

# PMA Real-Time PCR Bacterial Viability Kit – Staphylococcus aureus (nuc)

nietum market proportion of the state of th

A kit designed to perform viability PCR on *Staphylococcus aureus*. It contains the viability PCR dye PMA and primers to amplify the nuc gene.

#### **Product attributes**

Call us: 800-304-5357

Viability dye	PMA		
Storage Conditions	Store at -10 to -35 °C		
Assay type/options	Live/dead discrimination, Viability PCR		
Detection method/readout	PCR/qPCR		

## **Product Description**

This kit is designed to perform viability PCR on *Staphylococcus aureus*. It contains the viability PCR dye PMA and primers to amplify the nuc gene.

#### About this kit

Staphylococcus aureus is an anaerobic gram positive bacterium that can cause of a variety of infections in humans, and methicillin-resistant *S. aureus* is a rising cause of hospital-acquired infections. The primers included in this kit have been tested for amplification of the S. aureus-specific gene *nuc* gene from genomic DNA from both methicillin-susceptible and methicillin-resistant strains of *S. aureus* and have been validated at Biotium for real-time qPCR using Forget-Me-Not<sup>TM</sup> EvaGreen® qPCR Master Mix. Biotium also offers the <u>Staphylococcus aureus (mecA) PMA-PCR kit</u>, which includes primers against the methicillin-resistance gene *mecA*.

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™ References and a list of PMA and PMAxx™ Validated Bacterial Strains.

This datasheet was generated on November 23, 2025 at 12:30:38 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-staphylococcus-aureus-nuc/">https://biotium.com/product/pma-real-time-pcr-bacterial-viability-kit-staphylococcus-aureus-nuc/</a>