

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



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

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™ EvaGreen® qPCR Master Mix. Biotium also offers the [Staphylococcus aureus \(mecA\) PMA-PCR kit](#), which includes primers against the methicillin-resistance gene *mecA*.

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	<a href="#">31033</a>	<a href="#">31033-X</a>
<i>Mycobacterium tuberculosis</i>	groEL2	<a href="#">31034</a>	N/A
<i>Staphylococcus aureus</i>	nuc	<a href="#">31035</a>	N/A
MRSA	mecA	<a href="#">31036</a>	N/A
<i>E. coli</i> O157:H7	Z3276	<a href="#">31037</a>	<a href="#">31037-X</a>
<i>E. coli</i>	uidA	<a href="#">31050</a>	<a href="#">31050-X</a>
<i>Listeria monocytogenes</i>	hly	<a href="#">31051</a>	<a href="#">31051-X</a>
<i>Legionella pneumophila</i>	mip	<a href="#">31053</a>	N/A

Don't see your favorite strain? Let us know at [techsupport@biotium.com](mailto: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 June 13, 2026 at 08:57:46 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/pma-real-time-pcr-bacterial-viability-kit-staphylococcus-aureus-nuc/>