

## NanoDrop Ultra Release Notes

<b>About This Document.....</b>	<b>2</b>
<b>NanoDrop Ultra Local Control Software v1.0.1.....</b>	<b>2</b>
Release Date: January 2025 .....	2
Supported Operating Systems.....	2
New Features.....	2
Resolved Issues.....	2
Plant and Bacterial Deconvolution.....	2
Improved Fluorescence applications algorithms .....	2
Known Issues.....	2
<b>NanoDrop Ultra PC Control Software v1.0.1.....</b>	<b>2</b>
Release Date: January 2025 .....	2
Supported Operating Systems.....	2
New Features.....	2
Resolved Issues.....	2
Plant and Bacterial Deconvolution.....	2
Improved Fluorescence applications algorithms .....	2
Known Issues.....	2
<b>NanoDrop Ultra Local Control Software v1.0 .....</b>	<b>2</b>
Release Date: January 2025 .....	2
Supported Operating Systems.....	2
New Features.....	2
Resolved Issues.....	3
Known Issues.....	3
<b>NanoDrop Ultra PC Control Software v1.0.....</b>	<b>3</b>
Release Date: January 2025 .....	3
Supported Operating Systems.....	3
New Features.....	3
Resolved Issues.....	3
Known Issues.....	3
<b>How to Contact Us.....</b>	<b>3</b>
<b>Trademarks .....</b>	<b>3</b>

## About This Document

This document contains a revision history of Thermo Scientific NanoDrop™ Ultra/Ultra<sup>C</sup>/Ultra FL/Ultra<sup>C</sup> FL local control and PC control software which includes new features, resolved and known issues.

### **NanoDrop Ultra Local Control Software v1.0.1**

**Release Date:** January 2025

#### **Supported Operating Systems**

NanoDrop Ultra Instrument Linux OS

#### **New Features**

None

#### **Resolved Issues**

##### **Plant and Bacterial Deconvolution**

Improved algorithms of plant and bacterial in dsDNA/RNA applications

**Improved Fluorescence applications algorithms**

#### **Known Issues**

None

### **NanoDrop Ultra PC Control Software v1.0.1**

**Release Date:** January 2025

#### **Supported Operating Systems**

Windows 10 64-bit

Windows 11 64-bit

#### **New Features**

None

#### **Resolved Issues**

##### **Plant and Bacterial Deconvolution**

Improved algorithms of plant and bacterial in dsDNA/RNA applications

**Improved Fluorescence applications algorithms**

#### **Known Issues**

None

### **NanoDrop Ultra Local Control Software v1.0**

**Release Date:** January 2025

#### **Supported Operating Systems**

NanoDrop Ultra Instrument Linux OS

#### **New Features**

Initial Release, all features are new

**Resolved Issues**

Initial release, there are no resolved Issues

**Known Issues**

None

**NanoDrop Ultra PC Control Software v1.0**

**Release Date: January 2025**

**Supported Operating Systems**

Windows 10 64-bit

Windows 11 64-bit

**New Features**

Initial Release, all features are new

**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](http://www.thermofisher.com)

Select "Contact Us" at the bottom of the screen under the Support section

**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.