

## Workflows for Detection of *E. coli* in Food Samples

Pub. No. MAN0009419 Rev. C.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.

### NF VALIDATION™-certified workflows (AFNOR™) for PCR detection

Workflow overview			
<b>Species</b>	<i>E. coli</i> O157:H7	<i>E. coli</i> O157:H7	<i>E. coli</i> O157:H7
<b>NF VALIDATION™ categories</b>	Raw beef meat and raw produce	Raw beef meat and raw produce	Raw beef meat
<b>Nucleic acid (NA) isolation (sample prep)</b>	Applied Biosystems™ PrepSEQ™ Nucleic Acid Extraction Kit (Cat. Nos. 4480466, 4428176)	Applied Biosystems™ PrepSEQ™ Rapid Spin Sample Preparation Kit (Cat. Nos. 4407760, 4426714, 4413269, 4426715)	Thermo Scientific™ SureTect™ <i>E. coli</i> O157:H7 PCR Assay (Cat. No. PT0400A)
<b>NA isolation workflow</b>	Automated	Spin column	Direct lysis
<b>PCR detection</b>	Applied Biosystems™ MicroSEQ™ <i>E. coli</i> O157:H7 Detection Kit (Cat. No. 4427409)	Applied Biosystems™ MicroSEQ™ <i>E. coli</i> O157:H7 Detection Kit (Cat. No. 4427409)	Thermo Scientific™ SureTect™ <i>E. coli</i> O157:H7 PCR Assay (Cat. No. PT0400A)
<b>User guide (Pub. No.)</b>	4426513: Sample enrichment and NA isolation 4426511: PCR <sup>[1]</sup> and confirmation	4426519: Sample enrichment and NA isolation 4426511: PCR <sup>[1]</sup> and confirmation	D12162

<sup>[1]</sup> Refer to the specified user guide for the real-time PCR instruments and analysis software included in the NF VALIDATION™-certified workflow.

### AOAC™ Performance Tested Methods<sup>SM</sup> - certified workflows for PCR detection

Workflow overview					
<b>Species</b>	<i>E. coli</i> O157:H7	<i>E. coli</i> O157:H7	<i>E. coli</i> O157:H7 and non-O157 STECs <sup>[1]</sup>	<i>E. coli</i> O157:H7 and non-O157 STECs <sup>[1]</sup>	<i>E. coli</i> O157:H7
<b>Performance Tested Methods<sup>SM</sup> matrices</b>	Meat products, leafy green products, and juices	Meat products, leafy green products, and juices	Raw beef	Raw beef	Raw beef, leafy green products, and juices
<b>Nucleic acid (NA) isolation (sample prep)</b>	Applied Biosystems™ PrepSEQ™ Nucleic Acid Extraction Kit (Cat. Nos. 4480466, 4428176)	Applied Biosystems™ PrepSEQ™ Rapid Spin Sample Preparation Kit (Cat. Nos. 4407760, 4426714, 4413269, 4426715)	Applied Biosystems™ PrepSEQ™ Nucleic Acid Extraction Kit (Cat. Nos. 4480466, 4428176)	Applied Biosystems™ PrepSEQ™ Rapid Spin Sample Preparation Kit – Extra Clean with Proteinase K (Cat. Nos. 4426715)	Thermo Scientific™ SureTect™ <i>E. coli</i> O157:H7 PCR Assay (Cat. No. PT0400A)
<b>NA isolation workflow</b>	Automated	Spin column	Automated	Spin column	Direct lysis
<b>PCR detection</b>	Applied Biosystems™ MicroSEQ™ <i>E. coli</i> O157:H7 Detection Kit (Cat. No. 4427409)	Applied Biosystems™ MicroSEQ™ <i>E. coli</i> O157:H7 Detection Kit (Cat. No. 4427409)	Applied Biosystems™ RapidFinder™ STEC Screening Assay (Cat. No. 4476886) Applied Biosystems™ RapidFinder™ STEC Confirmation Assay (Cat. No. 4476901)	Applied Biosystems™ RapidFinder™ STEC Screening Assay (Cat. No. 4476886) Applied Biosystems™ RapidFinder™ STEC Confirmation Assay (Cat. No. 4476901)	Thermo Scientific™ SureTect™ <i>E. coli</i> O157:H7 PCR Assay (Cat. No. PT0400A)
<b>User guide (Pub. No.)</b>	4426513: Sample enrichment and NA isolation 4426511: PCR <sup>[2]</sup> and confirmation	4426519: Sample enrichment and NA isolation 4426511: PCR <sup>[2]</sup> and confirmation	4477013	MAN0015934	D12162

<sup>[1]</sup> Serogroups O26, O45, O103, O111, O121, and O145.

<sup>[2]</sup> Refer to the specified user guide for the real-time PCR instruments and analysis software included in the Performance Tested Methods<sup>SM</sup>-certified workflow.

## AOAC™ Performance Tested Methods<sup>SM</sup>-certified workflows for selective agar detection

Workflow overview			
<b>Species</b>	<i>E. coli</i> O157:H7	<i>E. coli</i> O157:H7	<i>E. coli</i> O157:H7
<b>Performance Tested Methods<sup>SM</sup> matrices</b>	Raw ground beef	Raw ground beef	Raw ground beef
<b>Sample preparation</b>	Applied Biosystems™ Pathatrix™ <i>E. coli</i> O157:H7 Kit (Cat. No. APE50SD; individual samples)	Applied Biosystems™ Pathatrix™ 5-Pooling <i>E. coli</i> O157:H7 Kit (Cat. No. APE250SDP; pooled samples)	Applied Biosystems™ Pathatrix™ 10-Pooling <i>E. coli</i> O157:H7 Kit (Cat. No. APE500SDP; pooled samples)
<b>Sample preparation workflow</b>	Immunomagnetic separation	Immunomagnetic separation	Immunomagnetic separation
<b>Detection</b>	Selective agar	Selective agar	Selective agar
<b>User guide (Pub. No.)</b>	MAN0007092	MAN0007090	MAN0007091

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.

**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. AOAC is a trademark and Performance Tested Methods is a service mark of AOAC International. AFNOR and NF VALIDATION are trademarks of Association Française de Normalisation (AFNOR).