

AccuSEQ™ Real-Time PCR Detection Software v2.1.2

Windows™ 10 Operating System

Catalog Numbers A45662

Publication Number MAN0018821 Revision A.0

Overview

The AccuSEQ™ Real-Time PCR Detection Software v2.1.2 (Windows™ 10 Operating System) is for use with the 7500 Fast Real-Time PCR Instrument.

Minimum computer requirements

IMPORTANT! We strongly recommend that you use computers that are provided by Thermo Fisher Scientific. All validation and verification studies were performed using Thermo Fisher Scientific computers. If you install the software on your own computer that meets the minimum specifications listed below, you must perform comprehensive testing to ensure proper operation of the software.

| Specification | Requirement |
|------------------|---|
| Processor | Intel™ Core™ i5 (Quad Core 2.6 GHz) Processor |
| Hard drive | ≥ 250 GB for software installation and data storage |
| Memory | ≥16 GB |
| Operating system | Windows™ 10, 64-bit operating system (IOT Enterprise) |

The following computers are compatible with AccuSEQ™ Real-Time PCR Detection Software v2.1.2 software and are available from Thermo Fisher Scientific:

- OptiPlex™ XE2 with Windows™ 10 IOT 2016-64Bit
- Dell™ Latitude™ E5570 Laptop with Windows™ 10 IOT 2016-64Bit
- Dell™ Latitude™ E5580 Laptop with Windows™ 10 IOT 2016-64Bit

Changes in the AccuSEQ™ software v2.1, v2.1.1, and v2.1.2

| Software version | Changes |
|------------------|---|
| 2.1 | <ul style="list-style-type: none"> • Automatic backup of .eds files function—After each run, a backup file is saved in the "<Install Dir>:\Applied Biosystems\AccuSEQ\experiments_temp" folder. Timestamps will be appended to backup file names. • Dilution Calculator—In Custom Mode > Setup > Plate Setup > Define Targets and Samples, the dilution factor used to prepare a sample can be entered. During analysis, the Dilution Factor is multiplied by the Quantity to automatically calculate the Dilution Adjusted concentration. • Efficiency/Recovery Calculator—In Custom Mode > Setup > Plate Setup > Define Targets and Samples, the spike concentration and reference information can be entered. During analysis, Spike Input and Reference are used to automatically calculate the value for Recovery/Efficiency. • Administrator MycoSEQ Call Defaults—In MycoSEQ Module > Analysis (top menu) > Analysis Settings..., a set of "Administrator Analysis Settings" mycoSEQ call settings can be set and saved. The "Factory Defaults" are always stored. "Administrator Defaults" have priority over the "Factory Defaults". |
| 2.1.1 | Defect fixes. |
| 2.1.2 | Windows™ 10 operating system. No software features have been changed or added. |

Related documentation

| Document | Publication number |
|--|--------------------|
| <i>AccuSEQ™ software Help</i> Access the help from within the software. | N/A |
| <i>AccuSEQ™ software: Mycoplasma Getting Started Guide</i> | 4425587 |
| <i>AccuSEQ™ software: Mycoplasma Quick Reference Card</i> | 4425586 |
| <i>AccuSEQ™ software: Custom Quick Reference Card</i> | 4425585 |

Customer and technical support

Visit thermofisher.com/support for the latest service and support information.

- Worldwide contact telephone numbers
- Product support information
 - Product FAQs
 - Software, patches, and updates
 - Training for many applications and instruments
- Order and web support

- Product documentation
 - User guides, manuals, and protocols
 - Certificates of Analysis
 - Safety Data Sheets (SDSs; also known as MSDSs)

Note: For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



Life Technologies Corporation | 200 Oyster Point Blvd | South San Francisco, CA 94080 | USA

For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Revision history: Pub. No. MAN0018821

| Revision | Date | Description |
|----------|-------------------|--|
| A.0 | 20 September 2019 | New document for the AccuSEQ™ Real-Time PCR Detection Software v2.1.2 on the Windows™ 10 Operating System. |

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

Trademarks: All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. Dell and Latitude are trademarks of Dell Computer Corporation.

©2019 Thermo Fisher Scientific Inc. All rights reserved.

thermofisher.com/support | thermofisher.com/askaquestion

thermofisher.com

20 September 2019

ThermoFisher
SCIENTIFIC

Draft-Comments