

NanoDrop Eight Release Notes

About This Document	2
NanoDrop Eight PC Software v1.0	2
<i>Release Date: September 2021.....</i>	<i>2</i>
<i>Supported Operating Systems.....</i>	<i>2</i>
<i>New Features.....</i>	<i>2</i>
<i>Nucleic Acid</i>	<i>2</i>
dsDNA application or RNA application.....	2
Custom Factor application.....	2
<i>Protein.....</i>	<i>2</i>
Protein Editor.....	2
Protein A280.....	2
<i>UV-Vis Application</i>	<i>2</i>
<i>Diagnostics.....</i>	<i>2</i>
Performance Verification	2
Intensity Check.....	3
<i>IQ/OQ.....</i>	<i>3</i>
<i>Security with 21CFR part 11 compliance Support.....</i>	<i>3</i>
<i>General</i>	<i>3</i>
History	3
Settings.....	3
Help	3
<i>Resolved Issues.....</i>	<i>3</i>
<i>Known Issues.....</i>	<i>3</i>
How to Contact Us.....	3
Trademarks	3

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 v1.0

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.