



Glowing products for science

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

PMA Real-Time PCR Bacterial Viability Kit – *Listeria monocytogenes* (hly)

A kit designed to perform viability PCR on *Listeria monocytogenes*. It contains a viability dye (PMAxx™ or PMA) and primers to amplify the hly 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 *Listeria monocytogenes*. It contains a viability PCR dye (PMA or PMAxx™) and primers to amplify the hly gene.

About this kit

Listeria monocytogenes is a food-borne pathogen that can cause digestive illness and fetal harm. The gene hly encodes the Listeria Lysin O (LLO) protein that is specific for *Listeria*. PMA-PCR using these primers for viability PCR in *Listeria* has been published. In addition, these primers 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 | 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 January 8, 2026 at 08:25:29 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/pma-real-time-pcr-bacterial-viability-kit-listeria-monocytogenes-hly/>