



Glowing products for science

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

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

A kit designed to perform viability PCR on MRSA. It contains the viability PCR dye PMA and primers to amplify the *mecA* 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 Methicillin-resistant *Staphylococcus aureus*. It contains the viability PCR dye PMA and primers to amplify the *mecA* gene.

About this kit

Staphylococcus aureus is an anaerobic gram negative bacterium that can cause a variety of infections in humans, and methicillin-resistant *S. aureus* is a rising cause of hospital-acquired infections. The *mecA* gene confers methicillin-resistance in *S. aureus* and other methicillin-resistant bacterial species. The *mecA* primers included in the kit have been used for detection of methicillin-resistant *S. aureus*, and have been validated for qPCR using Biotium's Forget-Me-Not™ EvaGreen® qPCR Master Mix. Note that the *mecA* primers are not specific to *S. aureus*, because *mecA* is expressed by other methicillin-resistant bacteria. Biotium also offers the [Staphylococcus aureus \(nuc\) PMA-PCR kit](#), which includes primers against the *S. aureus*-specific *nuc* gene, expressed by methicillin-susceptible and methicillin-resistant *S. aureus* strains.

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 December 20, 2025 at 02:20:57 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/pma-real-time-pcr-bacterial-viability-kit-staphylococcus-meca/>