



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

PMA Real-Time PCR Bacterial Viability Kit – *E. coli* O157:H7 (Z3276)

A kit designed to perform viability PCR on *E. coli* O157:H7. It contains a viability PCR dye (PMAxx™ or PMA) and primers to amplify the Z3276 gene.



Product attributes

Viability dye	PMA, PMAxx™
Storage Conditions	Store at -10 to -35 °C
Assay type/options	Live/dead discrimination, Viability PCR
Detection method/readout	PCR/qPCR

Product Description

This kit is designed to perform viability PCR on *E. coli* O157:H7. It contains a viability PCR dye (PMA or PMAxx™) and primers to amplify the Z3276 gene.

About this kit

E. coli O157:H7 is a pathogenic strain of *Escherichia coli* bacteria. Often a food-borne pathogen, it can cause digestive illness. PMA-based viability PCR to detect *E. coli* O157:H7 has been reported using the primers provided in the kit, and ORF Z3276 was shown to be specific for strain O157:H7. These primers have been validated at Biotium for real-time qPCR using Forget-Me-Not™ EvaGreen® qPCR Master Mix. For pan-*E. coli* primers to detect other strains of *E. coli* besides O157:H7, please see kit [31050](#).

To learn more about the advantages of determining bacterial viability using viability PCR, visit the [Viability PCR Technology Page](#).

Strain-Specific Bacterial Viability PCR Kits

Bacteria Strain	Gene name	Kit with PMA (catalog #)	Kit with PMAxx™ (catalog #)
<i>Salmonella enterica</i>	invA	31033	31033-X
<i>Mycobacterium tuberculosis</i>	groEL2	31034	N/A
<i>Staphylococcus aureus</i>	nuc	31035	N/A
MRSA	mecA	31036	N/A
<i>E. coli</i> O157:H7	Z3276	31037	31037-X
<i>E. coli</i>	uidA	31050	31050-X
<i>Listeria monocytogenes</i>	hly	31051	31051-X
<i>Legionella pneumophila</i>	mip	31053	N/A

Don't see your favorite strain? Let us know at techsupport@biotium.com.

PMAxx technology is covered by granted and/or pending US and international patents.

References

Download [PMA and PMAxx™ References](#) and a list of [PMA and PMAxx™ Validated Bacterial Strains](#).

This datasheet was generated on September 1, 2026 at 10:27:54 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/pma-real-time-pcr-bacterial-viability-kit-e-coli-0157h7-z3276/>