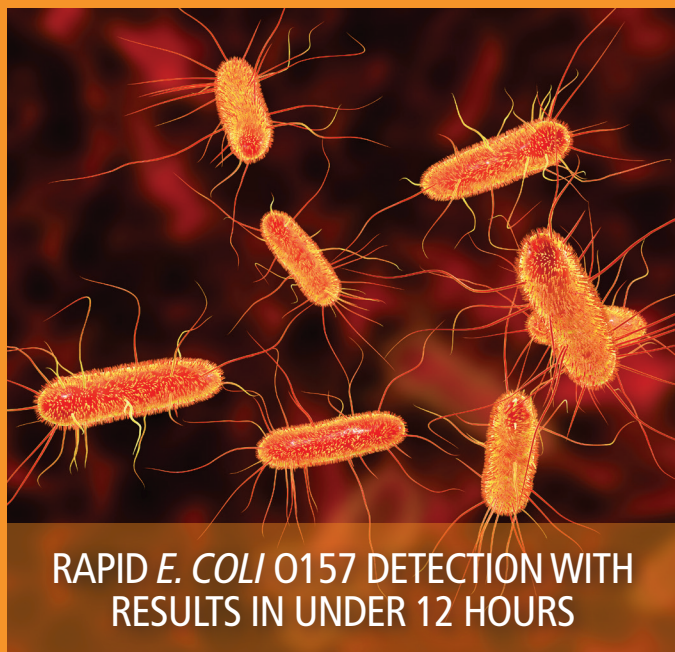


RAPID E. COLI 0157 TESTING



Solus One *E. coli* 0157
Assay and Automated Workflow Package



Rapid Testing and Enhanced Throughput for Raw Beef Samples

When Time to Results is Critical

In a fast-paced food testing environment, it's critical to process samples quickly and efficiently. Providing rapid, accurate results allows production facilities to act quickly if *E. coli* O157 is detected.

You also need the flexibility to cope with changeable sample volumes to fit with the natural fluctuations in supply and demand. Your lab must be able to maximize sample testing throughput without additional resources. That's the kind of capability our Solus One *E. coli* O157 detection system gives you.

A highly efficient assay for same-day detection of *E. coli* O157 in raw meat samples, our Solus One *E. coli* O157 provides results from a single enrichment step in less than 12 hours. The Solus One *E. coli* O157 assay extends our pathogen detection systems' range, which comprises specifically designed immunoassay kits, dedicated selective enrichment media, and automated liquid handling.

The Platform of Choice for *E. coli* O157 Detection

Our Solus pathogen systems have been specifically developed around the constraints of food testing environments. We employ proven, industry-standard immunoassay technologies enhanced to be as sensitive and robust as other pathogen screening methods. Solus assay kits are manufactured in our own production facility using tightly-controlled protocols within our ISO 9001-accredited quality system. Solus pathogen detection systems are validated by independent laboratories to the highest standards, are easy to use, and can be adapted for use by labs of all sizes.



Solus One *E. coli* O157 Benefits

With Solus One *E. coli* O157, you get a reliable and robust *E. coli* O157 detection method, yielding results within 12 hours. And you have the option to run assays manually or by automation. By choosing our highly efficient automated processing, you can achieve high-sample throughput on a single machine, with negative or presumptive positive results in around two hours, post enrichment. Plus, you'll benefit from:

- The ability to cope with fluctuating sample volumes and the capacity to grow
- Excellent sample traceability when using the onboard barcode facility
- A small instrument footprint that maximizes bench space
- Compact product packaging that reduces cold storage space and packaging waste
- An AOAC PTM-certified method



HIGH THROUGHPUT
INCREASES YOUR EFFICIENCY



EASY TO IMPLEMENT,
EASY TO USE

Solus One *E. coli* O157 Automation

Solus One *E. coli* O157 ELISA automated process is carried out on a Dynex DS2 instrument. A single DS2 processes two 96-well plates simultaneously, generating results for up to 558 samples in an eight-hour shift.

Once the instrument is loaded, the technician can walk away for up to two hours during each run, providing time to carry out other essential laboratory tasks.



Dynex DS2

Ordering information

Product Code	Product Description	Pack Size
EC1-0096	Solus One <i>E. coli</i> O157	96 wells
EC1-0480	Solus One <i>E. coli</i> O157	480 wells
MED017-10KG	Buffered Peptone Water	10 KG
SALSUPP22.5	Solus One Supplement	10 vials – for up to 45 liters media
SALSUPP112.5	Solus One Supplement	10 vials – for up to 225 liters media
SALSUPP200	Solus One Supplement	10 vials – for up to 400 liters media



Solus One *E. coli* O157 is certified by the AOAC Performance Tested Method (PTM) program certificate number 112001 for the detection of *E. coli* O157.

Test Methodology

Solus Automated Process



STEP 1

Add warm supplemented BPW to sample and mix

Incubation:

41.5 °C, 10-12 hrs.

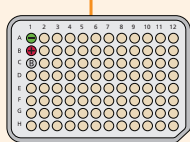


STEP 2

Transfer 1 mL culture to boiling tube

Heat:

85-100 °C, 15-20 min.

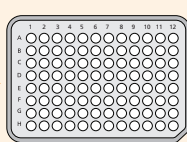


STEP 3

Pipette 100 µL sample into wells (plus controls)

Incubation:

37 °C, 30 min.

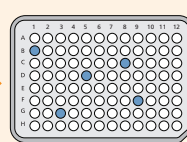


STEP 4

Wash plate and add 100 µL of conjugate

Incubation:

37 °C, 30 min.

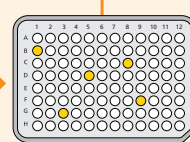


STEP 5

Wash plate and add 100 µL of substrate

Incubation:

18-25 °C, 30 min.



STEP 6

Add 100 µL of stop solution

Read Plate:

450 nm.

Negative Control
 Positive Control
 Blank Well
 Sample wells
 Positive (After Substrate Addition)
 Positive (After Stop Solution Addition)

For sales and support contacts, please visit www.solusscientific.com

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2020, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.