

# PMA Real-Time PCR Bacterial Viability Kit – E. coli (uidA)

 Viability dye
 PMA, PMAxx™

 Storage Conditions
 Store at -10 to -35 °C

 Assay type/options
 Live/dead discrimination, Viability PCR

PCR/qPCR, HRM®

Call us: 800-304-5357

Product attributes

Detection method/readout

A kit designed to perform viability PCR on E. coli. It contains a viability dye (PMAxx<sup>TM</sup> or PMA) and primers to amplify the uidA gene.

## **Product Description**

This kit is designed to perform viability PCR on *E. coli*. It contains a viability PCR dye (PMA or PMAxx<sup>TM</sup>) and primers to amplify the uidA gene.

#### About this kit

Escherichia coli is a species of bacteria that is commonly used in laboratories. Some strains of *E. coli* are food-borne pathogens and can cause digestive illness. For specific detection of the pathogenic *E. coli* strain O157:H7, please see kit number 31037. PCR to amplify the gene uidA has been published and shown to be highly specific for *E. coli*. The primers provided in the kit have been validated at Biotium for real-time qPCR using Forget-Me-Not<sup>TM</sup> EvaGreen® qPCR Master Mix.

To learn more about the advantages of determining bacterial viability using viability PCR, visit the <u>Viability PCR Technology Page</u>.

### Strain-Specific Bacterial Viability PCR Kits

Bacteria Strain	Gene name	Kit with PMA (catalog #)	Kit with PMAxx™ (catalog #)
Salmonella enterica	invA	<u>31033</u>	31033-X
Mycobacterium tuberculosis	groEL2	<u>31034</u>	N/A
Staphylococcus aureus	nuc	<u>31035</u>	N/A
MRSA	mecA	<u>31036</u>	N/A
E. coli O157:H7	Z3276	<u>31037</u>	31037-X
E. coli	uidA	<u>31050</u>	31050-X
Listeria monocytogenes	hly	<u>31051</u>	<u>31051-X</u>
Legionella pneumophila	mip	<u>31053</u>	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 November 23, 2025 at 12:36:43 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/pma-real-time-pcr-bacterial-viability-kit-e-coli-uida/">https://biotium.com/product/pma-real-time-pcr-bacterial-viability-kit-e-coli-uida/</a>