

PMA Real-Time PCR Bacterial Viability Kit – *Mycobacterium tuberculosis* (groEL2)



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

A kit designed to perform viability PCR on *Mycobacterium tuberculosis*. It contains the viability PCR dye PMA and primers to amplify the groEL2 gene.

Product Description

This kit is designed to perform viability PCR on *Mycobacterium tuberculosis*. It contains the viability PCR dye PMA and primers to amplify the groEL2 gene.

About this kit

Mycobacterium tuberculosis is a pathogenic bacteria that infects the lungs and causes the disease tuberculosis. PCR to detect *Mycobacterium tuberculosis* 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. Note: groEL2 primers also amplify other mycobacteria species, but products may be distinguishable by melt curve analysis.

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 June 18, 2026 at 09:17:39 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/pma-real-time-pcr-bacterial-viability-kit-mycobacterium-tuberculosis-groel2/>