

NanoDrop Lite Plus Release Notes

| | |
|---|----------|
| About This Document..... | 3 |
| NanoDrop Lite Plus Software v1.2.1 | 3 |
| <i>Release Date: November 2023.....</i> | <i>3</i> |
| <i>Supported Hardware Systems.....</i> | <i>3</i> |
| <i>New Features.....</i> | <i>3</i> |
| <i>Resolved Issues.....</i> | <i>3</i> |
| This version resolves an issue updating the wavelength offset calculation method | <i>3</i> |
| This version resolves an issue which was triggering an instrument initialization failure to occur | <i>3</i> |
| <i>Known Issues.....</i> | <i>3</i> |
| NanoDrop Lite Plus Software v1.2 | 3 |
| <i>Release Date: July 2023</i> | <i>3</i> |
| <i>Supported Hardware Systems.....</i> | <i>3</i> |
| <i>New Features.....</i> | <i>3</i> |
| <i>Support new hardware system MCU STM32H743.....</i> | <i>3</i> |
| <i>Resolved Issues.....</i> | <i>3</i> |
| <i>Known Issues.....</i> | <i>3</i> |
| NanoDrop Lite Plus Software v1.1 | 3 |
| <i>Release Date: July 2022</i> | <i>3</i> |
| <i>Supported Hardware Systems.....</i> | <i>3</i> |
| <i>New Features.....</i> | <i>3</i> |
| <i>Chinese language support in Settings menu.....</i> | <i>3</i> |
| <i>Resolved Issues.....</i> | <i>3</i> |
| <i>Known Issues.....</i> | <i>3</i> |
| NanoDrop Lite Plus Software v1.0 | 4 |
| <i>Release Date: May 2022.....</i> | <i>4</i> |
| <i>Supported Hardware Systems.....</i> | <i>4</i> |
| <i>New Features.....</i> | <i>4</i> |
| <i>Nucleic Acid.....</i> | <i>4</i> |
| dsDNA application, RNA application and ssDNA application | <i>4</i> |
| Custom Factor application..... | <i>4</i> |
| <i>Protein</i> | <i>4</i> |
| 1Abs=1mg/mL, BSA, IgG, Lysozyme..... | <i>4</i> |
| Other Protein E & MW..... | <i>4</i> |
| Other Protein E1%..... | <i>4</i> |
| <i>Maintenance</i> | <i>4</i> |
| Performance Verification..... | <i>4</i> |

Diagnostics 5
History 5
Resolved Issues 5
Known Issues 5
How to Contact Us 5

About This Document

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

NanoDrop Lite Plus Software v1.2.1

Release Date: November 2023

Supported Hardware Systems

STM32F469, STM32H743, FreeRTOS

New Features

None

Resolved Issues

This version resolves an issue updating the wavelength offset calculation method

This version resolves an issue which was triggering an instrument initialization failure to occur

Known Issues

None

NanoDrop Lite Plus Software v1.2

Release Date: July 2023

Supported Hardware Systems

STM32F469, STM32H743, FreeRTOS

New Features

Support new hardware system MCU STM32H743

Resolved Issues

None

Known Issues

None

NanoDrop Lite Plus Software v1.1

Release Date: July 2022

Supported Hardware Systems

STM32F469, FreeRTOS

New Features

Chinese language support in Settings menu

Resolved Issues

None

Known Issues

None

NanoDrop Lite Plus Software v1.0

Release Date: May 2022

Supported Hardware Systems

STM32F469, FreeRTOS

New Features

Nucleic Acid

dsDNA application, RNA application and ssDNA application

Run blank and measure samples, supports baseline correction at 340nm and customization of the sample name is available. dsDNA, ssDNA and RNA experiment data can be viewed in the History detail page. If the printer is connected, you can click the printer button on the History detail screen or the Measurement results screen to print the measurement results to the label. In the History page data can be exported to a USB flash driver in .csv file format.

Custom Factor application

Calculate nucleic acid concentration based on user defined factor. Run blank and measure samples, supports baseline correction at 340nm and customization of the sample name is available. Custom factor experiment data can be viewed in the History detail page. If the printer is connected, you can click the printer button on the History detail screen or the Measurement results screen to print the measurement results to the label. In the History page data can be exported to a USB flash driver in .csv file format.

Protein

1Abs=1mg/mL, BSA, IgG, Lysozyme

Each application uses a different E1% parameter. Run blank and measure samples, supports baseline correction on 340nm and customize the sample name. Protein application experiment data can be viewed in the History detail page. In the History page data can be exported to a USB flash driver in .csv file format.

Other Protein E & MW

This application allows the user to input Molar Extinction Coefficient and Molecular Weight parameters. Run blank and measure samples, supports baseline correction at 340nm and customization of the sample name is available. Other Protein E&MW experiment data can be viewed in the History detail page. In the History page data can be exported to a USB flash driver in .csv file format.

Other Protein E1%

This application allows the user to input E1% Extinction Coefficient parameter. Run blank and measure samples, supports baseline correction at 340nm and customization of the sample name is available. Other Protein E1% experiment data can be viewed in the History detail page. In the History page data can be exported to a USB flash driver in .csv file format.

Maintenance

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 View Previous PV detail page. Exportation of data to a USB flash driver is available in .txt format.

Diagnostics

Press Run button to start system diagnostics. The test results are immediately displayed on the measurement interface and show the PASS or FAIL criteria to the right of the results. The diagnostic results reflect the strength of the energy of the instrument's lamp and the wavelength accuracy of the CGS to pass or fail.

History

Measurement results will be displayed here in a list, 100 records will be displayed on each page, supporting up to 10 pages with a total of 1,000 history records; support for viewing and printing data and exporting data to a USB flash drive.

Resolved Issues

Initial release, there are no resolved issues

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