

**NanoDrop 2000- 2000c Release Notes**

August 2014

**About This Document** ..... 3

**About NanoDrop 2000 Software** ..... 3

**NanoDrop 2000 1.6.198 – August 2014** ..... 3

*New Features* ..... 3

        January 2017: Tested compatibility with Windows 10 Pro ..... 3

        Compatibility with Windows 8.1 ..... 3

        Region Control added to Kinetics ..... 3

*Resolved Issues* ..... 3

        Workbooks could be corrupted and only partially recoverable ..... 3

        User was not required to re-blank when switching application "modes" ..... 3

        Auto path selection not properly displayed ..... 3

        NanoDrop 2000 Simulator did not install on Windows 7 or Windows 8 ..... 3

        Removed the check box "Recalibrate Wavelengths upon measure" from Diagnostics ..... 3

        Calibration check diagnostic did not flag out of limit standard deviation value ..... 3

        No serial number in calibration check ..... 3

        Protein A280 Reprocess Button Issue ..... 4

**NanoDrop 2000 1.5 – May 2013** ..... 4

*New Features* ..... 4

        NanoDrop installs on Windows 8 but requires manual installation of .NET 3.5 ..... 4

*Resolved Issues* ..... 4

        No new issues resolved ..... 4

**NanoDrop 2000 1.4.2 – March 17, 2011** ..... 4

*New Features* ..... 4

        No new features ..... 4

*Resolved Issues* ..... 4

        "Cannot communicate" errors ..... 4

**NanoDrop 2000 1.4 – February 22, 2010** ..... 4

*New Features* ..... 4

        No new features ..... 4

*Resolved Issues* ..... 4

        Windows 7 (32-bit and 64-bit) compatibility ..... 4

**NanoDrop 2000 1.3.1 – June 17, 2009** ..... 4

*Resolved Issues* ..... 4

        "Low Bias" errors in Intensity Check/Diagnostics ..... 4

**NanoDrop 2000 1.2.2 – May 2009** ..... 5

*New Features* ..... 5

No new features ..... 5

*Resolved Issues* ..... 5

    “Low Bias” errors in Intensity Check/Diagnostics ..... 5

**NanoDrop 2000 1.1.9 – April 2009** ..... **5**

*New Features* ..... 5

        No new features ..... 5

*Resolved Issues* ..... 5

        Conflict with running non-English Regional and Language Options ..... 5

**Technical Support** ..... **5**

## About This Document

This document contains information about the NanoDrop 2000/2000c software features and resolved issues.

## About NanoDrop 2000 Software

Thermo Scientific NanoDrop 2000 software is used to perform measurements with the NanoDrop 2000/2000c instrument. The most recent version of NanoDrop 2000 software has been verified on Professional versions of Windows 7 (both 32-bit and 64-bit), Windows 8.1 (64-bit) and Windows 10 (64-bit).

Consult the computer requirements document for additional requirements for individual hardware or software components.

## NanoDrop 2000 1.6.198 – August 2014

### New Features

#### ***Compatibility with Windows 10***

January 2017: NanoDrop version 1.6 has been tested and approved for use on Windows 10 (64-bit) Pro.

#### ***Compatibility with Windows 8.1***

NanoDrop version 1.6 has been tested and approved for use on Windows 8.1 (64-bit) Professional.

#### ***Region Control added to Kinetics***

Region control is now available in the post acquisition screen. It can be used to establish end-points for the curve.

### Resolved Issues

#### ***Workbooks could be corrupted and only partially recoverable***

Large workbooks being saved on a slow system could be corrupted if the user closed the software before the save operation has completed. Preventative measures have been added to NanoDrop to minimize the possibility a user could prematurely terminate the save process.

#### ***User was not required to re-blank when switching application "modes"***

Fixed: If a user did not re-blank when entering an application, internal settings from the previous application could alter the measured data and give incorrect results.

#### ***Auto path selection not properly displayed***

Fixed: Although the software utilizes the optimal path length when the auto path option was selected for a custom method, the check box for the option was incorrectly displayed as not checked once the method is saved.

#### ***NanoDrop 2000 Simulator did not install on Windows 7 or Windows 8***

Fixed: The simulator installer blocked installation on all Windows 7 (and newer) operating systems.

#### ***Removed the check box "Recalibrate Wavelengths upon measure" from Diagnostics***

Inactive function: button removed.

#### ***Calibration check diagnostic did not flag out of limit standard deviation value***

When the calibration check procedure on the diagnostic page is run the standard deviation of the measurements for each pedestal position is tracked. If the value is outside the limits coded into the procedure, a calibration failure is now flagged.

#### ***No serial number in calibration check***

The instrument serial number is now displayed or recorded in the calibration check screen shot.

***Protein A280 Reprocess Button Issue***

Fixed: When either BSA, IGG or 1 ABS = 1mg per mil sample type was selected, a sample was measured and the user selected the Report tab, and then Reprocess, if they input E and MW (other protein) as the sample type, a concentration value is not generated.

**NanoDrop 2000 1.5 – May 2013****New Features*****NanoDrop installs on Windows 8 but requires manual installation of .NET 3.5***

Although the software has been approved for installation and operation on Windows 8, the user must first manually obtain and install Microsoft .Net version 3.5 as a required precursor.

**Resolved Issues**

*No new issues resolved*

**NanoDrop 2000 1.4.2 – March 17, 2011****New Features**

*No new features*

**Resolved Issues*****“Cannot communicate” errors***

Fixed: Various instrument communication errors and driver installation issues. Revision of the driver to correct for the “Cannot Communicate” and “Cannot Start Driver” errors.

**NanoDrop 2000 1.4 – February 22, 2010****New Features**

*No new features*

**Resolved Issues*****Windows 7 (32-bit and 64-bit) compatibility***

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

**NanoDrop 2000 1.3.1 – June 17, 2009****Resolved Issues*****“Low Bias” errors in Intensity Check/Diagnostics***

Corrected low and high bias limits for Intensity Check/Diagnostics. It now updates the bias value on the internal spectrometer when needed.

## **NanoDrop 2000 1.2.2 – May 2009**

### **New Features**

*No new features*

### **Resolved Issues**

#### ***“Low Bias” errors in Intensity Check/Diagnostics***

Corrected low and high bias limits for Intensity Check/Diagnostics.

## **NanoDrop 2000 1.1.9 – April 2009**

### **New Features**

*No new features*

### **Resolved Issues**

#### ***Conflict with running non-English Regional and Language Options***

Corrected issues which occurred upon application startup when using the software with Windows XP or Vista with regional languages that handle number strings differently than the English (United States) setting.

## **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: [nanodrop@thermofisher.com](mailto:nanodrop@thermofisher.com)  
[thermofisher.com/nanodrop](http://thermofisher.com/nanodrop)

For International Support, please contact your local distributor.

Microsoft, Windows is either a trademark or registered trademark 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.