

NanoDrop One Release Notes

About This Document	11
NanoDrop One PC Software v2.11.0	11
<i>Release Date: October 2023</i>	<i>11</i>
<i>Supported Operating Systems.....</i>	<i>11</i>
<i>New Features.....</i>	<i>11</i>
Increase to 16 unique colors when overlaying spectra in the software.....	11
Update extinction coefficient of Alexa Fluor 488, Alexa Fluor 546, Alexa Fluor 555, Alexa Fluor 594, and Alexa Fluor 647 based on new published sources.....	11
Add an option to save the spectrum as a .png file in the PC software.....	11
<i>Resolved Issues.....</i>	<i>11</i>
This version updates the correct NanoDrop One.xml in the NanoDrop One PC installation folder.....	11
This version resolves a crash issue when adding 13 U characters in the Oligo RNA application.....	11
<i>Known Issues.....</i>	<i>11</i>
NanoDrop One Local Control Software v2.11.0.....	11
<i>Release Date: October 2023</i>	<i>11</i>
<i>New Features.....</i>	<i>11</i>
Update extinction coefficient of Alexa Fluor 488, Alexa Fluor 546, Alexa Fluor 555, Alexa Fluor 594, and Alexa Fluor 647 based on new published sources.....	11
<i>Resolved Issues.....</i>	<i>11</i>
This version resolves a crash issue when adding 13 U characters in the Oligo RNA application.....	11
This version improves an issue seen with the SD and adds a dialog prompt to change the SD card if storage is full during the initialization check.....	11
This version updates Proteins & Labels baseline correction from 750nm to 850nm.....	11
<i>Known Issues.....</i>	<i>11</i>
NanoDrop One PC Software v2.10.0	11
<i>Release Date: March 2023</i>	<i>11</i>
<i>Supported Operating Systems.....</i>	<i>11</i>
<i>New Features.....</i>	<i>12</i>
Add absorbance and concentration data for standard curve samples in exported files	12
Add one policy 'Ability to back up and restore NanoDrop One measurement database' in Security Administration.....	12
Add the database backup and restore functionality of Settings database tab in the PC software.	12
Add the serial number of the instrument to the printable report of the Pedestal Image Check	12
<i>Resolved Issues.....</i>	<i>12</i>
This version improves the consistency of decimal places displayed in Protein A280	12
This version resolves an issue importing Kinetics experiments from the Local control when the measurement run time is over one hour.....	12
This version resolves an issue with Custom Methods where concentration column shows "out of range", not the correct value, in exported files.....	12
This version improves the printing of log oligo sequences	12
<i>Known Issues.....</i>	<i>12</i>

NanoDrop One Local Control Software v2.10.0.....12

Release Date: March 2023 12

New Features..... 12

 NanoDrop One with Security installation package adds the Security OQ documentation..... 12

Resolved Issues..... 12

 This version resolves a limitation in specifying a network path for networked data export using SMB2.0 12

 This version resolves an issue where data export to a shared network location from the Local Control could not be uploaded to the PC control software. 12

 This version resolves a formatting issue in which is shifted over one column in .tsv files exported from Local Control software 12

 This version resolves an issue with displayed wavelengths when exporting UV-Vis application data in .tsv file format. 12

 This version increases the number of absorbance value decimal places displayed in the UV-Vis application to 3 to keep consistent with the PC software 12

Known Issues..... 12

NanoDrop One PC Software v2.9.012

Release Date: September 2022..... 12

Supported Operating Systems..... 12

New Features..... 12

 Export data to add new Local control spreadsheet measurement data .csv file format 12

Resolved Issues..... 13

 This version resolves an issue with the auto export audit trail data creation and deletion records in a temporary folder 13

 This version resolves an issue with two samples displayed incompletely from printed reports with more than one page..... 13

Known Issues..... 13

NanoDrop One Local Control Software v2.9.0.....13

Release Date: Aug 2022..... 13

New Features..... 13

 The firmware package has been separated from NanoDrop One Local control software installation..... 13

Resolved Issues..... 13

 This version resolves an issue where two user accounts cannot simultaneously login to the instrument Cloud 13

 This version resolves an issue where the instrument cannot remember Cloud user account and password . 13

 This version resolves an issue where the Cloud user account cannot re-establish the connection with Thermo Cloud once the user disconnects 13

Known Issues..... 13

NanoDrop One PC Software v2.8.013

Release Date: June 2022 13

Supported Operating Systems..... 13

New Features..... 13

 Export data to add .csv file format..... 13

 Dye Editor wavelength update to 350 – 840 nm 13

 Add new access control: Ability to edit Auto Export..... 13

Add new system policies: Prevent export of .ned(experiment files) from the PC control software.....	13
Add new system policies: Prevent import of .ned(experiment files) into the PC control software	13
Intensity Check update fail information message.....	13
<i>Resolved Issues</i>	13
This version resolves an issue with Security secure data directory permission	13
This version resolves an issue with Protein A205 Extinction Coefficient incorrect display	13
<i>Known Issues</i>	13
NanoDrop One Local Control Software v2.8.0	14
<i>Release Date: June 2022</i>	14
<i>New Features</i>	14
LC software supports DYMO 550 Turbo printer.....	14
<i>Resolved Issues</i>	14
This version resolves an issue with Protein A205 Extinction Coefficient incorrect display	14
This version resolves an issue where Protein Bradford reuse standard curve did not display the standard curve data when imported into PC software	14
This version resolves an issue that .tsv files did not include the standard curve absorbance and concentration data when a standard curve was reused	14
<i>Known Issues</i>	14
Instrument software did not indicate the printer error.....	14
NanoDrop One PC Software v2.7.0	14
<i>Release Date: Mar 2022</i>	14
<i>Supported Operating Systems</i>	14
<i>New Features</i>	14
Protein Bradford application	14
Protein Lowry application	14
Protein Pierce 660	14
Add the instrument serial number of printed report Diagnostics application.....	14
Ability to turn on/off auto export feature	14
<i>Resolved Issues</i>	14
Intensity Check PC software and LC software results difference issue.....	14
This version resolved an issue where the .sql file from the LC software would display an incorrect result when imported to the PC software.....	14
<i>Known Issues</i>	14
NanoDrop One Local Control Software v2.7.0	14
<i>Release Date: Mar 2022</i>	14
<i>New Features</i>	15
LC software supports lock out feature.....	15
Protein Lowry supports standard curve import feature	15
Protein Pierce 660 supports standard curve import feature.....	15
<i>Resolved Issues</i>	15
<i>Known Issues</i>	15
NanoDrop One PC Software v2.6.0	15
<i>Release Date: Dec 2021</i>	15

Supported Operating Systems.....	15
New Features.....	15
None	15
Resolved Issues.....	15
Dutch and French system	15
Known Issues.....	15
NanoDrop One Local Control Software v2.6.0.....	15
Release Date: Dec 2021.....	15
New Features.....	15
LC software upgrades protocol from SMBv1.0 to SMBv2.0.....	15
Resolved Issues.....	15
Known Issues.....	15
NanoDrop One PC Software v2.5.0	15
Release Date: Oct 2021.....	15
Supported Operating Systems.....	15
New Features.....	15
OD600 application	15
Resolved Issues.....	16
Standard Deviation in PV	16
Print report.....	16
German system	16
Security lock inactive system policy	16
Known Issues.....	16
NanoDrop One Local Control Software v2.5.0.....	16
Release Date: Oct 2021.....	16
New Features.....	16
NA.....	16
Resolved Issues.....	16
Standard Deviation in PV	16
NanoDrop One PC Software v2.4.0	16
Release Date: July 2021.....	16
Supported Operating Systems.....	16
New Features.....	16
Kinetics application.....	16
Resolved Issues.....	16
Known Issues.....	16
NanoDrop One Local Control Software v2.4.0.....	17
Release Date: July 2021.....	17
New Features.....	17

Protein BCA supports standard curve import feature.....	17
Protein Bradford supports standard curve import feature	17
Remember default password when test connection for selected network location	17
<i>Resolved Issues</i>	17
NanoDrop One PC Software v2.3.0	17
<i>Release Date: Mar 2021</i>	17
<i>Supported Operating Systems</i>	17
<i>New Features</i>	17
Protein & Labels application.....	17
Protein BCA application	17
Print reports added the units with the concentration column	17
<i>Resolved Issues</i>	17
Negative concentration display	17
Export files.....	17
Analytical Wavelength in UV-Vis	17
<i>Known Issues</i>	17
NanoDrop One Local Control Software v2.3.0.....	17
<i>Release Date: Mar 2021</i>	17
<i>New Features</i>	18
Protein & Labels added A260/A280 result of sample details screen	18
<i>Resolved Issues</i>	18
Analytical Wavelength in UV-Vis	18
NanoDrop One PC Software v2.2.0	18
<i>Release Date: Dec 2020</i>	18
<i>Supported Operating Systems</i>	18
<i>New Features</i>	18
RNA/DNA Deconvolution	18
Microarray application	18
Released application support of exporting .sql instrument files from the PC software	18
Add policy in Security Administration	18
<i>Resolved Issues</i>	18
Print reports	18
Export files.....	18
<i>Known Issues</i>	18
NanoDrop One Local Control Software v2.2.0.....	18
<i>Release Date: Dec 2020</i>	18
<i>New Features</i>	18
RNA/DNA Deconvolution	18
<i>Resolved Issues</i>	18
History icon.....	18
NanoDrop One PC Software v2.1.0	19

<i>Release Date: Oct 2020</i>	19
<i>Supported Operating Systems</i>	19
<i>New Features</i>	19
RNA/DNA Deconvolution	19
Oligo DNA/Oligo RNA applications	19
Protein A205 application	19
Show monitored wavelength	19
<i>Resolved Issues</i>	19
Export files.....	19
PC viewer Custom method.....	19
<i>Known Issues</i>	19
NanoDrop One Local Control Software v2.1.0	19
<i>Release Date: Oct 2020</i>	19
<i>New Features</i>	19
RNA/DNA Deconvolution	19
Dye Editor wavelength	19
<i>Resolved Issues</i>	20
Launcher screen.....	20
Pedestal image sensor is not responding	20
NanoDrop One PC Software v2.0.1	20
<i>Release Date: Sep 2020</i>	20
<i>Supported Operating Systems</i>	20
<i>New Features</i>	20
<i>Resolved Issues</i>	20
Intensity Check.....	20
<i>Known Issues</i>	20
NanoDrop One PC Software v2.0.0	20
<i>Release Date: July 2020</i>	20
<i>Supported Operating Systems</i>	20
<i>New Features</i>	20
<i>Nucleic Acid</i>	20
dsDNA application	20
ssDNA application	20
RNA application.....	21
Custom Factor application.....	21
<i>Protein</i>	21
Protein Editor.....	21
Protein A280.....	21
<i>Custom</i>	21
Custom Methods	21
UV-Vis	21
<i>Diagnostics</i>	21
Performance Verification	21

Intensity Check.....	21
Pedestal Image Check	22
<i>IQ/OQ</i>	22
<i>Security with 21CFR part 11 compliance Support</i>	22
General	22
History	22
Settings.....	22
Help/Documentation.....	22
<i>Resolved Issues</i>	22
<i>Known Issues</i>	22
NanoDrop One Local Control Software v2.0.0	22
<i>Release Date: July 2020</i>	22
<i>New Features</i>	22
<i>Resolved Issues</i>	22
NanoDrop One Local Control Software v1.4.5	22
<i>Release Date: February 2020</i>	22
<i>Resolved Issues</i>	22
Fixed issue with running manufacturing and service utilities	22
NanoDrop One Local Control Software v1.4.4	23
<i>Release Date: November 2019</i>	23
<i>Resolved Issues</i>	23
California SB-327 Compliance	23
NanoDrop One Local Control Software v1.4.3	23
<i>Release Date: October 2019</i>	23
<i>Resolved Issues</i>	23
Customer Special Software.....	23
NanoDrop One Local Control Software v1.4.2	23
<i>Release Date: November 2017</i>	23
<i>Resolved Issues</i>	23
Blue Screen Lockup	23
NanoDrop One Local Control Software v1.4	23
<i>Release Date: May 2017</i>	23
<i>New Features</i>	23
Wi-Fi and Ethernet Networking and Export:.....	23
Cloud Connect:.....	23
Wi-Fi/USB Printing.....	24
16GB Data Storage	24
Export Logs	24
“All” Option in Search Time Range.....	24
Edit Experiment Name	24

Export .sql files for individual experiments	24
<i>Resolved Issues</i>	24
Custom Method Issues.....	24
Protein Bradford	24
<i>Known Issues</i>	24
NanoDrop One PC Viewer Software v1.4	25
Release Date: May 2017	25
Supported Operating Systems.....	25
New Features.....	25
Edit Experiment Name	25
Resolved Issues.....	25
Printing Intensity Check.....	25
Exporting to .tsv Format	25
Known Issues.....	25
NanoDrop One Local Control Software v1.3 Hot Fix 1	25
Release Date: November 2016.....	25
Supported Operating Systems.....	25
New Features.....	25
Export for Cloud.....	25
Resolved Issues.....	25
Known Issues.....	25
NanoDrop One Local Control Software v1.3.....	26
Release Date: September 2016.....	26
Supported Operating Systems.....	26
New Features.....	26
Translations: Software and Help	26
Resolved Issues.....	26
Blank PC Name	26
Kinetics Interval Error Message	26
GuSCN or GUHCI contamination:	26
Execution of Custom Methods:	26
Known Issues.....	26
NanoDrop One PC Viewer Software v1.3	26
Release Date: September 2016.....	26
Supported Operating Systems.....	26
New Features.....	27
Translations: Software and Help	27
Resolved Issues.....	27
Display Issues	27
Y-axis Label for Blank	27

<i>Known Issues</i>	27
NanoDrop One Local Control Software v1.2	27
<i>Release Date: February 2016</i>	27
<i>Supported Operating Systems</i>	27
<i>New Features</i>	27
UV-Vis Application.....	27
Kinetics Application.....	27
Help System Translation.....	28
IQ/OQ.....	28
Additional Custom Methods.....	28
Acclaro Blanking.....	28
Wi-Fi Data Transfer.....	28
<i>Resolved Issues</i>	28
<i>Known Issues</i>	28
Data Size Limit.....	28
NanoDrop One PC Viewer Software v1.2	28
<i>Release Date: February 2016</i>	28
<i>Supported Operating Systems</i>	28
<i>New Features</i>	28
Wi-Fi Data Transfer.....	28
Additional Custom Method.....	28
Export .sql File Format.....	29
<i>Resolved Issues</i>	29
<i>Known Issues</i>	29
NanoDrop One Local Control Software v1.1	29
<i>Release Date: December 2015</i>	29
<i>Supported Operating Systems</i>	29
<i>New Features</i>	29
Direct Connect (PC Tethering).....	29
Pedestal Image Check.....	29
Software Translation.....	29
<i>Resolved Issues</i>	29
<i>Known Issues</i>	29
NanoDrop One PC Viewer Software v1.1	29
<i>Release Date: December 2015</i>	29
<i>Supported Operating Systems</i>	29
<i>New Features</i>	30
Direct Connect (PC Tethering).....	30
<i>Resolved Issues</i>	30
<i>Known Issues</i>	30

NanoDrop One Local Control Software v1.0 **30**
 Release Date: November 2015..... 30
 Supported Operating Systems..... 30
 New Features..... 30
 Resolved Issues..... 30
 Known Issues..... 30

NanoDrop One PC Viewer Software v1.0 **30**
 Release Date: November 2015..... 30
 Supported Operating Systems..... 30
 New Features..... 30
 Resolved Issues..... 30
 Known Issues..... 30

How to Contact Us **30**

Trademarks **31**

About This Document

This document contains a revision history of Thermo Scientific NanoDrop™ One/One^C PC Software and Thermo Scientific NanoDrop™ One/ One^C Local Control Software which includes new features, resolved and known issues.

NanoDrop One PC Software v2.11.0

Release Date: October 2023

Supported Operating Systems

Windows 10 64-bit
Windows 11 64-bit

New Features

Increase to 16 unique colors when overlaying spectra in the software.
Update extinction coefficient of Alexa Fluor 488, Alexa Fluor 546, Alexa Fluor 555, Alexa Fluor 594, and Alexa Fluor 647 based on new published sources.
Add an option to save the spectrum as a .png file in the PC software.

Resolved Issues

This version updates the correct NanoDrop One.xml in the NanoDrop One PC installation folder.
This version resolves a crash issue when adding 13 U characters in the Oligo RNA application.

Known Issues

None

NanoDrop One Local Control Software v2.11.0

Release Date: October 2023

New Features

Update extinction coefficient of Alexa Fluor 488, Alexa Fluor 546, Alexa Fluor 555, Alexa Fluor 594, and Alexa Fluor 647 based on new published sources.

Resolved Issues

This version resolves a crash issue when adding 13 U characters in the Oligo RNA application.
This version improves an issue seen with the SD and adds a dialog prompt to change the SD card if storage is full during the initialization check.
This version updates Proteins & Labels baseline correction from 750nm to 850nm.

Known Issues

None

NanoDrop One PC Software v2.10.0

Release Date: March 2023

Supported Operating Systems

Windows 10 64-bit
Windows 11 64-bit

New Features

- Add absorbance and concentration data for standard curve samples in exported files
- Add one policy 'Ability to back up and restore NanoDrop One measurement database' in Security Administration
- Add the database backup and restore functionality of Settings database tab in the PC software.
- Add the serial number of the instrument to the printable report of the Pedestal Image Check

Resolved Issues

- This version improves the consistency of decimal places displayed in Protein A280
- This version resolves an issue importing Kinetics experiments from the Local control when the measurement run time is over one hour
- This version resolves an issue with Custom Methods where concentration column shows "out of range", not the correct value, in exported files
- This version improves the printing of log oligo sequences

Known Issues

None

NanoDrop One Local Control Software v2.10.0

Release Date: March 2023

New Features

- NanoDrop One with Security installation package adds the Security OQ documentation

Resolved Issues

- This version resolves a limitation in specifying a network path for networked data export using SMB2.0
- This version resolves an issue where data export to a shared network location from the Local Control could not be uploaded to the PC control software.
- This version resolves a formatting issue in which is shifted over one column in .tsv files exported from Local Control software
- This version resolves an issue with displayed wavelengths when exporting UV-Vis application data in .tsv file format.
- This version increases the number of absorbance value decimal places displayed in the UV-Vis application to 3 to keep consistent with the PC software

Known Issues

None

NanoDrop One PC Software v2.9.0

Release Date: September 2022

Supported Operating Systems

- Windows 10 64-bit
- Windows 11 64-bit

New Features

- Export data to add new Local control spreadsheet measurement data .csv file format

Resolved Issues

This version resolves an issue with the auto export audit trail data creation and deletion records in a temporary folder

This version resolves an issue with two samples displayed incompletely from printed reports with more than one page

Known Issues

None

NanoDrop One Local Control Software v2.9.0

Release Date: Aug 2022

New Features

The firmware package has been separated from NanoDrop One Local control software installation

Resolved Issues

This version resolves an issue where two user accounts cannot simultaneously login to the instrument Cloud

This version resolves an issue where the instrument cannot remember Cloud user account and password

This version resolves an issue where the Cloud user account cannot re-establish the connection with Thermo Cloud once the user disconnects

Known Issues

None

NanoDrop One PC Software v2.8.0

Release Date: June 2022

Supported Operating Systems

Windows 10 64-bit

Windows 11 64-bit

New Features

Export data to add .csv file format

Dye Editor wavelength update to 350 – 840 nm

Add new access control: Ability to edit Auto Export

Add new system policies: Prevent export of .ned(experiment files) from the PC control software

Add new system policies: Prevent import of .ned(experiment files) into the PC control software

Intensity Check update fail information message

Resolved Issues

This version resolves an issue with Security secure data directory permission

This version resolves an issue with Protein A205 Extinction Coefficient incorrect display

Known Issues

None

NanoDrop One Local Control Software v2.8.0

Release Date: June 2022

New Features

LC software supports DYMO 550 Turbo printer

Resolved Issues

This version resolves an issue with Protein A205 Extinction Coefficient incorrect display

This version resolves an issue where Protein Bradford reuse standard curve did not display the standard curve data when imported into PC software

This version resolves an issue that .tsv files did not include the standard curve absorbance and concentration data when a standard curve was reused

Known Issues

Instrument software did not indicate the printer error

This version does not include software alerts for printer errors

NanoDrop One PC Software v2.7.0

Release Date: Mar 2022

Supported Operating Systems

Windows 10 64-bit

Windows 11 64-bit

New Features

Protein Bradford application

Protein Lowry application

Protein Pierce 660

A standard curve is required for colorimetric protein analysis and support to import a standard curve. Run blank and measurement on pedestal and cuvette. Protein Bradford, Protein Lowry and Protein Pierce 660 experiment data can be viewed in the History page. Ability to customize which columns are displayed in History and in the measurement acquisition screen. Export data in multiple file formats (.tsv and .ned). Support import of .ned/.sql files and PC Viewer database files into History.

Add the instrument serial number of printed report Diagnostics application

Ability to turn on/off auto export feature

Resolved Issues

Intensity Check PC software and LC software results difference issue

This version resolved an issue where the .sql file from the LC software would display an incorrect result when imported to the PC software.

Known Issues

None

NanoDrop One Local Control Software v2.7.0

Release Date: Mar 2022

New Features

- LC software supports lock out feature
- Protein Lowry supports standard curve import feature
- Protein Pierce 660 supports standard curve import feature

Resolved Issues

None

Known Issues

None

NanoDrop One PC Software v2.6.0

Release Date: Dec 2021

Supported Operating Systems

- Windows 10 64-bit
- Windows 11 64-bit

New Features

None

Resolved Issues

Dutch and French system

This version resolved an issue where the pathlength choice is incorrect on Dutch and French operating systems

Known Issues

None

NanoDrop One Local Control Software v2.6.0

Release Date: Dec 2021

New Features

- LC software upgrades protocol from SMBv1.0 to SMBv2.0

Resolved Issues

None

Known Issues

None

NanoDrop One PC Software v2.5.0

Release Date: Oct 2021

Supported Operating Systems

- Windows 10 64-bit

New Features

OD600 application

Run blank and measurement to monitor the growth rate of bacterial or other microbial cell cultures in growth media at 600nm. Ability to customize which columns are displayed in History and in the measurement

acquisition screen. Export data in multiple file formats (. tsv and .ned). Support import of . ned/.sql files and PC Viewer database files into History.

Resolved Issues

Standard Deviation in PV

This version resolved an issue where the Standard Deviation calculation at 1.0 mm was not functioning properly in the Performance Verification application

Print report

This version resolved issues printing truncated sample names and supporting the printing of pathlength length data.

German system

This version resolved an issue where some Custom Methods could not run on German operating systems.

Security lock inactive system policy

This version resolved a bug in the “Automatically lock inactive system after(minutes)” policy in Security Administration

Known Issues

None

NanoDrop One Local Control Software v2.5.0

Release Date: Oct 2021

New Features

NA

Resolved Issues

Standard Deviation in PV

This version resolved an issue where the Standard Deviation calculation at 1.0 mm was not functioning properly in the Performance Verification application

NanoDrop One PC Software v2.4.0

Release Date: July 2021

Supported Operating Systems

Windows 10 64-bit

New Features

Kinetics application

Run blank and measurement to make time-based experiments in cuvettes, up to 3 wavelengths between 190 nm and 850 nm can be designated for continuous absorbance monitoring at user-defined intervals in up to 5 stages. Ability to customize which columns are displayed in History and in the measurement acquisition screen. Export data in multiple file formats (. tsv and .ned). Support import of . ned/.sql files and PC Viewer database files into History.

Resolved Issues

None

Known Issues

None

NanoDrop One Local Control Software v2.4.0

Release Date: July 2021

New Features

- Protein BCA supports standard curve import feature**
- Protein Bradford supports standard curve import feature**
- Remember default password when test connection for selected network location**

Resolved Issues

None

NanoDrop One PC Software v2.3.0

Release Date: Mar 2021

Supported Operating Systems

Windows 10 64-bit

New Features

Protein & Labels application

Run blank and measurement on pedestal and cuvette, supports Analysis Correction. Protein & Labels experiment data can be viewed in the History page. Ability to customize which columns are displayed in History and in the measurement acquisition screen. Export data in multiple file formats (.tsv and .ned). Support import of .ned/.sql files and PC Viewer database files into History.

Protein BCA application

A standard curve is required for colorimetric protein analysis and support to import a standard curve. Run blank and measurement on pedestal and cuvette. Protein BCA experiment data can be viewed in the History page. Ability to customize which columns are displayed in History and in the measurement acquisition screen. Export data in multiple file formats (.tsv and .ned). Support import of .ned/.sql files and PC Viewer database files into History.

Print reports added the units with the concentration column

Resolved Issues

Negative concentration display

Released applications display the negative concentration value instead of showing 'NA'.

Export files

Protein A280 concentration results exported in .tsv files supporting three decimal points.

Analytical Wavelength in UV-Vis

This version resolved an issue where the analytical wavelength was not worked in UV-Vis application

Known Issues

None

NanoDrop One Local Control Software v2.3.0

Release Date: Mar 2021

New Features

Protein & Labels added A260/A280 result of sample details screen

Resolved Issues

Analytical Wavelength in UV-Vis

This version resolved an issue where the analytical wavelength was not worked in UV-Vis application.

NanoDrop One PC Software v2.2.0

Release Date: Dec 2020

Supported Operating Systems

Windows 10 64-bit

New Features

RNA/DNA Deconvolution

Improved algorithms for differentiation of mammalian DNA from RNA and vice versa in dsDNA/RNA applications

Microarray application

Run, blank, and measurement on pedestal and cuvette, supports Analysis Correction. Microarray experiment data can be viewed in the History page. Ability to customize which columns are displayed in History and in the measurement acquisition screen. Export data in multiple file formats (.tsv and .ned). Support import of .ned/.sql files and PC Viewer database files into History.

Released application support of exporting .sql instrument files from the PC software

Add policy in Security Administration

Prevent export of .sql (instrument files) from the PC software

Resolved Issues

Print reports

Signature Log did not print out with the Intensity Check and Performance Verification

Export files

Cuvette Intensity Check results exported in .tsv format misidentified the sample location

Known Issues

None

NanoDrop One Local Control Software v2.2.0

Release Date: Dec 2020

New Features

RNA/DNA Deconvolution

Improved algorithms for differentiation of mammalian DNA from RNA and vice versa in dsDNA/RNA applications

Resolved Issues

History icon

The word "Local" in certain languages would cause the History icon to disappear

NanoDrop One PC Software v2.1.0

Release Date: Oct 2020

Supported Operating Systems

Windows 10 64-bit

New Features

RNA/DNA Deconvolution

New analyses that differentiate mammalian DNA from RNA and vice versa in dsDNA/RNA applications

Oligo DNA/Oligo RNA applications

Run blank and measurement on pedestal and cuvette, supports baseline correction. Oligo DNA/Oligo RNA experiment data can be viewed in the History page. Ability to customize what columns are displayed in History and in the measurement acquisition screen. Export data in multiple file formats (.tsv and .ned). Support import of .ned/.sql files and PC viewer database files into History.

Protein A205 application

Calculate peptide concentration based on user defined factor, run blank and measurement on pedestal and cuvette, supports baseline correction. Protein A205 experiment data can be viewed in the History page. Ability to customize what columns are displayed in History and in the measurement acquisition screen. Acclaro bubble detection includes a check for blank and measurement integrity. Export data in multiple file formats (.tsv and .ned). Support import of .ned/.sql file and PC viewer database files into History.

Show monitored wavelength

Show and Print extra monitored wavelength in Nucleic Acids and Protein applications.

Resolved Issues

Export files

Fixed a formatting issue when exporting .tsv data on French and Danish operation systems.

PC viewer Custom method

Fixed Custom Method from PC viewer cannot import to NanoDrop One PC software

Known Issues

None

NanoDrop One Local Control Software v2.1.0

Release Date: Oct 2020

New Features

RNA/DNA Deconvolution

New analyses that differentiate mammalian DNA from RNA and vice versa in dsDNA/RNA applications

Dye Editor wavelength

Wider Dye Editor wavelength update to 350-850nm

Resolved Issues**Launcher screen**

The version resolves an issue related to the software launch interface.

Pedestal image sensor is not responding

Added message prompting users to restart the instrument when pedestal image sensor is not responding.

NanoDrop One PC Software v2.0.1

Release Date: Sep 2020

Supported Operating Systems

Windows 10 64-bit

New Features

NA

Resolved Issues**Intensity Check**

This version resolves an issue in how the experiment timestamp was displayed in History.

Known Issues

None

NanoDrop One PC Software v2.0.0

Release Date: July 2020

Supported Operating Systems

Windows 10 64-bit

New Features

Initial Release of software from which users can control the instrument

- NanoDrop One PC Viewer Software is retired.
- Integrate features required by 21 CFR Part 11 regulations
- Chinese language support
- All new Features list as below:

Nucleic Acid**dsDNA application**

Run blank and measurement on pedestal and cuvette, supports baseline correction. dsDNA experiment data can be viewed in the History detail page. Ability to customize what columns are displayed in History and in the measurement acquisition screen. Acclaro contaminant analysis/ bubble detection includes a check for blank and measurement. Export data in multiple file formats (.tsv and .ned). Support to import .ned/.sql file and PC viewer database files in History.

ssDNA application

Run blank and measurement on pedestal and cuvette, supports baseline correction. ssDNA experiment data can be viewed in the History detail page. Ability to customize what columns are displayed in History and in the measurement acquisition screen. Acclaro includes a bubble detection check for blank and measurement. Export data in multiple file formats (.tsv and .ned). Support to import .ned/.sql file and PC viewer database files in History.

RNA application

Run blank and measurement on pedestal and cuvette, supports baseline correction. RNA experiment data can be viewed in the History detail page. Ability to customize what columns are displayed in History and in the measurement acquisition screen. Acclaro contaminant analysis/ bubble detection includes a check for blank and measurement. Export data in multiple file formats (.tsv and .ned). Support to import .ned/.sql file and PC viewer database files in History.

Custom Factor application

Calculate nucleic acid concentration based on user defined factor. Run blank and measurement on pedestal and cuvette, supports baseline correction. Custom factor experiment data can be viewed in the History detail page. Ability to customize what columns are displayed in History and in the measurement acquisition screen. Acclaro includes a bubble detection check for blank and measurement. Export data in multiple file formats (.tsv and .ned). Support to import .ned/ .sql file and PC viewer database files in History.

Protein**Protein Editor**

User-entered protein extinction coefficients in the Protein Editor can be selected from the Protein A280 drop-down menu

Protein A280

Calculate protein concentration based on user defined factor, run blank and measurement on pedestal and cuvette, supports baseline correction. Protein A280 experiment data can be viewed in the History detail page. Ability to customize what columns are displayed in History and in the measurement acquisition screen. Acclaro contaminant analysis/ bubble detection includes a check for blank and measurement. Export data in multiple file formats (.tsv and .ned). Support to import .ned/ .sql file and PC viewer database files in History.

Custom**Custom Methods**

Formula and standard curve custom methods can be created in the Custom method application. Run blank and measurement on pedestal and cuvette, supports analysis wavelength and quantification step. Export data in multiple file formats (.tsv and .ned). Support to import .ned/ .sql file and NanoDrop One PC viewer database file in History.

UV-Vis

The UV-Vis application allows the NanoDrop One/NanoDrop One PC to function as a conventional spectrophotometer. Wavelengths can be entered for absorbance monitoring and inclusion in the results. Sample absorbance is displayed on the screen from 190nm to 840nm.

Diagnostics**Performance Verification**

Users enter the target absorbance value on the PV-1 ampoule in the Performance Verification setup page, then blank and measure aliquots of the standard. Pass or fail instructions show at the top of the screen. Experiment data can be viewed in the history detail page. Export data in multiple file formats (.tsv and .ned). Support to import .ned/ .sql file and PC viewer database files in History.

Intensity Check

Run an Intensity Check through the pedestal or cuvette light paths. Pass or fail instructions show at the top. Experiment data can be viewed in the History detail page. Export data in multiple file formats (.tsv and .ned). Support to import .ned/ .sql file and PC viewer database files in History.

Pedestal Image Check

Run a Pedestal Image Check. Supports Image optimization. Pass or fail instructions show at the top. Experiment data can be viewed in the History detail page. Export data in multiple file formats (. tsv and. ned). Support to import .ned/ .sql file and PC viewer database files in History

IQ/OQ

Qualification procedures are now available for the NanoDrop One PC Control software

Security with 21CFR part 11 compliance Support

NanoDrop One Security software, together with Thermo Security Administration software, lets the system administrator set system policies and control user access to NanoDrop One software feature. Supports access control, system policies, and signature meanings

General

History

Data can be viewed and updated from the History page

Settings

Security server location

Automatic .ned file export to specific location (Support local and network location)

Protein Editor

Language settings, support Chinese language

Help/Documentation

Resolved Issues

Initial release, there are no resolved Issues

Known Issues

None

NanoDrop One Local Control Software v2.0.0

Release Date: July 2020

New Features

NanoDrop One instrument directly connects to the NanoDrop One PC Control software with an Ethernet cable. Instrument can be controlled from the PC and all data are stored on the computer.

Change "Data Viewer" to "History"

Resolved Issues

None

NanoDrop One Local Control Software v1.4.5

Release Date: February 2020

Resolved Issues

Fixed issue with running manufacturing and service utilities

NanoDrop One Local Control Software is updated to be compatible with manufacturing and service utilities.

NanoDrop One Local Control Software v1.4.4**Release Date: November 2019****Resolved Issues****California SB-327 Compliance**

NanoDrop One Local Control Software is updated to comply with California SB-327: unique password on each instrument to turn off kiosk mode.

NanoDrop One Local Control Software v1.4.3**Release Date: October 2019****Resolved Issues****Customer Special Software**

NanoDrop One Local Control Software and PC Data Viewer are updated to support two customer special orders.

NanoDrop One Local Control Software v1.4.2**Release Date: November 2017****Resolved Issues****Blue Screen Lockup**

Some NanoDrop One/NanoDrop One^C instruments experienced a lockup where the instrument software did not get past the blue droplet screen. This version fixes this issue. It is recommended that all users upgrade to this version of the local control software.

NanoDrop One Local Control Software v1.4**Release Date: May 2017****New Features****Wi-Fi and Ethernet Networking and Export:**

Users can connect the NanoDrop One/NanoDrop One^C to a Wi-Fi network by selecting **System → Networking → Wi-Fi**.

Users can select **System → Networking → Ethernet** to connect to a wired network or directly to a PC. A list of network locations can be pre-defined under the **System → Export** tab. This list is presented when the user selects Export in the End Experiment dialog or in the Data Viewer.

Cloud Connect:

The NanoDrop One/NanoDrop One^C can now connect to the Thermo Fisher Cloud and export data to the user's cloud account. The exported data may be viewed or manipulated using the NanoDrop Cloud application.

Users can setup their account details on the **Settings → Network** tab. Once an account is setup, a new option “Thermo Fisher Cloud” is added to the Export dropdown. When “Thermo Fisher Cloud” is selected as the export location, a user account list is presented. Data can be exported by selecting the appropriate user account name.

Wi-Fi/USB Printing

All NanoDrop One/NanoDrop One^C instruments with v1.4 of Local Control Software can print reports to a standard USB or Wi-Fi connected printer. To configure a printer for the first time, select the Print option from the main menu when in an experiment and follow the instructions.

If you have upgraded the NanoDrop One/NanoDrop One^C Local Control Software from an earlier version to v1.4, your NanoDrop One/NanoDrop One^C instrument will need to be connected to the Internet over Wi-Fi or Ethernet. Once connected, select the Print option from the main menu of an experiment and follow the instructions.

When a Dymo printer is connected, the instrument will honor the Dymo printer exclusively. In order to use regular printer, the Dymo printer needs to be disconnected.

Note:

A variety of printer manufacturers and models are supported. Many printers from the same printer manufacturer behave the same way. If your printer model is not found in the list, try selecting the closest printer model and test it.

16GB Data Storage

Data storage limit on the NanoDrop One/NanoDrop One^C is extended to 16GB. This extended memory will also be available to NanoDrop One/NanoDrop One^C instruments upgrading to v1.4.

Export Logs

Users can export logs to an USB flash drive by selecting **Diagnostics → Export Logs** option. Logs can only be exported to a connected USB flash drive.

“All” Option in Search Time Range

It is now possible to search all the experiments by selecting the “All” option in the Search Dialog’s time range dropdown.

Edit Experiment Name

In Experiment Details dialog the experiment name field is editable and the experiment name can be changed and saved.

Export .sql files for individual experiments

Users can now export .sql files for individual experiments, rather than a single .sql file for all experiments.

Resolved Issues

Custom Method Issues

This version resolves an issue where the entered wavelength was not used during analysis.

Protein Bradford

When saving standards under the Protein Bradford method, all numbers after the decimal are now included.

Known Issues

None

NanoDrop One PC Viewer Software v1.4

Release Date: May 2017

Supported Operating Systems

Windows 7 32-Bit
Windows 7 64-bit
Windows 10

New Features

Edit Experiment Name

In Experiment Details dialog the experiment name field is editable and the experiment name can be changed and saved.

Resolved Issues

Printing Intensity Check

An issue where the Xe monitored peaks were printed incorrectly when printing an Intensity Check experiment is now resolved.

Exporting to .tsv Format

Fixed an issue where an incorrect delimiter was used when exporting in .TSV format has now been resolved.

Known Issues

None

NanoDrop One Local Control Software v1.3 Hot Fix 1

Release Date: November 2016

Supported Operating Systems

NanoDrop One Instrument Android OS

New Features

Export for Cloud

A new feature has been added to Export to export the currently selected experiment to Cloud format to read into the NanoDrop Cloud application

Resolved Issues

None

Known Issues

None

NanoDrop One Local Control Software v1.3

Release Date: September 2016

Supported Operating Systems

NanoDrop One Instrument Android OS

New Features

Translations: Software and Help

Software and Help translations are available in the following languages:

- Chinese
- Japanese
- Polish
- Korean
- German
- French
- Spanish

The instrument is delivered with English software and Help. Language packs can be downloaded from our website: www.thermoscientific.com/nanodrop

Resolved Issues

Blank PC Name

PC Name in PC Viewer cannot be blank to connect to Wi-Fi, if it is blank the instrument will recognize that it is connected to a PC via Wi-Fi but it will not be available in the drop down menu.

Kinetics Interval Error Message

A dialog was displayed in error indicating that the kinetics interval was less than two seconds when the interval was acceptable.

GuSCN or GUHCl contamination:

When a sample is contaminated with GuSCN or GUHCl the Acclaro results displayed the incorrect result under the contaminate row-A260 column. This issue has been fixed.

Execution of Custom Methods:

Custom methods failed to run correctly, this issue has been fixed.

Known Issues

None

NanoDrop One PC Viewer Software v1.3

Release Date: September 2016

Supported Operating Systems

Windows 7 32-Bit
Windows 7 64-bit
Windows 10

New Features

Translations: Software and Help

Software and Help translations are available in the following languages:

- Chinese
- Japanese
- Polish
- Korean
- German
- French
- Spanish

The instrument is delivered with English software and Help. Language packs can be downloaded from our website: www.thermoscientific.com/nanodrop

Resolved Issues

Display Issues

When Overlay mode has been turned off, there were some issues with selection of sample. These issues have been fixed

Y-axis Label for Blank

The Y-axis label for blank measurements in the PC Data Viewer software was displaying 10mm Absorbance. This is incorrect and has been fixed

Known Issues

None

NanoDrop One Local Control Software v1.2

Release Date: February 2016

Supported Operating Systems

Windows 7 32-Bit
Windows 7 64-bit
Windows 10

New Features

UV-Vis Application

The UV-Vis application allows the NanoDrop One/NanoDrop One^C to function as a conventional spectrophotometer. Up to 40 wavelengths can be selected for absorbance monitoring and inclusion in the final results. Sample absorbance is displayed on the screen from 190nm to 840nm. UV-Vis experiment data can be viewed in the PC Viewer software.

Kinetics Application

The NanoDrop One^C can be used to make time-based kinetic measurements. Setting up kinetics experiments is easy to do and a visual display shows the stage information such as delay, measurement intervals and stage duration. Up to 5 stages can be created and up to 3 wavelengths can be monitored during the experiment.

The user also has the option to Pause or Stop the kinetics experiment. Kinetics data can be displayed as Time vs ABS or Time vs Delta ABS. Kinetics data can also be viewed in the NanoDrop One PC Viewer software.

Help System Translation

Translation for the help system and Acclaro is now included in the following languages; Chinese, French, German, Japanese, Korean, Polish and Spanish.

IQ/OQ

Qualification procedures are now available for the NanoDrop One and NanoDrop One^C. New part numbers have been released for full qualification kits in both US binder size and A4 binder size.

Additional Custom Methods

New custom method options are now available in the NanoDrop One PC Viewer software. These methods are easily created on the PC and executed on the NanoDrop One system. Standard curve custom method options are now available along with the formula and equation types that were available in an earlier version.

Acclaro Blanking

Acclaro now includes a check for blank. The software will notify the user if residue remains on the pedestal during blanking operations. If Acclaro has a concern you will be notified with a message and have the option to find our more information from our help system.

Wi-Fi Data Transfer

WiFi is now available for transferring data to a computer within your WiFi network. Any NanoDrop One instrument can be configured to send data or execute Custom Methods from a specified PC on the WiFi network. The data can then be reviewed and analyzed in more detail at your workstation away from the NanoDrop One instrument.

Resolved Issues**Known Issues****Data Size Limit**

There is a 4 GB limit on the android file system, which means that our database will be at the maximum storage capacity at 4GB. When the data storage is reaching the maximum data size, the user will be prompted to delete or export data to make room for additional data storage. Data can be imported to store and view using the NanoDrop One PC Viewer software that will not have this storage limitation.

NanoDrop One PC Viewer Software v1.2

Release Date: February 2016

Supported Operating Systems

Windows 7 32-Bit
Windows 7 64-bit
Windows 10

New Features**Wi-Fi Data Transfer**

WiFi is now possible using version 1.2 NanoDrop One PC Viewer software. Data can be seamlessly transferred from the NanoDrop One instrument to any WiFi connected PC. All data will be saved on the attached PC rather than on the local control instrument. Data can then be easily viewed using this software during the acquisition and after the experiment has completed.

Additional Custom Method

New custom method options are now available in the NanoDrop One PC Viewer software. These methods are easily created on the PC and executed on the NanoDrop One system. Standard curve

custom method options are now available along with the formula and equation types that were available in an earlier version.

Export .sql File Format

The database can now be exported from the NanoDrop One PC Viewer software. This was previously only an option from the local control but now it is available from the PC software as well.

Resolved Issues

None

Known Issues

None

NanoDrop One Local Control Software v1.1

Release Date: December 2015

Supported Operating Systems

NanoDrop One Instrument Android OS

New Features

Direct Connect (PC Tethering)

Ethernet direct connect is now possible using version 1.1 NanoDrop One software. Data can be seamlessly transferred from the NanoDrop One instrument to an Ethernet direct connected PC. All data will be saved on the attached PC rather than on the local control system when this options is selected. Data can be viewed from the NanoDrop One PC Viewer software.

Pedestal Image Check

Enhancements have been completed for the Pedestal Image Check diagnostic. Additional pass/fail information is now displayed along with more detailed information about any issues that may be encountered.

Software Translation

Translation for the local control software is now included in the following languages; Chinese, French, German, Japanese, Korean, Polish and Spanish

Resolved Issues

None

Known Issues

None

NanoDrop One PC Viewer Software v1.1

Release Date: December 2015

Supported Operating Systems

Windows 7 32-Bit
Windows 7 64-bit
Windows 10

New Features

Direct Connect (PC Tethering)

Ethernet direct connect is now possible using version 1.1 NanoDrop One software. Data can be seamlessly transferred from the NanoDrop One instrument to an Ethernet direct connected PC. All data will be saved on the attached PC rather than on the local control system when this options is selected. Data can be viewed from the NanoDrop One PC Viewer software.

Resolved Issues

None

Known Issues

None

NanoDrop One Local Control Software v1.0

Release Date: November 2015

Supported Operating Systems

NanoDrop One Instrument Android OS

New Features

Initial Release, all features are new

Resolved Issues

Initial release, there are no resolved Issues

Known Issues

None

NanoDrop One PC Viewer Software v1.0

Release Date: November 2015

Supported Operating Systems

Windows 7 32bit

Windows 7 64bit

Windows 10

New Features

Initial Release, all features are new

Resolved Issues

Initial release, there are no resolved Issues

Known Issues

None

How to Contact Us

Current contact information is located at www.thermofisher.com
Select the "Contact Us" icon at the top of the screen

Trademarks

Microsoft and Windows are either trademarks or registered trademarks of Microsoft Corporation in the United States and/or other countries.

All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.