

NanoDrop 3300 Release Notes

August 2020

1. About This Document	2
2. About the NanoDrop 3300	2
3. Released Versions	2
NanoDrop 3300- version 2.8.1 – August 2020	2
NanoDrop 3300- version 2.8.0.1 – December 2016	2
Older versions	2
4. Technical Support	2

1. About This Document

This document contains information about the NanoDrop™ 3300 software features and resolved issues.

2. About the NanoDrop 3300

Thermo Scientific™ NanoDrop 3300 software is used to perform measurements with the NanoDrop 3300 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 3300- version 2.8.1 – August 2020

New Features

- No new software capabilities

Resolved Issues

- Added compatibility with Windows 10, 64-bit for build versions 1903 or greater

NanoDrop 3300- version 2.8.0.1 – 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

Older versions

- No release notes are available for previous versions

4. Technical Support

Thermo Fisher Scientific
3411 Silverside Road
Tatnall Building, Suite 100
Wilmington, DE 19810 U.S.A.
Telephone: 302-479-7707
Fax: 302-792-7155
Email: nanodrop@thermofisher.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.