



Glowing products for science

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

## PMA Real-Time PCR Bacterial Viability Kit – *Salmonella enterica* (invA)

A kit designed to perform viability PCR on *Salmonella enterica*. It contains a viability PCR dye (PMAxx™ or PMA) and primers to amplify the invA gene.



### Product attributes

Viability dye	PMA, PMAxx™
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 *Salmonella enterica*. It contains a viability PCR dye (PMA or PMAxx™) and primers to amplify the invA gene.

### About this kit

*Salmonella enterica* is a gram-negative bacteria that causes the food-borne illness salmonellosis. PMA-based viability PCR for *S. enterica* has been reported using the primers provided in the kit. 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 [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 January 7, 2026 at 10:44:39 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/pma-real-time-pcr-bacterial-viability-kit-salmonella-enterica-inva/>