

# Microscope Companion™ 1.17

## User Guide

PN 1553273

Revision A • April 2025

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 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, April 2025

# Introduction

Thermo Fisher Scientific Microscope Companion™ software is a microscope or hardware-oriented application to prepare Thermo Fisher Scientific Transmission Electron Microscopes for data acquisition in other dedicated software applications like Thermo Fisher Scientific Velox, EPU, etc. In the remainder of this document, and wherever practical, we use MiCo as an abbreviation of Microscope Companion™.

MiCo software is aimed at supervisors (thus not at all users) of Thermo Fisher Scientific Transmission Electron Microscopes. It is intended as a helpful assistant in identifying and correcting issues. Additionally, MiCo allows you to adjust basic microscope settings, do some sample navigation and perform hardware management activities like corrections or calibrations. The supervisor should be able to successfully perform the following:

- Evaluate the quality of acquired data, possibly in other data acquisition software.
- Identify the cause of accuracy and quality issues with the acquired data.
- Perform the necessary corrective actions to resolve the identified issues, provided that these actions are available to the user. For higher level corrective actions, the assistance of a Thermo Fisher Scientific engineer may be required.

The available functionality of the Microscope Companion software depends on the microscope type, configuration, and software versions. Therefore, some functionality described in this user manual may not be available, and the appearance of the User Interface (UI) may differ from those shown.

# Release Notes of MiCo 1.17

This release includes the following features to be introduced on applicable systems:

- Field emission source and High-Tension state display with options to change the state on the home page
- Several additions for Autoloader (requiring a software update of Autoloader):
  - Autoloader Docking/Undocking on the actions page
  - Autoloader Cassette page with inventory for cassette and individual slots
  - Vibration Prevention for the automated alignments
- Scheduling through a new planning page
  - With a cryo-cycle or temperature mode change for Autoloader
- Eco mode to save power when the microscope is not in use



# Microscope Compatibility of MiCo 1.17

MiCo can be used with any Thermo Fisher Scientific and FEI Transmission Electron Microscope (TEM) with compatible software.

The best user experience of MiCo and access to the latest features is achieved in combination with the latest Microscope Software version. The table below details the range of compatible Microscope Software versions for this version of MiCo. It is recommended to upgrade MiCo to this version for the listed Microscope Software versions.

MiCo	Titan	Talos
MiCo 1.17	3.24 - 3.19	2.24 - 2.19

MiCo is (backward) compatible with six Microscope Software versions and this enables an independent and more frequent software upgrade of MiCo to deliver improvements on a quarterly basis without the requirement of a Microscope Software version upgrade. Of course, it may be that features are only available from certain Microscope Software version onwards and may not be found on your Microscope (just yet). Software version combinations outside of the range specified in the table are not supported.

# Installation and Start-Up

MiCo can be installed or updated in the following ways:

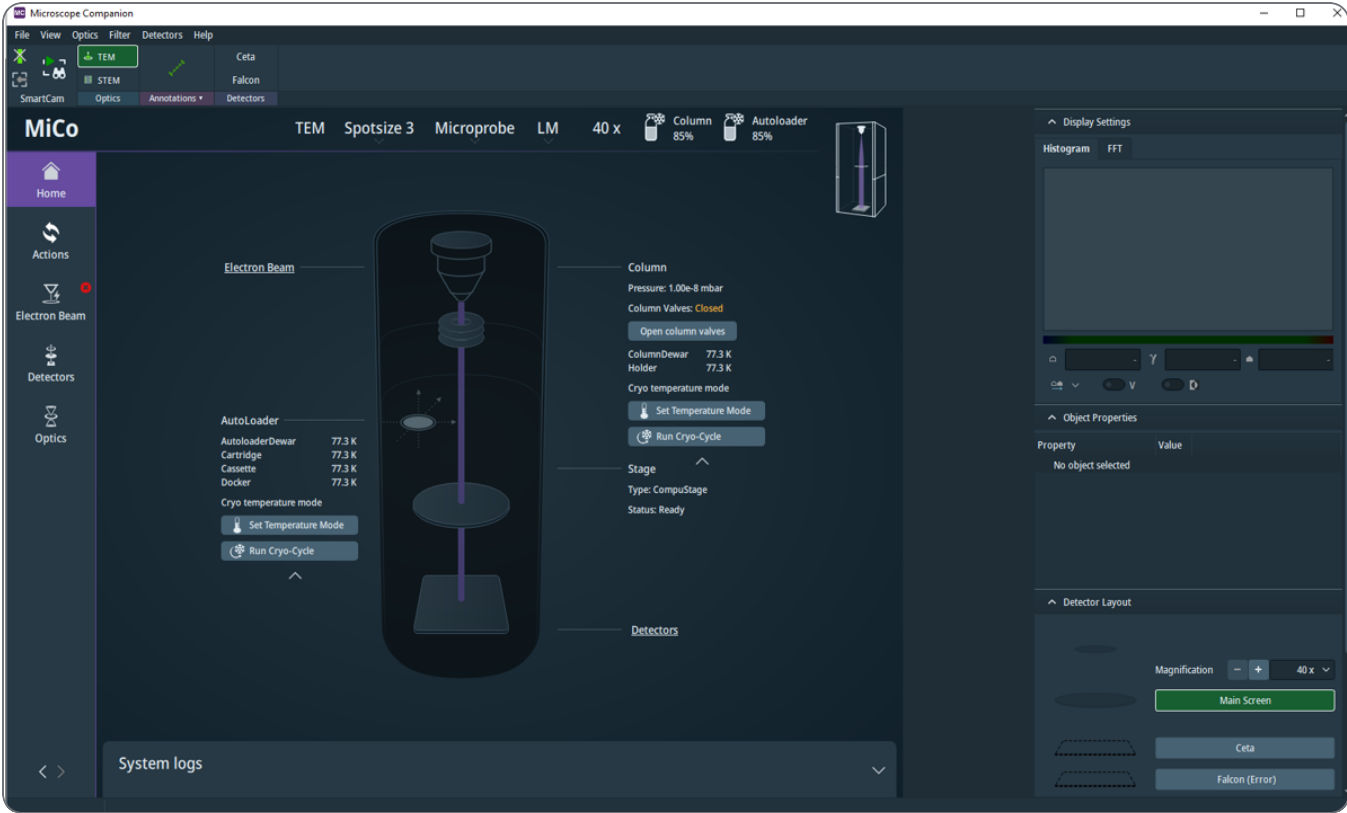
- Using an installer provided through the service, typically available quarterly
- Through the MyMicroscope service, if the microscope supports it
- As part of a Microscope Software installation by a qualified Thermo Fisher Scientific service engineer

The MiCo software is included in the TEM Server software installation and should be installed by a qualified Thermo Fisher Scientific service engineer during a microscope software upgrade. Additionally, the software can be installed or updated using an installer provided by a Thermo Fisher Scientific service engineer. To ensure compatibility with your microscope software versions, it is highly recommended that you upgrade to the latest version of MiCo. For compatibility specifics, please consult the Release Notes for your particular MiCo software version.

Before initiating the MiCo software, ensure that the TEM server is operational. To start MiCo, you can either use the Windows Start menu (navigate to All Programs > Thermo Scientific MiCo > MiCo) or the microscope software launcher. Upon starting the software, the main window will appear, as shown in the image below.



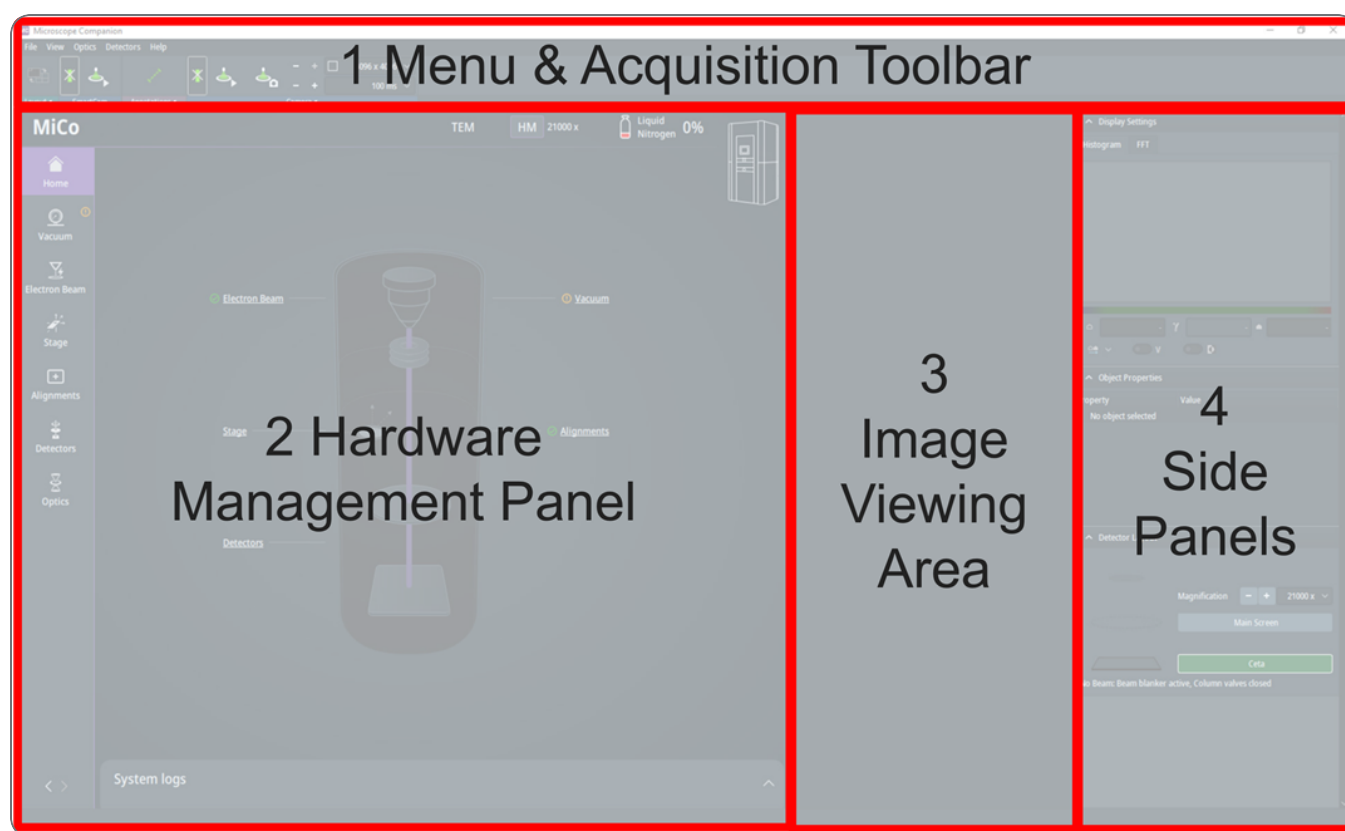
**Note:** The images in this document depend on the microscope and may look slightly different.



# User Interface Layout and Features

The Microscope Companion user interface is divided into different parts as shown in the following image. These parts are:

1. Menu and Acquisition toolbar.
2. Hardware Management panel.
3. Image Viewing Area is the location where the live-view windows of cameras and detectors appear in the UI.
4. Side panels.



## Menu and Acquisition Toolbar

The main menu area contains the following items.

Menu item	Description
<b>File</b>	Close the application.
<b>View</b>	Layout of the image windows.
<b>Optics</b>	Optical settings like magnification, optical mode (STEM/TEM), projection mode (Imaging/Diffraction) and probe mode (microprobe/nanoprobe).
<b>Filter</b>	Set the system to EFTEM mode and set up or modify settings of the energy filter on the system.
<b>Detectors</b>	Available cameras on the system.
<b>Help</b>	User manual and release notes.

The acquisition toolbar primarily focuses on sample navigation using live-view windows from cameras and detectors, rather than data acquisition, as seen in other applications. The toolbar consists of the items outlined in the table below.

Toolbar	Description / Main purpose
<b>SmartCam</b>	Start SmartCam live-view and beam blank button.
<b>Optics</b>	Change optical mode (TEM/STEM).
<b>Annotations</b>	Measure the dimension of a feature on a sample in the image window.
<b>Detectors</b>	Select a Camera or detector for live-view.
<b>Camera</b>	Set up properties of live-view of various cameras.
<b>Scan Rotation</b>	Change scanning direction of STEM live-view.
<b>STEM Imaging</b>	Set up properties of live-view of various detectors.
<b>Energy Filter</b>	Set up properties of the energy filter.
<b>Dose rate</b>	See and measure dose on the sample.

## Side Panels

The right side of the UI grants access to the following panels:

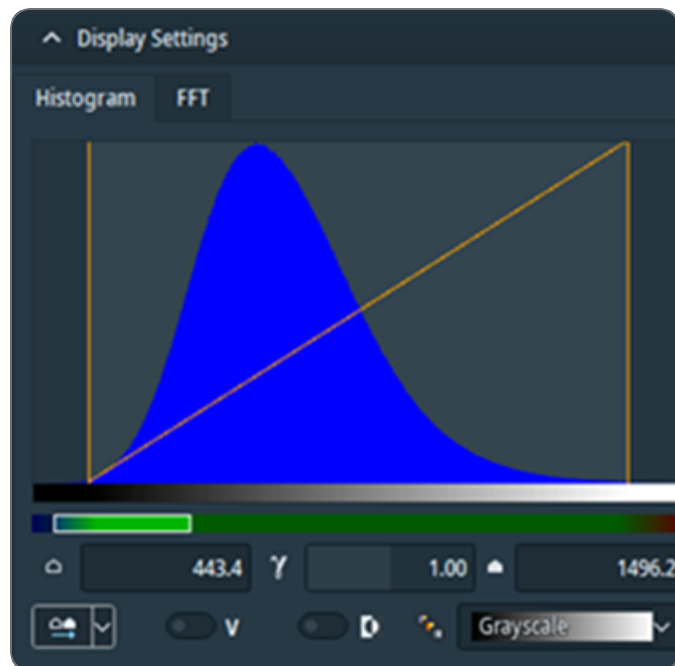
1. Histogram.
2. Fast Fourier Transform (FFT).
3. Object Properties, which displays parameters for a selected live-view or annotation, such as image size and conditions under which a live-view is being recorded.
4. Detector Layout.
5. Detector Settings, which offers parameters and functions to fine-tune the STEM detectors.

The Histogram, FFT, and Detector Layout panels are discussed in the subsequent sections. The Detector Settings panel is addressed in conjunction with live-views of STEM detectors.

## Histogram



The Histogram is shown for a live-view window as the first tab of the display settings panel and as shown in the figure below. It should be noted that it is possible to have multiple live-view windows open and there is only one histogram at a time. The Histogram is displayed for the selected image window where the title of the image window is high-lighted.

The purpose of the Histogram is to adjust the live-viewing of cameras without changing the controls of the camera or detector itself.



## Options of a Live-View Window



The appearance of the live-view is determined by a number of parameters:

- Black-level is the histogram intensity value represented by black. Any intensity value below the black-level will be black. In the figure above, the black-level of 443.4 is applied.
- White-level is the histogram intensity value represented by white. Any intensity value above the white-level will be white. In the figure above, the white-level of 1496.2 is applied.
- Gamma value is the gradient applied to the histogram intensity value. In the figure above, a linear gradient is applied with a gamma value of 1.00. The linear gradient is represented by the diagonal line in the graph. Select the **Gamma** button () to reset the Gamma value to 0.50 for diffraction pattern images, or to 1.00 for all other image types.
- Look-up-Table (LUT) describes the relation between the histogram intensity value and screen color for the range of black-level to white-level. There are a number of predefined LUTs available. In the figure above, the Grayscale LUT is used. It is also possible to invert the LUT (.

For a reference of Cubehelix LUT please see: [Green, D. A., 2011, 'A colour scheme for the display of astronomical intensity images', Bulletin of the Astronomical Society of India, 39, 289.](#)

## Adjust a Live-View Window

There are several ways in which to adjust the live-viewing window:

- Automatically adjust Black and White level () or White level () only. The automatic adjustment of Black and White level is the default option for all images, except diffraction pattern images.
- Manually interact with the graph by
  - Moving the vertical lines using the mouse to adjust the minimum intensity value (black-level) and maximum intensity value (white-level).
  - Changing the shape of the curve by moving the line using the mouse.
- Manually use the controls underneath the graph.

## Optimize a Live-View Window

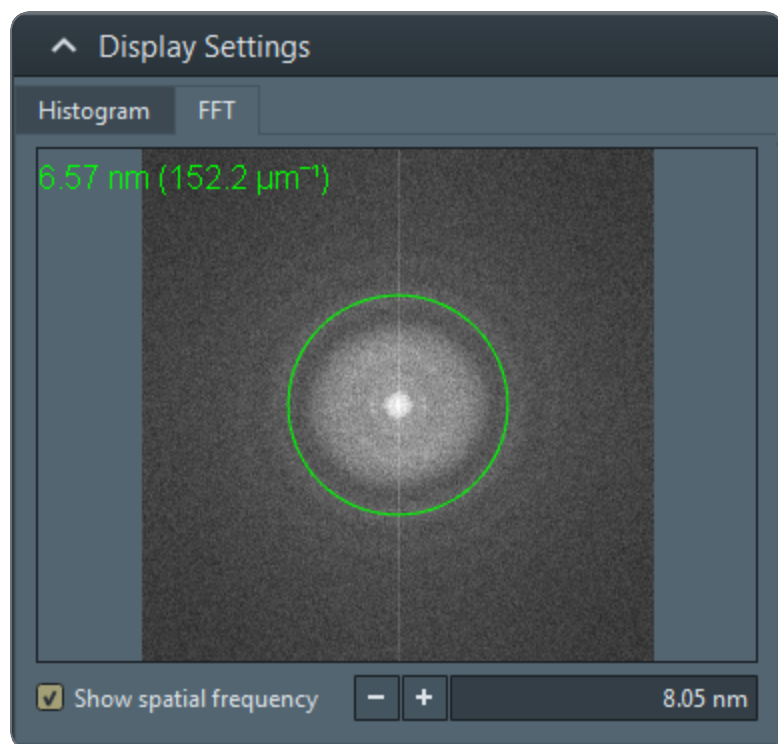
The Detector Saturation Indicator shows the saturation of the detector as shown below. The Detector Saturation Indicator can be used to optimize the imaging conditions, so the detector

captures the full range of intensities without overexposure.



## Fast Fourier Transform

The Fast Fourier Transform (FFT) can be found in the second tab of the Display settings side panel, as illustrated in the figure below. It displays a high-speed live FFT of the selected image window. The interactive window allows for zooming in or zooming out using the mouse wheel when the cursor hovers over the FFT.



You can draw a circle at a specified spatial frequency by selecting the **Show spatial frequency** check box and entering the value and unit (pm, nm, mm). Alternatively, double-clicking within the FFT sets the spatial frequency at the cursor's location.

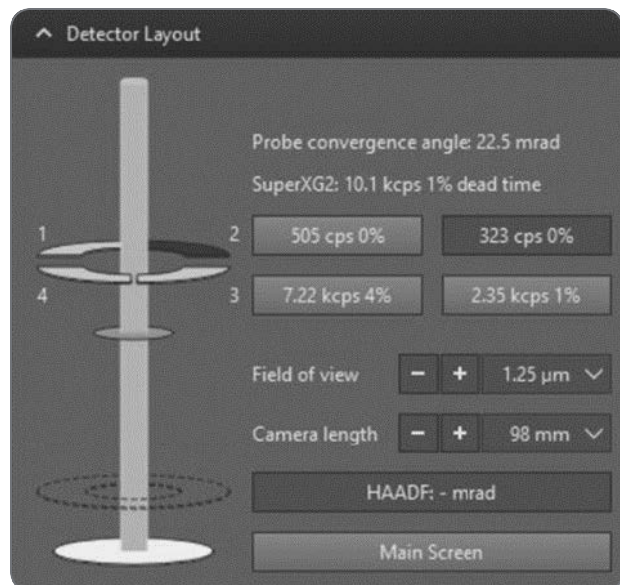
A right-click in the FFT reveals a menu with several additional options, including:

- Resetting the zoom.
- Zooming to one FFT pixel per display pixel (zoom 1:1).
- Utilizing interpolation (anti-aliasing) to enhance the smoothness of the displayed image.

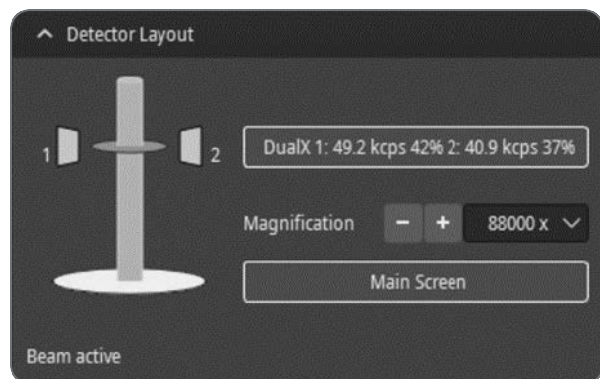


- Applying a Hamming window (windowed FFT) to reduce horizontal and vertical lines caused by image borders.

## Detector Layout



*Detector Layout panel with an HAADF and a Super-X G2 detector. For the Super-X G2 detector, segments 1, 3 and 4 are enabled.*



*Detector Layout panel with a Dual-X detector.*

Depending on the system configuration and Optics mode, the Detector Layout panel shows:

- Detector status or Detector Segment status**  
The inserted/retracted or enabled/disabled status for each detector or detector segment. Select the detector-specific button to toggle its status. Depending on the type of detector, additional properties can be displayed.

- Super-X and Dual-X sensors: the count rate and dead time percentage.
- STEM detectors: the collection angle.

- **Main Screen**

Insert or retract the FluScreen.

- **Beam status**

No Beam: Column valves closed, FEG Standby, High Tension off

If no beam is present, the reason is shown here.

- **Magnification**

Magnification   -   +   5.5 kx ▼

The available magnification values are nominal.

Select [+] or [-] or select a value from the drop-down list.

- **Camera length**

Camera length   -   +   44 mm ▼

Select [+] or [-] or select a value from the drop-down list.

- **Field of View**

Field of view   -   +   1234 nm ▼

The available magnification values are calibrated.

Select [+] or [-] or select a value from the drop-down list, or enter the desired value in the list box.

- **Proj. Mag.**

Proj. Mag.   -   +   6500 x ▼







Select [+] or [-] or select a value from the drop-down list.

# Hardware Management Panel

The hardware management panel is divided into four distinct parts: the Toolbar, the Header, the management pages, and the log window. The following sections describe each part.






## Hardware Toolbar

The toolbar grants access to pages as outlined in the table below. The selected page will be highlighted on the toolbar, with the button appearing in a purple color. Please note that pages marked with an asterisk (\*) are only available on supported systems.

Page	Icon	Brief description of main functionalities
Home		Status overview of main system functionality, Column valve status and open/close, Column pressure, Stage type and status
Actions		Cryo-cycle, Evacuate Column, Recover Source and High-Tension, Recover Stage, Recover Optics boards
Alignments*		Overview of automated alignments Executing automated alignments with progress indication, information and diagnostic images
Detectors		Camera status EDS status and controls
Optics		Change beam diameter in TEM imaging*
Planning		Schedule Cryo-cycle and temperature mode change for Autoloader*. View eco mode status*.

## Status Indications

The hardware management toolbar and panels provide feedback on the status of various subsystems, modules and functionality of the microscope. Status symbols and coloring are used to indicate status and it should be obvious where issues occur. The symbols and coloring are shown in the following table.

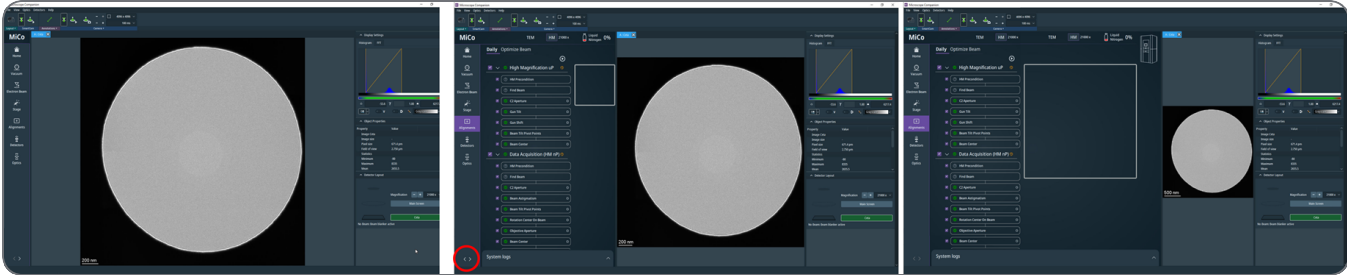
Symbol	Status	Remarks
	Good, Ready, On	Good, Ready, On
	Warning	Indicates something may require attention. For example, it indicates that the column valves are closed on the vacuum page. These valves are opened automatically when required by some other actions on the microscope, like viewing an live-view.
	Expired	Indicates something is expired. Typically, the status of something like an alignment may expire after not being executed for a specific time period.
	Unknown	Not sure what the status is at this moment. Status will be overwritten after execution. However, it should be noted that some status may not be persistent and/or stored and the indication may return after restarting the application.
	Bad, Not Ready, Off	Something is not in an operational state and requires attention and/or recovery actions. For example, alignments may fail when they were executed under non-intended conditions.

## Resizing the Panel

The hardware management panel is fully open by default and the home page is shown when starting the application. The button for the currently displayed page becomes purple.

It is possible to collapse the panel completely by clicking on the icon of the selected page in the toolbar.

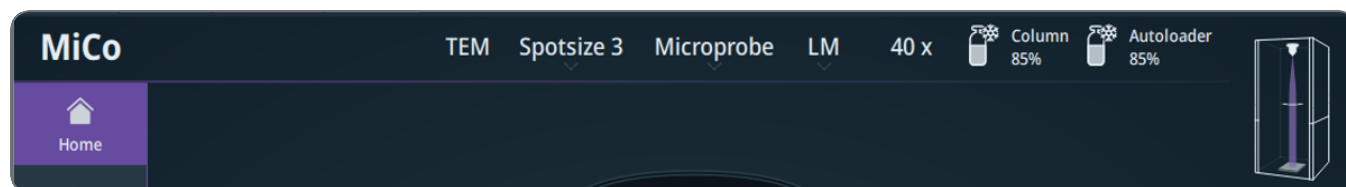
Additionally, the toolbar provides a way to re-size the Hardware Management Panel at the bottom as indicated in the figure below. The panel can be open (left), half-way open (middle) or closed (right).



## Hardware Management Header

The header of the hardware management panel, located above the pages, displays the system status as depicted in the figure below. The status elements in the header include:

- The optical mode, which may comprise TEM or STEM.
- The EF represents the switch between TEM and EFTEM (in case of energy filter)
- Spotsize (1-11).
- Probe mode (microprobe, nanoprobe).
- The Magnification range, which can be either High-Magnification (HM) or Low-Magnification (LM). Furthermore, you can toggle between the HM and LM by clicking the button.
- The Liquid Nitrogen level in the Column and Autoloader or Cryo Loading Station (CLS). The liquid nitrogen level will be shown on Autoloader systems only when in Cryo temperature mode. In all other cases, the highest temperature will be displayed to indicate the system is (not yet) in Cryo temperature mode. It is recommended to look at the home page for all temperatures available for column or Autoloader.



## Hardware Management Pages

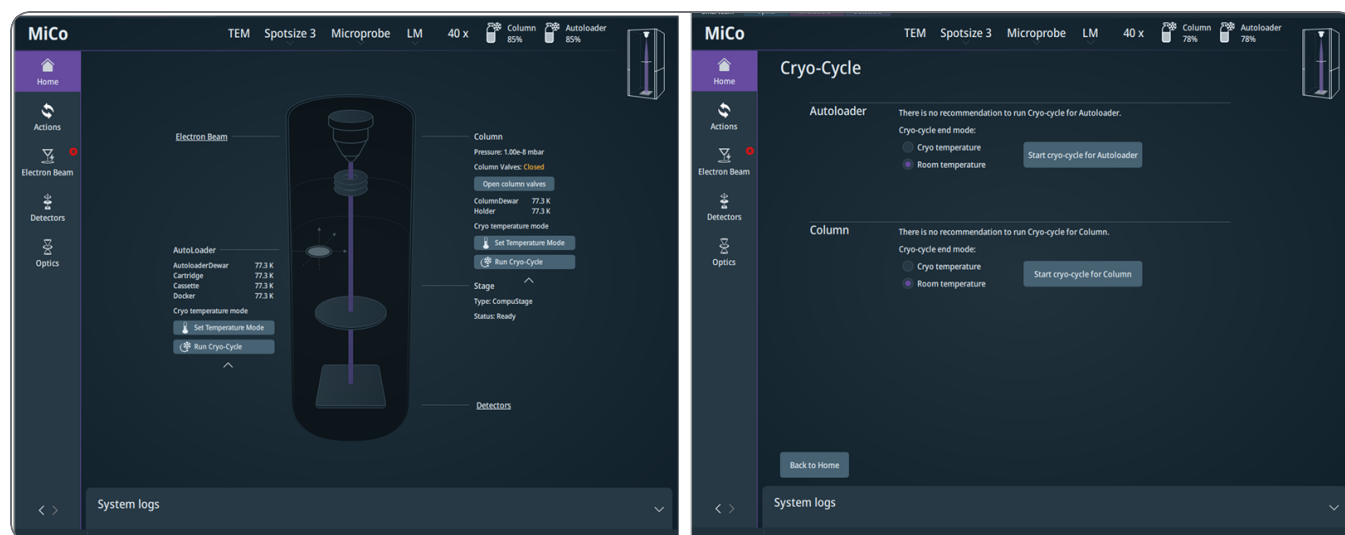
The toolbar provides access to various pages, which are described in the sections below.

### Home Page

The Home Page offers a status overview of the main system functionality, along with a generalized microscope image. It does not display the status of every component within the microscope.

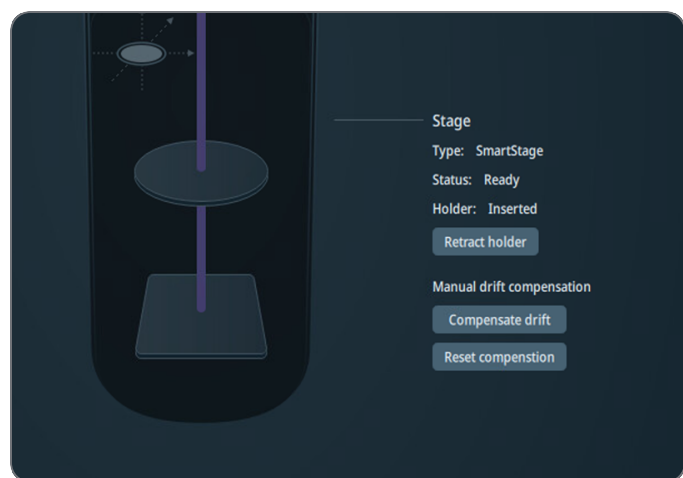
The page contains control over the column valves and shows its status. Also, it will show something when in Eco Mode.

It shows stage type and status. Optionally, holder status and insert/retract with progress information is available on supported systems.



## SmartStage Functionality

Additional functionality will be available on the Home page on microscopes equipped with a SmartStage (Metrios only). The microscope will have an extended Stage section with status information as shown below.



The page provides access to the automatic insert and retracts of the holder with progress reporting.

Additionally, the SmartStage supports a manual drift compensation that can be applied directly after a holder is inserted using a live-view of a camera or STEM detector to observe the drift. The user is left to determine the drift compensation vector. The drift compensation allows to capture of an image under quasi-static conditions. However, the drift will change over time as the holder will (thermally) stabilize inside the microscope. Therefore, the manual compensation should or

could be reduced to zero after some point in time. It should be reduced to zero (reset) when the (automated and) active drift compensation takes over. The compensate drift button provides access to the drift compensation graph. Here, a three-dimensional drift compensation vector can be set. The two-dimensional graph corresponds roughly to the imaging plane. The vertical direction can be compensated if the image is moving out of focus. The actual compensation vector is shown below the graph.



**Note:** The compensation is not automatically reset and is left up to the user. It is recommended to reset after holder stabilization within the microscope or after holder retraction.

**Note:** The scan rotation in STEM mode or TEM image rotation is not taken into account.

## Actions Page

The Actions page displays multiple buttons and allows scrolling when there are too many to fit on the screen.

Evacuate Column (previously known as "evacuate all")

Eco Mode (see, [Eco Mode](#) section)

Dock/Undock Cassette (see, [Auto Loader](#) section)

Recover Source and High-Tension, the following should be taken into account:

- The Recover Source and High-Tension button enables to start source from an off or power-on state to emitting.
- Set the High-Tension to a default value, which may not match the intended (maximum) value for the microscope. Adjust the High-Tension value on the home page.
- The system reports the progress of the source, which may take more than 10 minutes to reach an emitting state.
- It is compatible with all FEGs, but not with Thermionic sources.

Recover Stage, previously known as Enable stage.

Recovery of Optics Board degradation.

Recover Autoloader, Autoloader temperature system, and column temperature system with recommendations if available. The temperature system recovery will only be enabled when needed, but the Recover Autoloader will remain enabled. It is recommended to try this functionality in case of problems experienced with the Autoloader. If it does not resolve your issues, then it is recommended to contact your local service engineer.

Start Cryo-cycle for supported systems.

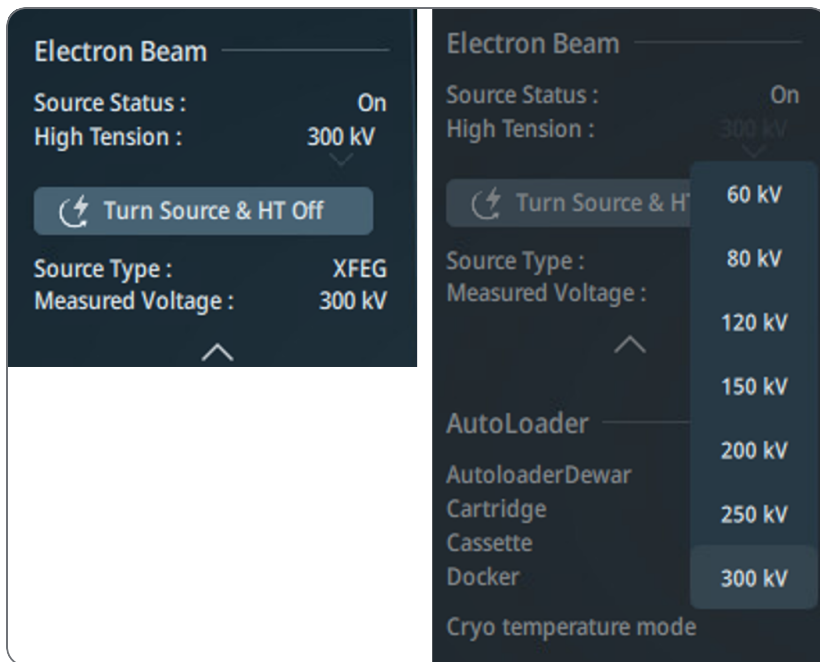
Start Heating Cycle for supported system, documentation for this feature will be provided elsewhere.

## Source and High-Tension

The Home page provides access to Source and High-Tension (HT) functionality. Users can view and switch the Source and HT states. This feature applies only to Field-Emission-Gun (FEG) sources and is not supported for thermionic sources.

For most FEG sources, the Source and High Tension (HT) are coupled and can only be turned on or off together. This process is automated. The HT value should be selected after they are turned on, as it will go to default values and not automatically go to the highest HT value. The HT value is shown when the HT is on, otherwise it will show the Off state. It is possible to change the HT value using the dropdown when hovering over the value.





Only for Cold-FEG it is possible to switch the state of source independent of the HT and it has the option to go to Power On state in addition to either On or Off. See screenshot below.




**Note:** The Source status is not supported for Thermionic sources

## Alignments Page

The Alignments Page is only available on systems with automated alignments, such as Tundra. The page displays various sequences with automated alignment modules. These modules are tailored to specific systems and workflows for other applications, like EPU.

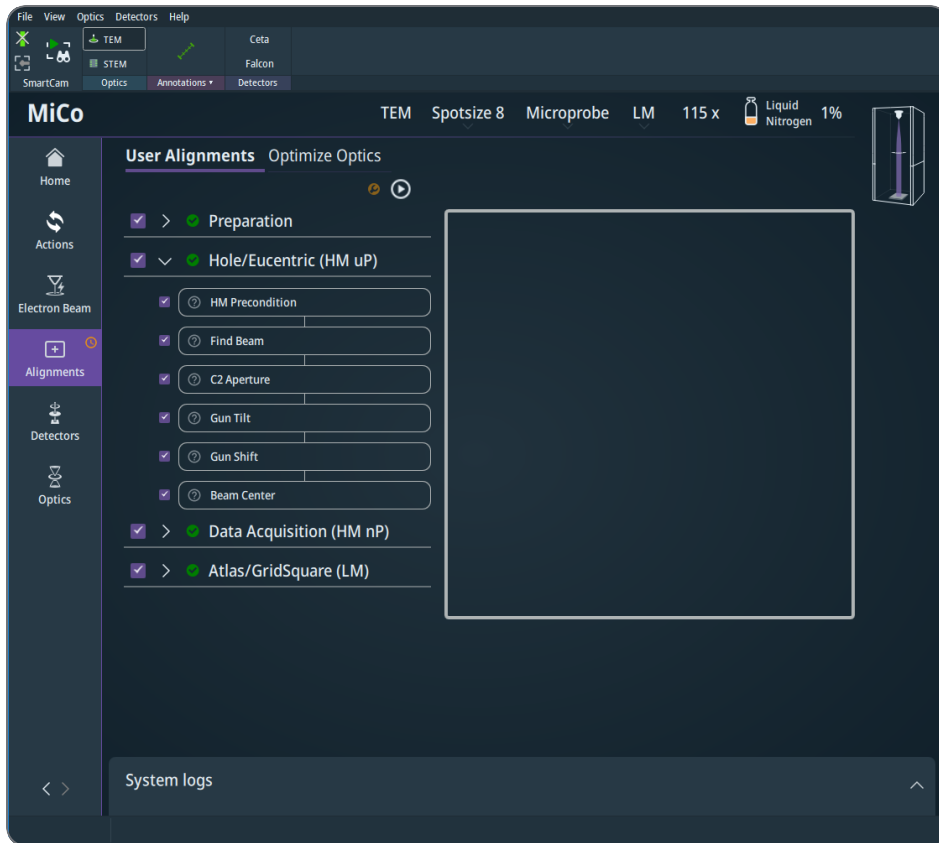
Consider the following points when interacting with the Alignments Page:

- Automated alignment modules offer an alternative to manual alignments using the TEM user interface. However, it should be assessed (by the supervisor) whether the alignments should be performed on the microscope.
- These automated processes are designed to operate under the default microscope settings defined in the sequences. Deviations from these defaults may result in module failure.
- Most (but not all) alignment modules have default expiration times. These should be used as guidelines rather than strict timelines.
- Most (but not all) alignment modules store the status. Some precondition and fast or necessary modules do not store the result and may show unknown status after restart.
- The execution of any automated alignment module can be initiated on-demand from MiCo, regardless of its expiration time.
- Adjusting microscope settings while automated alignments are running is not advised, even though it may be technically possible.

 **Note:** The SmartCam live-view is not automatically paused and its continuation may hinder successful execution of alignments (Known Issues).

This page includes:

- Alignment status, error messages, and expiration times.
- Alignment sequence execution with progress messages and diagnostic images.
- Start/Stop: Begins the automated alignment procedure for the selected Alignments set.
- Other features: Individual alignment selection, progress indication per running alignment, and messages indicating the progress and results of the automated alignments procedure.



It should be noted that the microscope state is automatically saved upon successful completion of each complete alignment sequence. In case of a failure for one or more modules, it is possible to manually perform these alignments using TEM User Interface. In such case, it is recommended to save the state (wrench icon) of the microscope manually after finishing.

In the case of Tundra microscopes, it's possible to manually save/load the state to perform alignments using the TEM user interface as well. This procedure typically involves saving the state, opening the TEM UI, loading the state, executing alignments, saving the state again, closing the TEM UI, and finally, loading the state.

## Alignments Sequence

### User Alignments Sequence

This sequence performs the basic alignment of the microscope. It is meant to run without a sample, and aligns appropriate alignments in LM, HM  $\mu$ P, and HM nP modes.

It consists of a Preparation part, and three sub-sequences. The sub-sequences and their purpose are as follows:

- **Vibration Prevention**

Makesure that the automated alignments are not interrupted by vibration sources such as automated liquid nitrogen filling. This also applies to the Optimize Optics Sequence.

- **Hole/Eucentric (HM  $\mu$ P)**

Roughly aligns the system in HM  $\mu$ P mode at a magnification defined in HM Precondition block (about 45 kx).

- **Data Acquisition (HM nP)**

Aligns the system in HM nP mode at a magnification defined in HM Precondition block (above 100 kx) to prepare it for the Data Acquisition.

- **Atlas (LM)**

Aligns the system in LM mode at its highest magnification (around 2 kx) to prepare it for the Atlas screening.

### Optimize Optics Sequence

This sequence performs fine tuning of the microscope. It is meant to run on a sample and aligns appropriate alignments in HM nP mode.

It is executed on the current optical settings. In particular, the following parameters are used:

- Gun lens (does not apply for Tundra)
- Spot size
- C2 Aperture
- Magnification
- Objective Aperture

### Alignments and Blocks Description

This blocks can be divided into two groups system actions, and Alignments.

#### System Actions

##### Load Aligned State

- Makes sure that the initial state of the microscope is the last aligned state.
- It loads the \*APM\*.alg file located in C:\Tecnai\Alg folder (as it comes aligned from factory and service).
- It also loads the last aligned state of all apertures (currently, this file is not available for user

access).

### HM $\mu$ P Preconditions

- Sets the default optical settings for the HM  $\mu$ P sequence. These settings are:
  - Sets Eucentric focus
  - Probe mode:  $\mu$ P
  - Magnification:
    - Tundra, F200C: 43 kx
    - Glacios, Arctica: TEM 45kx, EFTEM 540 kx
  - Spot size: 4
    - Glacios with C-FEG: variable (based on Gun lens value)
    - C2 aperture: 150  $\mu$ m

### HM nP Preconditions

- Sets the default optical settings for the HM nP sequence. These settings are:
  - Sets Eucentric focus
  - Probe mode: nP
  - Magnification:
    - Tundra, F200C: 110 kx
    - Glacios non-FFI non-EFTEM, Arctica: 120 kx
    - Glacios FFI non-EFTEM: 45 kx
    - Glacios EFTEM (both non-FFI and FFI): 540kx
  - Spot size: 4
    - Glacios with C-FEG: variable (based on Gun lens value)
  - C2 aperture:
    - Tundra: 70  $\mu$ m
    - Glacios and Glacios FFI: 50  $\mu$ m

### LM Preconditions

- Sets the default optical settings for the LM sequence. These settings are:

## Hardware Management Panel

- Sets Eucentric focus
- Probe mode:  $\mu\text{P}$
- Magnification: Highest LM magnification
- Spot size: 4
  - Glacios with C-FEG: variable (based on Gun lens value)
- C2 aperture: 150  $\mu\text{m}$

### Optimize Optics Preconditions

- Sets the default optical settings for the optimize optics sequence. These settings are:
  - Sets Eucentric focus
  - Magnification: Glacios FFI non-EFTEM: 17500 x
  - C2 aperture: Glacios FFI: 50  $\mu\text{m}$

### Move Stage To Vacuum

- Moves stage to a vacuum position defined within the EPU.
- It is not possible to define the position in MiCo (this block is skipped when run from MiCo).

### Move Stage To Site

- Moves stage to a sample position defined within the EPU.
- It is not possible to define the position in MiCo (this block is skipped when run from MiCo).

### Restore C2 Aperture

- Inserts the C2 aperture defined by the user.

### Set Magnification For C2

- Decreases the magnification to see the whole beam in the FluCam Viewer.
- The target magnification depends on the size of the C2 aperture. The larger the C2 aperture, the lower the magnification.

### Restore Magnification

- Restores the magnification to the one defined by the user.

## Alignments

### Find Beam

- Assures the electron beam is visible on the FluScreen.
- It maximizes the screen current value by changing the beam position and its size.
- Find beam may also change the magnification to find the beam faster.
- Other alignments can internally call the Find Beam procedure as well, if they need help with finding the electron beam.

### C2 Aperture

- Aligns the important C2 apertures.
- It minimizes the beam movement while changing the beam size.

### Gun Tilt

- Centers the image of the extractor aperture onto the C2 aperture.
- It maximizes the screen current value by changing the position of the extractor's image with respect to the C2 aperture.

### Gun Shift

- Minimizes the beam movement while changing the Spot size value.
- It minimizes the beam movement while changing the Spot size values between 3 and 9.
- To do that, it centers the beam using a beam shift parameter on Spot size 9, while on Spot size 3, it centers the beam using a gun shift parameter.

### Beam Center

- Moves the electron beam to the center of the FluCam.
- It minimizes the distance from the center by changing the beam position.

### Beam Astigmatism

- Makes the electron beam circular.
- It minimizes the eccentricity of the electron beam's shape using the condenser stigmators.

### Objective Aperture

- Centers the objective apertures with respect to the electron beam in diffraction.

## Hardware Management Panel

- First, it finds the target position by focusing the beam to a spot in diffraction.
- Then, it spreads the beam to see the objective aperture edge and minimizes the objective aperture distance from the target position using the objective aperture movement.

### Parallel Illumination

- Sets the parallel illumination on the sample.
- It switches to diffraction mode and minimizes the beam size.

### Image Astigmatism

- Minimizes the astigmatism within an image.
- It acquires the image of an amorphous area, performs a FFT of the image, fits the Thon rings, calculates the image astigmatism, and minimizes it using the objective stigmators.

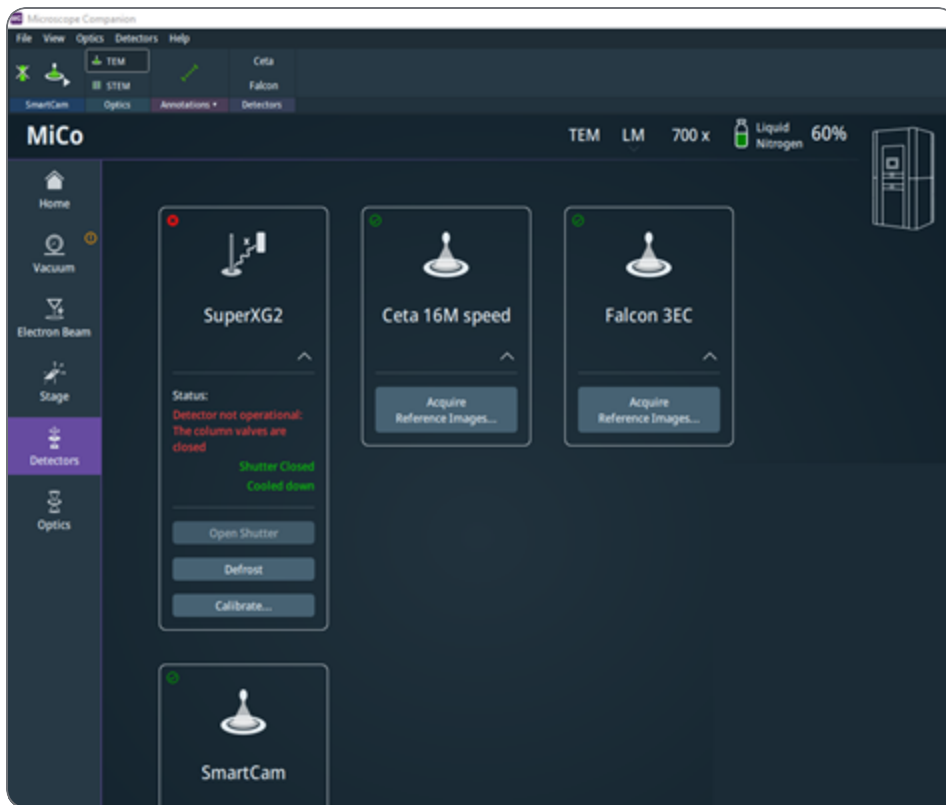
### Image Coma

- Minimizes the coma aberration within an image.
- It acquires more images for different electron beam angles, calculates the coma value, and minimizes it using the beam tilt parameter.

## Detectors Page

The Detectors Page contains tiles related to the detectors available on the system, such as the SuperXG2, Ceta 16M speed, Falcon 3EC, and SmartCam in the following image.





## EDS detector section

Depending on the type of EDS detector, the following functions may be offered:

- **Shutter / Position**

Depending on the type of detector, either:

- Shutter status: *Open* or *Closed*
- Detector position: *Inserted* or *Retracted*

The microscope software has automatic safety functions. Depending on the status of the microscope and its subsystems, these safety functions can deny the execution of actions that could damage the EDS detector.

- **Temperature**

The cooling status of the EDS detector.

- **Status toggle**

Depending on the type of detector, either:

- Shutter status: *Open* or *Close*

- Detector position: *Insert* or *Retract*
- **Defrost**  
Warm up the EDS detector.  
While warming up, a timer is displayed below the Defrost button. The defrost procedure can be aborted at any time.
- **Calibrate**  
Opens an external calibration tool.
- **Status**  
Displays status and progress messages for the EDS detector.
- **External Tools**  
Opens external tools such as EDX Calibration.

### Energy Filter section

The Energy Filter section shows the name of the filter and information on its status.

### Camera section

The Camera's on the system are displayed.

## Optics Page

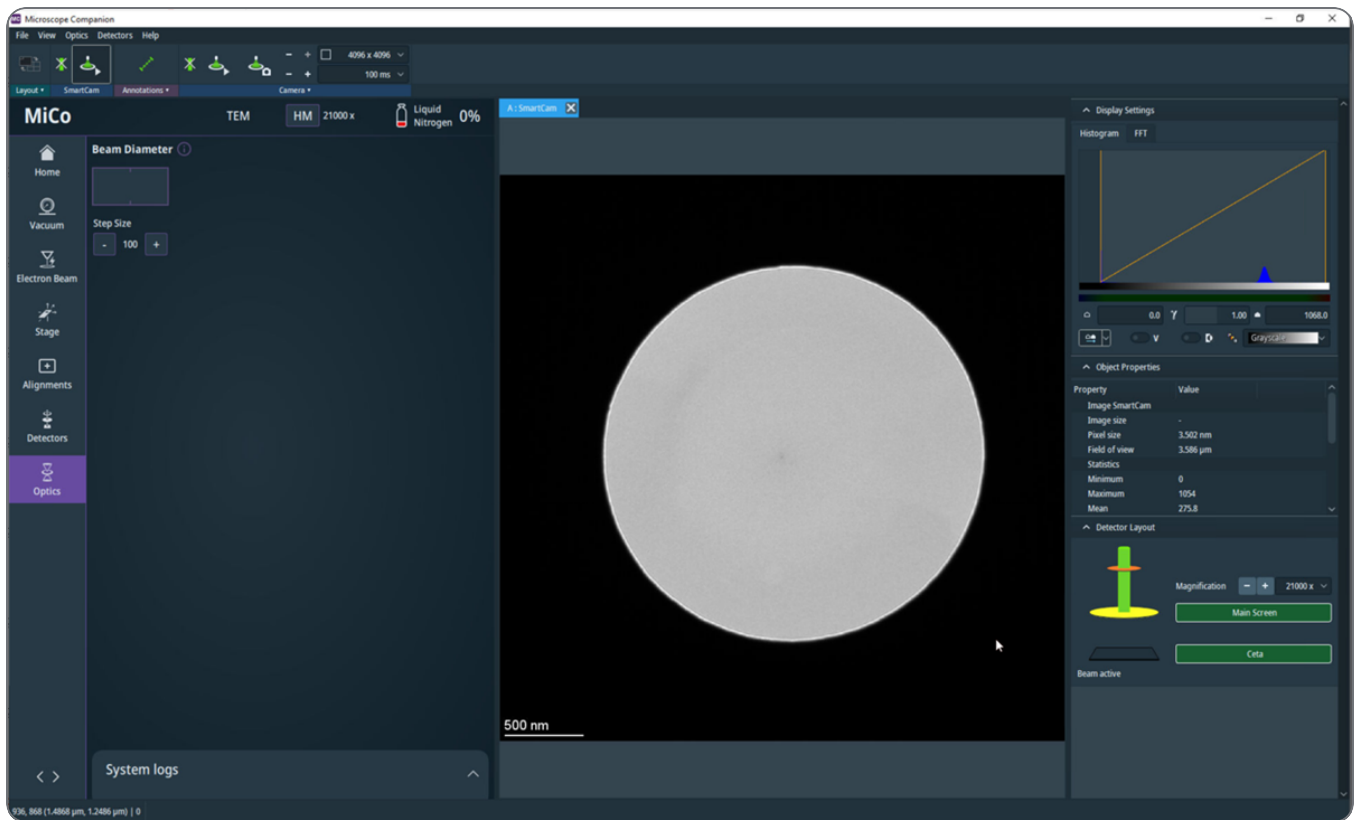
This page includes control of the beam diameter, only available for TEM imaging on supported systems.

### Changing beam diameter

It is recommended to use a SmartCam or camera live-view and half-way open hardware management panel, as shown in the figure below. The beam diameter control consists of a purple-boxed area with two options for manipulating the beam diameter:

- Use the mouse-wheel when the cursor is over the boxed area.
- Use mouse-clicks on the two halves of the purple-boxed area, where the left and right halves represent the mouse buttons.

Currently, no keyboard shortcuts are available for adjusting the beam diameter.



## Autoloader Functionality

**Note:** The Autoloader functionality in MiCo requires an upgrade of Autoloader Software through a Microscope Version upgrade.

## Introduction

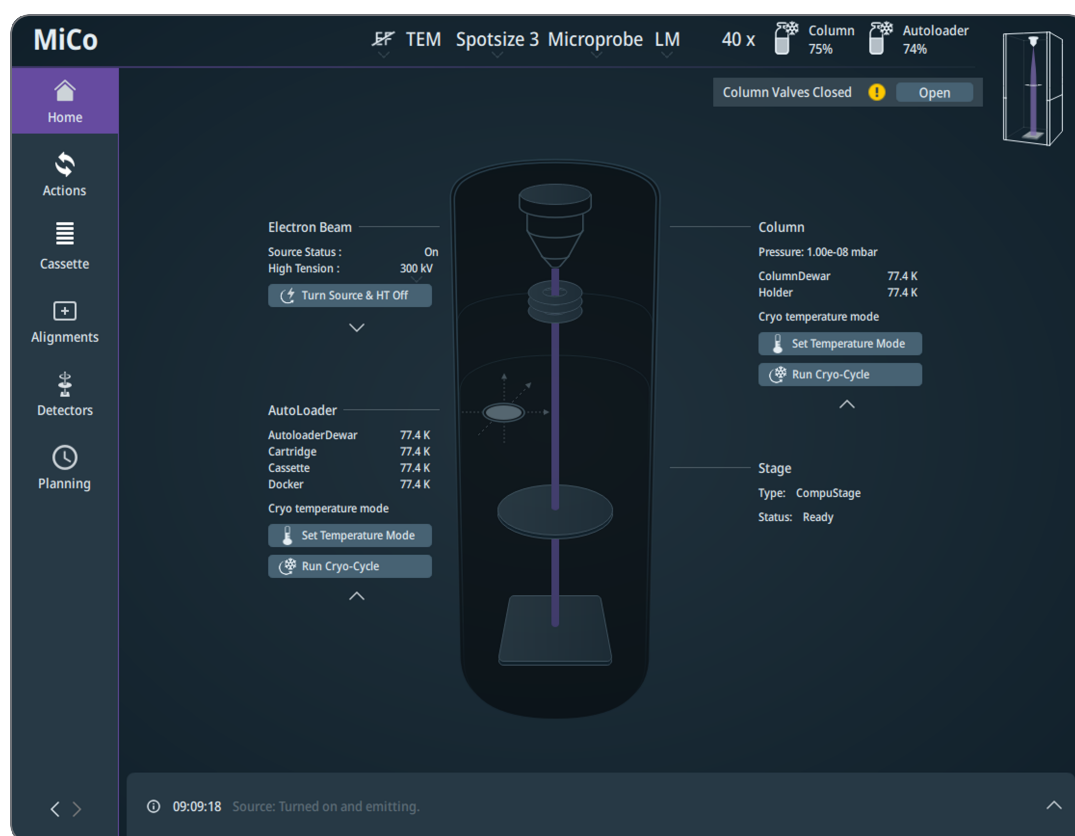
The autoloader functionality can be found on several locations within MiCo:

- The Header toolbar displays the liquid nitrogen levels of Autoloader and Column when in cryo-mode only. Otherwise, it will display temperatures.
- The Home page displays the temperature mode, temperatures and pressure information. Additionally, there is the ability to change the temperature mode and initiate a cryo-cycle instantly.
- The Actions page contains the Cassette docking/undocking and recovery functionality.

- The Cassette page contains the Cassette overview and the ability to run an inventory on the cassette. This can be done for the whole cassette at once, or by selecting individual slots.
- The Planning page contains the ability to schedule a cryo-cycle or temperature mode change.

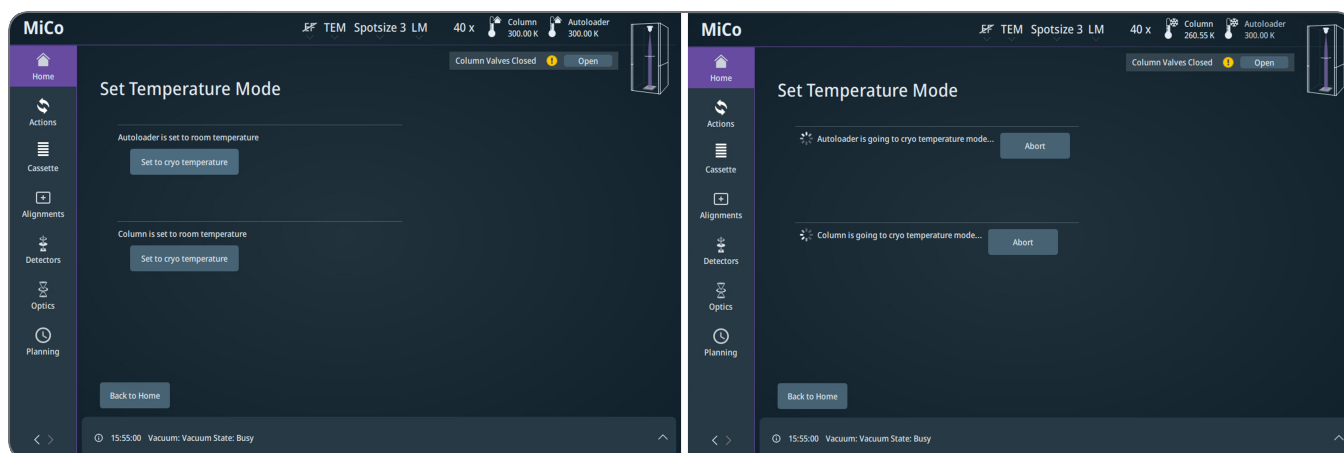
## Temperature handling

The Autoloader and Column can be set to room-temperature or cryo-temperature mode on the Home page. It should be noted that the Autoloader and the Column are treated as two separate compartments and the temperature mode and cryo-cycle are taken care of individually. It is possible to start them at the same time or independently.



## Set temperature mode

The temperature mode can be changed through the Set Temperature Mode button on the Home page. The button will lead to a page where the mode can be set or changed, see screenshots below.



The cryo-mode will be achieved by automatically filling the liquid nitrogen dewars in the system and waiting until the measured temperatures are below fixed thresholds. Liquid nitrogen will be filled automatically when the levels are running low while in cryo mode.

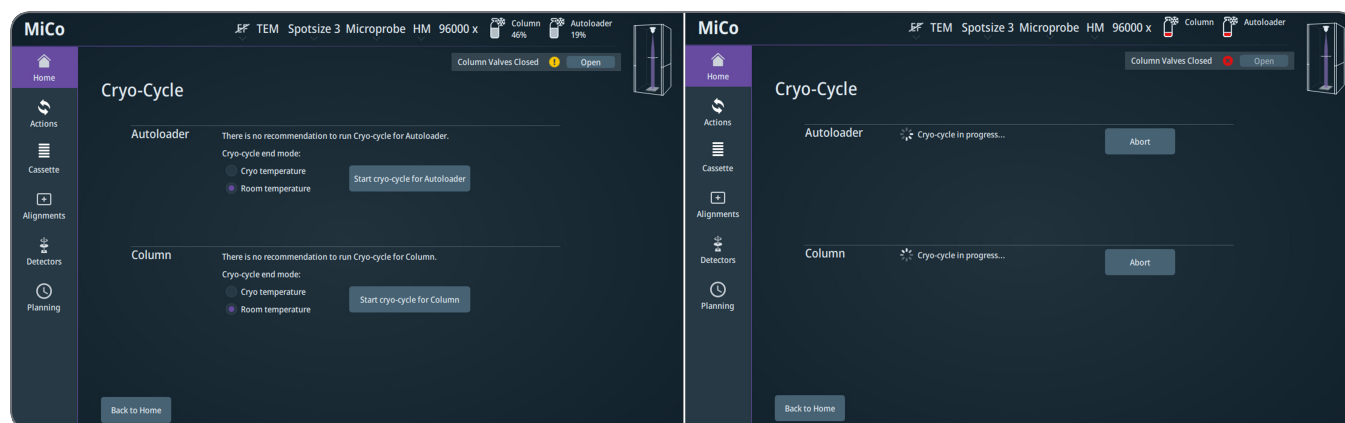
There are multiple temperature sensors in the Column and Autoloader compartments. Their temperature values are visible on the home page. There is currently no indication of whether the temperatures are above or below the thresholds for the cryo-mode.

Temperatures will be displayed in the header toolbar when transitioning to cryo mode and in room temperature mode. The change in temperature mode will show a blinking Autoloader or Column temperature in the header toolbar. The header will display the liquid nitrogen levels only when in cryo mode.

In case there are problems with the filling of liquid nitrogen, an error dialog will be presented. For example, it may happen when the liquid nitrogen in the vessel has run out and no liquid nitrogen level increase is detected during the filling.

## Run a Cryo-cycle

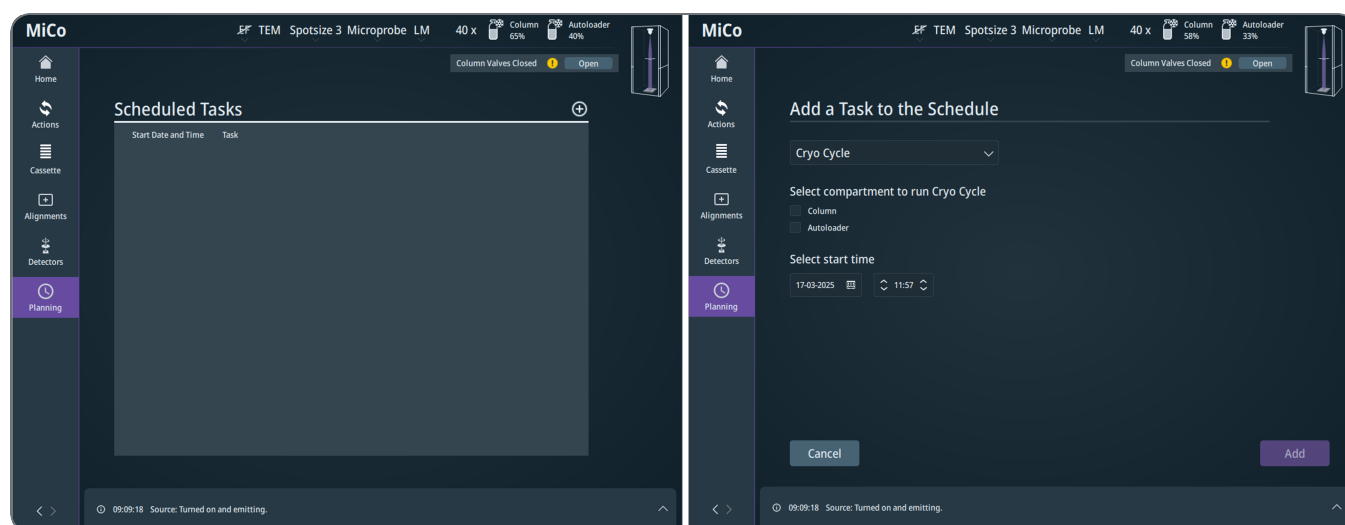
The cryo-cycle can be started from the Home page and the Run Cryo-cycle button will lead to the screenshots below.



The temperature mode can be selected after the cryo-cycle and the state will be set right after the cryo-cycle is finished. It should be noted that the temperature mode will only be set as long as MiCo is not closed as there is no scheduling of the temperature mode change in this process.

## Scheduling

The cryo-cycle and temperature mode change can be scheduled on the planning page. These two actions could then be scheduled independently. For example, it would be possible to schedule the cryo-cycle at a Friday night and do a temperature mode change early Monday morning. The planning of a cryo-cycle is shown in the image below. It is also possible to remove the scheduled actions from the planning page. It should be noted that the scheduled actions are not automatically cancelled or removed when the actions are triggered manually elsewhere, so they may fail to run or cause conflicts, and they are best removed manually as well after manually performing the operation.

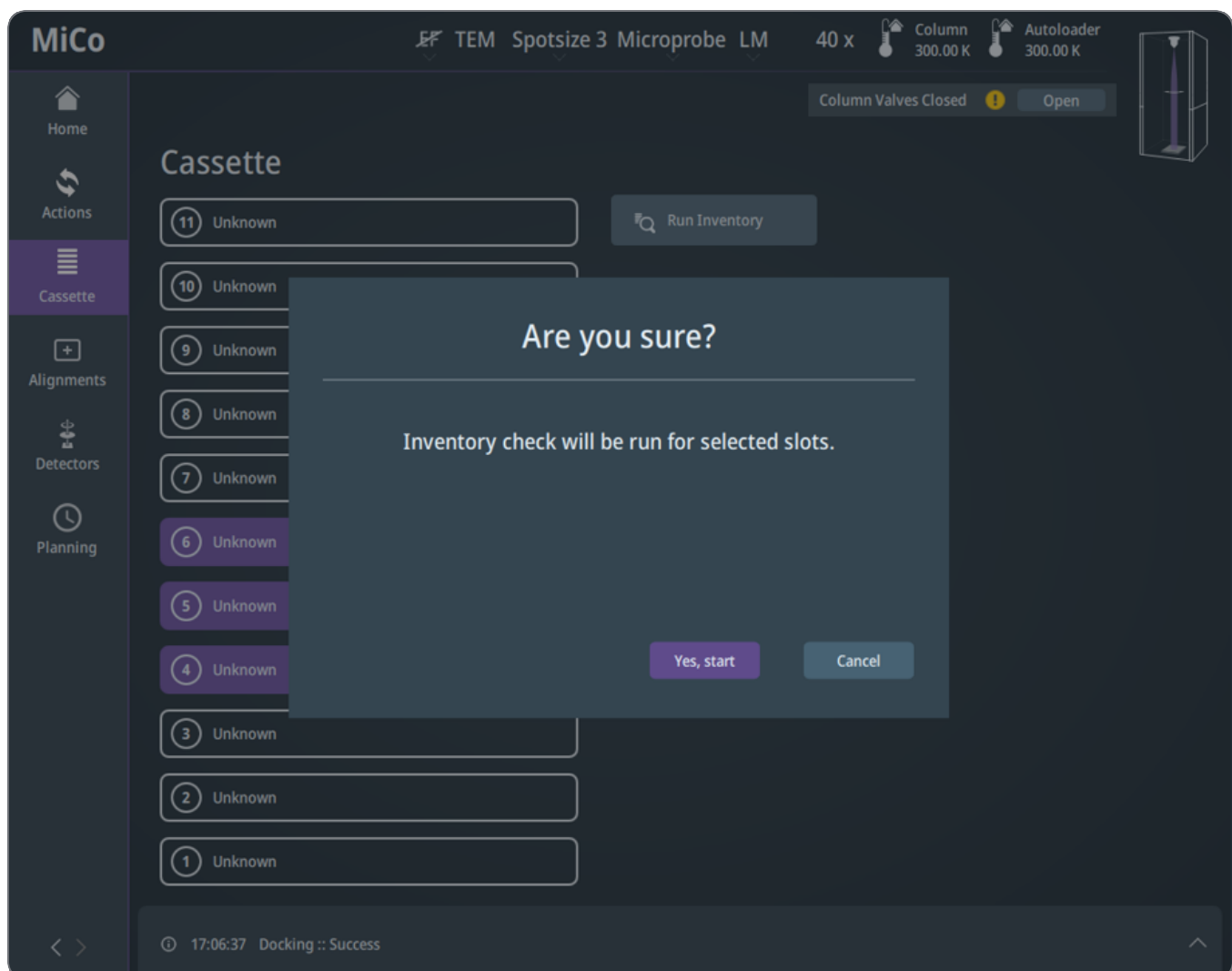


## Sample handling

The Cassette can be docked or undocked on the Actions page within MiCo, or using the On-System-Display at the microscope.

Once the cassette is docked, it will be possible to run the inventory on the Cassette. Otherwise, the button will be disabled. By default, the inventory will run on the whole cassette when no slots are selected. The button can then be used to abort the inventory, but it will finish the slot that is in progress.

It is possible to run the inventory on specific slots that can be selected by clicking on them. Multiple slots can be selected as the screenshot below shows.



### Recovery

There are three recovery actions available on the Actions page. One is a general Autoloader recovery, which is recommended when Autoloader is not behaving as expected. The functionality will assess the state of Autoloader and try to automatically resolve any issue.

The other two are for the temperature system of Autoloader and Column and will only be enabled in case of an issue encountered.

### Known issues

1. The Autoloader and Column will be filled automatically when in cryo-mode. There is no information or notification available within MiCo when the filling will happen.
2. There is no option to load/unload a sample.
3. The cassette page always shows the 12 slots of the cassette, even though there may not be a Cassette docked in the Autoloader.
4. Running a cryo-cycle or changing the temperature mode will show progress, but no time information is currently available.
5. There is an "undefined state" in the planning of a cryo-cycle, which is not recommended to use and will be removed. It is advised to use either room or cryo temperature selection.

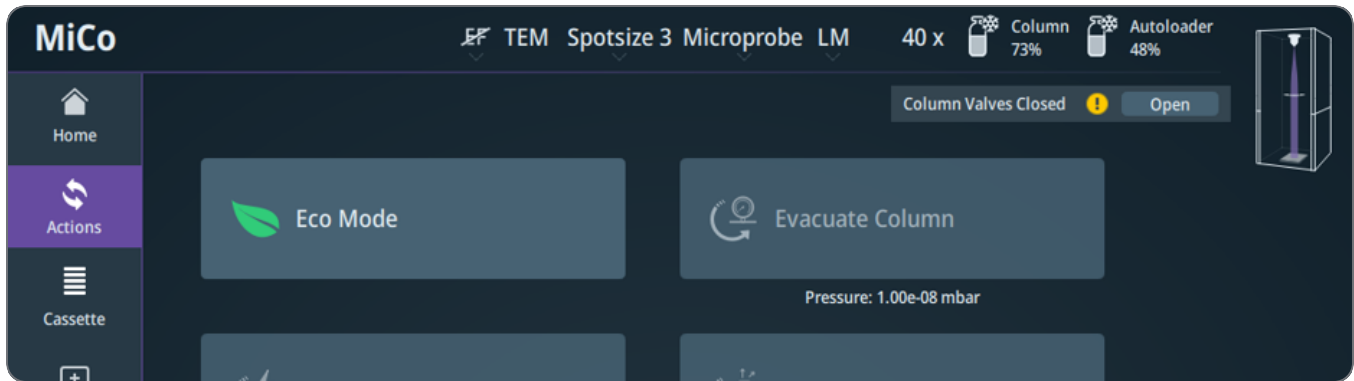
### Eco Mode

Designed for sustainability, low-power Eco modes deactivate parts of the microscope when not in use, reducing power consumption by up to 80%, lowering the carbon footprint, and cutting running costs. There are two mode available, standby saving 5-10% of power and sleep saving more than 80% of power. The process of going through the eco mode takes time and the following should be considered:

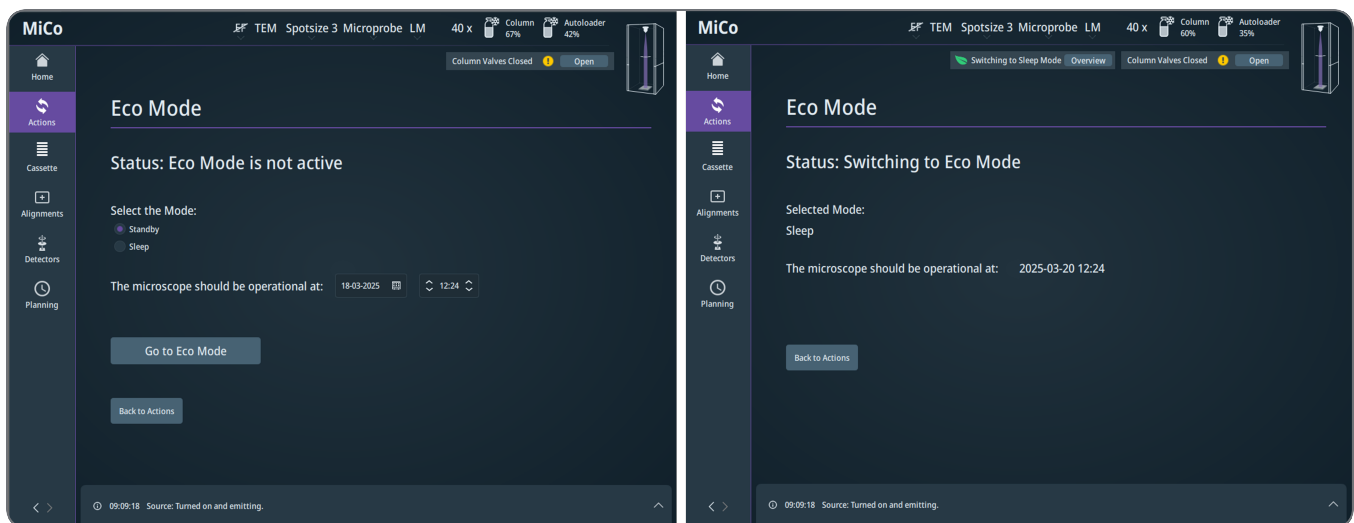
- The standby mode is best used when there is 24 hours or more idle time.
- The sleep mode is best used when there is more than 48 hours of idle time and is recommended over the weekend or holidays. It is not recommended to schedule more than a week of sleep for the vacuum may deteriorate over such time.
- The microscope will not be operational during this time and making changes to the state of server or components through MiCo and TEM UI is not advised while it can lead to issues with the automated process.
- It is recommended to remove any sample before initiating the sleep mode.

The Eco mode can be found on the Actions page.



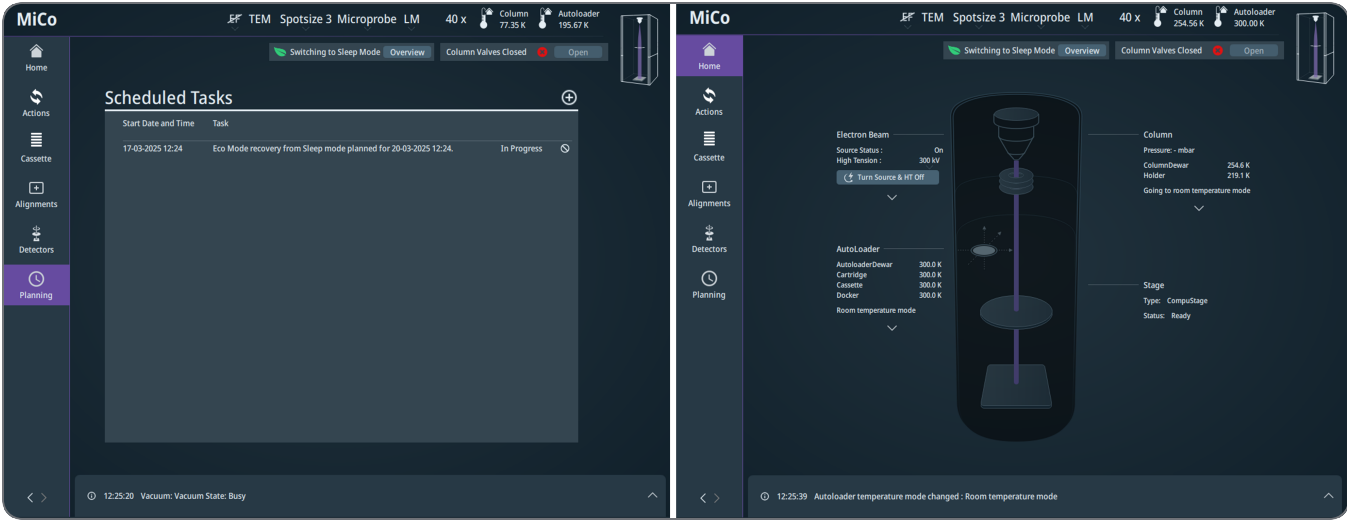


The button opens a page where the standby or sleep mode can be selected and the operational time of the microscope can be scheduled, see the screenshot below. The microscope will switch to the selected mode, stay in the mode for as long as the time allows, and it will be recovered at the requested operational time.



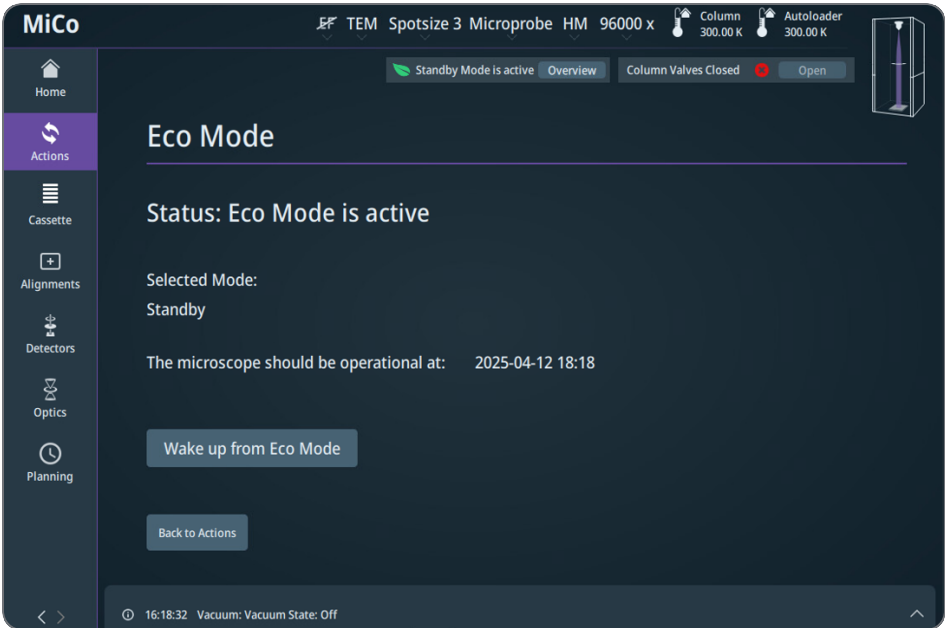
There will be an entry on the planning page when the eco mode is in progress and it is possible to see when the microscope will be operational again from there as well. On the top of the pages, there is also an eco mode status reported with the option to go to the eco mode overview page (similar to the screenshot on the right side as shown above).

# Hardware Management Panel



In case of Autoloader, the sleep mode will transition the microscope to room temperature. It is possible to schedule the transition to cryo-mode some time before the microscope is operational again. In such case, it is necessary to make sure sufficient liquid nitrogen is available.

It is possible to stop the Eco Mode early by using the Wakeup button on the overview page as shown below. This will immediately trigger the recovery, and the microscope will be operational as soon as possible. It should be noted it will still take considerable time from this point when in sleep mode. Also, the wakeup will invalidate the scheduled task on the planning page and it will report a failure.



# Live-View of Cameras and Detectors

The live-views of the cameras and detectors can be initiated from the Acquisition Toolbar. The toolbar also includes a beam blank button for manually blanking/unblanking the beam. Typically, the required camera settings include image size and exposure time. The live-view of the cameras and detectors will be displayed in the Image Viewing area of the User Interface.

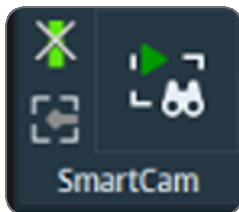
The following topics are discussed in separate sections:

- SmartCam (previously known as FluCam).
- STEM detectors.

## SmartCam

The SmartCam is a low-resolution and low-sensitivity camera that can be used to set up the microscope or for sample navigation, for example. The camera looks at the retractable Main Screen (a fluorescent plate, previously known as FluScreen) inside the microscope as indicated in the detector layout on the right side.

The live-view is started from the Acquisition Toolbar at the top left of the application. The small SmartCam toolbar contains a button for blanking and unblanking the beam. Also, it has a button to insert the beam stop, which will alternate between retracted, half-inserted, and fully inserted.



## Automatic Start of SmartCam Live-View

The SmartCam toolbar initiates the automatic start-up of a SmartCam live view. The start-up consists of the following steps:

- A Message Box is presented to ask whether the Column valves should be open to see a beam on the camera.
- The Main Screen is inserted. Additionally, the Main Screen can be manually inserted and retracted from the detector layout.
- The beam is unblanked and is visible in the detector layout. The beam blank button is

available next to the SmartCam start button in the Acquisition Toolbar.

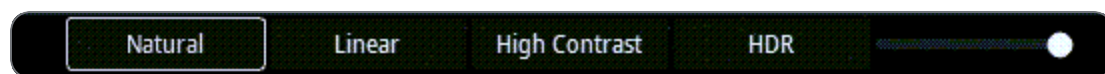
- A SmartCam live-view is shown in the Image and Data Viewing area in the middle of the application.

## Manual Actions for SmartCam Live-View

There are several possible manual actions:

- It is possible to run SmartCam with a blanked beam or with column valves closed, although the image will be black. The beam can be blanked by clicking the button next to the SmartCam start button. The column valves can be closed via the Home page.
- It is not possible to run SmartCam with a retracted Main Screen.
- The appearance of the live-view can be adjusted using the histogram (Black-level, White-level, Gamma, Color Lookup Table).
- Right-clicking with the computer mouse on the SmartCam display provides access to additional features such as the data bar and screen markings.

## Modes



The SmartCam display can be set in different modes to optimize the imaging conditions.

- **Natural:** Natural mode displays the image as it would appear on the flu-screen. The camera sensitivity is adapted to the highest intensity in the image. This mode is most suitable for general use.
- **Linear:** In Linear mode gamma is set to 1, this mode gives more contrast image compared to *Natural*.
- **High Contrast:** Features with very low contrast can be difficult to observe in *Natural* mode. *High Contrast* adapts the gamma level to produce more contrast compared to Linear and Natural.
- **High Dynamic Range:** Some images (mostly diffraction patterns) have an extreme intrascene dynamic range, e.g. there are both bright and dark areas with information. Adapting the exposure time is not enough, the camera lacks dynamic range to image both the bright spots and the patterns in the background. To enhance the dynamic range the exposure time is alternated between a short and long exposure and an image is

reconstructed from these images. To fit the dynamic range in the display a gamma curve is chosen that enhances the dark areas and suppresses contrast in the highlights.

Several exposure time pairs can be chosen in the camera control panel or by using the mouse wheel in the image.

- **Manual (using the slider):** Automatic adjustment is paused, the camera parameters can be adjusted by hand by using the slider on the SmartCam display.

## Notes on SmartCam

The reference images for SmartCam (dark and gain) are applied automatically and should be recorded elsewhere (FlucamViewer or the new Reference Image Manager from the Microscope Software Launcher).

## STEM Detectors

### Use the STEM Live-View to Prepare and Acquire a Single STEM Image

The basic steps to acquire a STEM image with MiCo are as follows:

1. Select the **Optics > STEM** mode.



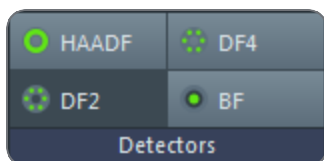
2. Insert the **Detector Layout > Main Screen** and open the column valves.
3. Select the **SmartCam** to start a live view of the Main Screen.



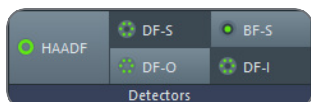
A Ronchigram should be visible in the SmartCam live view.

4. Use the **handpanels** to optimize the Ronchigram.
5. Depending on the configuration, insert one or more detectors.  
Either:

- In the **Detectors** toolbar, activate one or more detectors or detector segments.

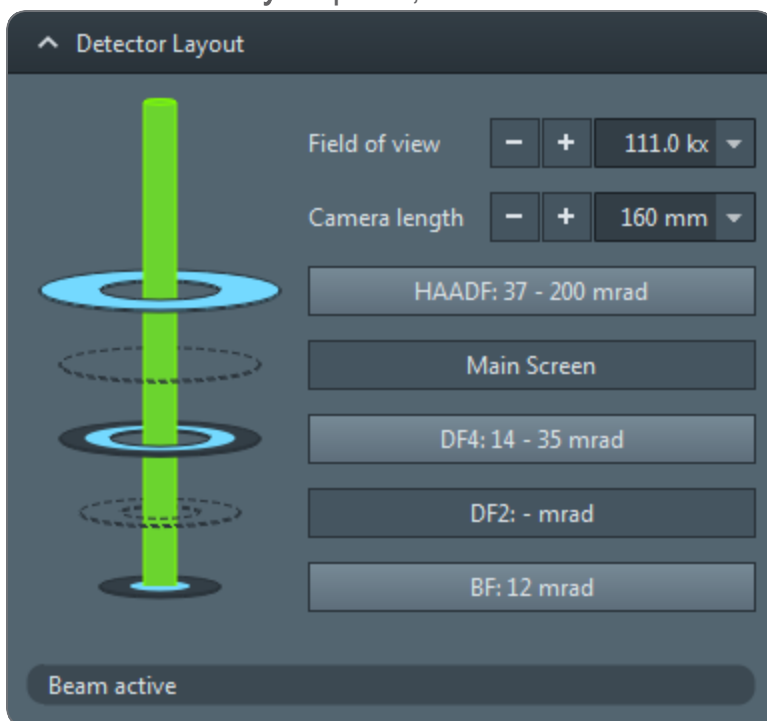


*Detectors ribbon bar for a system with HAADF and a BF/DF detector.*



*Detectors ribbon bar for a system with HAADF and a Panther STEM BF-S/DF-S detector. The DF-S option combines the DF-I (inner) and DF-O (outer) signals.*

- In the **Detector Layout** panel, activate one or more detectors or detector segments.



6. Select **STEM Imaging > Start Scanning** to start a STEM live view.

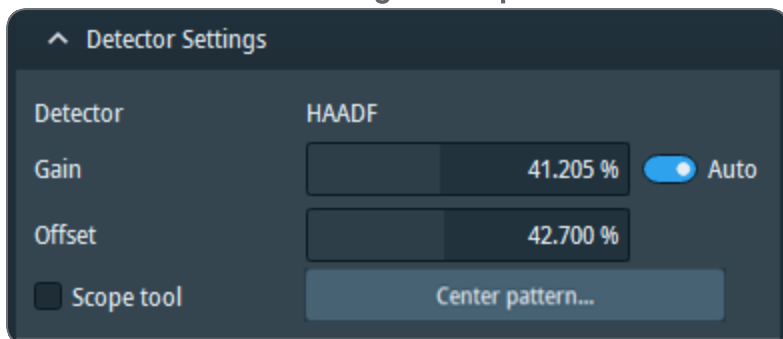


For each selected detector, MiCo opens an Image Display tab and starts scanning in View mode. While scanning, the current scan line position is indicated by an orange line that runs along the side of the image(s).

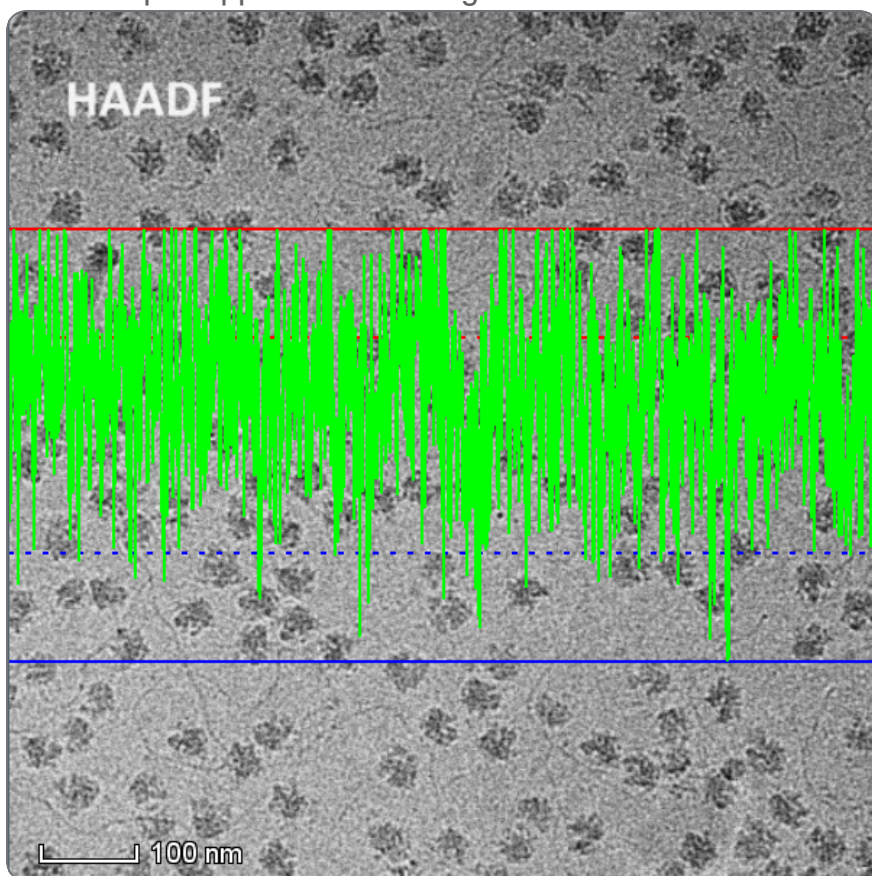
If a detector is retracted or deactivated, then the related image is removed again.

7. If necessary, use the **handpanels** to optimize the STEM live view.

8. (Optional) Maximize the dynamic range for the selected detector image.  
For the HAADF detector and for the Panther STEM BF-S and DF-S detectors it is also possible to use the Auto Gain function for an automatically optimized dynamic range. For instructions, see [How to use Auto Gain](#).
  - a. Select the **STEM image** of the detector for which Offset and/or Gain need to be adjusted.
  - b. Select the **Detector Settings > Scope tool**.



A *counts* plot appears in the image of each selected detector.



- c. Select **STEM Imaging > Beam Blank**.



- d. Adjust **Detector Settings > Offset**, so that the entire *counts* plot stays just above the lower blue line (not the dotted blue line).
- e. Select **STEM Imaging > Beam Blank** again to unblank the beam.
- f. Adjust **Detector Settings > Gain**, so that the entire *counts* plot stays just below the upper red line.
- g. Verify that the **Display Settings > Histogram > Detector Saturation Indicator** is entirely in the green, but close to the blue and red areas. This means that the Offset and Gain are set properly.



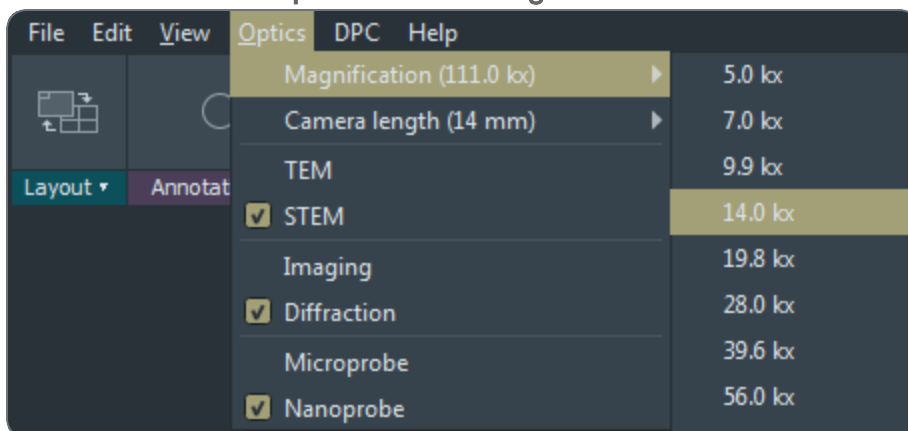
9. Select or specify the desired **Field of view** or **Magnification**.

Either:

- Use the **Handpanels > Magnification** knob.
- Select or specify the desired **Detector Layout** side panel > **Field of view**.
  - Select **[+]** or **[-]**
  - Select a **value** from the drop list.
  - Specify a **custom value**



- Select the desired **Optics** menu > **Magnification**.



In STEM Imaging mode, the Optics menu does not offer the *Field of View* parameter. Instead it offers a set of *Magnification* values, where  $\text{Magnification} = 0.1/\text{Field}$



of view with the assumption that the width of the field of view is 10 cm.

10. Select the **STEM Imaging > Scan Size**.



The preset Scan Size values include square and non-square areas of various aspect ratios and dimensions. Non-square scan sizes are not supported for DPC and iDPC acquisitions.

11. Select the **STEM Imaging > Dwell Time** or specify a custom value.

- For most acquisitions, a dwell time between 5  $\mu$ s and 40  $\mu$ s results in good quality images. A very short dwell time may cause noisy images.
- The overhead time is an estimation. For very short dwell times, the overhead time may be significantly larger than estimated.
- The frame time is calculated from the dwell time, the scan size, and the overhead time.

12. Select **STEM Imaging > Start Scanning** again to stop the STEM live view.

13. (Optional) Select **STEM Imaging > Start Acquiring** to acquire a single image.



It is also possible to start a single image acquisition without first stopping the STEM live view.

## How to Use Auto Gain

The Auto Gain function is available for the HAADF detector and for the Panther STEM BF-S and DF-S detectors. The Auto Gain function automatically adjusts the Gain value, so that the optimal dynamic range of the STEM detector is used.

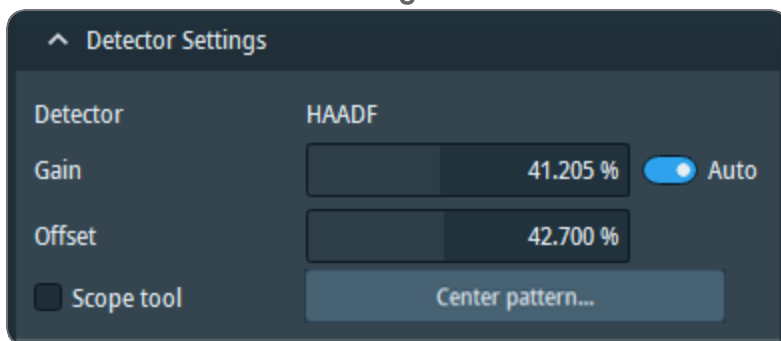
- In live viewing mode, the Auto Gain function checks the Gain value after each frame and adjusts it if necessary.
- The Gain value is not updated during the acquisition of an image or series.

The Auto Gain function also sets the Offset value to a pre-defined default value, so that the count is (nearly) zero in areas where no current is detected.

1. In the **Image Display**, select the image of the detector for which the Gain and Offset value need to be adjusted.



2. Activate the **Detector Settings > Auto** function.



To deactivate the Auto Gain function, either:

- Select **Detector Settings > Auto** again,
- Manually adjust the **Gain** or **Offset** value.

## The Live Scan Rotation Functionality

The Scan Rotation functionality adjusts the orientation of the scanning pattern.



- **Rotate tool**



Specify the Image Rotation angle by drawing an artificial horizon in the Image Display.

## Rotation angle



Either: Select **[+]** and **[-]** to fine tune the current orientation. Select the input field and enter the desired rotation angle. Select the input field and drag the mouse to the left or to the right. The *Rotation angle* value must be in the  $[-180, 180]$  degrees range. If the entered value is outside the valid range, then MiCo corrects the angle to the  $[-180, 180]$  degrees range.

- Rotate left / right 90 degrees



- Reset rotation to 0 degrees



Navigation with the joystick takes the Rotation Angle into account, so that the image on screen moves in the same direction as the joystick.

If the specimen is moved via the *TEM User Interface* > *Stage* control panel or another software application, then the Rotation Angle is *not* accounted for.

## Sample Navigation

Any camera or detector window (live-view or paused) can be used for Sample Navigation by using the Click-Center functionality. This feature makes it easier and faster to center a feature or region of interest in the Field of View by double-clicking on it.

Depending on the preceding stage moves, the accuracy of the first Click-Center move may vary. The Click-Center functionality does not apply backlash correction to absorb any mechanical tolerances of the stage. If the selected feature is not accurately centered the first time, then center it once more.

Alternatively, it is possible to use the **Arrow** keys on the keyboard to move the stage, which will move the stage approximately 80% of the Field of View in the selected direction. Using the **Shift+Arrow** keys reduces the shift to approximately 40% of the Field of View.

## Known Issues

The following are known issues that are being released with this version of the software.

ID	Description	Workaround
1	The SmartCam in MiCo may give an error when FlucamViewer is running.	It is recommended to close FlucamViewer before starting SmartCam in MiCo.
2	The startup time of MiCo may be longer than usual for some systems during the first time after TEM server start, taking from about 10 to 60 seconds.	None available for now.
3	Grey progress images are shown for few alignments using a Falcon (3) camera.	None available for now.
4	Dose measurement for Falcon camera's is too sensitive and may prevent live-viewing of the camera.	Use other applications, like EPU.
5	Running SmartCam in the application blocks other software from using the same (like automated alignments) and it may cause alignments to fail quickly.	Close SmartCam manually.  Note that MiCo does not need to remain open when working in other applications.
6	Gun Shift alignment requires an inserted holder for X-Ray safety on systems without Autoloader. However, it requires vacuum/no sample in beam.	Holder can be navigated to hole or holder can be slightly retracted from microscope.
7	Stage recovery error is thrown with recovery of stage.	Ignore the error, the stage will be enabled.