maps now supports TEM server versions 3.17-3.22 for Titan and 2.17-2.22 for Talos.

Bug Fixes

The following bugs were fixed in this release of the Maps software.

Bug ID	Description
MAPS-11821	The Job Control button located in the status bar now supports a toggle pattern with a chevron icon. When you click the Job Control button repeatedly, the Job Control screen toggles between expanded and collapsed states. Previously, you could not collapse the screen by clicking the Job Control button.
MAPS-12455	A resolution of 512 x 512 pixels is now available on Phenom microscopes.
MAPS-12655	Maps now removes temporary acquisition files on Phenom microscopes when the files are no longer needed. This previously caused problems during long acquisitions of large tile sets, including errors and crashes when a system's temporary space filled up.

Maps 3.30.2 Release Notes

This document specifies changes made to the Thermo Scientific[™] Maps 3.30.2 software application

Known Issues

See Known issues for known issues introduced with this version of Maps software.

Maps 3.30.1 Release Notes

This document specifies changes made to the Thermo Scientific[™] Maps 3.30.1 software application.

New Features

Phenom: Default license includes basic features

Maps on your Phenom microscope now includes basic features without a need for product registration and activation. The basic features allow you to control the Phenom microscopes; create and edit projects; create and acquire tile sets; and perform stitching of acquired data.

TEM: Support for TEM 3.21 and TEM 2.21

Maps now supports TEM server versions 3.16-3.21 for Titan and 2.16-2.21 for Talos.

Known Issues

See Known issues for known issues introduced with this version of Maps software.

Bug Fixes

The following bugs were fixed in this release of the Maps software.

Bug ID	Description
MAPS- 11391	TEM: Autofunctions now run correctly in EFTEM mode on all supported cameras. Previously, the autofunctions failed to run on some cameras.
MAPS- 11407	Maps no longer freezes when exporting a group layer to a new or existing project.
MAPS- 11418	TEM: The first tile is now acquired in a more precise position after eucentric height autofunction runs on the first tile. Previously, it could be visibly shifted relatively to the rest of the tile set in some use cases.
MAPS- 11478	Leica XLEF import: Confocal stacks are now oriented consistently with overview images.

Bug ID	Description
MAPS- 11479	The color indicator for fluorescence channels now shows a color gradient based on a lookup table. Previously, the color indicator showed as white if the lookup table was used.
MAPS- 11793	TEM: The Data Exchange Service runs correctly again, thus allowing data exchange between the latest versions of Maps and Tomography. This service was previously broken by a library update conflict, which caused the data exchange to fail.

Maps 3.29.1 Release Notes

This document specifies changes made to the Thermo Scientific[™] Maps 3.29.1 software application.

Product News

Mineralogy: Revised product name

The product name of the Maps Mineralogy plugin application has been revised to "Maps Min". The new product name is effective as of this release.

New Features

Mineralogy: Bruker XFlash® 7 EDS Detector support

Maps Min now supports the Bruker XFlash 7 EDS detector. This new option enables you to configure and utilize the advanced capabilities of the Bruker 7 Series detector for your elemental mapping analysis. You can select this detector when you configure the Maps software for mineralogy acquisition.

See Configure Mineralogy Acquisition for more information.

Mineralogy: Maps Min Reporting plugin application

- You now have the ability to import a reporting list from another project and integrate it into your current project.
- You can now easily delete project and associated data. Select a project, click the delete button, confirm the deletion, and then project will be permanently removed.
- A new report named 'Mineral Association' is available and is visible in tabular form. This report provides an analysis of the minerals grains relationships and interactions through the perimeter proportion between grains. This report also includes particle filter; data set selection; reporting list selection and combination by replicates; and can it can be exported in .csv format.
- Previously, you were limited to publishing particles with BSE (Backscattered Electron) data less than 20 MB. Now, you can publish and view large particles without any size limitations. This enhancement ensures compatibility with image sizes up to 50-82K pixels, allowing you to work with high-resolution particles.

• In the Reporting user interface, you can open the dataset and select the Particle View report. All particles (including those exceeding 20 MB), will be displayed, thus, providing a comprehensive view of the dataset.

Phenom: Advanced tab control enhancements

The following enhancements have been added to the advanced tile acquisition controls.

- High Voltage: The preset options now include 20 kV.
- Detector: The options now include the SED detector.

See Phenom Tile Set Tab for more information.

Phenom: New advanced tab controls

There are new advanced tile acquisition controls in the user interface.

- Vacuum: Use this control to specify the required vacuum pressure during acquisition.
- Acquisition Order: Use this control to determine the order in which the contained tiles are acquired.

See Phenom Tile Set Tab for more information.

Phenom: Post-acquistion actions

Post-acquisiton actions are now available for Phenom. After all jobs are completed, you can return the microscope to HiVac mode, move the sample to the NavCam, and put the microscope in the standby operational state.

See Phenom Post-Acquisition Actions for more information.

SEM: Support for Quattro 35.0

Maps now fully supports Quattro 35.0. The corresponding xT server version is automatically recognized during Maps installation.

TEM: Support for TEM 3.20 and TEM 2.20

Maps now supports TEM server versions 3.15-3.20 for Titan and 2.15-2.20 for Talos.

Updated system requirements for installation

- Windows compatibility: The Maps application can now run on Windows 11.
- .Net Framework: The required version of .Net Framework has been updated to 4.8, which will be now installed during the Maps installation process if it is not already installed on the system.

Known Issues

See Known issues for known issues introduced with this version of Maps software.

Bug Fixes

The following bugs were fixed in this release of the Maps software.

Bug ID	Description
MAPS-10804	Stitching of tile sets with some tiles deactivated no longer causes an exception to occur.
MAPS-11136	Importing TIF files with invalid metadata no longer causes Maps to crash.
MAPS-7863	Mineralogy: Detector names in the xT quadrants are now being retained correctly.
MAPS-10158	Mineralogy: Exporting size distribution to CSV now produces correct ranges for the X-axis depending on the dynamic or fixed bins.
MAPS-10680	Mineralogy: The Mineralogy Reporting server graphs are now accurate.
MAPS-10901	Mineralogy: You can now export images with Porosity overlay on BSE in Nanomin.
MAPS-11034	Mineralogy: The Sort by Mineral or Element proportion are now functional within the Particle View report.
MAPS-10967	STEM: The beam convergence angle now displays in a more user friendly manner, rounded to two decimal places and with units displayed after the angle value.

Maps 3.28.1 Release Notes

This document specifies changes made to the Thermo Scientific[™] Maps 3.28.1 software application.

New Features

Bright Phase Search acquisition

The Maps application now supports Bright Phase Search acquisition, which produces faster acquisition and results within the web-based Maps Mineralogy Reporting application. Bright Phase Search acquisition is available in either Grid mode or Centroid mode (beta).

See Optional: Bright Phase Search Acquisition Workflow for more information.

Maps improvement program

Maps has introduced an improvement program that will send your anonymous usage data back to the Maps team for product improvement purposes. If desired, you can opt out of this program after you first launch the Maps application and accept the End User License Agreement (EULA). Go to the application settings and deselect the "Allow usage logging" setting under the Improvement Program section.

See Options menu for more information.

Mineralogy: Bruker and Nanomin connection states

The Analytics connection indicator has been expanded to separate the Bruker and Nanomin status icons, which indicate the online connection states between the Maps application and the microscope. This change provides more status support by communicating if a connection is in use, if a connection is attempting to connect, or if a connection has timed out.

See Microscope connection states for more information.

Mineralogy: Improved Nanomin performance for large datasets

Pyramid rendering has been updated in the Nanomin application to improve panning, zooming, and overlay rendering speeds while you navigate large datasets.

Mineralogy: Maps Mineralogy Reporting application

- A desktop shortcut is now available on your desktop soon after you successfully install the Maps Mineralogy Reporting application. Double-click the desktop shortcut and the application will launch the user interface within your browser.
- You can now generate Locking and Liberation reports by Area % within the Maps Mineralogy Reporting application. This means that you can choose between Weight % or Area % as the output type for increased flexibility and precision.
- When you select either a locking or liberation report while using the Maps Mineralogy Reporting application, the initial option within the Liberation Type field has been changed from "Area %" to "Composition" to clearly identify particle composition.
- You can now view product information within the Maps Mineralogy Reporting application. This information includes the version number of the application, server, and client for you to share with Thermo Fisher Scientific while troubleshooting a software issue.

Mineralogy: Plugin enhancements

- New and improved data tracking of Mineralogy tile set metadata.
- You can now view Mineralogy tile set parameters offline.
- Mining tags have moved to a new tab.
- Analytics tab for mineralogy has been renamed to "Mineralogy Acquisition" tab.

Phenom: Simplified connectivity

The microscope connection process for the Phenom has been simplified, so that the connection dialog no longer requires you to enter your user name and password. The Maps application contains a security tool that allows it to connect to any available Phenom within your network.

See Phenom Connectivity for more information.

TEM: Added support for TEM 3.19 and TEM 2.19

Maps 3.28.1 supports TEM server versions 3.14-3.19 for Titan and 2.14-2.19 for Talos.

TEM: Standalone Velox plugin

Acquisition of EDS tile sets through Velox is now available as a standalone plugin, so connectivity to Velox and the acquisition process is more robust.

See Activate Velox Software for more information.

TEM: Tiling correction

You can use the Tiling Correction beta feature to achieve better tile alignment during acquisition.

See TEM Tiling Corrections and Tiling Corrections Quality for more information.

Known Issues

See Known issues for known issues introduced with this version of Maps software.

Bug Fixes

The following bugs were fixed in this release of the Maps software.

Bug ID	Description
MAPS-7863	Mineralogy: Detector names in the xT quadrants are now being retained correctly.
MAPS-10158	Mineralogy: Exporting size distribution to CSV now produces correct ranges for the X-axis depending on the dynamic or fixed bins.
MAPS-10774	The beam convergence angle on STEM is now displayed in mrad units instead of degrees.
MAPS-10781	Taking a STEM snapshot on a system without the HAADF detector no longer causes an error. Previously, the HAADF detector was used as a default detector for snapshots on all systems, regardless of the actual configuration.

Maps 3.26.1 Release Notes

This document specifies changes made to the Thermo Scientific[™] Maps 3.26.1 software application.

New Features

Array Tomography workflow controls

The Array Tomography workflow has been streamlined with new and improved controls including:

- New Line Tool to order sections
- New keyboard shortcuts for specific steps to expedite workflow
- Increased flexibility for section previews creation
- Revised refinement step to improve viewing and editing of results
- New "stitch all" option during tile set array acquisitions

See the Array Tomography Workflow for more details.

CRYO: Arctis data features

This bundle of new features allow you to import and view Arctis Cryo-Plasma-FIB data in the Maps software application on TEM systems.

See Import and view Arctis data for more information.

TEM: OptiSTEM autofunction automatically restores last known good state

When the OptiSTEM autofunction fails to converge or the quality of the result is poor, the Maps software application automatically restores the focus and stigmator values to the last known good state.

TEM: Added support for TEM 3.17 and TEM 2.17

Maps now supports the latest TEM server versions for Titan (TEM 3.17) and Talos (TEM 2.17).

Documentation Update

A comprehensive Phenom section has been added to the Maps software application user guide. This new section describes the Phenom connectivity, settings, operational states, and user interface elements.

See Phenom for more information.

Bug Fixes

The following bugs were fixed in this release of the Maps software.

Bug ID	Description
MAPS-9687	TEM: The position of the crosshair and cross-correlation peak annotations is now correct. Previously, the annotations in the intermediate cross-correlation images during the execution of autofunctions were not positioned correctly when a camera with a rectangular image resolution was used (for exmaple, a Gatan K3 camera).
MAPS- 10276	TEM: Camera settings are now properly applied when used in the autofunction templates on TEM and EFTEM. Previously, autofunctions used the default camera settings and ignored settings like Frames Summed for Ceta-F camera or Electron Counting Mode for Falcon camera.

Maps 3.25.1 Release Notes

This document specifies changes made to the Thermo Scientific[™] Maps 3.25.1 software application.

New Features

Animation for viewer functions

The "Center View on Sample," "Center View on Selection," and "Go to Microscope FOV" viewer functions in the toolbar now animate pan and zoom to help keep you oriented in the sample space. If desired, you can deactivate this feature under the "Animate Pan & Zoom" View menu item.

See Toolbar and View menu for more information.

EFTEM slit control

Slit insertion state and energy-selecting slit width can now be set directly on a tile set. Slit parameters are also now part of EFTEM tile set templates, and are propagated from Tomography presets when importing the presets to Maps.

See Advanced group, EFTEM mode for more information.

Phenom SEM support

This version of Maps now supports the Thermo Scientific[™] Phenom[™] desktop SEM series. As a Phenom user, you will be able to access all the Maps features that are currently available on your other Thermo Fisher SEM, SDB, and TEM systems. This integrated support also allows you to expand the correlative workflows within Maps from desktop SEMs to high-end TEM systems for a variety of scientific experiments. A Maps Phenom license is required, so please contact your Thermo Fisher account manager for additional information.

Note: The Maps software application user guide will be updated in a future release to include the Phenom functionality.

Project History search bar

The Project History window now has a search bar available. When you use this search bar, it immediately filters the project history and provides results that match either the Project Name or Project Description fields.

See Open a Project for more information.

TEM/EFTEM: Automated image filtering for autofunctions

TEM and EFTEM autofunctions support a new setting called "Use Image Filtering". You can use this new setting to improve the reliability and accuracy of these autofunctions.

See Automated Image Filtering for more information.

TEM: Added support for TEM 3.16 and TEM 2.16

Maps now supports the latest TEM server versions for Titan (TEM 3.16) and Talos (TEM 2.16).

Bug Fixes

The following bugs were fixed in this release of the Maps software.

Bug ID	Description
MAPS-9271	TEM: Correct milling angle of a lamella is now displayed in projects imported from SDB to TEM. Previously, the sample pre-tilt angle used on SDB was either ignored, or assumed to be 35 degrees (Maps 3.24), causing the incorrect milling angle to be computed on TEM. Maps on TEM systems now reads the sample pre-tilt angle that was used on the SDB system and adjusts the computed milling angle accordingly.
MAPS-9274	TEM: Falcon 4 camera settings now work properly in EFTEM mode. Previously, the camera mode was not properly initialized based on microscope state, and Linear mode was used for acquisition even if Counted mode was specified on the tile set.
MAPS-9442	TEM: Templates created from a new tile set can now be used immediately as an autofunction template within the new tile set. Previously, there was a synchronization issue that caused the new template to appear only after switching to another layer in the layer tree and then back to the tile set.

License Information

The following licensed content is included in Maps 3.32.1.

Accord.NET Framework

The Accord.NET Framework Copyright © 2009-2017, Accord.NET Authors <authors@accord-framework.net>

This library is free software; you can redistribute it and/or modify it under the terms of the GNU Lesser General Public License as published by the Free Software Foundation; either version 2.1 of the License, or (at your option) any later version.

The copyright holders provide no reassurances that the source code provided does not infringe any patent, copyright, or any other intellectual property rights of third parties. The copyright holders disclaim any liability to any recipient for claims brought against recipient by any third party for infringement of that parties intellectual property rights.

This library is distributed in the hope that it will be useful, but WITHOUT ANY WARRANTY; without even the implied warranty of MERCHANTABILITY or FITNESS FOR A PARTICULAR PURPOSE.

See the GNU Lesser General Public License for more details.

You should have received a copy of the GNU Lesser General Public License along with this program. If not, see http://www.gnu.org/licenses/>.

Some modules of this library are distributed under the GPL, and not the LGPL. Those modules have been marked with .GPL.dll in their name. All relevant licensing information should also be accompanying each respective source code. Other extra modules may carry other restrictions.

Please see relevant source headers for details.

The documentation sections contains references to Wikipedia and is thus entirely available under the Creative Commons Attribution/Share-Alike License. All referenced sections include a hyperlink to the original article page. If you think you have found something which has not been properly linked, please let me know so it can be corrected.

Overall framework architecture and style was greatly inspired by AForge.NET. In May 2015, this project has been merged with the AForge.NET framework since public support for AForge.NET has ended. The original AForge.NET Framework is a copyrighted work by Andrew Kirillov, altough developed and shared under the same LGPL license.

This library also contains routines adapted from Lutz Roeder's Mapack for .NET, LAPACK and Jama routines. Jama has been released under the public domain. LAPACK has been released under a compatible 3-clause BSD license, listed below. Original ArrayDataView code is copyright © 2004 Mihail Stefanov, licensed under the LGPL with permission of the original author. Original SortableBindingList source code is copyright © 2008 Tim Van Wassenhove, licensed under the LGPL with permission of the original author. Statistics module contains code for computing the Kolmogorov-Smirnov distribution based on original code by Richard Simard, redistributed under the LGPL with permission of the original author. Imaging module contains code for Gabor filter based on the original code by Max Bügler, redistributed under the LGPL with permission of the original author.

The Augmented Lagrangian Solver implementation is based on the NLopt Numerical Optimization Library, copyright © 2008 Steven G. Johnson, distributed under the MIT license. Some image filters have been contributed by Diego Catalano, distributed under the LGPL license and included with permission of the original author. Mathematics module includes the Kinematic joint models by Rémy Dispagne, licensed under the LGPL with permission of the original author. The Haar Object detector contains contributions from Darko Jurić, added with permission of the original author.

The Accord.Vision module includes definitions for Haar Cascades created by Modesto Castrillon-Santana, distributed under the LGPL with permission of the original author. The SURF interest points detector implementation is based on code by Christopher Evans, licensed under the LGPL with permission of the original author. The FAST corners detector has been created using a code generation tool created by Edward Rosten, distributed under a compatible 3-clause BSD license, listed below. The Camshift tracker implementation is based on the FaceIt ActionScript Camshift tracker, copyright © 2009 Benjamin Jung, distributed under the MIT license, listed below. The Project Marilena is copyright © 2008 Masakazu Ohtsuka, distributed under a compatible 2-clause BSD license, listed below. See actual sources for details.

Below are the licenses applying to specific sections of the code. Those are not applicable to the work as a whole (the Framework), but to the specifics of each portion of the framework. See actual sources for details.

Circular Statistics Toolbox (Directional Statistics)

Copyright © Philipp Berens, 2006-2012, All rights reserved.

Redistribution and use in source and binary forms, with or without modification, are permitted provided that the following conditions are met:

- 1. Redistributions of source code must retain the above copyright notice, this list of conditions and the following disclaimer.
- Redistributions in binary form must reproduce the above copyright notice, this list of conditions and the following disclaimer in the documentation and/or other materials provided with the distribution

THIS SOFTWARE IS PROVIDED BY THE COPYRIGHT HOLDERS AND CONTRIBUTORS "AS IS" AND ANY EXPRESS OR IMPLIED WARRANTIES, INCLUDING, BUT NOT LIMITED TO, THE IMPLIED WARRANTIES OF MERCHANTABILITY AND FITNESS FOR A PARTICULAR PURPOSE ARE DISCLAIMED. IN NO EVENT SHALL THE COPYRIGHT OWNER OR CONTRIBUTORS BE LIABLE FOR ANY DIRECT, INDIRECT, INCIDENTAL, SPECIAL, EXEMPLARY, OR CONSEQUENTIAL DAMAGES (INCLUDING, BUT NOT LIMITED TO, PROCUREMENT OF SUBSTITUTE GOODS OR SERVICES; LOSS OF USE, DATA, OR PROFITS; OR BUSINESS INTERRUPTION) HOWEVER CAUSED AND ON ANY THEORY OF LIABILITY, WHETHER IN CONTRACT, STRICT LIABILITY, OR TORT (INCLUDING NEGLIGENCE OR OTHERWISE) ARISING IN ANY WAY OUT OF THE USE OF THIS SOFTWARE, EVEN IF ADVISED OF THE POSSIBILITY OF SUCH DAMAGE.

LIBLINEAR - A library for large linear classification

Copyright © 2007-2011 The LIBLINEAR Project. All rights reserved.

Redistribution and use in source and binary forms, with or without modification, are permitted provided that the following conditions are met:

- 1. Redistributions of source code must retain the above copyright notice, this list of conditions and the following disclaimer.
- 2. Redistributions in binary form must reproduce the above copyright notice, this list of conditions and the following disclaimer in the documentation and/or other materials provided with the distribution.
- 3. Neither name of copyright holders nor the names of its contributors may be used to endorse or promote products derived from this software without specific prior written permission.

THIS SOFTWARE IS PROVIDED BY THE COPYRIGHT HOLDERS AND CONTRIBUTORS "AS IS" AND ANY EXPRESS OR IMPLIED WARRANTIES, INCLUDING, BUT NOT LIMITED TO, THE IMPLIED WARRANTIES OF MERCHANTABILITY AND FITNESS FOR A PARTICULAR PURPOSE ARE DISCLAIMED. IN NO EVENT SHALL THE REGENTS OR CONTRIBUTORS BE LIABLE FOR ANY DIRECT, INDIRECT, INCIDENTAL, SPECIAL, EXEMPLARY, OR CONSEQUENTIAL DAMAGES (INCLUDING, BUT NOT LIMITED TO, PROCUREMENT OF SUBSTITUTE GOODS OR SERVICES; LOSS OF USE, DATA, OR PROFITS; OR BUSINESS INTERRUPTION) HOWEVER CAUSED AND ON ANY THEORY OF LIABILITY, WHETHER IN CONTRACT, STRICT LIABILITY, OR TORT (INCLUDING NEGLIGENCE OR OTHERWISE) ARISING IN ANY WAY OUT OF THE USE OF THIS SOFTWARE, EVEN IF ADVISED OF THE POSSIBILITY OF SUCH DAMAGE.

LibSVM & liblinear library for Support Vector Machines

Copyright © 2000-2014 Chih-Chung Chang and Chih-Jen Lin, All rights reserved.

Redistribution and use in source and binary forms, with or without modification, are permitted provided that the following conditions are met:

- 1. Redistributions of source code must retain the above copyright notice, this list of conditions and the following disclaimer.
- Redistributions in binary form must reproduce the above copyright notice, this list of conditions and the following disclaimer in the documentation and/or other materials provided with the distribution.
- 3. Neither name of copyright holders nor the names of its contributors may be used to endorse or promote products derived from this software without specific prior written permission.

THIS SOFTWARE IS PROVIDED BY THE COPYRIGHT HOLDERS AND CONTRIBUTORS "AS IS" AND ANY EXPRESS OR IMPLIED WARRANTIES, INCLUDING, BUT NOT LIMITED TO, THE IMPLIED WARRANTIES OF MERCHANTABILITY AND FITNESS FOR A PARTICULAR PURPOSE ARE DISCLAIMED. IN NO EVENT SHALL THE REGENTS OR CONTRIBUTORS BE LIABLE FOR ANY DIRECT, INDIRECT, INCIDENTAL, SPECIAL, EXEMPLARY, OR CONSEQUENTIAL DAMAGES (INCLUDING, BUT NOT LIMITED TO, PROCUREMENT OF SUBSTITUTE GOODS OR SERVICES; LOSS OF USE, DATA, OR PROFITS; OR BUSINESS INTERRUPTION) HOWEVER CAUSED AND ON ANY THEORY OF LIABILITY, WHETHER IN CONTRACT, STRICT LIABILITY, OR TORT (INCLUDING NEGLIGENCE OR OTHERWISE) ARISING IN ANY WAY OUT OF THE USE OF THIS SOFTWARE, EVEN IF ADVISED OF THE POSSIBILITY OF SUCH DAMAGE.

LumenWorks.Framework.IO.CSV.CsvReader - A Fast CSV Reader

 $Copyright @ Sebastien \ Lorion, \ 2005-2011, \ All \ rights \ reserved.$

MIT license (http://en.wikipedia.org/wiki/MIT_License)

Permission is hereby granted, free of charge, to any person obtaining a copy of this software and associated documentation files (the "Software"), to deal in the Software without restriction, including without limitation the rights to use, copy, modify, merge, publish, distribute, sublicense, and/or sell copies of the Software, and to permit persons to whom the Software is furnished to do so, subject to the following conditions:

The above copyright notice and this permission notice shall be included in all copies or substantial portions of the Software.

THE SOFTWARE IS PROVIDED "AS IS", WITHOUT WARRANTY OF ANY KIND, EXPRESS OR IMPLIED, INCLUDING BUT NOT LIMITED TO THE WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE AND NONINFRINGEMENT. IN NO EVENT SHALL THE AUTHORS OR COPYRIGHT HOLDERS BE LIABLE FOR ANY CLAIM, DAMAGES OR OTHER LIABILITY, WHETHER IN AN ACTION OF CONTRACT, TORT OR OTHERWISE, ARISING FROM, OUT OF OR IN CONNECTION WITH THE SOFTWARE OR THE USE OR OTHER DEALINGS IN THE SOFTWARE.

MathExtension: Mathematical extensions and structures for the .NET framework

MIT License

Copyright © 2015 AJ Richardson

Permission is hereby granted, free of charge, to any person obtaining a copy of this software and associated documentation files (the "Software"), to deal in the Software without restriction, including without limitation the rights to use, copy, modify, merge, publish, distribute, sublicense, and/or sell copies of the Software, and to permit persons to whom the Software is furnished to do so, subject to the following conditions:

The above copyright notice and this permission notice shall be included in all copies or substantial portions of the Software.

THE SOFTWARE IS PROVIDED "AS IS", WITHOUT WARRANTY OF ANY KIND, EXPRESS OR IMPLIED, INCLUDING BUT NOT LIMITED TO THE WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE AND NONINFRINGEMENT. IN NO EVENT SHALL THE AUTHORS OR COPYRIGHT HOLDERS BE LIABLE FOR ANY CLAIM, DAMAGES OR OTHER LIABILITY, WHETHER IN AN ACTION OF CONTRACT, TORT OR OTHERWISE, ARISING FROM, OUT OF OR IN CONNECTION WITH THE SOFTWARE OR THE USE OR OTHER DEALINGS IN THE SOFTWARE.

Munkres algorithm - The Hungarian method for solving the assignment problem

MIT License

Copyright © 2000 Robert A. Pilgrim, Murray State University, Dept. of Computer Science & Information Systems, Murray, Kentucky

Permission is hereby granted, free of charge, to any person obtaining a copy of this software and associated documentation files (the "Software"), to deal in the Software without restriction, including without limitation the rights to use, copy, modify, merge, publish, distribute, sublicense, and/or sell copies of the Software, and to permit persons to whom the Software is furnished to do so, subject to the following conditions:

The above copyright notice and this permission notice shall be included in all copies or substantial portions of the Software.

THE SOFTWARE IS PROVIDED "AS IS", WITHOUT WARRANTY OF ANY KIND, EXPRESS OR IMPLIED, INCLUDING BUT NOT LIMITED TO THE WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE AND NONINFRINGEMENT. IN NO EVENT SHALL THE AUTHORS OR COPYRIGHT HOLDERS BE LIABLE FOR ANY CLAIM, DAMAGES OR OTHER LIABILITY, WHETHER IN AN ACTION OF CONTRACT, TORT OR OTHERWISE, ARISING FROM, OUT OF OR IN CONNECTION WITH THE SOFTWARE OR THE USE OR OTHER DEALINGS IN THE SOFTWARE.

Partitioning around medoids and Voronoi iteration algorithm for the K-Medoids clustering: algorithm implementation and unit tests

MIT License

Copyright © 2017 Ivan Pizhenko

Permission is hereby granted, free of charge, to any person obtaining a copy of this software and associated documentation files (the "Software"), to deal in the Software without restriction, including without limitation the rights to use, copy, modify, merge, publish, distribute, sublicense, and/or sell copies of the Software, and to permit persons to whom the Software is furnished to do so, subject to the following conditions:

The above copyright notice and this permission notice shall be included in all copies or substantial portions of the Software.

THE SOFTWARE IS PROVIDED "AS IS", WITHOUT WARRANTY OF ANY KIND, EXPRESS OR IMPLIED, INCLUDING BUT NOT LIMITED TO THE WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE AND NONINFRINGEMENT. IN NO EVENT SHALL THE AUTHORS OR COPYRIGHT HOLDERS BE LIABLE FOR ANY CLAIM, DAMAGES OR OTHER LIABILITY, WHETHER IN AN ACTION OF CONTRACT, TORT OR OTHERWISE, ARISING FROM, OUT OF OR IN CONNECTION WITH THE SOFTWARE OR THE USE OR OTHER DEALINGS IN THE SOFTWARE.

NLopt Numerical Optimization Library - Copyright © 2008-2014 Steven G. Johnson

Copyright © 2007-2011 Massachusetts Institute of Technology

Permission is hereby granted, free of charge, to any person obtaining a copy of this software and associated documentation files (the "Software"), to deal in the Software without restriction, including without limitation the rights to use, copy, modify, merge, publish, distribute, sublicense, and/or sell copies of the Software, and to permit persons to whom the Software is furnished to do so, subject to the following conditions:

The above copyright notice and this permission notice shall be included in all copies or substantial portions of the Software.

THE SOFTWARE IS PROVIDED "AS IS", WITHOUT WARRANTY OF ANY KIND, EXPRESS OR IMPLIED, INCLUDING BUT NOT LIMITED TO THE WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE AND NONINFRINGEMENT. IN NO EVENT SHALL THE AUTHORS OR COPYRIGHT HOLDERS BE LIABLE FOR ANY CLAIM, DAMAGES OR OTHER LIABILITY, WHETHER IN AN ACTION OF CONTRACT, TORT OR OTHERWISE, ARISING FROM, OUT OF OR IN CONNECTION WITH THE SOFTWARE OR THE USE OR OTHER DEALINGS IN THE SOFTWARE.

Peter Kovesi's Computer Vision functions - Copyright © 1995-2010 Peter Kovesi

Copyright © 1995-2010 Peter Kovesi, Centre for Exploration Targeting, School of Earth and Environment, The University of Western Australia

Permission is hereby granted, free of charge, to any person obtaining a copy of this software and associated documentation files (the "Software"), to deal in the Software without restriction, including without limitation the rights to use, copy, modify, merge, publish, distribute, sublicense, and/or sell copies of the Software, and to permit persons to whom the Software is furnished to do so, subject to the following conditions:

The above copyright notice and this permission notice shall be included in all copies or substantial portions of the Software.

The software is provided "as is", without warranty of any kind, express or implied, including but not limited to the warranties of merchantability, fitness for a particular purpose and noninfringement. In no event shall the authors or copyright holders be liable for any claim, damages or other liability, whether in an action of contract, tort or otherwise, arising from, out of or in connection with the software or the use or other dealings in the software.

Snowball: A language for stemming algorithms

Copyright © 2001, Dr Martin Porter, Copyright © 2004,2005, Richard Boulton, All rights reserved.

Redistribution and use in source and binary forms, with or without modification, are permitted provided that the following conditions are met:

- 1. Redistributions of source code must retain the above copyright notice, this list of conditions and the following disclaimer.
- Redistributions in binary form must reproduce the above copyright notice, this list of conditions and the following disclaimer in the documentation and/or other materials provided with the distribution.
- 3. Neither the name of the Snowball project nor the names of its contributors may be used to endorse or promote products derived from this software without specific prior written permission.

THIS SOFTWARE IS PROVIDED BY THE COPYRIGHT HOLDERS AND CONTRIBUTORS "AS IS" AND ANY EXPRESS OR IMPLIED WARRANTIES, INCLUDING, BUT NOT LIMITED TO, THE IMPLIED WARRANTIES OF MERCHANTABILITY AND FITNESS FOR A PARTICULAR PURPOSE ARE DISCLAIMED. IN NO EVENT SHALL THE COPYRIGHT OWNER OR CONTRIBUTORS BE LIABLE FOR ANY DIRECT, INDIRECT, INCIDENTAL, SPECIAL, EXEMPLARY, OR CONSEQUENTIAL DAMAGES (INCLUDING, BUT NOT LIMITED TO, PROCUREMENT OF SUBSTITUTE GOODS OR SERVICES; LOSS OF USE, DATA, OR PROFITS; OR BUSINESS INTERRUPTION) HOWEVER CAUSED AND ON ANY THEORY OF LIABILITY, WHETHER IN CONTRACT, STRICT LIABILITY, OR TORT (INCLUDING NEGLIGENCE OR OTHERWISE) ARISING IN ANY WAY OUT OF THE USE OF THIS SOFTWARE, EVEN IF ADVISED OF THE POSSIBILITY OF SUCH DAMAGE.

Statistical Quantiles: implementation and unit tests

MIT License

Copyright © 2017 Ivan Pizhenko

Permission is hereby granted, free of charge, to any person obtaining a copy of this software and associated documentation files (the "Software"), to deal in the Software without restriction, including without limitation the rights to use, copy, modify, merge, publish, distribute, sublicense, and/or sell copies of the Software, and to permit persons to whom the Software is furnished to do so, subject to the following conditions:

The above copyright notice and this permission notice shall be included in all copies or substantial portions of the Software.

THE SOFTWARE IS PROVIDED "AS IS", WITHOUT WARRANTY OF ANY KIND, EXPRESS OR IMPLIED, INCLUDING BUT NOT LIMITED TO THE WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE AND NONINFRINGEMENT. IN NO EVENT SHALL THE AUTHORS OR COPYRIGHT HOLDERS BE LIABLE FOR ANY CLAIM, DAMAGES OR OTHER LIABILITY, WHETHER IN AN ACTION OF CONTRACT, TORT OR OTHERWISE, ARISING FROM, OUT OF OR IN CONNECTION WITH THE SOFTWARE OR THE USE OR OTHER DEALINGS IN THE SOFTWARE.

The Cephes Math Library - Copyright © 1984-2000

Release 2.8 - June, 2000 Copyright 1984, 1987, 1988, 2000 by Stephen L. Moshier.

Some software in this archive may be from the book _Methods and Programs for Mathematical Functions_ (Prentice-Hall or Simon & Schuster International, 1989) or from the Cephes Mathematical Library, a commercial product. In either event, it is copyrighted by the author. What you see here may be used freely but it comes with no support or guarantee.

The two known misprints in the book are repaired here in the source listings for the gamma function and the incomplete beta integral.

Stephen L. Moshier, moshier@na-net.ornl.gov.

The Facelt Library - Copyright © 2009 Benjamin Jung

Permission is hereby granted, free of charge, to any person obtaining a copy of this software and associated documentation files (the "Software"), to deal in the Software without restriction, including without limitation the rights to use, copy, modify, merge, publish, distribute, sublicense, and/or sell copies of the Software, and to permit persons to whom the Software is furnished to do so, subject to the following conditions:

The above copyright notice and this permission notice shall be included in all copies or substantial portions of the Software.

THE SOFTWARE IS PROVIDED "AS IS", WITHOUT WARRANTY OF ANY KIND, EXPRESS OR IMPLIED, INCLUDING BUT NOT LIMITED TO THE WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE AND NONINFRINGEMENT. IN NO EVENT SHALL THE AUTHORS OR COPYRIGHT HOLDERS BE LIABLE FOR ANY CLAIM, DAMAGES OR OTHER LIABILITY, WHETHER IN AN ACTION OF CONTRACT, TORT OR OTHERWISE, ARISING FROM, OUT OF OR IN CONNECTION WITH THE SOFTWARE OR THE USE OR OTHER DEALINGS IN THE SOFTWARE.

The FAST corners detector - Copyright © 2006-2010 Edward Rosten

Redistribution and use in source and binary forms, with or without modification, are permitted provided that the following conditions are met:

- 1. Redistributions of source code must retain the above copyright notice, this list of conditions and the following disclaimer.
- 2. Redistributions in binary form must reproduce the above copyright notice, this list of conditions and the following disclaimer in the documentation and/or other materials provided with the distribution.
- Neither the name of the University of Cambridge nor the names of its contributors may be used to endorse or promote products derived from this software without specific prior written permission.

THIS SOFTWARE IS PROVIDED BY THE COPYRIGHT HOLDERS AND CONTRIBUTORS "AS IS" AND ANY EXPRESS OR IMPLIED WARRANTIES, INCLUDING, BUT NOT LIMITED TO, THE IMPLIED WARRANTIES OF MERCHANTABILITY AND FITNESS FOR A PARTICULAR PURPOSE ARE DISCLAIMED. IN NO EVENT SHALL THE COPYRIGHT OWNER OR CONTRIBUTORS BE LIABLE FOR ANY DIRECT, INDIRECT, INCIDENTAL, SPECIAL, EXEMPLARY, OR CONSEQUENTIAL DAMAGES (INCLUDING, BUT NOT LIMITED TO, PROCUREMENT OF SUBSTITUTE GOODS OR SERVICES; LOSS OF USE, DATA, OR PROFITS; OR BUSINESS INTERRUPTION) HOWEVER CAUSED AND ON ANY THEORY OF LIABILITY, WHETHER IN CONTRACT, STRICT LIABILITY, OR TORT (INCLUDING NEGLIGENCE OR OTHERWISE) ARISING IN ANY WAY OUT OF THE USE OF THIS SOFTWARE, EVEN IF ADVISED OF THE POSSIBILITY OF SUCH DAMAGE.

The LAPACK Linear Algebra Package - Copyright © 1992-2010

Copyright © 1992-2010 The University of Tennessee and The University of Tennessee Research Foundation. All rights reserved.

Copyright © 2000-2010 The University of California Berkeley. All rights reserved.

Copyright © 2006-2010 The University of Colorado Denver. All rights reserved.

Redistribution and use in source and binary forms, with or without modification, are permitted provided that the following conditions are met:

- 1. Redistributions of source code must retain the above copyright notice, this list of conditions and the following disclaimer.
- 2. Redistributions in binary form must reproduce the above copyright notice, this list of conditions and the following disclaimer listed in this license in the documentation and/or other materials provided with the distribution.
- 3. Neither the name of the copyright holders nor the names of its contributors may be used to endorse or promote products derived from this software without specific prior written permission.

The copyright holders provide no reassurances that the source code provided does not infringe any patent, copyright, or any other intellectual property rights of third parties. The copyright holders disclaim any liability to any recipient for claims brought against recipient by any third party for infringement of that parties intellectual property rights.

THIS SOFTWARE IS PROVIDED BY THE COPYRIGHT HOLDERS AND CONTRIBUTORS "AS IS" AND ANY EXPRESS OR IMPLIED WARRANTIES, INCLUDING, BUT NOT LIMITED TO, THE IMPLIED WARRANTIES OF MERCHANTABILITY AND FITNESS FOR A PARTICULAR PURPOSE ARE DISCLAIMED. IN NO EVENT SHALL THE COPYRIGHT OWNER OR CONTRIBUTORS BE LIABLE FOR ANY DIRECT, INDIRECT, INCIDENTAL, SPECIAL, EXEMPLARY, OR CONSEQUENTIAL DAMAGES (INCLUDING, BUT NOT LIMITED TO, PROCUREMENT OF SUBSTITUTE GOODS OR SERVICES; LOSS OF USE, DATA, OR PROFITS; OR BUSINESS INTERRUPTION) HOWEVER CAUSED AND ON ANY THEORY OF LIABILITY, WHETHER IN CONTRACT, STRICT LIABILITY, OR TORT (INCLUDING NEGLIGENCE OR OTHERWISE) ARISING IN ANY WAY OUT OF THE USE OF THIS SOFTWARE, EVEN IF ADVISED OF THE POSSIBILITY OF SUCH DAMAGE.

The Marilena Project - Copyright © 2008 Masakazu Ohtsuka

Redistribution and use in source and binary forms, with or without modification, are permitted provided that the following conditions are met:

- 1. Redistributions of source code must retain the above copyright notice, this list of conditions and the following disclaimer.
- 2. Redistributions in binary form must reproduce the above copyright notice, this list of conditions and the following disclaimer in the documentation and/or other materials provided with the distribution.

This software is provided by the copyright holders and contributors "as is" and any express or implied warranties, including, but not limited to, the implied warranties of merchantability and fitness for a particular purpose are disclaimed. In no event shall the Intel Corporation or contributors be liable for any direct, indirect, incidental, special, exemplary, or consequential damages (including, but not limited to, procurement of substitute goods or services; loss of use, data, or profits; or business interruption) however caused and on any theory of liability, whether in contract, strict liability, or tort (including negligence or otherwise) arising in any way out of the use of this software, even if advised of the possibility of such damage.

Autofac

MIT License

Copyright © 2014 Autofac Project

Permission is hereby granted, free of charge, to any person obtaining a copy of this software and associated documentation files (the "Software"), to deal in the Software without restriction, including without limitation the rights to use, copy, modify, merge, publish, distribute, sublicense, and/or sell copies of the Software, and to permit persons to whom the Software is furnished to do so, subject to the following conditions:

The above copyright notice and this permission notice shall be included in all copies or substantial portions of the Software.

THE SOFTWARE IS PROVIDED "AS IS", WITHOUT WARRANTY OF ANY KIND, EXPRESS OR IMPLIED, INCLUDING BUT NOT LIMITED TO THE WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE AND NONINFRINGEMENT. IN NO EVENT SHALL THE AUTHORS OR COPYRIGHT HOLDERS BE LIABLE FOR ANY CLAIM, DAMAGES OR OTHER LIABILITY, WHETHER IN AN ACTION OF CONTRACT, TORT OR OTHERWISE, ARISING FROM, OUT OF OR IN CONNECTION WITH THE SOFTWARE OR THE USE OR OTHER DEALINGS IN THE SOFTWARE.

AutoMapper

MIT License

Copyright © 2010 Jimmy Bogard

Permission is hereby granted, free of charge, to any person obtaining a copy of this software and associated documentation files (the "Software"), to deal in the Software without restriction, including without limitation the rights to use, copy, modify, merge, publish, distribute, sublicense, and/or sell copies of the Software, and to permit persons to whom the Software is furnished to do so, subject to the following conditions:

The above copyright notice and this permission notice shall be included in all copies or substantial portions of the Software.

THE SOFTWARE IS PROVIDED "AS IS", WITHOUT WARRANTY OF ANY KIND, EXPRESS OR IMPLIED, INCLUDING BUT NOT LIMITED TO THE WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE AND NONINFRINGEMENT. IN NO EVENT SHALL THE AUTHORS OR COPYRIGHT HOLDERS BE LIABLE FOR ANY CLAIM, DAMAGES OR OTHER LIABILITY, WHETHER IN AN ACTION OF CONTRACT, TORT OR OTHERWISE, ARISING FROM, OUT OF OR IN CONNECTION WITH THE SOFTWARE OR THE USE OR OTHER DEALINGS IN THE SOFTWARE.

Bio-Formats

Copyright © 2005-2018 Open Microscopy Environment, Board of Regents of the University of Wisconsin-Madison

- Glencoe Software, Inc.
- University of Dundee

BitMiracle.LibTiff.NET

Copyright © 2008-2019, Bit Miracle, All rights reserved.

Redistribution and use in source and binary forms, with or without modification, are permitted provided that the following conditions are met:

- 1. Redistributions of source code must retain the above copyright notice, this list of conditions and the following disclaimer.
- 2. Redistributions in binary form must reproduce the above copyright notice, this list of conditions and the following disclaimer in the documentation and/or other materials provided with the distribution.
- 3. Neither the name of the Bit Miracle nor the names of its contributors may be used to endorse or promote products derived from this software without specific prior written permission.

THIS SOFTWARE IS PROVIDED BY THE COPYRIGHT HOLDERS AND CONTRIBUTORS "AS IS" AND ANY EXPRESS OR IMPLIED WARRANTIES, INCLUDING, BUT NOT LIMITED TO, THE IMPLIED WARRANTIES OF MERCHANTABILITY AND FITNESS FOR A PARTICULAR PURPOSE ARE DISCLAIMED. IN NO EVENT SHALL BIT MIRACLE BE LIABLE FOR ANY DIRECT, INDIRECT, INCIDENTAL, SPECIAL, EXEMPLARY, OR CONSEQUENTIAL DAMAGES (INCLUDING, BUT NOT LIMITED TO, PROCUREMENT OF SUBSTITUTE GOODS OR SERVICES; LOSS OF USE, DATA, OR PROFITS; OR BUSINESS INTERRUPTION) HOWEVER CAUSED AND ON ANY THEORY OF LIABILITY, WHETHER IN CONTRACT, STRICT LIABILITY, OR TORT (INCLUDING NEGLIGENCE OR OTHERWISE) ARISING IN ANY WAY OUT OF THE USE OF THIS SOFTWARE, EVEN IF ADVISED OF THE POSSIBILITY OF SUCH DAMAGE. NOTE: This software is based in part on the The Independent JPEG Group's JPEG software, http://www.ijg.org/, Copyright © 1991-1998, Thomas G. Lane.

NOTE: This software is based in part on the LibTIFF - TIFF Library and Utilities http://www.remotesensing.org/libtiff/, Copyright © 1988-1997 Sam Leffler

Copyright © 1991-1997 Silicon Graphics, Inc., Portions Copyright © 2006, ComponentAce, http://www.componentace.com, All rights reserved.

Redistribution and use in source and binary forms, with or without modification, are permitted provided that the following conditions are met:

- 1. Redistributions of source code must retain the above copyright notice, this list of conditions and the following disclaimer.
- 2. Redistributions in binary form must reproduce the above copyright notice, this list of conditions and the following disclaimer in the documentation and/or other materials provided with the distribution.
- 3. Neither the name of ComponentAce nor the names of its contributors may be used to endorse or promote products derived from this software without specific prior written permission.

THIS SOFTWARE IS PROVIDED BY THE COPYRIGHT HOLDERS AND CONTRIBUTORS "AS IS" AND ANY EXPRESS OR IMPLIED WARRANTIES, INCLUDING, BUT NOT LIMITED TO, THE IMPLIED WARRANTIES OF MERCHANTABILITY AND FITNESS FOR A PARTICULAR PURPOSE ARE DISCLAIMED. IN NO EVENT SHALL THE COPYRIGHT OWNER OR CONTRIBUTORS BE LIABLE FOR ANY DIRECT, INDIRECT, INCIDENTAL, SPECIAL, EXEMPLARY, OR CONSEQUENTIAL DAMAGES (INCLUDING, BUT NOT LIMITED TO, PROCUREMENT OF SUBSTITUTE GOODS OR SERVICES; LOSS OF USE, DATA, OR PROFITS; OR BUSINESS INTERRUPTION) HOWEVER CAUSED AND ON ANY THEORY OF LIABILITY, WHETHER IN CONTRACT, STRICT LIABILITY, OR TORT (INCLUDING NEGLIGENCE OR OTHERWISE) ARISING IN ANY WAY OUT OF THE USE OF THIS SOFTWARE, EVEN IF ADVISED OF THE POSSIBILITY OF SUCH DAMAGE.

Boost

Boost Software License - Version 1.0 - August 17th, 2003

Permission is hereby granted, free of charge, to any person or organization obtaining a copy of the software and accompanying documentation covered by this license (the "Software") to use, reproduce, display, distribute, execute, and transmit the Software, and to prepare derivative works of the Software, and to permit third-parties to whom the Software is furnished to do so, all subject to the following:

The copyright notices in the Software and this entire statement, including the above license grant, this restriction and the following disclaimer, must be included in all copies of the Software, in whole or in part, and all derivative works of the Software, unless such copies or derivative works are solely in the form of machine-executable object code generated by a source language processor.

THE SOFTWARE IS PROVIDED "AS IS", WITHOUT WARRANTY OF ANY KIND, EXPRESS OR IMPLIED, INCLUDING BUT NOT LIMITED TO THE WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, TITLE AND NON-INFRINGEMENT. IN NO EVENT SHALL THE COPYRIGHT HOLDERS OR ANYONE DISTRIBUTING THE SOFTWARE BE LIABLE FOR ANY DAMAGES OR OTHER LIABILITY, WHETHER IN CONTRACT, TORT OR OTHERWISE, ARISING FROM, OUT OF OR IN CONNECTION WITH THE SOFTWARE OR THE USE OR OTHER DEALINGS IN THE SOFTWARE.
Castle Core

Copyright (c) 2004-2019 Castle Project - http://www.castleproject.org/

CefSharp

 $Copyright @ 2010-2017 \ The \ CefSharp \ Authors. \ All \ rights \ reserved. \ Redistribution \ and \ use \ in \ source \ and \ binary \ forms, \ with \ or \ without \ modification, \ are \ permitted \ provided \ that \ the \ following \ conditions \ are \ met:$

- 1. Redistributions of source code must retain the above copyright notice, this list of conditions and the following disclaimer.
- 2. Redistributions in binary form must reproduce the above copyright notice, this list of conditions and the following disclaimer in the documentation and/or other materials provided with the distribution.
- 3. Neither the name of Google Inc. nor the name Chromium Embedded Framework nor the name CefSharp nor the names of its contributors may be used to endorse or promote products derived from this software without specific prior written permission.

THIS SOFTWARE IS PROVIDED BY THE COPYRIGHT HOLDERS AND CONTRIBUTORS "AS IS" AND ANY EXPRESS OR IMPLIED WARRANTIES, INCLUDING, BUT NOT LIMITED TO, THE IMPLIED WARRANTIES OF MERCHANTABILITY AND FITNESS FOR A PARTICULAR PURPOSE ARE DISCLAIMED. IN NO EVENT SHALL THE COPYRIGHT OWNER OR CONTRIBUTORS BE LIABLE FOR ANY DIRECT, INDIRECT, INCIDENTAL, SPECIAL, EXEMPLARY, OR CONSEQUENTIAL DAMAGES (INCLUDING, BUT NOT LIMITED TO, PROCUREMENT OF SUBSTITUTE GOODS OR SERVICES; LOSS OF USE, DATA, OR PROFITS; OR BUSINESS INTERRUPTION) HOWEVER CAUSED AND ON ANY THEORY OF LIABILITY, WHETHER IN CONTRACT, STRICT LIABILITY, OR TORT (INCLUDING NEGLIGENCE OR OTHERWISE) ARISING IN ANY WAY OUT OF THE USE OF THIS SOFTWARE, EVEN IF ADVISED OF THE POSSIBILITY OF SUCH DAMAGE.

Chromium

Copyright © 2015 The Chromium Authors. All rights reserved. Redistribution and use in source and binary forms, with or without modification, are permitted provided that the following conditions are met:

- 1. Redistributions of source code must retain the above copyright notice, this list of conditions and the following disclaimer.
- 2. Redistributions in binary form must reproduce the above copyright notice, this list of conditions and the following disclaimer in the documentation and/or other materials provided with the distribution.
- 3. Neither the name of Google Inc. nor the names of its contributors may be used to endorse or promote products derived from this software without specific prior written permission.

THIS SOFTWARE IS PROVIDED BY THE COPYRIGHT HOLDERS AND CONTRIBUTORS "AS IS" AND ANY EXPRESS OR IMPLIED WARRANTIES, INCLUDING, BUT NOT LIMITED TO, THE IMPLIED WARRANTIES OF MERCHANTABILITY AND FITNESS FOR A PARTICULAR PURPOSE ARE DISCLAIMED. IN NO EVENT SHALL THE COPYRIGHT OWNER OR CONTRIBUTORS BE LIABLE FOR ANY DIRECT, INDIRECT, INCIDENTAL, SPECIAL, EXEMPLARY, OR CONSEQUENTIAL DAMAGES (INCLUDING, BUT NOT LIMITED TO, PROCUREMENT OF SUBSTITUTE GOODS OR SERVICES; LOSS OF USE, DATA, OR PROFITS; OR BUSINESS INTERRUPTION) HOWEVER CAUSED AND ON ANY THEORY OF LIABILITY, WHETHER IN CONTRACT, STRICT LIABILITY, OR TORT (INCLUDING NEGLIGENCE OR OTHERWISE) ARISING IN ANY WAY OUT OF THE USE OF THIS SOFTWARE, EVEN IF ADVISED OF THE POSSIBILITY OF SUCH DAMAGE.

Common Service Locator

Copyright © Microsoft 2008

CsvHelper

Copyright © 2009-2017 Josh Close and Contributors

DotNetZip Library

Copyright © 2006-2011 Dino Chiesa

Copyright © 2006, 2007, 2008, 2009 Dino Chiesa and Microsoft Corporation.

The managed ZLIB code included in Ionic.Zlib.dll and Ionic.Zip.dll is derived from jzlib.

jzlib (https://github.com/ymnk/jzlib) is provided under a BSD-style (3 clause)

Copyright © 2000,2001,2002,2003 ymnk, JCraft, Inc.

Redistribution and use in source and binary forms, with or without modification, are permitted provided that the following conditions are met:

- 1. Redistributions of source code must retain the above copyright notice, this list of conditions and the following disclaimer.
- 2. Redistributions in binary form must reproduce the above copyright notice, this list of conditions and the following disclaimer in the documentation and/or other materials provided with the distribution.
- 3. The names of the authors may not be used to endorse or promote products derived from this software without specific prior written permission.

THIS SOFTWARE IS PROVIDED "AS IS" AND ANY EXPRESSED OR IMPLIED WARRANTIES, INCLUDING, BUT NOT LIMITED TO, THE IMPLIED WARRANTIES OF MERCHANTABILITY AND FITNESS FOR A PARTICULAR PURPOSE ARE DISCLAIMED. IN NO EVENT SHALL JCRAFT, INC. OR ANY CONTRIBUTORS TO THIS SOFTWARE BE LIABLE FOR ANY DIRECT, INDIRECT, INCIDENTAL, SPECIAL, EXEMPLARY, OR CONSEQUENTIAL DAMAGES (INCLUDING, BUT NOT LIMITED TO, PROCUREMENT OF SUBSTITUTE GOODS OR SERVICES; LOSS OF USE, DATA, OR PROFITS; OR BUSINESS INTERRUPTION) HOWEVER CAUSED AND ON ANY THEORY OF LIABILITY, WHETHER IN CONTRACT, STRICT LIABILITY, OR TORT (INCLUDING NEGLIGENCE OR OTHERWISE) ARISING IN ANY WAY OUT OF THE USE OF THIS SOFTWARE, EVEN IF ADVISED OF THE POSSIBILITY OF SUCH DAMAGE.

The jzlib library, itself, is a re-implementation of ZLIB v1.1.3 in pure Java.

zlib is provided under the zlib license:

Copyright © 1995-2004 Jean-loup Gailly and Mark Adler

The ZLIB software is provided 'as-is', without any express or implied warranty. In no event will the authors be held liable for any damages arising from the use of this software.

Permission is granted to anyone to use this software for any purpose, including commercial applications, and to alter it and redistribute it freely, subject to the following restrictions:

1. The origin of this software must not be misrepresented; you must not claim that you wrote the original software. If you use this software in a product, an acknowledgment in the product documentation would be appreciated but is not required.

2. Altered source versions must be plainly marked as such, and must not be misrepresented as being the original software.

3. This notice may not be removed or altered from any source distribution.

Jean-loup Gailly jloup@gzip.org

Mark Adler madler@alumni.caltech.edu

The managed BZIP2 code included in Ionic.BZip2.dll and Ionic.Zip.dll is modified code, based on Java code in the Apache commons compress library.

Apache Commons Compress (http://commons.apache.org/proper/commons-compress/) is provided under the Apache 2 license:

Apache Commons Compress

Copyright 2002-2014 The Apache Software Foundation

Licensed to the Apache Software Foundation (ASF) under one or more contributor license agreements. See the NOTICE file distributed with this work for additional information regarding copyright ownership. The ASF licenses this file to you under the Apache License, Version 2.0 (the "License"); you may not use this file except in compliance with the License. You may obtain a copy of the License at

http://www.apache.org/licenses/LICENSE-2.0

Unless required by applicable law or agreed to in writing, software distributed under the License is distributed on an "AS IS" BASIS, WITHOUT WARRANTIES OR CONDITIONS OF ANY KIND, either express or implied. See the License for the specific language governing permissions and limitations under the License.

Many thanks to Julian Seward for the original C implementation of BZip2 (http://www.bzip.org/).

FluentMigrator

Copyright - Sean Chambers 2008-2014

Fody

MIT License

Copyright (c) 2012 Simon Cropp and contributors

Permission is hereby granted, free of charge, to any person obtaining a copy of this software and associated documentation files (the "Software"), to deal in the Software without restriction, including without limitation the rights to use, copy, modify, merge, publish, distribute, sublicense, and/or sell copies of the Software, and to permit persons to whom the Software is furnished to do so, subject to the following conditions:

The above copyright notice and this permission notice shall be included in all copies or substantial portions of the Software.

THE SOFTWARE IS PROVIDED "AS IS", WITHOUT WARRANTY OF ANY KIND, EXPRESS OR IMPLIED, INCLUDING BUT NOT LIMITED TO THE WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE AND NONINFRINGEMENT. IN NO EVENT SHALL THE AUTHORS OR COPYRIGHT HOLDERS BE LIABLE FOR ANY CLAIM, DAMAGES OR OTHER LIABILITY, WHETHER IN AN ACTION OF CONTRACT, TORT OR OTHERWISE, ARISING FROM, OUT OF OR IN CONNECTION WITH THE SOFTWARE OR THE USE OR OTHER DEALINGS IN THE SOFTWARE.

Framework Microsoft SqlClient Data Provider

MIT License

Copyright (c) Microsoft Corporation

Permission is hereby granted, free of charge, to any person obtaining a copy of this software and associated documentation files (the "Software"), to deal in the Software without restriction, including without limitation the rights to use, copy, modify, merge, publish, distribute, sublicense, and/or sell copies of the Software, and to permit persons to whom the Software is furnished to do so, subject to the following conditions: The above copyright notice and this permission notice shall be included in all copies or substantial portions of the Software.

THE SOFTWARE IS PROVIDED "AS IS", WITHOUT WARRANTY OF ANY KIND, EXPRESS OR IMPLIED, INCLUDING BUT NOT LIMITED TO THE WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE AND NONINFRINGEMENT. IN NO EVENT SHALL THE AUTHORS OR COPYRIGHT HOLDERS BE LIABLE FOR ANY CLAIM, DAMAGES OR OTHER LIABILITY, WHETHER IN AN ACTION OF CONTRACT, TORT OR OTHERWISE, ARISING FROM, OUT OF OR IN CONNECTION WITH THE SOFTWARE OR THE USE OR OTHER DEALINGS IN THE SOFTWARE.

GongSolutions.WPF.DragDrop 1.1.0.0

MIT License

Copyright © 2013-2019, Jan Karger, Steven Kirk. All rights reserved.

Redistribution and use in source and binary forms, with or without modification, are permitted provided that the following conditions are met:

- 1. Redistributions of source code must retain the above copyright notice, this list of conditions and the following disclaimer.
- 2. Redistributions in binary form must reproduce the above copyright notice, this list of conditions and the following disclaimer in the documentation and/or other materials provided with the distribution.
- 3. Neither the name of gong-wpf-dragdrop nor the names of its contributors may be used to endorse or promote products derived from this software without specific prior written permission.

THIS SOFTWARE IS PROVIDED BY THE COPYRIGHT HOLDERS AND CONTRIBUTORS "AS IS" AND ANY EXPRESS OR IMPLIED WARRANTIES, INCLUDING, BUT NOT LIMITED TO, THE IMPLIED WARRANTIES OF MERCHANTABILITY AND FITNESS FOR A PARTICULAR PURPOSE ARE DISCLAIMED. IN NO EVENT SHALL THE COPYRIGHT OWNER OR CONTRIBUTORS BE LIABLE FOR ANY DIRECT, INDIRECT, INCIDENTAL, SPECIAL, EXEMPLARY, OR CONSEQUENTIAL DAMAGES (INCLUDING, BUT NOT LIMITED TO, PROCUREMENT OF SUBSTITUTE GOODS OR SERVICES; LOSS OF USE, DATA, OR PROFITS; OR BUSINESS INTERRUPTION) HOWEVER CAUSED AND ON ANY THEORY OF LIABILITY, WHETHER IN CONTRACT, STRICT LIABILITY, OR TORT (INCLUDING NEGLIGENCE OR OTHERWISE) ARISING IN ANY WAY OUT OF THE USE OF THIS SOFTWARE, EVEN IF ADVISED OF THE POSSIBILITY OF SUCH DAMAGE.

GoogleAnalyticsTracker

Copyright © GoogleAnalyticsTracker 2012-2017

gRPC

Apache License

Version 2.0, January 2004

http://www.apache.org/licenses/

TERMS AND CONDITIONS FOR USE, REPRODUCTION, AND DISTRIBUTION

1. Definitions.

"License" shall mean the terms and conditions for use, reproduction, and distribution as defined by Sections 1 through 9 of this document.

"Licensor" shall mean the copyright owner or entity authorized by the copyright owner that is granting the License.

"Legal Entity" shall mean the union of the acting entity and all other entities that control, are controlled by, or are under common control with that entity. For the purposes of this definition, "control" means (i) the power, direct or indirect, to cause the direction or management of such entity, whether by contract or otherwise, or (ii) ownership of fifty percent (50%) or more of the outstanding shares, or (iii) beneficial ownership of such entity.

"You" (or "Your") shall mean an individual or Legal Entity exercising permissions granted by this License.

"Source" form shall mean the preferred form for making modifications, including but not limited to software source code, documentation source, and configuration files.

"Object" form shall mean any form resulting from mechanical transformation or translation of a Source form, including but not limited to compiled object code, generated documentation, and conversions to other media types.

"Work" shall mean the work of authorship, whether in Source or Object form, made available under the License, as indicated by a copyright notice that is included in or attached to the work (an example is provided in the Appendix below).

"Derivative Works" shall mean any work, whether in Source or Object form, that is based on (or derived from) the Work and for which the editorial revisions, annotations, elaborations, or other modifications represent, as a whole, an original work of authorship. For the purposes of this License, Derivative Works shall not include works that remain separable from, or merely link (or bind by name) to the interfaces of, the Work and Derivative Works thereof.

"Contribution" shall mean any work of authorship, including the original version of the Work and any modifications or additions to that Work or Derivative Works thereof, that is intentionally submitted to Licensor for inclusion in the Work by the copyright owner or by an individual or Legal Entity authorized to submit on behalf of the copyright owner. For the purposes of this definition, "submitted" means any form of electronic, verbal, or written communication sent to the Licensor or its representatives, including but not limited to communication on electronic mailing lists, source code control systems, and issue tracking systems that are managed by, or on behalf of, the Licensor for the purpose of discussing and improving the Work, but excluding communication that is conspicuously marked or otherwise designated in writing by the copyright owner as "Not a Contribution."

"Contributor" shall mean Licensor and any individual or Legal Entity on behalf of whom a Contribution has been received by Licensor and subsequently incorporated within the Work.

2. Grant of Copyright License.

Subject to the terms and conditions of this License, each Contributor hereby grants to You a perpetual, worldwide, nonexclusive, no-charge, royalty-free, irrevocable copyright license to reproduce, prepare Derivative Works of, publicly display, publicly perform, sublicense, and distribute the Work and such Derivative Works in Source or Object form. 3. Grant of Patent License. Subject to the terms and conditions of this License, each Contributor hereby grants to You a perpetual, worldwide, non-exclusive, no-charge, royalty-free, irrevocable (except as stated in this section) patent license to make, have made, use, offer to sell, sell, import, and otherwise transfer the Work, where such license applies only to those patent claims licensable by such Contributor that are necessarily infringed by their Contribution(s) alone or by combination of their Contribution(s) with the Work to which such Contribution(s) was submitted. If You institute patent litigation against any entity (including a cross-claim or counterclaim in a lawsuit) alleging that the Work or a Contribution incorporated within the Work constitutes direct or contributory patent infringement, then any patent licenses granted to You under this License for that Work shall terminate as of the date such litigation is filed.

4. Redistribution.

You may reproduce and distribute copies of the Work or Derivative Works thereof in any medium, with or without modifications, and in Source or Object form, provided that You meet the following conditions:

(a) You must give any other recipients of the Work or Derivative Works a copy of this License; and

(b) You must cause any modified files to carry prominent notices stating that You changed the files; and

(c) You must retain, in the Source form of any Derivative Works that You distribute, all copyright, patent, trademark, and attribution notices from the Source form of the Work, excluding those notices that do not pertain to any part of the Derivative Works; and

(d) If the Work includes a "NOTICE" text file as part of its distribution, then any Derivative Works that You distribute must include a readable copy of the attribution notices contained within such NOTICE file, excluding those notices that do not pertain to any part of the Derivative Works, in at least one of the following places: within a NOTICE text file distributed as part of the Derivative Works; within the Source form or documentation, if provided along with the Derivative Works; or, within a display generated by the Derivative Works, if and wherever such third-party notices normally appear. The contents of the NOTICE file are for informational purposes only and do not modify the License. You may add Your own attribution notices within Derivative Works that You distribute, alongside or as an addendum to the NOTICE text from the Work, provided that such additional attribution notices cannot be construed as modifying the License. You may add Your own copyright statement to Your modifications and may provide additional or different license terms and conditions for use, reproduction, or distribution of Your modifications, or for any such Derivative Works as a whole, provided Your use, reproduction, and distribution of the Work otherwise complies with the conditions stated in this License.

5. Submission of Contributions.

Unless You explicitly state otherwise, any Contribution intentionally submitted for inclusion in the Work by You to the Licensor shall be under the terms and conditions of this License, without any additional terms or conditions. Notwithstanding the above, nothing herein shall supersede or modify the terms of any separate license agreement you may have executed with Licensor regarding such Contributions.

6. Trademarks.

This License does not grant permission to use the trade names, trademarks, service marks, or product names of the Licensor, except as required for reasonable and customary use in describing the origin of the Work and reproducing the content of the NOTICE file.

7. Disclaimer of Warranty.

Unless required by applicable law or agreed to in writing, Licensor provides the Work (and each Contributor provides its Contributions) on an "AS IS" BASIS, WITHOUT WARRANTIES OR CONDITIONS OF ANY KIND, either express or implied, including, without limitation, any warranties or conditions of TITLE, NON-INFRINGEMENT, MERCHANTABILITY, or FITNESS FOR A PARTICULAR PURPOSE. You are solely responsible for determining the appropriateness of using or redistributing the Work and assume any risks associated with Your exercise of permissions under this License.

8. Limitation of Liability.

In no event and under no legal theory, whether in tort (including negligence), contract, or otherwise, unless required by applicable law (such as deliberate and grossly negligent acts) or agreed to in writing, shall any Contributor be liable to You for damages, including any direct, indirect, special, incidental, or consequential damages of any character arising as a result of this License or out of the use or inability to use the Work (including but not limited to damages for loss of goodwill, work stoppage, computer failure or malfunction, or any and all other commercial damages or losses), even if such Contributor has been advised of the possibility of such damages.

9. Accepting Warranty or Additional Liability.

While redistributing the Work or Derivative Works thereof, You may choose to offer, and charge a fee for, acceptance of support, warranty, indemnity, or other liability obligations and/or rights consistent with this License. However, in accepting such obligations, You may act only on Your own behalf and on Your sole responsibility, not on behalf of any other Contributor, and only if You agree to indemnify, defend, and hold each Contributor harmless for any liability incurred by, or claims asserted against, such Contributor by reason of your accepting any such warranty or additional liability.

END OF TERMS AND CONDITIONS

APPENDIX: How to apply the Apache License to your work.

To apply the Apache License to your work, attach the following boilerplate notice, with the fields enclosed by brackets "[]" replaced with your own identifying information. (Don't include the brackets!) The text should be enclosed in the appropriate comment syntax for the file format. We also recommend that a file or class name and description of purpose be included on the same "printed page" as the copyright notice for easier identification within third-party archives.

Copyright [yyyy] [name of copyright owner]

Licensed under the Apache License, Version 2.0 (the "License"); you may not use this file except in compliance with the License.

You may obtain a copy of the License at http://www.apache.org/licenses/LICENSE-2.0

Unless required by applicable law or agreed to in writing, software distributed under the License is distributed on an "AS IS" BASIS, WITHOUT WARRANTIES OR CONDITIONS OF ANY KIND, either express or implied.

See the License for the specific language governing permissions and limitations under the License.

BSD 3-Clause License

Copyright 2016, Google Inc.

Redistribution and use in source and binary forms, with or without modification, are permitted provided that the following conditions are met:

1. Redistributions of source code must retain the above copyright notice, this list of conditions and the following disclaimer.

2. Redistributions in binary form must reproduce the above copyright notice, this list of conditions and the following disclaimer in the documentation and/or other materials provided with the distribution.

3. Neither the name of the copyright holder nor the names of its contributors may be used to endorse or promote products derived from this software without specific prior written permission.

THIS SOFTWARE IS PROVIDED BY THE COPYRIGHT HOLDERS AND CONTRIBUTORS "AS IS" AND ANY EXPRESS OR IMPLIED WARRANTIES, INCLUDING, BUT NOT LIMITED TO, THE IMPLIED WARRANTIES OF MERCHANTABILITY AND FITNESS FOR A PARTICULAR PURPOSE ARE DISCLAIMED. IN NO EVENT SHALL THE COPYRIGHT HOLDER OR CONTRIBUTORS BE LIABLE FOR ANY DIRECT, INDIRECT, INCIDENTAL, SPECIAL, EXEMPLARY, OR CONSEQUENTIAL DAMAGES (INCLUDING, BUT NOT LIMITED TO, PROCUREMENT OF SUBSTITUTE GOODS OR SERVICES; LOSS OF USE, DATA, OR PROFITS; OR BUSINESS INTERRUPTION) HOWEVER CAUSED AND ON ANY THEORY OF LIABILITY, WHETHER IN CONTRACT, STRICT LIABILITY, OR TORT (INCLUDING NEGLIGENCE OR OTHERWISE) ARISING IN ANY WAY OUT OF THE USE OF THIS SOFTWARE, EVEN IF ADVISED OF THE POSSIBILITY OF SUCH DAMAGE.

Mozilla Public License, v. 2.0

This Source Code Form is subject to the terms of the Mozilla Public License, v. 2.0.

If a copy of the MPL was not distributed with this file, You can

obtain one at https://mozilla.org/MPL/2.0/.

Java(TM) Platform SE 8

Oracle Binary Code License Agreement for the Java SE Platform Products and JavaFX ORACLE AMERICA, INC. ("ORACLE"), FOR AND ON BEHALF OF ITSELF AND ITS SUBSIDIARIES AND AFFILIATES UNDER COMMON CONTROL, IS WILLING TO LICENSE THE SOFTWARE TO YOU ONLY UPON THE CONDITION THAT YOU ACCEPT ALL OF THE TERMS CONTAINED IN THIS BINARY CODE LICENSE AGREEMENT AND SUPPLEMENTAL LICENSE TERMS (COLLECTIVELY "AGREEMENT"). PLEASE READ THE AGREEMENT CAREFULLY. BY SELECTING THE "ACCEPT LICENSE AGREEMENT" (OR THE EQUIVALENT) BUTTON AND/OR BY USING THE SOFTWARE YOU ACKNOWLEDGE THAT YOU HAVE READ THE TERMS AND AGREE TO THEM. IF YOU ARE AGREEING TO THESE TERMS ON BEHALF OF A COMPANY OR OTHER LEGAL ENTITY, YOU REPRESENT THAT YOU HAVE THE LEGAL AUTHORITY TO BIND THE LEGAL ENTITY TO THESE TERMS. IF YOU DO NOT HAVE SUCH AUTHORITY, OR IF YOU DO NOT WISH TO BE BOUND BY THE TERMS, THEN SELECT THE "DECLINE LICENSE AGREEMENT" (OR THE EQUIVALENT) BUTTON AND YOU MUST NOT USE THE SOFTWARE ON THIS SITE OR ANY OTHER MEDIA ON WHICH THE SOFTWARE IS CONTAINED. DEFINITIONS. "Software" means the software identified above in binary form that you selected for download, install or use (in the version You selected for download, install or use) from Oracle or its authorized licensees and/or those portions of such software produced by jlink as output using a Program's code, when such output is in unmodified form in combination, and for sole use with, that Program, as well as any other machine readable materials (including, but not limited to, libraries, source files, header files, and data files), any updates or error corrections provided by Oracle, and any user manuals, programming guides and other documentation provided to you by Oracle under this Agreement. The Java Linker (jlink) is available with Java 9 and later versions. "General Purpose Desktop Computers and Servers" means computers, including desktop and laptop computers, or servers, used for general computing functions under end user control (such as but not specifically limited to email, general purpose Internet browsing, and office suite productivity tools). The use of Software in systems and solutions that provide dedicated functionality (other than as mentioned above) or designed for use in embedded or functionspecific software applications, for example but not limited to: Software embedded in or bundled with industrial control systems, wireless mobile telephones, wireless handheld devices, kiosks, TV/STB, Blu-ray Disc devices, telematics and network control switching equipment, printers and storage management systems, and other related systems are excluded from this definition and not licensed under this Agreement. "Programs" means (a) Java technology applets and applications intended to run on the Java Platform. Standard Edition platform on Java-enabled General Purpose Desktop Computers and Servers; and (b) JavaFX technology applications intended to run on the JavaFX Runtime on JavaFX-enabled General Purpose Desktop Computers and Servers. "Java SE LIUM" means the Licensing Information User Manual - Oracle Java SE and Oracle Java Embedded Products Document accessible at http://www.oracle.com/java/technologies/java-se-doc.html. "Commercial Features" means those features that are identified as such in the Java SE LIUM under the "Description of Product Editions and Permitted Features" section.

LICENSE TO USE. Subject to the terms and conditions of this Agreement including, but not limited to, the Java Technology Restrictions of the Supplemental License Terms, Oracle grants you a non-exclusive, non-transferable, limited license without license fees to reproduce and use internally the Software complete and unmodified for the sole purpose of running Programs. THE LICENSE SET FORTH IN THIS SECTION 2 DOES NOT EXTEND TO THE COMMERCIAL FEATURES. YOUR RIGHTS AND OBLIGATIONS RELATED TO THE COMMERCIAL FEATURES ARE AS SET FORTH IN THE SUPPLEMENTAL TERMS ALONG WITH ADDITIONAL LICENSES FOR DEVELOPERS AND PUBLISHERS.

RESTRICTIONS. Software is copyrighted. Title to Software and all associated intellectual property rights is retained by Oracle and/or its licensors. Unless enforcement is prohibited by applicable law, you may not modify, decompile, or reverse engineer Software. You acknowledge that the Software is developed for general use in a variety of information management applications; it is not developed or intended for use in any inherently dangerous applications, including applications that may create a risk of personal injury. If you use the Software in dangerous applications, then you shall be responsible to take all appropriate fail-safe, backup, redundancy, and other measures to ensure its safe use. Oracle disclaims any express or implied warranty of fitness for such uses. No right, title or interest in or to any trademark, service mark, logo or trade name of Oracle or its licensors is granted under this Agreement. Additional restrictions for developers and/or publishers licenses are set forth in the Supplemental License Terms.

DISCLAIMER OF WARRANTY. THE SOFTWARE IS PROVIDED "AS IS" WITHOUT WARRANTY OF ANY KIND. ORACLE FURTHER DISCLAIMS ALL WARRANTIES, EXPRESS AND IMPLIED, INCLUDING WITHOUT LIMITATION, ANY IMPLIED WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE OR NONINFRINGEMENT.

LIMITATION OF LIABILITY. IN NO EVENT SHALL ORACLE BE LIABLE FOR ANY INDIRECT, INCIDENTAL, SPECIAL, PUNITIVE OR CONSEQUENTIAL DAMAGES, OR DAMAGES FOR LOSS OF PROFITS, REVENUE, DATA OR DATA USE, INCURRED BY YOU OR ANY THIRD PARTY, WHETHER IN AN ACTION IN CONTRACT OR TORT, EVEN IF ORACLE HAS BEEN ADVISED OF THE POSSIBILITY OF SUCH DAMAGES. ORACLE'S ENTIRE LIABILITY FOR DAMAGES HEREUNDER SHALL IN NO EVENT EXCEED ONE THOUSAND DOLLARS (U.S. \$1,000).

TERMINATION. This Agreement is effective until terminated. You may terminate this Agreement at any time by destroying all copies of Software. This Agreement will terminate immediately without notice from Oracle if you fail to comply with any provision of this Agreement. Either party may terminate this Agreement immediately should any Software become, or in either party's opinion be likely to become, the subject of a claim of infringement of any intellectual property right. Upon termination, you must destroy all copies of Software.

EXPORT REGULATIONS. You agree that U.S. export control laws and other applicable export and import laws govern your use of the Software, including technical data; additional information can be found on Oracle's Global Trade Compliance web site (http://www.oracle.com/us/products/export). You agree that neither the Software nor any direct product thereof will be exported, directly, or indirectly, in violation of these laws, or will be used for any purpose prohibited by these laws including, without limitation, nuclear, chemical, or biological weapons proliferation.

TRADEMARKS AND LOGOS. You acknowledge and agree as between you and Oracle that Oracle owns the ORACLE and JAVA trademarks and all ORACLE- and JAVA-related trademarks, service marks, logos and other brand designations ("Oracle Marks"), and you agree to comply with the Third Party Usage Guidelines for Oracle Trademarks currently located at http://www.oracle.com/legal/trademarks.html. Any use you make of the Oracle Marks inures to Oracle's benefit.

U.S. GOVERNMENT LICENSE RIGHTS. If Software is being acquired by or on behalf of the U.S. Government or by a U.S. Government prime contractor or subcontractor (at any tier), then the Government's rights in Software and accompanying documentation shall be only those set forth in this Agreement.

GOVERNING LAW. This agreement is governed by the substantive and procedural laws of California. You and Oracle agree to submit to the exclusive jurisdiction of, and venue in, the courts of San Francisco, or Santa Clara counties in California in any dispute arising out of or relating to this agreement.

SEVERABILITY. If any provision of this Agreement is held to be unenforceable, this Agreement will remain in effect with the provision omitted, unless omission would frustrate the intent of the parties, in which case this Agreement will immediately terminate.

INTEGRATION. This Agreement is the entire agreement between you and Oracle relating to its subject matter. It supersedes all prior or contemporaneous oral or written communications, proposals, representations and warranties and prevails over any conflicting or additional terms of any quote, order, acknowledgment, or other communication between the parties relating to its subject matter during the term of this Agreement. No modification of this Agreement will be binding, unless in writing and signed by an authorized representative of each party.

SUPPLEMENTAL LICENSE TERMS

These Supplemental License Terms add to or modify the terms of the Binary Code License Agreement. Capitalized terms not defined in these Supplemental Terms shall have the same meanings ascribed to them in the Binary Code License Agreement. These Supplemental Terms shall supersede any inconsistent or conflicting terms in the Binary Code License Agreement, or in any license contained within the Software.

A. COMMERCIAL FEATURES. You may not use the Commercial Features for running Programs, Java applets or applications in your internal business operations or for any commercial or production purpose, or for any purpose other than as set forth in Sections B, C, D and E of these Supplemental Terms. If You want to use the Commercial Features for any purpose other than as permitted in this Agreement, You must obtain a separate license from Oracle.

B. SOFTWARE INTERNAL USE FOR DEVELOPMENT LICENSE GRANT. Subject to the terms and conditions of this Agreement and restrictions and exceptions set forth in the Java SE LIUM incorporated herein by reference, including, but not limited to the Java Technology Restrictions of these Supplemental Terms, Oracle grants you a non-exclusive, non-transferable, limited license without fees to reproduce internally and use internally the Software complete and unmodified for the purpose of designing, developing, and testing your Programs.

C. LICENSE TO DISTRIBUTE SOFTWARE. Subject to the terms and conditions of this Agreement and restrictions and exceptions set forth in the Java SE LIUM, including, but not limited to the Java Technology Restrictions and Limitations on Redistribution of these Supplemental Terms, Oracle grants you a non-exclusive, non-transferable, limited license without fees to reproduce and distribute the Software, provided that (i) you distribute the Software complete and unmodified and only bundled as part of, and for the sole purpose of running, your Programs, (ii) the Programs add significant and primary functionality to the Software, (iii) you do not distribute additional software intended to replace any component(s) of the Software, (iv) you do not remove or alter any proprietary legends or notices contained in the Software, (v) you only distribute the Software subject to a license agreement that: (a) is a complete, unmodified reproduction of this Agreement; or (b) protects Oracle's interests consistent with the terms contained in this Agreement and that includes the notice set forth in Section H, and (vi) you agree to defend and indemnify Oracle and its licensors from and against any damages, costs, liabilities, settlement amounts and/or expenses (including attorneys' fees) incurred in connection with any claim, lawsuit or action by any third party that arises or results from the use or distribution of any and all Programs and/or Software. The license set forth in this Section C does not extend to the Software identified in Section G.

D. LICENSE TO DISTRIBUTE REDISTRIBUTABLES. Subject to the terms and conditions of this Agreement and restrictions and exceptions set forth in the Java SE LIUM, including but not limited to the Java Technology Restrictions and Limitations on Redistribution of these Supplemental Terms, Oracle grants you a non-exclusive, non-transferable, limited license without fees to reproduce and distribute those files specifically identified as redistributable in the Java SE LIUM ("Redistributables") provided that: (i) you distribute the Redistributables complete and unmodified, and only bundled as part of Programs, (ii) the Programs add significant and primary functionality to the Redistributables, (iii) you do not distribute additional software intended to supersede any component(s) of the Redistributables (unless otherwise specified in the applicable Java SE LIUM), (iv) you do not remove or alter any proprietary legends or notices contained in or on the Redistributables, (v) you only distribute the Redistributables pursuant to a license agreement that: (a) is a complete, unmodified reproduction of this Agreement; or (b) protects Oracle's interests consistent with the terms contained in the Agreement and includes the notice set forth in Section H, (vi) you agree to defend and indemnify Oracle and its licensors from and against any damages, costs, liabilities, settlement amounts and/or expenses (including attorneys' fees) incurred in connection with any claim, lawsuit or action by any third party that arises or results from the use or distribution of any and all Programs and/or Software. The license set forth in this Section D does not extend to the Software identified in Section G.

E. DISTRIBUTION BY PUBLISHERS. This section pertains to your distribution of the JavaTM SE Development Kit Software ("JDK") with your printed book or magazine (as those terms are commonly used in the industry) relating to Java technology ("Publication"). Subject to and conditioned upon your compliance with the restrictions and obligations contained in the Agreement, Oracle hereby grants to you a non-exclusive, nontransferable limited right to reproduce complete and unmodified copies of the JDK on electronic media (the "Media") for the sole purpose of inclusion and distribution with your Publication(s), subject to the following terms: (i) You may not distribute the JDK on a stand-alone basis; it must be distributed with your Publication(s): (ii) You are responsible for downloading the JDK from the applicable Oracle web site; (iii) You must refer to the JDK as JavaTM SE Development Kit; (iv) The JDK must be reproduced in its entirety and without any modification whatsoever (including with respect to all proprietary notices) and distributed with your Publication subject to a license agreement that is a complete, unmodified reproduction of this Agreement; (v) The Media label shall include the following information: "Copyright [YEAR], Oracle America, Inc. All rights reserved. Use is subject to license terms. ORACLE and JAVA trademarks and all ORACLE- and JAVA-related trademarks, service marks, logos and other brand designations are trademarks or registered trademarks of Oracle in the U.S. and other countries." [YEAR] is the year of Oracle's release of the Software; the year information can typically be found in the Software's "About" box or screen. This information must be placed on the Media label in such a manner as to only apply to the JDK; (vi) You must clearly identify the JDK as Oracle's product on the Media holder or Media label, and you may not state or imply that Oracle is responsible for any third-party software contained on the Media; (vii) You may not include any third party software on the Media which is intended to be a replacement or substitute for the JDK; (viii) You agree to defend and indemnify Oracle and its licensors from and against any damages, costs, liabilities, settlement amounts and/or expenses (including attorneys' fees) incurred in connection with any claim, lawsuit or action by any third party that arises or results from the use or distribution of the JDK and/or the Publication; and (ix) You shall provide Oracle with a written notice for each Publication; such notice shall include the following information: (1) title of Publication, (2) author(s), (3) date of Publication, and (4) ISBN or ISSN numbers. Such notice shall be sent to Oracle America, Inc., 500 Oracle Parkway, Redwood Shores, California 94065 U.S.A , Attention: General Counsel.

F. JAVA TECHNOLOGY RESTRICTIONS. You may not create, modify, or change the behavior of, or authorize your licensees to create, modify, or change the behavior of, classes, interfaces, or subpackages that are in any way identified as "java", "javax", "sun", "oracle" or similar convention as specified by Oracle in any naming convention designation.

G. LIMITATIONS ON REDISTRIBUTION. You may not redistribute or otherwise transfer patches, bug fixes or updates made available by Oracle through Oracle Premier Support, including those made available under Oracle's Java SE Support program.

H. COMMERCIAL FEATURES NOTICE. For purpose of complying with Supplemental Term Section C.(v)(b) and D.(v)(b), your license agreement shall include the following notice, where the notice is displayed in a manner that anyone using the Software will see the notice:

Use of the Commercial Features for any commercial or production purpose requires a separate license from Oracle. "Commercial Features" means those features that are identified as such in the Licensing Information User Manual - Oracle Java SE and Oracle Java Embedded Products Document, accessible at http://www.oracle.com/java/technologies/java-sedoc.html, under the "Description of Product Editions and Permitted Features" section.

I. SOURCE CODE. Software may contain source code that, unless expressly licensed for other purposes, is provided solely for reference purposes pursuant to the terms of this Agreement. Source code may not be redistributed unless expressly provided for in this Agreement.

J. THIRD PARTY CODE. Additional copyright notices and license terms applicable to portions of the Software are set forth in the Java SE LIUM accessible at http://www.oracle.com/java/technologies/java-se-doc.html. In addition to any terms and conditions of any third party opensource/freeware license identified in the Java SE LIUM, the disclaimer of warranty and limitation of liability provisions in paragraphs 4 and 5 of the Binary Code License Agreement shall apply to all Software in this distribution.

K. TERMINATION FOR INFRINGEMENT. Either party may terminate this Agreement immediately should any Software become, or in either party's opinion be likely to become, the subject of a claim of infringement of any intellectual property right.

L. INSTALLATION AND AUTO-UPDATE. The Software's installation and auto-update processes transmit a limited amount of data to Oracle (or its service provider) about those specific processes to help Oracle understand and optimize them. Oracle does not associate the data with personally identifiable information. You can find more information about the data Oracle collects as a result of your Software download at http://www.oracle.com/java/technologies/java-se-doc.html.

For inquiries please contact: Oracle America, Inc., 500 Oracle Parkway,

Redwood Shores, California 94065, USA.

Last updated 21 September 2017

JsonSubTypes

MIT License

Permission is hereby granted, free of charge, to any person obtaining a copy of ______ (the "Software"), to deal in the Software without restriction, including without limitation the rights to use, copy, modify, merge, publish, distribute, sublicense, and/or sell copies of the Software, and to permit persons to whom the Software is furnished to do so, subject to the following conditions:

The above copyright notice and this permission notice (including the next paragraph) shall be included in all copies or substantial portions of the Software.

THE SOFTWARE IS PROVIDED "AS IS", WITHOUT WARRANTY OF ANY KIND, EXPRESS OR IMPLIED, INCLUDING BUT NOT LIMITED TO THE WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE AND NONINFRINGEMENT. IN NO EVENT SHALL ______ BE LIABLE FOR ANY CLAIM, DAMAGES OR OTHER LIABILITY, WHETHER IN AN ACTION OF CONTRACT, TORT OR OTHERWISE, ARISING FROM, OUT OF OR IN CONNECTION WITH THE SOFTWARE OR THE USE OR OTHER DEALINGS IN THE SOFTWARE.

LibJpeg

This software is based in part on the work of the Independent JPEG Group.

Copyright © 1991-2016, Thomas G. Lane, Guido Vollbeding.

LibTiff

Copyright © 1988-1997 Sam Leffler, Copyright © 1991-1997 Silicon Graphics, Inc.

Permission to use, copy, modify, distribute, and sell this software and its documentation for any purpose is hereby granted without fee, provided that (i) the above copyright notices and this permission notice appear in all copies of the software and related documentation, and (ii) the names of Sam Leffler and Silicon Graphics may not be used in any advertising or publicity relating to the software without the specific, prior written permission of Sam Leffler and Silicon Graphics.

THE SOFTWARE IS PROVIDED "AS-IS" AND WITHOUT WARRANTY OF ANY KIND, EXPRESS, IMPLIED OR OTHERWISE, INCLUDING WITHOUT LIMITATION, ANY WARRANTY OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE.

IN NO EVENT SHALL SAM LEFFLER OR SILICON GRAPHICS BE LIABLE FOR ANY SPECIAL, INCIDENTAL, INDIRECT OR CONSEQUENTIAL DAMAGES OF ANY KIND, OR ANY DAMAGES WHATSOEVER RESULTING FROM LOSS OF USE, DATA OR PROFITS, WHETHER OR NOT ADVISED OF THE POSSIBILITY OF DAMAGE, AND ON ANY THEORY OF LIABILITY, ARISING OUT OF OR IN CONNECTION WITH THE USE OR PERFORMANCE OF THIS SOFTWARE.

LibTiff.net

Copyright 2008-2011, Bit Miracle. All rights reserved.

This software is provided by the copyright holders and contributors "as is" and any express or implied warranties, including, but not limited to, the implied warranties of merchantability and fitness for a particular purpose are disclaimed. In no event shall Bit Miracle be liable for any direct, indirect, incidental, special, exemplary, or consequential damages (including, but not limited to, procurement of substitute goods or services; loss of use, data, or profits; or business interruption) however caused and on any theory of liability, whether in contract, strict liability, or tort (including negligence or otherwise) arising in any way out of the use of this software, even if advised of the possibility of such damage.

log4net

Copyright 2004-2013 The Apache Software Foundation.

LZ4 Library

Copyright (c) 2011-2020, Yann Collet

All rights reserved.

Redistribution and use in source and binary forms, with or without modification, are permitted provided that the following conditions are met:

* Redistributions of source code must retain the above copyright notice, this list of conditions and the following disclaimer.

* Redistributions in binary form must reproduce the above copyright notice, this list of conditions and the following disclaimer in the documentation and/or other materials provided with the distribution.

THIS SOFTWARE IS PROVIDED BY THE COPYRIGHT HOLDERS AND CONTRIBUTORS "AS IS" AND ANY EXPRESS OR IMPLIED WARRANTIES, INCLUDING, BUT NOT LIMITED TO, THE IMPLIED WARRANTIES OF MERCHANTABILITY AND FITNESS FOR A PARTICULAR PURPOSE ARE DISCLAIMED. IN NO EVENT SHALL THE COPYRIGHT HOLDER OR CONTRIBUTORS BE LIABLE FOR ANY DIRECT, INDIRECT, INCIDENTAL, SPECIAL, EXEMPLARY, OR CONSEQUENTIAL DAMAGES (INCLUDING, BUT NOT LIMITED TO, PROCUREMENT OF SUBSTITUTE GOODS OR SERVICES; LOSS OF USE, DATA, OR PROFITS; OR BUSINESS INTERRUPTION) HOWEVER CAUSED AND ON ANY THEORY OF LIABILITY, WHETHER IN CONTRACT, STRICT LIABILITY, OR TORT (INCLUDING NEGLIGENCE OR OTHERWISE) ARISING IN ANY WAY OUT OF THE USE OF THIS SOFTWARE, EVEN IF ADVISED OF THE POSSIBILITY OF SUCH DAMAGE.

Math.NET Numerics

Copyright © 2002-2020 Math.NET

Permission is hereby granted, free of charge, to any person obtaining a copy of this software and associated documentation files (the "Software"), to deal in the Software without restriction, including without limitation the rights to use, copy, modify, merge, publish, distribute, sublicense, and/or sell copies of the Software, and to permit persons to whom the Software is furnished to do so, subject to the following conditions:

The above copyright notice and this permission notice shall be included in all copies or substantial portions of the Software.

THE SOFTWARE IS PROVIDED "AS IS", WITHOUT WARRANTY OF ANY KIND, EXPRESS OR IMPLIED, INCLUDING BUT NOT LIMITED TO THE WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE AND NONINFRINGEMENT. IN NO EVENT SHALL THE AUTHORS OR COPYRIGHT HOLDERS BE LIABLE FOR ANY CLAIM, DAMAGES OR OTHER LIABILITY, WHETHER IN AN ACTION OF CONTRACT, TORT OR OTHERWISE, ARISING FROM, OUT OF OR IN CONNECTION WITH THE SOFTWARE OR THE USE OR OTHER DEALINGS IN THE SOFTWARE.

Microsoft Enterprise Library for .NET

Copyright © Microsoft Corporation. All rights reserved.

Microsoft Entity Framework

© Microsoft Corporation. All rights reserved.

Nito Libraries

MIT License

Permission is hereby granted, free of charge, to any person obtaining a copy of ______ (the "Software"), to deal in the Software without restriction, including without limitation the rights to use, copy, modify, merge, publish, distribute, sublicense, and/or sell copies of the Software, and to permit persons to whom the Software is furnished to do so, subject to the following conditions:

The above copyright notice and this permission notice (including the next paragraph) shall be included in all copies or substantial portions of the Software.

THE SOFTWARE IS PROVIDED "AS IS", WITHOUT WARRANTY OF ANY KIND, EXPRESS OR IMPLIED, INCLUDING BUT NOT LIMITED TO THE WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE AND NONINFRINGEMENT. IN NO EVENT SHALL _____ BE LIABLE FOR ANY CLAIM, DAMAGES OR OTHER LIABILITY, WHETHER IN AN ACTION OF CONTRACT, TORT OR OTHERWISE, ARISING FROM, OUT OF OR IN CONNECTION WITH THE SOFTWARE OR THE USE OR OTHER DEALINGS IN THE SOFTWARE.

NLog v3.2.0.0 (PrivateBuild)

Copyright © 2004-2020 Jaroslaw Kowalski <jaak@jkowalski.net>, Kim Christensen, Julian Verdurmen

All rights reserved.

Redistribution and use in source and binary forms, with or without modification, are permitted provided that the following conditions are met:

- 1. Redistributions of source code must retain the above copyright notice, this list of conditions and the following disclaimer.
- 2. Redistributions in binary form must reproduce the above copyright notice, this list of conditions and the following disclaimer in the documentation and/or other materials provided with the distribution.
- 3. Neither the name of Jaroslaw Kowalski nor the names of its contributors may be used to endorse or promote products derived from this software without specific prior written permission.

THIS SOFTWARE IS PROVIDED BY THE COPYRIGHT HOLDERS AND CONTRIBUTORS "AS IS" AND ANY EXPRESS OR IMPLIED WARRANTIES, INCLUDING, BUT NOT LIMITED TO, THE IMPLIED WARRANTIES OF MERCHANTABILITY AND FITNESS FOR A PARTICULAR PURPOSE ARE DISCLAIMED. IN NO EVENT SHALL THE COPYRIGHT OWNER OR CONTRIBUTORS BE LIABLE FOR ANY DIRECT, INDIRECT, INCIDENTAL, SPECIAL, EXEMPLARY, OR CONSEQUENTIAL DAMAGES (INCLUDING, BUT NOT LIMITED TO, PROCUREMENT OF SUBSTITUTE GOODS OR SERVICES; LOSS OF USE, DATA, OR PROFITS; OR BUSINESS INTERRUPTION) HOWEVER CAUSED AND ON ANY THEORY OF LIABILITY, WHETHER IN CONTRACT, STRICT LIABILITY, OR TORT (INCLUDING NEGLIGENCE OR OTHERWISE) ARISING IN ANY WAY OUT OF THE USE OF THIS SOFTWARE, EVEN IF ADVISED OF THE POSSIBILITY OF SUCH DAMAGE.

OpenCvSharp3

BSD 3-Clause License

Copyright © 2019, shimat

All rights reserved.

Redistribution and use in source and binary forms, with or without modification, are permitted provided that the following conditions are met:

- 1. Redistributions of source code must retain the above copyright notice, this list of conditions and the following disclaimer.
- 2. Redistributions in binary form must reproduce the above copyright notice, this list of conditions and the following disclaimer in the documentation and/or other materials provided with the distribution.
- 3. Neither the name of the copyright holder nor the names of its contributors may be used to endorse or promote products derived from this software without specific prior written permission.

THIS SOFTWARE IS PROVIDED BY THE COPYRIGHT HOLDERS AND CONTRIBUTORS "AS IS" AND ANY EXPRESS OR IMPLIED WARRANTIES, INCLUDING, BUT NOT LIMITED TO, THE IMPLIED WARRANTIES OF MERCHANTABILITY AND FITNESS FOR A PARTICULAR PURPOSE ARE DISCLAIMED. IN NO EVENT SHALL THE COPYRIGHT HOLDER OR CONTRIBUTORS BE LIABLE FOR ANY DIRECT, INDIRECT, INCIDENTAL, SPECIAL, EXEMPLARY, OR CONSEQUENTIAL DAMAGES (INCLUDING, BUT NOT LIMITED TO, PROCUREMENT OF SUBSTITUTE GOODS OR SERVICES; LOSS OF USE, DATA, OR PROFITS; OR BUSINESS INTERRUPTION) HOWEVER CAUSED AND ON ANY THEORY OF LIABILITY, WHETHER IN CONTRACT, STRICT LIABILITY, OR TORT (INCLUDING NEGLIGENCE OR OTHERWISE) ARISING IN ANY WAY OUT OF THE USE OF THIS SOFTWARE, EVEN IF ADVISED OF THE POSSIBILITY OF SUCH DAMAGE.

OxyPlot

MIT License

Copyright © 2014 OxyPlot contributors

Permission is hereby granted, free of charge, to any person obtaining a copy of this software and associated documentation files (the "Software"), to deal in the Software without restriction, including without limitation the rights to use, copy, modify, merge, publish, distribute, sublicense, and/or sell copies of the Software, and to permit persons to whom the Software is furnished to do so, subject to the following conditions:

The above copyright notice and this permission notice shall be included in all copies or substantial portions of the Software.

THE SOFTWARE IS PROVIDED "AS IS", WITHOUT WARRANTY OF ANY KIND, EXPRESS OR IMPLIED, INCLUDING BUT NOT LIMITED TO THE WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE AND NONINFRINGEMENT. IN NO EVENT SHALL THE AUTHORS OR COPYRIGHT HOLDERS BE LIABLE FOR ANY CLAIM, DAMAGES OR OTHER LIABILITY, WHETHER IN AN ACTION OF CONTRACT, TORT OR OTHERWISE, ARISING FROM, OUT OF OR IN CONNECTION WITH THE SOFTWARE OR THE USE OR OTHER DEALINGS IN THE SOFTWARE.

Polly

Redistribution and use in source and binary forms, with or without modification, are permitted provided that the following conditions are met:

1. Redistributions of source code must retain the above copyright notice, this list of conditions and the following disclaimer.

2. Redistributions in binary form must reproduce the above copyright notice, this list of conditions and the following disclaimer in the documentation and/or other materials provided with the distribution.

3. _____ be used to endorse or promote products derived from this software without specific prior written permission.

THIS SOFTWARE IS PROVIDED BY _____ "AS IS" AND ANY _____ OR IMPLIED WARRANTIES, INCLUDING, BUT NOT LIMITED TO, THE IMPLIED WARRANTIES OF MERCHANTABILITY AND FITNESS FOR A PARTICULAR PURPOSE ARE DISCLAIMED. IN NO EVENT SHALL _____ BE LIABLE FOR ANY DIRECT, INDIRECT, INCIDENTAL, SPECIAL, EXEMPLARY, OR CONSEQUENTIAL DAMAGES (INCLUDING, BUT NOT LIMITED TO, PROCUREMENT OF SUBSTITUTE GOODS OR SERVICES; LOSS OF USE, DATA, OR PROFITS; OR BUSINESS INTERRUPTION) HOWEVER CAUSED AND ON ANY THEORY OF LIABILITY, WHETHER IN CONTRACT, STRICT LIABILITY, OR TORT (INCLUDING NEGLIGENCE OR OTHERWISE) ARISING IN ANY WAY OUT OF THE USE OF THIS SOFTWARE, EVEN IF ADVISED OF THE POSSIBILITY OF SUCH DAMAGE.

Protocol Buffers .NET

The core Protocol Buffers technology is provided courtesy of Google. At the time of writing, this is released under the BSD license.

Full details can be found here: http://code.google.com/p/protobuf/

This .NET implementation is Copyright 2008 Marc Gravell Licensed under the Apache License, Version 2.0 (the "License"); you may not use this file except in compliance with the License. You may obtain a copy of the License at http://www.apache.org/licenses/LICENSE-2.0

Unless required by applicable law or agreed to in writing, software distributed under the License is distributed on an "AS IS" BASIS, WITHOUT WARRANTIES OR CONDITIONS OF ANY KIND, either express or implied.

See the License for the specific language governing permissions and limitations under the License.

RBush.dll

MIT License

Copyright © 2017 viceroypenguin

Permission is hereby granted, free of charge, to any person obtaining a copy of this software and associated documentation files (the "Software"), to deal in the Software without restriction, including without limitation the rights to use, copy, modify, merge, publish, distribute, sublicense, and/or sell copies of the Software, and to permit persons to whom the Software is furnished to do so, subject to the following conditions:

The above copyright notice and this permission notice shall be included in all copies or substantial portions of the Software.

THE SOFTWARE IS PROVIDED "AS IS", WITHOUT WARRANTY OF ANY KIND, EXPRESS OR IMPLIED, INCLUDING BUT NOT LIMITED TO THE WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE AND NONINFRINGEMENT. IN NO EVENT SHALL THE AUTHORS OR COPYRIGHT HOLDERS BE LIABLE FOR ANY CLAIM, DAMAGES OR OTHER LIABILITY, WHETHER IN AN ACTION OF CONTRACT, TORT OR OTHERWISE, ARISING FROM, OUT OF OR IN CONNECTION WITH THE SOFTWARE OR THE USE OR OTHER DEALINGS IN THE SOFTWARE.

Reactive Extensions (Rx) for .NET

MIT License

Copyright © .NET Foundation and Contributors

All rights reserved.

Permission is hereby granted, free of charge, to any person obtaining a copy of this software and associated documentation files (the "Software"), to deal in the Software without restriction, including without limitation the rights to use, copy, modify, merge, publish, distribute, sublicense, and/or sell copies of the Software, and to permit persons to whom the Software is furnished to do so, subject to the following conditions:

The above copyright notice and this permission notice shall be included in all copies or substantial portions of the Software.

THE SOFTWARE IS PROVIDED "AS IS", WITHOUT WARRANTY OF ANY KIND, EXPRESS OR IMPLIED, INCLUDING BUT NOT LIMITED TO THE WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE AND NONINFRINGEMENT. IN NO EVENT SHALL THE AUTHORS OR COPYRIGHT HOLDERS BE LIABLE FOR ANY CLAIM, DAMAGES OR OTHER LIABILITY, WHETHER IN AN ACTION OF CONTRACT, TORT OR OTHERWISE, ARISING FROM, OUT OF OR IN CONNECTION WITH THE SOFTWARE OR THE USE OR OTHER DEALINGS IN THE SOFTWARE.

RestSharp

Apache License

Version 2.0, January 2004

http://www.apache.org/licenses/

TERMS AND CONDITIONS FOR USE, REPRODUCTION, AND DISTRIBUTION

1. Definitions.

"License" shall mean the terms and conditions for use, reproduction, and distribution as defined by Sections 1 through 9 of this document.

"Licensor" shall mean the copyright owner or entity authorized by the copyright owner that is granting the License.

"Legal Entity" shall mean the union of the acting entity and all other entities that control, are controlled by, or are under common control with that entity. For the purposes of this definition, "control" means (i) the power, direct or indirect, to cause the direction or management of such entity, whether by contract or otherwise, or (ii) ownership of fifty percent (50%) or more of the outstanding shares, or (iii) beneficial ownership of such entity.

"You" (or "Your") shall mean an individual or Legal Entity exercising permissions granted by this License.

"Source" form shall mean the preferred form for making modifications, including but not limited to software source code, documentation source, and configuration files.

"Object" form shall mean any form resulting from mechanical transformation or translation of a Source form, including but not limited to compiled object code, generated documentation, and conversions to other media types.

"Work" shall mean the work of authorship, whether in Source or Object form, made available under the License, as indicated by a copyright notice that is included in or attached to the work (an example is provided in the Appendix below).

"Derivative Works" shall mean any work, whether in Source or Object form, that is based on (or derived from) the Work and for which the editorial revisions, annotations, elaborations, or other modifications represent, as a whole, an original work of authorship. For the purposes of this License, Derivative Works shall not include works that remain separable from, or merely link (or bind by name) to the interfaces of, the Work and Derivative Works thereof.

"Contribution" shall mean any work of authorship, including the original version of the Work and any modifications or additions to that Work or Derivative Works thereof, that is intentionally submitted to Licensor for inclusion in the Work by the copyright owner or by an individual or Legal Entity authorized to submit on behalf of the copyright owner. For the purposes of this definition, "submitted" means any form of electronic, verbal, or written communication sent to the Licensor or its representatives, including but not limited to communication on electronic mailing lists, source code control systems, and issue tracking systems that are managed by, or on behalf of, the Licensor for the purpose of discussing and improving the Work, but excluding communication that is conspicuously marked or otherwise designated in writing by the copyright owner as "Not a Contribution."

"Contributor" shall mean Licensor and any individual or Legal Entity on behalf of whom a Contribution has been received by Licensor and subsequently incorporated within the Work.

2. Grant of Copyright License.

Subject to the terms and conditions of this License, each Contributor hereby grants to You a perpetual, worldwide, nonexclusive, no-charge, royalty-free, irrevocable copyright license to reproduce, prepare Derivative Works of, publicly display, publicly perform, sublicense, and distribute the Work and such Derivative Works in Source or Object form.

3. Grant of Patent License.

Subject to the terms and conditions of this License, each Contributor hereby grants to You a perpetual, worldwide, nonexclusive, no-charge, royalty-free, irrevocable (except as stated in this section) patent license to make, have made, use, offer to sell, sell, import, and otherwise transfer the Work, where such license applies only to those patent claims licensable by such Contributor that are necessarily infringed by their Contribution(s) alone or by combination of their Contribution(s) with the Work to which such Contribution(s) was submitted. If You institute patent litigation against any entity (including a crossclaim or counterclaim in a lawsuit) alleging that the Work or a Contribution incorporated within the Work constitutes direct or contributory patent infringement, then any patent licenses granted to You under this License for that Work shall terminate as of the date such litigation is filed.

4. Redistribution.

You may reproduce and distribute copies of the Work or Derivative Works thereof in any medium, with or without modifications, and in Source or Object form, provided that You meet the following conditions:

You must give any other recipients of the Work or Derivative Works a copy of this License; and

You must cause any modified files to carry prominent notices stating that You changed the files; and

You must retain, in the Source form of any Derivative Works that You distribute, all copyright, patent, trademark, and attribution notices from the Source form of the Work, excluding those notices that do not pertain to any part of the Derivative Works; and

If the Work includes a "NOTICE" text file as part of its distribution, then any Derivative Works that You distribute must include a readable copy of the attribution notices contained within such NOTICE file, excluding those notices that do not pertain to any part of the Derivative Works, in at least one of the following places: within a NOTICE text file distributed as part of the Derivative Works; within the Source form or documentation, if provided along with the Derivative Works; or, within a display generated by the Derivative Works, if and wherever such third-party notices normally appear. The contents of the NOTICE file are for informational purposes only and do not modify the License. You may add Your own attribution notices within Derivative Works that You distribute, alongside or as an addendum to the NOTICE text from the Work, provided that such additional attribution notices cannot be construed as modifying the License.

You may add Your own copyright statement to Your modifications and may provide additional or different license terms and conditions for use, reproduction, or distribution of Your modifications, or for any such Derivative Works as a whole, provided Your use, reproduction, and distribution of the Work otherwise complies with the conditions stated in this License.

5. Submission of Contributions.

Unless You explicitly state otherwise, any Contribution intentionally submitted for inclusion in the Work by You to the Licensor shall be under the terms and conditions of this License, without any additional terms or conditions. Notwithstanding the above, nothing herein shall supersede or modify the terms of any separate license agreement you may have executed with Licensor regarding such Contributions.

6. Trademarks.

This License does not grant permission to use the trade names, trademarks, service marks, or product names of the Licensor, except as required for reasonable and customary use in describing the origin of the Work and reproducing the content of the NOTICE file.

7. Disclaimer of Warranty.

Unless required by applicable law or agreed to in writing, Licensor provides the Work (and each Contributor provides its Contributions) on an "AS IS" BASIS, WITHOUT WARRANTIES OR CONDITIONS OF ANY KIND, either express or implied, including, without limitation, any warranties or conditions of TITLE, NON-INFRINGEMENT, MERCHANTABILITY, or FITNESS FOR A PARTICULAR PURPOSE. You are solely responsible for determining the appropriateness of using or redistributing the Work and assume any risks associated with Your exercise of permissions under this License.

8. Limitation of Liability.

In no event and under no legal theory, whether in tort (including negligence), contract, or otherwise, unless required by applicable law (such as deliberate and grossly negligent acts) or agreed to in writing, shall any Contributor be liable to You for damages, including any direct, indirect, special, incidental, or consequential damages of any character arising as a result of this License or out of the use or inability to use the Work (including but not limited to damages for loss of goodwill, work stoppage, computer failure or malfunction, or any and all other commercial damages or losses), even if such Contributor has been advised of the possibility of such damages.

9. Accepting Warranty or Additional Liability.

While redistributing the Work or Derivative Works thereof, You may choose to offer, and charge a fee for, acceptance of support, warranty, indemnity, or other liability obligations and/or rights consistent with this License. However, in accepting such obligations, You may act only on Your own behalf and on Your sole responsibility, not on behalf of any other Contributor, and only if You agree to indemnify, defend, and hold each Contributor harmless for any liability incurred by, or claims asserted against, such Contributor by reason of your accepting any such warranty or additional liability.

END OF TERMS AND CONDITIONS

Serilog

Apache License Version 2.0, January 2004

http://www.apache.org/licenses/

TERMS AND CONDITIONS FOR USE, REPRODUCTION, AND DISTRIBUTION

1. Definitions.

"License" shall mean the terms and conditions for use, reproduction, and distribution as defined by Sections 1 through 9 of this document. "Licensor" shall mean the copyright owner or entity authorized by the copyright owner that is granting the License. "Legal Entity" shall mean the union of the acting entity and all other entities that control, are controlled by, or are under common control with that entity. For the purposes of this definition, "control" means (i) the power, direct or indirect, to cause the direction or management of such entity, whether by contract or otherwise, or (ii) ownership of fifty percent (50%) or more of the outstanding shares, or (iii) beneficial ownership of such entity. "You" (or "Your") shall mean an individual or Legal Entity exercising permissions granted by this License. "Source" form shall mean the preferred form for making modifications, including but not limited to software source code, documentation source, and configuration files.

"Object" form shall mean any form resulting from mechanical transformation or translation of a Source form, including but not limited to compiled object code, generated documentation, and conversions to other media types. "Work" shall mean the work of authorship, whether in Source or Object form, made available under the License, as indicated by a copyright notice that is included in or attached to the work (an example is provided in the Appendix below).

"Derivative Works" shall mean any work, whether in Source or Object form, that is based on (or derived from) the Work and for which the editorial revisions, annotations, elaborations, or other modifications represent, as a whole, an original work of authorship. For the purposes of this License, Derivative Works shall not include works that remain separable from, or merely link (or bind by name) to the interfaces of, the Work and Derivative Works thereof. "Contribution" shall mean any work of authorship, including the original version of the Work and any modifications or additions to that Work or Derivative Works thereof, that is intentionally submitted to Licensor for inclusion in the Work by the copyright owner or by an individual or Legal Entity authorized to submit on behalf of the copyright owner. For the purposes of this definition, "submitted" means any form of electronic, verbal, or written communication sent to the Licensor or its representatives, including but not limited to communication on electronic mailing lists, source code control systems, and issue tracking systems that are managed by, or on behalf of, the Licensor for the purpose of discussing and improving the Work, but excluding communication that is conspicuously marked or otherwise designated in writing by the copyright owner as "Not a Contribution."

"Contributor" shall mean Licensor and any individual or Legal Entity on behalf of whom a Contribution has been received by Licensor and subsequently incorporated within the Work.

2. Grant of Copyright License.

Subject to the terms and conditions of this License, each Contributor hereby grants to You a perpetual, worldwide, nonexclusive, no-charge, royalty-free, irrevocable copyright license to reproduce, prepare Derivative Works of, publicly display, publicly perform, sublicense, and distribute the Work and such Derivative Works in Source or Object form.

3. Grant of Patent License.

Subject to the terms and conditions of this License, each Contributor hereby grants to You a perpetual, worldwide, non-exclusive, no-charge, royalty-free, irrevocable (except as stated in this section) patent license to make, have

made, use, offer to sell, sell, import, and otherwise transfer the Work, where such license applies only to those patent claims licensable by such Contributor that are necessarily infringed by their Contribution(s) alone or by combination of their Contribution(s) with the Work to which such Contribution(s) was submitted. If You institute patent litigation against any entity (including a cross-claim or counterclaim in a lawsuit) alleging that the Work or a Contribution incorporated within the Work constitutes direct or contributory patent infringement, then any patent licenses granted to You under this License for that Work shall terminate as of the date such litigation is filed.

4. Redistribution.

You may reproduce and distribute copies of the Work or Derivative Works thereof in any medium, with or without modifications, and in Source or Object form, provided that You meet the following conditions:

You must give any other recipients of the Work or Derivative Works a copy of this License; and You must cause any modified files to carry prominent notices stating that You changed the files; and You must retain, in the Source form of any Derivative Works that You distribute, all copyright, patent, trademark, and attribution notices from the Source form of the Work, excluding those notices that do not pertain to any part of the Derivative Works; and If the Work includes a "NOTICE" text file as part of its distribution, then any Derivative Works that You distribute must include a readable copy of the attribution notices contained within such NOTICE file, excluding those notices that do not pertain to any part of the Derivative Works; within the Source form or documentation, if provided along with the Derivative Works; or, within a display generated by the Derivative Works, if and wherever such third-party notices normally appear. The contents of the NOTICE file are for informational purposes only and do not modify the License. You may add Your own attribution notices within Derivative Works that You distribute, alongside or as an addendum to the NOTICE text from the Work, provided that such additional attribution notices cannot be construed as modifying the License. You may add Your own copyright statement to Your modifications, or for any such Derivative Works as a whole, provided Your use, reproduction, and distribution of the Work otherwise complies with the conditions stated in this License.

5. Submission of Contributions.

Unless You explicitly state otherwise, any Contribution intentionally submitted for inclusion in the Work by You to the Licensor shall be under the terms and conditions of this License, without any additional terms or conditions. Notwithstanding the above, nothing herein shall supersede or modify the terms of any separate license agreement you may have executed with Licensor regarding such Contributions.

6. Trademarks.

This License does not grant permission to use the trade names, trademarks, service marks, or product names of the Licensor, except as required for reasonable and customary use in describing the origin of the Work and reproducing the content of the NOTICE file.

7. Disclaimer of Warranty.

Unless required by applicable law or agreed to in writing, Licensor provides the Work (and each Contributor provides its Contributions) on an "AS IS" BASIS, WITHOUT WARRANTIES OR CONDITIONS OF ANY KIND, either express or implied, including, without limitation, any warranties or conditions of TITLE, NON-INFRINGEMENT, MERCHANTABILITY, or FITNESS FOR A PARTICULAR PURPOSE. You are solely responsible for determining the appropriateness of using or redistributing the Work and assume any risks associated with Your exercise of permissions under this License.

8. Limitation of Liability.

In no event and under no legal theory, whether in tort (including negligence), contract, or otherwise, unless required by applicable law (such as deliberate and grossly negligent acts) or agreed to in writing, shall any Contributor be liable to You for damages, including any direct, indirect, special, incidental, or consequential damages of any character arising as a result of this License or out of the use or inability to use the Work (including but not limited to damages for loss of goodwill, work stoppage, computer failure or malfunction, or any and all other commercial damages or losses), even if such Contributor has been advised of the possibility of such damages.

9. Accepting Warranty or Additional Liability.

While redistributing the Work or Derivative Works thereof, You may choose to offer, and charge a fee for, acceptance of support, warranty, indemnity, or other liability obligations and/or rights consistent with this License. However, in accepting such obligations, You may act only on Your own behalf and on Your sole responsibility, not on behalf of any other Contributor, and only if You agree to indemnify, defend, and hold each Contributor harmless for any liability incurred by, or claims asserted against, such Contributor by reason of your accepting any such warranty or additional liability.

END OF TERMS AND CONDITIONS

APPENDIX: How to apply the Apache License to your work

To apply the Apache License to your work, attach the following boilerplate notice, with the fields enclosed by brackets "[]" replaced with your own identifying information. (Don't include the brackets!) The text should be enclosed in the appropriate comment syntax for the file format. We also recommend that a file or class name and description of purpose be included on the same "printed page" as the copyright notice for easier identification within third-party archives.

Copyright [yyyy] [name of copyright owner]

Licensed under the Apache License, Version 2.0 (the "License"); you may not use this file except in compliance with the License. You may obtain a copy of the License at http://www.apache.org/licenses/LICENSE-2.0

Unless required by applicable law or agreed to in writing, software distributed under the License is distributed on an "AS IS" BASIS, WITHOUT WARRANTIES OR CONDITIONS OF ANY KIND, either express or implied. See the License for the specific language governing permissions and limitations under the License.

SQLitePCLRaw

Apache License

Version 2.0, January 2004

http://www.apache.org/licenses/

TERMS AND CONDITIONS FOR USE, REPRODUCTION, AND DISTRIBUTION

1. Definitions.

"License" shall mean the terms and conditions for use, reproduction, and distribution as defined by Sections 1 through 9 of this document.

"Licensor" shall mean the copyright owner or entity authorized by the copyright owner that is granting the License.

"Legal Entity" shall mean the union of the acting entity and all other entities that control, are controlled by, or are under common control with that entity. For the purposes of this definition, "control" means (i) the power, direct or indirect, to cause the direction or management of such entity, whether by contract or otherwise, or (ii) ownership of fifty percent (50%) or more of the outstanding shares, or (iii) beneficial ownership of such entity. "You" (or "Your") shall mean an individual or Legal Entity exercising permissions granted by this License.

"Source" form shall mean the preferred form for making modifications, including but not limited to software source code, documentation source, and configuration files.

"Object" form shall mean any form resulting from mechanical transformation or translation of a Source form, including but not limited to compiled object code, generated documentation, and conversions to other media types.

"Work" shall mean the work of authorship, whether in Source or Object form, made available under the License, as indicated by a copyright notice that is included in or attached to the work (an example is provided in the Appendix below).

"Derivative Works" shall mean any work, whether in Source or Object form, that is based on (or derived from) the Work and for which the editorial revisions, annotations, elaborations, or other modifications represent, as a whole, an original work of authorship. For the purposes of this License, Derivative Works shall not include works that remain separable from, or merely link (or bind by name) to the interfaces of, the Work and Derivative Works thereof.

"Contribution" shall mean any work of authorship, including the original version of the Work and any modifications or additions to that Work or Derivative Works thereof, that is intentionally submitted to Licensor for inclusion in the Work by the copyright owner or by an individual or Legal Entity authorized to submit on behalf of the copyright owner. For the purposes of this definition, "submitted" means any form of electronic, verbal, or written communication sent to the Licensor or its representatives, including but not limited to communication on electronic mailing lists, source code control systems, and issue tracking systems that are managed by, or on behalf of, the Licensor for the purpose of discussing and improving the Work, but excluding communication that is conspicuously marked or otherwise designated in writing by the copyright owner as "Not a Contribution."

"Contributor" shall mean Licensor and any individual or Legal Entity on behalf of whom a Contribution has been received by Licensor and subsequently incorporated within the Work.

2. Grant of Copyright License. Subject to the terms and conditions of this License, each Contributor hereby grants to You a perpetual, worldwide, non-exclusive, no-charge, royalty-free, irrevocable copyright license to reproduce, prepare Derivative Works of, publicly display, publicly perform, sublicense, and distribute the Work and such Derivative Works in Source or Object form.

3. Grant of Patent License. Subject to the terms and conditions of this License, each Contributor hereby grants to You a perpetual, worldwide, non-exclusive, no-charge, royalty-free, irrevocable (except as stated in this section) patent license to make, have made, use, offer to sell, sell, import, and otherwise transfer the Work, where such license applies only to those patent claims licensable by such Contributor that are necessarily infringed by their Contribution(s) alone or by combination of their Contribution(s) with the Work to which such Contribution(s) was submitted. If You institute patent litigation against any entity (including a cross-claim or counterclaim in a lawsuit) alleging that the Work or a Contribution incorporated within the Work constitutes direct or contributory patent infringement, then any patent licenses granted to You under this License for that Work shall terminate as of the date such litigation is filed.

4. Redistribution. You may reproduce and distribute copies of the Work or Derivative Works thereof in any medium, with or without modifications, and in Source or Object form, provided that You meet the following conditions:

(a) You must give any other recipients of the Work or Derivative Works a copy of this License; and

(b) You must cause any modified files to carry prominent notices stating that You changed the files; and

(c) You must retain, in the Source form of any Derivative Works that You distribute, all copyright, patent, trademark, and attribution notices from the Source form of the Work, excluding those notices that do not pertain to any part of the Derivative Works; and

(d) If the Work includes a "NOTICE" text file as part of its distribution, then any Derivative Works that You distribute must include a readable copy of the attribution notices contained within such NOTICE file, excluding those notices that do not pertain to any part of the Derivative Works, in at least one of the following places: within a NOTICE text file distributed as part of the Derivative Works; within the Source form or documentation, if provided along with the Derivative Works; or, within a display generated by the Derivative Works, if and wherever such third-party notices normally appear. The contents of the NOTICE file are for informational purposes only and do not modify the License. You may add Your own attribution notices within Derivative Works that You distribute, alongside or as an addendum to the NOTICE text from the Work, provided that such additional attribution notices cannot be construed as modifying the License.

You may add Your own copyright statement to Your modifications and may provide additional or different license terms and conditions for use, reproduction, or distribution of Your modifications, or for any such Derivative Works as a whole, provided Your use, reproduction, and distribution of the Work otherwise complies with the conditions stated in this License.

5. Submission of Contributions. Unless You explicitly state otherwise, any Contribution intentionally submitted for inclusion in the Work by You to the Licensor shall be under the terms and conditions of this License, without any additional terms or conditions. Notwithstanding the above, nothing herein shall supersede or modify the terms of any separate license agreement you may have executed with Licensor regarding such Contributions.

6. Trademarks. This License does not grant permission to use the trade names, trademarks, service marks, or product names of the Licensor, except as required for reasonable and customary use in describing the origin of the Work and reproducing the content of the NOTICE file.

7. Disclaimer of Warranty. Unless required by applicable law or agreed to in writing, Licensor provides the Work (and each Contributor provides its Contributions) on an "AS IS" BASIS, WITHOUT WARRANTIES OR CONDITIONS OF ANY KIND, either express or implied, including, without limitation, any warranties or conditions of TITLE, NON-INFRINGEMENT, MERCHANTABILITY, or FITNESS FOR A PARTICULAR PURPOSE. You are solely responsible for determining the appropriateness of using or redistributing the Work and assume any risks associated with Your exercise of permissions under this License.

8. Limitation of Liability. In no event and under no legal theory, whether in tort (including negligence), contract, or otherwise, unless required by applicable law (such as deliberate and grossly negligent acts) or agreed to in writing, shall any Contributor be liable to You for damages, including any direct, indirect, special, incidental, or consequential damages of any character arising as a result of this License or out of the use or inability to use the Work (including but not limited to damages for loss of goodwill, work stoppage, computer failure or malfunction, or any and all other commercial damages or losses), even if such Contributor has been advised of the possibility of such damages.

9. Accepting Warranty or Additional Liability. While redistributing the Work or Derivative Works thereof, You may choose to offer, and charge a fee for, acceptance of support, warranty, indemnity, or other liability obligations and/or rights consistent with this License. However, in accepting such obligations, You may act only on Your own behalf and on Your sole responsibility, not on behalf of any other Contributor, and only if You agree to indemnify, defend, and hold each Contributor harmless for any liability incurred by, or claims asserted against, such Contributor by reason of your accepting any such warranty or additional liability.

END OF TERMS AND CONDITIONS

APPENDIX: How to apply the Apache License to your work.

To apply the Apache License to your work, attach the following boilerplate notice, with the fields enclosed by brackets "[]" replaced with your own identifying information. (Don't include the brackets!) The text should be enclosed in the appropriate comment syntax for the file format. We also recommend that a file or class name and description of purpose be included on the same "printed page" as the copyright notice for easier identification within third-party archives.

Copyright [yyyy] [name of copyright owner]

Licensed under the Apache License, Version 2.0 (the "License"); you may not use this file except in compliance with the License.

You may obtain a copy of the License at http://www.apache.org/licenses/LICENSE-2.0

Unless required by applicable law or agreed to in writing, software distributed under the License is distributed on an "AS IS" BASIS, WITHOUT WARRANTIES OR CONDITIONS OF ANY KIND, either express or implied.

See the License for the specific language governing permissions and limitations under the License.

SystemWrapper

Copyright © 2008 Jared Barneck, Portions Copyright © 2014-2018 Jozef Izso

VCL Styles InnoSetup plugin

Copyright © Rodrigo Ruz

YamlDotNet

Copyright (c) 2008, 2009, 2010, 2011, 2012, 2013, 2014 Antoine Aubry and contributors Permission is hereby granted, free of charge, to any person obtaining a copy of this software and associated documentation files (the "Software"), to deal in the Software without restriction, including without limitation the rights to use, copy, modify, merge, publish, distribute, sublicense, and/or sell copies of the Software, and to permit persons to whom the Software is furnished to do so, subject to the following conditions:

The above copyright notice and this permission notice shall be included in all copies or substantial portions of the Software.

THE SOFTWARE IS PROVIDED "AS IS", WITHOUT WARRANTY OF ANY KIND, EXPRESS OR IMPLIED, INCLUDING BUT NOT LIMITED TO THE WARRANTIES OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE AND NONINFRINGEMENT. IN NO EVENT SHALL THE AUTHORS OR COPYRIGHT HOLDERS BE LIABLE FOR ANY CLAIM, DAMAGES OR OTHER LIABILITY, WHETHER IN AN ACTION OF CONTRACT, TORT OR OTHERWISE, ARISING FROM, OUT OF OR IN CONNECTION WITH THE SOFTWARE OR THE USE OR OTHER DEALINGS IN THE SOFTWARE.

ZedGraph Library

GNU LESSER GENERAL PUBLIC LICENSE, Version 2.1, February 1999, Copyright © 1991, 1999 Free Software Foundation, Inc., 51 Franklin Street, Fifth Floor, Boston, MA 02110-1301 USA

Everyone is permitted to copy and distribute verbatim copies of this license document, but changing it is not allowed.

[This is the first released version of the Lesser GPL. It also counts as the successor of the GNU Library Public License, version 2, hence the version number 2.1.]

Preamble

The licenses for most software are designed to take away your freedom to share and change it. By contrast, the GNU General Public Licenses are intended to guarantee your freedom to share and change free software--to make sure the software is free for all its users.

This license, the Lesser General Public License, applies to some specially designated software packages--typically librariesof the Free Software Foundation and other authors who decide to use it. You can use it too, but we suggest you first think carefully about whether this license or the ordinary General Public License is the better strategy to use in any particular case, based on the explanations below.

When we speak of free software, we are referring to freedom of use, not price. Our General Public Licenses are designed to make sure that you have the freedom to distribute copies of free software (and charge for this service if you wish); that you receive source code or can get it if you want it; that you can change the software and use pieces of it in new free programs; and that you are informed that you can do these things.

To protect your rights, we need to make restrictions that forbid distributors to deny you these rights or to ask you to surrender these rights. These restrictions translate to certain responsibilities for you if you distribute copies of the library or if you modify it.

For example, if you distribute copies of the library, whether gratis or for a fee, you must give the recipients all the rights that we gave you. You must make sure that they, too, receive or can get the source code. If you link other code with the library, you must provide complete object files to the recipients, so that they can relink them with the library after making changes to the library and recompiling it. And you must show them these terms so they know their rights.

We protect your rights with a two-step method: (1) we copyright the library, and (2) we offer you this license, which gives you legal permission to copy, distribute and/or modify the library.

To protect each distributor, we want to make it very clear that there is no warranty for the free library. Also, if the library is modified by someone else and passed on, the recipients should know that what they have is not the original version, so that the original author's reputation will not be affected by problems that might be introduced by others.

Finally, software patents pose a constant threat to the existence of any free program. We wish to make sure that a company cannot effectively restrict the users of a free program by obtaining a restrictive license from a patent holder. Therefore, we insist that any patent license obtained for a version of the library must be consistent with the full freedom of use specified in this license.

Most GNU software, including some libraries, is covered by the ordinary GNU General Public License. This license, the GNU Lesser General Public License, applies to certain designated libraries, and is quite different from the ordinary General Public License. We use this license for certain libraries in order to permit linking those libraries into non-free programs.

When a program is linked with a library, whether statically or using a shared library, the combination of the two is legally speaking a combined work, a derivative of the original library. The ordinary General Public License therefore permits such linking only if the entire combination fits its criteria of freedom. The Lesser General Public License permits more lax criteria for linking other code with the library.

We call this license the "Lesser" General Public License because it does Less to protect the user's freedom than the ordinary General Public License. It also provides other free software developers Less of an advantage over competing non-free programs. These disadvantages are the reason we use the ordinary General Public License for many libraries. However, the Lesser license provides advantages in certain special circumstances.

For example, on rare occasions, there may be a special need to encourage the widest possible use of a certain library, so that it becomes a de-facto standard. To achieve this, non-free programs must be allowed to use the library. A more frequent case is that a free library does the same job as widely used non-free libraries. In this case, there is little to gain by limiting the free library to free software only, so we use the Lesser General Public License.

In other cases, permission to use a particular library in non-free programs enables a greater number of people to use a large body of free software. For example, permission to use the GNU C Library in non-free programs enables many more people to use the whole GNU operating system, as well as its variant, the GNU/Linux operating system.

Although the Lesser General Public License is Less protective of the users' freedom, it does ensure that the user of a program that is linked with the Library has the freedom and the wherewithal to run that program using a modified version of the Library.

The precise terms and conditions for copying, distribution and modification follow. Pay close attention to the difference between a "work based on the library" and a "work that uses the library". The former contains code derived from the library, whereas the latter must be combined with the library in order to run.

GNU LESSER GENERAL PUBLIC LICENSE

TERMS AND CONDITIONS FOR COPYING, DISTRIBUTION AND MODIFICATION

0. This License Agreement applies to any software library or other program which contains a notice placed by the copyright holder or other authorized party saying it may be distributed under the terms of this Lesser General Public License (also called "this License"). Each licensee is addressed as "you".

A "library" means a collection of software functions and/or data prepared so as to be conveniently linked with application programs (which use some of those functions and data) to form executables.

The "Library", below, refers to any such software library or work which has been distributed under these terms. A "work based on the Library" means either the Library or any derivative work under copyright law: that is to say, a work containing the Library or a portion of it, either verbatim or with modifications and/or translated straightforwardly into another language. (Hereinafter, translation is included without limitation in the term "modification".)

"Source code" for a work means the preferred form of the work for making modifications to it. For a library, complete source code means all the source code for all modules it contains, plus any associated interface definition files, plus the scripts used to control compilation and installation of the library.

Activities other than copying, distribution and modification are not covered by this License; they are outside its scope. The act of running a program using the Library is not restricted, and output from such a program is covered only if its contents constitute a work based on the Library (independent of the use of the Library in a tool for writing it). Whether that is true depends on what the Library does and what the program that uses the Library does.

1. You may copy and distribute verbatim copies of the Library's complete source code as you receive it, in any medium, provided that you conspicuously and appropriately publish on each copy an appropriate copyright notice and disclaimer of warranty; keep intact all the notices that refer to this License and to the absence of any warranty; and distribute a copy of this License along with the Library.

You may charge a fee for the physical act of transferring a copy, and you may at your option offer warranty protection in exchange for a fee.

2. You may modify your copy or copies of the Library or any portion of it, thus forming a work based on the Library, and copy and distribute such modifications or work under the terms of Section 1 above, provided that you also meet all of these conditions:

a) The modified work must itself be a software library.

b) You must cause the files modified to carry prominent notices stating that you changed the files and the date of any change.

c) You must cause the whole of the work to be licensed at no charge to all third parties under the terms of this License.

d) If a facility in the modified Library refers to a function or a table of data to be supplied by an application program that uses the facility, other than as an argument passed when the facility is invoked, then you must make a good faith effort to ensure that, in the event an application does not supply such function or table, the facility still operates, and performs whatever part of its purpose remains meaningful.

(For example, a function in a library to compute square roots has a purpose that is entirely well-defined independent of the application. Therefore, Subsection 2d requires that any application-supplied function or table used by this function must be optional: if the application does not supply it, the square root function must still compute square roots.)

These requirements apply to the modified work as a whole. If identifiable sections of that work are not derived from the Library, and can be reasonably considered independent and separate works in themselves, then this License, and its terms, do not apply to those sections when you distribute them as separate works. But when you distribute the same sections as part of a whole which is a work based on the Library, the distribution of the whole must be on the terms of this License, whose permissions for other licensees extend to the entire whole, and thus to each and every part regardless of who wrote it.

Thus, it is not the intent of this section to claim rights or contest your rights to work written entirely by you; rather, the intent is to exercise the right to control the distribution of derivative or collective works based on the Library.

In addition, mere aggregation of another work not based on the Library with the Library (or with a work based on the Library) on a volume of a storage or distribution medium does not bring the other work under the scope of this License.

3. You may opt to apply the terms of the ordinary GNU General Public License instead of this License to a given copy of the Library. To do this, you must alter all the notices that refer to this License, so that they refer to the ordinary GNU General Public License, version 2, instead of to this License. (If a newer version than version 2 of the ordinary GNU General Public License has appeared, then you can specify that version instead if you wish.) Do not make any other change in these notices.

Once this change is made in a given copy, it is irreversible for that copy, so the ordinary GNU General Public License applies to all subsequent copies and derivative works made from that copy.

This option is useful when you wish to copy part of the code of the Library into a program that is not a library.

4. You may copy and distribute the Library (or a portion or derivative of it, under Section 2) in object code or executable form under the terms of Sections 1 and 2 above provided that you accompany it with the complete corresponding machine-readable source code, which must be distributed under the terms of Sections 1 and 2 above on a medium customarily used for software interchange.

If distribution of object code is made by offering access to copy from a designated place, then offering equivalent access to copy the source code from the same place satisfies the requirement to distribute the source code, even though third parties are not compelled to copy the source along with the object code.

5. A program that contains no derivative of any portion of the Library, but is designed to work with the Library by being compiled or linked with it, is called a "work that uses the Library". Such a work, in isolation, is not a derivative work of the Library, and therefore falls outside the scope of this License.

However, linking a "work that uses the Library" with the Library creates an executable that is a derivative of the Library (because it contains portions of the Library), rather than a "work that uses the library". The executable is therefore covered by this License. Section 6 states terms for distribution of such executables.

When a "work that uses the Library" uses material from a header file that is part of the Library, the object code for the work may be a derivative work of the Library even though the source code is not. Whether this is true is especially significant if the work can be linked without the Library, or if the work is itself a library. The threshold for this to be true is not precisely defined by law. If such an object file uses only numerical parameters, data structure layouts and accessors, and small macros and small inline functions (ten lines or less in length), then the use of the object file is unrestricted, regardless of whether it is legally a derivative work. (Executables containing this object code plus portions of the Library will still fall under Section 6.)

Otherwise, if the work is a derivative of the Library, you may distribute the object code for the work under the terms of Section 6. Any executables containing that work also fall under Section 6, whether or not they are linked directly with the Library itself.

6. As an exception to the Sections above, you may also combine or link a "work that uses the Library" with the Library to produce a work containing portions of the Library, and distribute that work under terms of your choice, provided that the terms permit modification of the work for the customer's own use and reverse engineering for debugging such modifications.

You must give prominent notice with each copy of the work that the Library is used in it and that the Library and its use are covered by this License. You must supply a copy of this License. If the work during execution displays copyright notices, you must include the copyright notice for the Library among them, as well as a reference directing the user to the copy of this License. Also, you must do one of these things:

a) Accompany the work with the complete corresponding machine-readable source code for the Library including whatever changes were used in the work (which must be distributed under Sections 1 and 2 above); and, if the work is an executable linked with the Library, with the complete machine-readable "work that uses the Library", as object code and/or source code, so that the user can modify the Library and then relink to produce a modified executable containing the modified Library. (It is understood that the user who changes the contents of definitions files in the Library will not necessarily be able to recompile the application to use the modified definitions.)

b) Use a suitable shared library mechanism for linking with the Library. A suitable mechanism is one that (1) uses at run time a copy of the library already present on the user's computer system, rather than copying library functions into the executable, and (2) will operate properly with a modified version of the library, if the user installs one, as long as the modified version is interface-compatible with the version that the work was made with.

c) Accompany the work with a written offer, valid for at least three years, to give the same user the materials specified in Subsection 6a, above, for a charge no more than the cost of performing this distribution.

d) If distribution of the work is made by offering access to copy from a designated place, offer equivalent access to copy the above specified materials from the same place.

e) Verify that the user has already received a copy of these materials or that you have already sent this user a copy.

For an executable, the required form of the "work that uses the Library" must include any data and utility programs needed for reproducing the executable from it. However, as a special exception, the materials to be distributed need not include anything that is normally distributed (in either source or binary form) with the major components (compiler, kernel, and so on) of the operating system on which the executable runs, unless that component itself accompanies the executable.

It may happen that this requirement contradicts the license restrictions of other proprietary libraries that do not normally accompany the operating system. Such a contradiction means you cannot use both them and the Library together in an executable that you distribute.

7. You may place library facilities that are a work based on the Library side-by-side in a single library together with other library facilities not covered by this License, and distribute such a combined library, provided that the separate distribution of the work based on the Library and of the other library facilities is otherwise permitted, and provided that you do these two things:

a) Accompany the combined library with a copy of the same work based on the Library, uncombined with any other library facilities. This must be distributed under the terms of the Sections above.

b) Give prominent notice with the combined library of the fact that part of it is a work based on the Library, and explaining where to find the accompanying uncombined form of the same work.

8. You may not copy, modify, sublicense, link with, or distribute the Library except as expressly provided under this License. Any attempt otherwise to copy, modify, sublicense, link with, or distribute the Library is void, and will automatically terminate your rights under this License. However, parties who have received copies, or rights, from you under this License will not have their licenses terminated so long as such parties remain in full compliance.

9. You are not required to accept this License, since you have not signed it. However, nothing else grants you permission to modify or distribute the Library or its derivative works. These actions are prohibited by law if you do not accept this License. Therefore, by modifying or distributing the Library (or any work based on the Library), you indicate your acceptance of this License to do so, and all its terms and conditions for copying, distributing or modifying the Library or works based on it.

10. Each time you redistribute the Library (or any work based on the Library), the recipient automatically receives a license from the original licensor to copy, distribute, link with or modify the Library subject to these terms and conditions. You may not impose any further restrictions on the recipients' exercise of the rights granted herein. You are not responsible for enforcing compliance by third parties with this License.

11. If, as a consequence of a court judgment or allegation of patent infringement or for any other reason (not limited to patent issues), conditions are imposed on you (whether by court order, agreement or otherwise) that contradict the conditions of this License, they do not excuse you from the conditions of this License. If you cannot distribute so as to satisfy simultaneously your obligations under this License and any other pertinent obligations, then as a consequence you may not distribute the Library at all. For example, if a patent license would not permit royalty-free redistribution of the Library by all those who receive copies directly or indirectly through you, then the only way you could satisfy both it and this License would be to refrain entirely from distribution of the Library.

If any portion of this section is held invalid or unenforceable under any particular circumstance, the balance of the section is intended to apply, and the section as a whole is intended to apply in other circumstances.

It is not the purpose of this section to induce you to infringe any patents or other property right claims or to contest validity of any such claims; this section has the sole purpose of protecting the integrity of the free software distribution system which is implemented by public license practices. Many people have made generous contributions to the wide range of software distributed through that system in reliance on consistent application of that system; it is up to the author/donor to decide if he or she is willing to distribute software through any other system and a licensee cannot impose that choice.

This section is intended to make thoroughly clear what is believed to be a consequence of the rest of this License.

12. If the distribution and/or use of the Library is restricted in certain countries either by patents or by copyrighted interfaces, the original copyright holder who places the Library under this License may add an explicit geographical distribution limitation excluding those countries, so that distribution is permitted only in or among countries not thus excluded. In such case, this License incorporates the limitation as if written in the body of this License.

13. The Free Software Foundation may publish revised and/or new versions of the Lesser General Public License from time to time. Such new versions will be similar in spirit to the present version, but may differ in detail to address new problems or concerns.

Each version is given a distinguishing version number. If the Library specifies a version number of this License which applies to it and "any later version", you have the option of following the terms and conditions either of that version or of any later version published by the Free Software Foundation. If the Library does not specify a license version number, you may choose any version ever published by the Free Software Foundation.

14. If you wish to incorporate parts of the Library into other free programs whose distribution conditions are incompatible with these, write to the author to ask for permission. For software which is copyrighted by the Free Software Foundation, write to the Free Software Foundation; we sometimes make exceptions for this. Our decision will be guided by the two goals of preserving the free status of all derivatives of our free software and of promoting the sharing and reuse of software generally.

NO WARRANTY

15. BECAUSE THE LIBRARY IS LICENSED FREE OF CHARGE, THERE IS NO WARRANTY FOR THE LIBRARY, TO THE EXTENT PERMITTED BY APPLICABLE LAW. EXCEPT WHEN OTHERWISE STATED IN WRITING THE COPYRIGHT HOLDERS AND/OR OTHER PARTIES PROVIDE THE LIBRARY "AS IS" WITHOUT WARRANTY OF ANY KIND, EITHER EXPRESSED OR IMPLIED, INCLUDING, BUT NOT LIMITED TO, THE IMPLIED WARRANTIES OF MERCHANTABILITY AND FITNESS FOR A PARTICULAR PURPOSE. THE ENTIRE RISK AS TO THE QUALITY AND PERFORMANCE OF THE LIBRARY IS WITH YOU. SHOULD THE LIBRARY PROVE DEFECTIVE, YOU ASSUME THE COST OF ALL NECESSARY SERVICING, REPAIR OR CORRECTION.

16. IN NO EVENT UNLESS REQUIRED BY APPLICABLE LAW OR AGREED TO IN WRITING WILL ANY COPYRIGHT HOLDER, OR ANY OTHER PARTY WHO MAY MODIFY AND/OR REDISTRIBUTE THE LIBRARY AS PERMITTED ABOVE, BE LIABLE TO YOU FOR DAMAGES, INCLUDING ANY GENERAL, SPECIAL, INCIDENTAL OR CONSEQUENTIAL DAMAGES ARISING OUT OF THE USE OR INABILITY TO USE THE LIBRARY (INCLUDING BUT NOT LIMITED TO LOSS OF DATA OR DATA BEING RENDERED INACCURATE OR LOSSES SUSTAINED BY YOU OR THIRD PARTIES OR A FAILURE OF THE LIBRARY TO OPERATE WITH ANY OTHER SOFTWARE), EVEN IF SUCH HOLDER OR OTHER PARTY HAS BEEN ADVISED OF THE POSSIBILITY OF SUCH DAMAGES.

END OF TERMS AND CONDITIONS

Zstandard

BSD License

For Zstandard software

Copyright (c) 2016-present, Facebook, Inc. All rights reserved.

Redistribution and use in source and binary forms, with or without modification, are permitted provided that the following conditions are met:

* Redistributions of source code must retain the above copyright notice, this list of conditions and the following disclaimer.

* Redistributions in binary form must reproduce the above copyright notice, this list of conditions and the following disclaimer in the documentation and/or other materials provided with the distribution.

* Neither the name Facebook nor the names of its contributors may be used to endorse or promote products derived from this software without specific prior written permission.

THIS SOFTWARE IS PROVIDED BY THE COPYRIGHT HOLDERS AND CONTRIBUTORS "AS IS" AND ANY EXPRESS OR IMPLIED WARRANTIES, INCLUDING, BUT NOT LIMITED TO, THE IMPLIED WARRANTIES OF MERCHANTABILITY AND FITNESS FOR A PARTICULAR PURPOSE ARE DISCLAIMED. IN NO EVENT SHALL THE COPYRIGHT HOLDER OR CONTRIBUTORS BE LIABLE FOR ANY DIRECT, INDIRECT, INCIDENTAL, SPECIAL, EXEMPLARY, OR CONSEQUENTIAL DAMAGES (INCLUDING, BUT NOT LIMITED TO, PROCUREMENT OF SUBSTITUTE GOODS OR SERVICES; LOSS OF USE, DATA, OR PROFITS; OR BUSINESS INTERRUPTION) HOWEVER CAUSED AND ON ANY THEORY OF LIABILITY, WHETHER IN CONTRACT, STRICT LIABILITY, OR TORT (INCLUDING NEGLIGENCE OR OTHERWISE) ARISING IN ANY WAY OUT OF THE USE OF THIS SOFTWARE, EVEN IF ADVISED OF THE POSSIBILITY OF SUCH DAMAGE.

ZstdNet

BSD License

For ZstdNet software

Copyright (c) 2016-present, SKB Kontur. All rights reserved.

Redistribution and use in source and binary forms, with or without modification, are permitted provided that the following conditions are met:

* Redistributions of source code must retain the above copyright notice, this list of conditions and the following disclaimer.

* Redistributions in binary form must reproduce the above copyright notice, this list of conditions and the following disclaimer in the documentation and/or other materials provided with the distribution.

* Neither the name SKB Kontur nor the names of its contributors may be used to endorse or promote products derived from this software without specific prior written permission.

THIS SOFTWARE IS PROVIDED BY THE COPYRIGHT HOLDERS AND CONTRIBUTORS "AS IS" AND ANY EXPRESS OR IMPLIED WARRANTIES, INCLUDING, BUT NOT LIMITED TO, THE IMPLIED WARRANTIES OF MERCHANTABILITY AND FITNESS FOR A PARTICULAR PURPOSE ARE DISCLAIMED. IN NO EVENT SHALL THE COPYRIGHT HOLDER OR CONTRIBUTORS BE LIABLE FOR ANY DIRECT, INDIRECT, INCIDENTAL, SPECIAL, EXEMPLARY, OR CONSEQUENTIAL DAMAGES (INCLUDING, BUT NOT LIMITED TO, PROCUREMENT OF SUBSTITUTE GOODS OR SERVICES; LOSS OF USE, DATA, OR PROFITS; OR BUSINESS INTERRUPTION) HOWEVER CAUSED AND ON ANY THEORY OF LIABILITY, WHETHER IN CONTRACT, STRICT LIABILITY, OR TORT (INCLUDING NEGLIGENCE OR OTHERWISE) ARISING IN ANY WAY OUT OF THE USE OF THIS SOFTWARE, EVEN IF ADVISED OF THE POSSIBILITY OF SUCH DAMAGE.