


MicroSEQ™ *Listeria monocytogenes* Detection Kit

Catalog Number 4403874

Pub. No. 4403929 Rev. E

 **WARNING!** Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from thermofisher.com/support.

Product description

The MicroSEQ™ *Listeria monocytogenes* Detection Kit rapidly assesses pre-enriched food and environmental samples for the presence of *Listeria monocytogenes*. The kit includes a lyophilized assay that provides all reagents (except for template DNA) required for real-time PCR detection of *Listeria monocytogenes* DNA. In addition, each assay bead includes an Internal Positive Control (IPC) to identify possible PCR inhibition.

Contents and storage

Component	Description	Amount	Cap color	Storage ^[1]
MicroSEQ™ <i>Listeria monocytogenes</i> Detection Kit	<i>Listeria monocytogenes</i> Assay Beads, 8-tube strips in rack	12 strips (96 tubes) 1 rack	Blue (rack)	5±3°C Protect from light and moisture. ^[2]
	MicroAmp™ Optical 8-Cap Strips	12 strips (96 caps)	N/A	
Pathogen Detection Negative Control ^[3]	Pathogen Detection Negative Control	1.5 mL	Red	5±3°C

^[1] See the product label for the expiration date.

^[2] Excessive exposure to light may affect the fluorescent probes. To protect the beads from moisture, do not remove the desiccant from the pouch, and seal the pouch tightly each time you remove assay bead strips.

^[3] The Pathogen Detection Negative Control is included in a separate box and may be shipped separately.

Procedural guidelines

- For detailed procedures, see the product web page at thermofisher.com; search for Cat. No. 4403874.
- For a list of workflows for detection of *Listeria*, see Pub. No. MAN0009418 at: tools.thermofisher.com/content/sfs/manuals/MAN0009418.pdf

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. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.

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

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES 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. 4403929

Revision	Date	Description
E	11 January 2017	<ul style="list-style-type: none">Remove expiration date; refer to the product labelUpdate format to current style
D	June 2009	Baseline for the revision

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

Limited Use Label License No. 492: Environmental Testing, Quality Control/Quality Assurance Testing, Food and Agricultural Testing: Notice to Purchaser: The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product (a) to perform internal research for the sole benefit of the purchaser; and (b) for environmental testing, quality control/quality assurance testing, food and agricultural testing, including reporting results of purchaser's activities in environmental testing, quality control/quality assurance testing, food and agricultural testing for a fee or other commercial consideration. No other right is hereby granted expressly, by implication, or by estoppel. This product is for environmental testing, quality control/ quality assurance testing, food and agricultural testing and research purposes only. The purchase of this product does not grant the purchaser any additional rights, including (without limitation) the right to transfer or resell the product in any form or the right to use the product as a therapeutic agent or diagnostics test component. For information on obtaining additional rights, please contact outlicensing@thermofisher.com or Out Licensing, Life Technologies Corporation (part of Thermo Fisher Scientific), 5823 Newton Drive, Carlsbad, California 92008.

Corporate entity: Life Technologies Corporation | Carlsbad, CA 92008 USA | Toll Free in USA 1 800 955 6288

©2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.