



Glowing products for science

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

PMA Real-Time PCR Bacterial Viability Kit – *Legionella pneumophila* (mip)

A kit designed to perform viability PCR on *Legionella pneumophila*. It contains the viability PCR dye PMA and primers to amplify the mip gene.



Product attributes

| | |
|--------------------------|---|
| Viability dye | PMA |
| 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 *Legionella pneumophila*. It contains the viability PCR dye PMA and primers to amplify the mip gene.

About this kit

Legionella pneumophila is a pathogenic species of gram-negative bacteria. It can infect the lungs and cause Legionnaire's Disease. *Legionella pneumophila* is sometimes detected in water sources such as cooling towers and swimming pools. PCR to amplify the gene mip has been published and is highly specific for Legionella. 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 | 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 12:36:46 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/pma-real-time-pcr-bacterial-viability-kit-legionella-pneumophila-mip/>