

NanoDrop Eight Release Notes

About This Document	3
NanoDrop Eight PC Software v1.5	3
<i>Release Date: April 2023</i>	3
<i>Supported Operating Systems</i>	3
<i>New Features</i>	3
Add database auto back up and restore functionality	3
Add the ability to edit oligo sequences on each channel of Oligo DNA and Oligo RNA applications	3
<i>Resolved Issues</i>	3
This version resolves an issue related to experiment data loss when two Windows users log into the software at the same time.	3
This version resolves an issues where molecular weight is not editable in Custom Methods when the Result concentration is in a molar unit and the extinction coefficient in a mass unit.	3
This version resolves an issue in software version 1.4 Performance Verification where Channel letters were not displayed properly during measurement. in the NanoDrop Eight software version 1.4 PV channel incorrectly displaying the data during measuring.	3
<i>Known Issues</i>	3
NanoDrop Eight PC Software v1.4	4
<i>Release Date: September 2022</i>	4
<i>Supported Operating Systems</i>	4
<i>New Features</i>	4
Protein Bradford application	4
Protein Lowry application	4
Protein Pierce 660 application	4
<i>Resolved Issues</i>	4
This version resolves an issue where the concentration is not correctly calculated when selecting a custom protein from the Protein Editor as the sample type	4
This version updates the export files to display “NA” as the negative corrected concentration	4
<i>Known Issues</i>	4
NanoDrop Eight PC Software v1.3	4
<i>Release Date: June 2022</i>	4
<i>Supported Operating Systems</i>	4
<i>New Features</i>	4
OD600 application	4
Protein BCA application	4
Performance Verification adds PV-1 Lot Number tracking	4
Performance Verification Avg %Error value is calculated based on the Average column	4
Dye Editor wavelength update to 350 – 840 nm	4
<i>Resolved Issues</i>	5
This version resolves an issue with Protein A205 Extinction Coefficient incorrect display	5
This version resolves an issue with PV application results status not matching Print report results	5
<i>Known Issues</i>	5

NanoDrop Eight PC Software v1.2.1HF	5
<i>Release Date: May 2022</i>	5
<i>Supported Operating Systems</i>	5
<i>New Features</i>	5
<i>Resolved Issues</i>	5
This version updates Performance Verification tolerances	5
<i>Known Issues</i>	5
NanoDrop Eight PC Software v1.2	5
<i>Release Date: March 2022</i>	5
<i>Supported Operating Systems</i>	5
<i>New Features</i>	5
Protein A205.....	5
Microarray application	5
Proteins & Labels application	5
Dye Editor.....	6
Diagnostic add instrument serial number of the printed report	6
<i>Resolved Issues</i>	6
This version resolves an issue with displaying incorrect result in Intensity Check print report.....	6
This version resolves an issue with experiment date and measurement date not matching when import experiment data between different time zone	6
Dutch and French system	6
<i>Known Issues</i>	6
NanoDrop Eight PC Software v1.1	6
<i>Release Date: December 2021</i>	6
<i>Supported Operating Systems</i>	6
<i>New Features</i>	6
<i>Nucleic Acid</i>	6
ssDNA application	6
Oligo DNA/Oligo RNA applications	6
<i>Custom</i>	6
Custom Methods	6
<i>Resolved Issues</i>	6
This version resolves an issue with displaying data in History when the experiment timestamp was in the future compared to the local computer date & time.	6
This version resolved an issue where the E1% and E & MW fields were not user editable in Protein A280	6
<i>Known Issues</i>	7
NanoDrop Eight PC Software v1.0	7
<i>Release Date: September 2021</i>	7
<i>Supported Operating Systems</i>	7
<i>New Features</i>	7
<i>Nucleic Acid</i>	7
dsDNA application or RNA application.....	7

Custom Factor application..... 7

Protein..... 7

 Protein Editor..... 7

 Protein A280..... 7

UV-Vis Application 7

Diagnostics..... 7

 Performance Verification 7

 Intensity Check..... 8

IQ/OQ..... 8

Security with 21CFR part 11 compliance Support..... 8

General 8

 History 8

 Settings..... 8

 Help 8

Resolved Issues..... 8

Known Issues..... 8

How to Contact Us.....8

Trademarks8

About This Document

This document contains a revision history of Thermo Scientific NanoDrop™ Eight PC Software which includes new features, resolved and known issues.

NanoDrop Eight PC Software v1.5

Release Date: April 2023

Supported Operating Systems

- Windows 10
- Windows 11

New Features

- Add database auto back up and restore functionality**
- Add the ability to edit oligo sequences on each channel of Oligo DNA and Oligo RNA applications**

Resolved Issues

- This version resolves an issue related to experiment data loss when two Windows users log into the software at the same time.**
- This version resolves an issues where molecular weight is not editable in Custom Methods when the Result concentration is in a molar unit and the extinction coefficient in a mass unit.**
- This version resolves an issue in software version 1.4 Performance Verification where Channel letters were not displayed properly during measurement. in the NanoDrop Eight software version 1.4 PV channel incorrectly displaying the data during measuring.**

Known Issues

None

NanoDrop Eight PC Software v1.4

Release Date: September 2022

Supported Operating Systems

Windows 10
Windows 11

New Features

Protein Bradford application

Protein Lowry application

Protein Pierce 660 application

A standard curve is required for colorimetric protein analysis and support to import a standard curve. Run blank and measure samples using either the full 8-channel mode or a convenient single sample mode. Protein Bradford, Lowry, and 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/ .csv and .n8db).

Resolved Issues

This version resolves an issue where the concentration is not correctly calculated when selecting a custom protein from the Protein Editor as the sample type

This version updates the export files to display “NA” as the negative corrected concentration

Known Issues

None

NanoDrop Eight PC Software v1.3

Release Date: June 2022

Supported Operating Systems

Windows 10
Windows 11

New Features

OD600 application

Run blank and measure samples using either the full 8-channel mode or a convenient single sample mode 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/ .csv and .n8db).

Protein BCA application

A standard curve is required for colorimetric protein analysis and support to import a standard curve. Run blank and measure samples using either the full 8-channel mode or a convenient single sample mode. 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/ .csv and .n8db).

Performance Verification adds PV-1 Lot Number tracking

Performance Verification Avg %Error value is calculated based on the Average column

Dye Editor wavelength update to 350 – 840 nm

Resolved Issues

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

This version resolves an issue with PV application results status not matching Print report results

Known Issues

None

NanoDrop Eight PC Software v1.2.1HF

Release Date: May 2022

Supported Operating Systems

Windows 10

Windows 11

New Features

None

Resolved Issues

This version updates Performance Verification tolerances

Known Issues

None

NanoDrop Eight PC Software v1.2

Release Date: March 2022

Supported Operating Systems

Windows 10

Windows 11

New Features

Protein A205

Calculate peptide concentration based on user defined factor, run blank and measure samples using either the full 8-channel mode or a convenient single sample mode, 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. Export data in multiple file formats (.tsv/ .csv and .n8db).

Microarray application

Run blank and measure samples using either the full 8-channel mode or a convenient single sample mode, supports Analysis Correction. Microarray 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/ .csv and .n8db).

Proteins & Labels application

Run blank and measure samples using either the full 8-channel mode or a convenient single sample mode, supports Analysis Correction. Proteins & Labels 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/ .csv and .n8db).

Dye Editor

Provide 11 pre-defined dyes used to label sample material or add custom dye.

Diagnostic add instrument serial number of the printed report

Resolved Issues

This version resolves an issue with displaying incorrect result in Intensity Check print report.
This version resolves an issue with experiment date and measurement date not matching when import experiment data between different time zone
Dutch and French system

This version resolves pathlength selection issue in the Dutch and French system

Known Issues

None

NanoDrop Eight PC Software v1.1

Release Date: December 2021

Supported Operating Systems

Windows 10
Windows 11

New Features

Nucleic Acid

ssDNA application

Run blank and measure samples using either the full 8-channel mode or a convenient single sample mode, 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. Export data in multiple file formats (.tsv/ .csv and .n8db).

Oligo DNA/Oligo RNA applications

Run blank and measure samples using either the full 8-channel mode or a convenient single sample mode, 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/ .csv and .n8db).

Custom

Custom Methods

Formula and standard curve custom methods can be created in the Custom method application. Run blank and measure samples using either the full 8-channel mode or a convenient single sample mode, supports analysis wavelength and quantification step. Export data in multiple file formats (.tsv/ .csv and .n8db).

Resolved Issues

This version resolves an issue with displaying data in History when the experiment timestamp was in the future compared to the local computer date & time.
This version resolved an issue where the E1% and E & MW fields were not user editable in Protein A280

Known Issues

None

NanoDrop Eight PC Software v1.0**Release Date: September 2021****Supported Operating Systems**

Windows 10

New Features***Nucleic Acid*****dsDNA application or RNA application**

Run blank and measure samples using either the full 8-channel mode or a convenient single sample mode, supports baseline correction. dsDNA and RNA experiment data can be viewed in the History detail page. In experiment details page the experiment name and sample name fields are editable. Ability to customize what columns are displayed in History and in the measurement acquisition screen. Analyses that differentiate mammalian DNA from RNA and vice versa in dsDNA/RNA applications. Acclaro contaminant analysis includes a check for measurement. Export data in multiple file formats (.tsv/ .csv and .n8db).

Custom Factor application

Calculate nucleic acid concentration based on user defined factor. Run blank and measure samples using either the full 8-channel mode or a convenient single sample mode, 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. Export data in multiple file formats (.tsv/ .csv and .n8db).

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 measure samples using either the full 8-channel mode or a convenient single sample mode, 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 includes a check for measurement. Export data in multiple file formats (.tsv/ .csv and .n8db).

UV-Vis Application

The UV-Vis application allows the NanoDrop Eight to function as a conventional spectrophotometer. Up to 40 wavelengths can be selected for absorbance monitoring and inclusion in the results. Sample absorbance is displayed on the screen from 190nm to 850nm. UV-Vis experiment data can be viewed in the history.

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, conditional pass or fail results show at the status column. Experiment data can be viewed in the history detail page. Export data in multiple file formats (.tsv/ .csv and .n8db).

Intensity Check

Run an Intensity Check through the pedestal light paths. Pass or fail result shows at the top. Experiment data can be viewed in the History detail page. Export data in multiple file formats (.tsv/ .csv and .n8db).

IQ/OQ

Qualification procedures are available for the NanoDrop Eight PC Control software

Security with 21CFR part 11 compliance Support

NanoDrop Eight SciVault software, includes the User Privileges feature to set system policies and control user access to NanoDrop Eight software feature. Supports access control, system policies, and signature meanings. Audit Logs records all SciVault and NanoDrop Eight software audit trails.

General

History

Data can be viewed and updated from the History page

Settings

System Settings allow user to change SciVault Server Address. Turn on or off automatic multiple files export to specific location feature (Support local and network location)

Protein Editor

Preferences

Help

NanoDrop Eight User Guide

E-mail question to technical support

Navigate to software logs

Resolved Issues

None

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.