

NanoDrop 8000 Release Notes

April 2020

1. About This Document	2
2. About the NanoDrop 8000	2
3. Released Versions	2
NanoDrop 8000- v.2.3.3 – April 2020	2
NanoDrop 8000- v.2.3.2a – December 2016.....	2
NanoDrop 8000- v.2.3.2 – June 2013.....	2
NanoDrop 8000- v.2.2.1 – June 2011.....	2
NanoDrop 8000- v.2.2.0 – January 2011.....	2
NanoDrop 8000- v.2.1.0 – September 2009.....	3
4. Technical Support	3

1. About This Document

This document contains information about the NanoDrop™ 8000 software features and resolved issues.

2. About the NanoDrop 8000

Thermo Scientific™ NanoDrop 8000 software is used to perform measurements with the NanoDrop 8000 instrument. See Windows Compatibility list for each version below.

Consult the computer requirements section of the User manual for additional requirements for individual hardware or software components.

3. Released Versions

NanoDrop 8000- v.2.3.3 – April 2020

New Features

- No new software capabilities

Resolved Issues

- Added improved compatibility with Windows 10 64-bit
- No longer compatible with Windows 2000, Windows XP or Windows Vista

NanoDrop 8000- v.2.3.2a – December 2016

New Features

- No new software capabilities

Resolved Issues

- Added compatibility with Windows 8.1 64 bit
- Added compatibility with Windows 10 64-bit
- No longer compatible with Windows 2000, Windows XP or Windows Vista

NanoDrop 8000- v.2.3.2 – June 2013

New Features

- No new software capabilities

Resolved Issues

- Improved linearity correction implementation
- Compatible with Windows™ XP (32-bit)
- Compatible with Windows Vista™ (32-bit) or
- Compatible with Windows 7 (32-bit or 64-bit)

NanoDrop 8000- v.2.2.1 – June 2011

New Features

- No new software capabilities

Resolved Issues

- Corrected setting that disallowed .nd8 files to open directly in Data Viewer.

NanoDrop 8000- v.2.2.0 – January 2011

New Features

- Revision of drivers and installer to support Windows 7 (32-bit and 64-bit) compatibility.

- Calibration Check results now save as .TXT files in addition to .JPG files.
- Revised Calibration Check User interface now displays absorbance profile and absorbance value for every measurement replicate.
- Historical Calibration Check Files may be imported into the Calibration Check application for review.

Resolved Issues

- No newly resolved issues

NanoDrop 8000- v.2.1.0 – September 2009

New Features

- Includes capability to run operate the instrument in a single position mode (colorimetric applications not included.) The new function includes sample ID import options and plot displays optimized for the single sample measurements.
- Added data export capability by plate row sorting as an alternative to plate column sorting.
- Expanded wavelength range for User Methods utilizing Standards.
- Included an autopath selection feature for User Methods.
- Included new prompt to Blank before loading a sample ID plate file for more efficient operation.

Resolved Issues

- Corrected methods list loading issue associated with opening the software via an archived file.
- Removed non-English LV support to prevent potential errors in the menus of some non-English installations.
- Modified graph ranges displayed on diagnostic screens.
- Improved application result units recalculation and display behavior.
- Corrected feature to allow for molecular weight based oligo nucleic acid concentration results upon switching back from other sample type selections.
- Corrected algorithms to treat dye correction factors as fractions rather than percent.
- Corrected report column title display error.
- Linked column break message to absorbance values at the wavelength associated with the cursor position for custom methods that do not specify a quantification mode.

4. Technical Support

Thermo Fisher Scientific
3411 Silverside Road
Bancroft Building, Suite 100
Wilmington, DE 19810 U.S.A.
Telephone: 302-479-7707
Fax: 302-792-7155
E-mail: info@nanodrop.com
www.nanodrop.com

For International Support, please contact your local distributor.

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.