

Workflows for Detection of *Listeria* in Food and Environmental Samples

Pub. No. MAN0009418 Rev. B.0

Note: Validated workflows include enrichment, sample preparation, detection, and confirmation methods described in the specified user guides. Refer to the appropriate publication for the specific matrices included in each validated workflow.

Table 1 NF Validation™-certified workflows (AFNOR) for PCR detection

Workflow overview				
Species	<i>Listeria</i> spp.	<i>Listeria</i> spp.	<i>Listeria monocytogenes</i>	<i>Listeria monocytogenes</i>
NF Validation categories	All food and feed categories and environmental samples	All food and feed categories and environmental samples	All food and feed categories and environmental samples	All food and feed categories and environmental samples
Nucleic acid (NA) isolation (sample prep)	PrepSEQ® Nucleic Acid Extraction Kit (Cat. nos. 4480466, 4428176)	PrepSEQ® Rapid Spin Sample Preparation Kits—Bead Beating [Cat. nos. 4464654, 4468304 (Extra Clean; for meat samples)]	PrepSEQ® Nucleic Acid Extraction Kit (Cat. nos. 4480466, 4428176)	PrepSEQ® Rapid Spin Sample Preparation Kits—Bead Beating [Cat. nos. 4464654, 4468304 (Extra Clean; for meat samples)]
NA isolation workflow	Automated	Spin column + mechanical/heat lysis	Automated	Spin column + mechanical/heat lysis
PCR detection	MicroSEQ® <i>Listeria</i> spp. Detection Kit (Cat. no. 4427410)	MicroSEQ® <i>Listeria</i> spp. Detection Kit (Cat. no. 4427410)	MicroSEQ® <i>Listeria monocytogenes</i> Detection Kit (Cat. no. 4403874)	MicroSEQ® <i>Listeria monocytogenes</i> Detection Kit (Cat. no. 4403874)
User guide (Pub. no.)	4470047: Sample enrichment and NA isolation 4489329 ^[1] : PCR ^[2] and confirmation	4467823: Sample enrichment and NA isolation 4489329 ^[1] : PCR ^[2] and confirmation	4470047: Sample enrichment and NA isolation 4489329 ^[1] : PCR ^[2] and confirmation	4467823: Sample enrichment and NA isolation 4489329 ^[1] : PCR ^[2] and confirmation

^[1] Replaces Pub. nos. 4405962 and 4426510.^[2] Refer to the specified user guide for the real-time PCR instruments and analysis software included in the NF Validation™-certified workflow.Table 2 AOAC® Performance Tested MethodsSM-certified workflows for PCR detection

Workflow overview				
Species	<i>Listeria</i> spp.	<i>Listeria</i> spp.	<i>Listeria monocytogenes</i>	<i>Listeria monocytogenes</i>
Performance Tested MethodsSM matrices	Food and environmental surfaces	Food and environmental surfaces	Food matrices	Food matrices
Nucleic acid (NA) isolation (sample prep)	PrepSEQ® Nucleic Acid Extraction Kit (Cat. nos. 4480466, 4428176)	PrepSEQ® Rapid Spin Sample Preparation Kits with Proteinase K [Cat. nos. 4426714, 4426715 (Extra Clean)]	PrepSEQ® Nucleic Acid Extraction Kit (Cat. nos. 4480466, 4428176)	PrepSEQ® Rapid Spin Sample Preparation Kits with Proteinase K [Cat. nos. 4426714, 4426715 (Extra Clean)]
NA isolation workflow	Automated	Spin column	Automated	Spin column
PCR detection	MicroSEQ® <i>Listeria</i> spp. Detection Kit (Cat. no. 4427410)	MicroSEQ® <i>Listeria</i> spp. Detection Kit (Cat. no. 4427410)	MicroSEQ® <i>Listeria monocytogenes</i> Detection Kit (Cat. no. 4403874)	MicroSEQ® <i>Listeria monocytogenes</i> Detection Kit (Cat. no. 4403874)
User guide (Pub. no.)	4426514: Sample enrichment and NA isolation 4489329 ^[1] : PCR ^[2] and confirmation	4426506: Sample enrichment and NA isolation 4489329 ^[1] : PCR ^[2] and confirmation	4405966: Sample enrichment and NA isolation 4489329 ^[1] : PCR ^[2] and confirmation	4412851: Sample enrichment and NA isolation 4489329 ^[1] : PCR ^[2] and confirmation

^[1] Replaces Pub. nos. 4405962 and 4426510.^[2] Refer to the specified user guide for the real-time PCR instruments and analysis software included in the Performance Tested MethodsSM-certified workflow.

Table 3 AOAC® Performance Tested MethodsSM-certified workflows for selective agar detection

Workflow overview		
Species	<i>Listeria</i> spp.	<i>Listeria</i> spp.
Performance Tested Methods SM matrices	Food matrices	Food matrices
Sample preparation	Pathatrix® <i>Listeria</i> spp. Kit (Cat. no. APL50; individual samples)	Pathatrix® 5-Pooling <i>Listeria</i> spp. Kit (Cat. no. APL250P; pooled samples)
Sample preparation workflow	Immunomagnetic separation	Immunomagnetic separation
Detection	Selective agar	Selective agar
User guide (Pub. no.)	MAN0007093	MAN0006909

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

For testing of Food and Environmental samples only.

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

DISCLAIMER

LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

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.

©2014 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. AOAC is a registered trademark and Performance Tested Methods is a service mark of AOAC International. AFNOR and NF Validation are trademarks of Association Française de Normalisation (AFNOR).

For support visit lifetechnologies.com/support or email techsupport@lifetech.com

lifetechnologies.com