



Glowing products for science

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

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

A kit designed to perform viability PCR on *E. coli*. It contains a viability dye (PMAxx™ or PMA) and primers to amplify the uidA 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, HRM®

### Product Description

This kit is designed to perform viability PCR on *E. coli*. It contains a viability PCR dye (PMA or PMAxx™) 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™ EvaGreen® qPCR Master Mix.

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	<a href="#">31033</a>	<a href="#">31033-X</a>
<i>Mycobacterium tuberculosis</i>	groEL2	<a href="#">31034</a>	N/A
<i>Staphylococcus aureus</i>	nuc	<a href="#">31035</a>	N/A
MRSA	mecA	<a href="#">31036</a>	N/A
<i>E. coli</i> O157:H7	Z3276	<a href="#">31037</a>	<a href="#">31037-X</a>
<i>E. coli</i>	uidA	<a href="#">31050</a>	<a href="#">31050-X</a>
<i>Listeria monocytogenes</i>	hly	<a href="#">31051</a>	<a href="#">31051-X</a>
<i>Legionella pneumophila</i>	mip	<a href="#">31053</a>	N/A

Don't see your favorite strain? Let us know at [techsupport@biotium.com](mailto: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 January 8, 2026 at 08:30:02 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-uidA/>