

# MycoSEQ™ Mycoplasma Real-Time PCR Detection Kit

**Catalog Numbers** 4460623, 4460626





**Pub. No.** 4482248 **Rev.** B

**Note:** For safety and biohazard guidelines, see the “Safety” appendix in the *MycoSEQ™ Mycoplasma Real-Time PCR Detection Kit User Guide* (Pub. No. 4465874). Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

## Product description

The MycoSEQ™ Mycoplasma Real-Time PCR Detection Kit detects *Mycoplasma* species simply, reliably, and rapidly.

## Contents and storage conditions

| Contents <sup>[1]</sup>   | Cap color   | Amount <sup>[2]</sup> | Storage   |
|---|---|-----------------------|---|
| <b>Box 1: MycoSEQ™ Mycoplasma Real-Time PCR Detection Kit</b>         |   |                       |   |
| <i>Mycoplasma</i> Real-Time PCR Primer Mix (10×)                      | <br>blue     | 325 µL                | –25°C to –15°C on receipt<br>2–8°C after first use  |
| Negative Control  | <br>white    | 1,000 µL              |   |
| <i>Power</i> SYBR™ Green PCR Master Mix (2×)                          | <br>white  | 2 × 1,000 µL          | –25°C to –15°C on receipt, protected from light.<br>2–8°C after first use, protected from light |
| <b>Box 2: MycoSEQ™ Discriminatory Positive/Extraction Control</b>     |   |                       |   |
| MycoSEQ™ Discriminatory Positive/ Extraction Control, 1,000 copies/µL | <br>yellow | 700 µL                | –25°C to –15°C  |

<sup>[1]</sup> To purchase the MycoSEQ™ Mycoplasma Real-Time PCR Detection Kit that includes the PrepSEQ™ Mycoplasma Sample Preparation Kit, use Catalog Number 4460626.

<sup>[2]</sup> The kit contains reagents for 100 reactions.

## 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 found on Life Technologies' website at [www.thermofisher.com/us/en/home/global/terms-and-conditions.html](http://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](http://www.thermofisher.com/support).



**Manufacturer:** Thermo Fisher Scientific | 7 Kingsland Grange | Warrington, Cheshire WA1 4SR | United Kingdom

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. 4482248

| Revision | Date        | Description                               |
|----------|-------------|---|
| B        | 24 May 2018 | Updated to the current document template. |
| A        | Dec 2013    | New document.                             |

**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.

©2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.