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	DCD/aDCD HDM®		

Call us: 800-304-5357 Email: btinfo@biotium.com

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-NotTM 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	31034	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>	<u>31037-X</u>
E. coli	uidA	<u>31050</u>	31050-X
Listeria monocytogenes	hly	<u>31051</u>	31051-X
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 5, 2025 at 09:14:25 PM. 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/