

# 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<sup>TM</sup> or PMA) and primers to amplify the hly gene.



#### **Product attributes**

Call us: 800-304-5357

Viability dye	PMA, PMAxx <sup>TM</sup>	
Storage Conditions	Store at -10 to -35 °C	
Assay type/options	Live/dead discrimination, Viability PCR	
Detection method/readout	PCR/cPCR HRM®	

## **Product Description**

This kit is designed to perform viability PCR on *Listeria monocytogenes*. It contains a viability PCR dye (PMA or PMAxx<sup>TM</sup>) 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 <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<sup>TM</sup> References and a list of PMA and PMAxx<sup>TM</sup> Validated Bacterial Strains.

This datasheet was generated on November 23, 2025 at 01:57:25 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-listeria-monocytogenes-hly/">https://biotium.com/product/pma-real-time-pcr-bacterial-viability-kit-listeria-monocytogenes-hly/</a>