

Viability PCR Starter Kits

Kits designed to make it simple to start performing viability PCR with your choice of viability dye (PMA or PMAxxTM).



Product attributes

Call us: 800-304-5357

Viability dye	PMA, PMAxx TM
Bacteria type	All species, Gram negative bacteria
Storage Conditions	Store at -10 to -35 °C
Assay type/options	Live/dead discrimination, Viability PCR
Detection method/readout	PCR/qPCR, HRM®

Product Description

Viability PCR starter kits contain the materials that you need for selective detection of viable cells using either PMA or PMAxx™ viability dye and real-time PCR. The kits contain Forget-Me-Not™ EvaGreen® qPCR Master Mix, and are available with or without Enhancer for gram-negative bacteria. The user will need to supply primers to amplify DNA from their species of interest (see this blog post for primer design considerations).

About this kit

This kit contains reagents sufficient to treat 80 cultures with PMA or PMAxxTM, and perform 200 PCR reactions. There is an option to choose either PMA or PMAxxTM for the viability dye. Original PMA dye has been validated for vPCR in hundreds of publications and a wide variety of bacteria, yeast, and viruses. PMAxxTM is a next-generation vPCR dye designed by Biotium scientists to provide superior live/dead discrimination compared to PMA.

Kit options are available with of PMA Enhnacer for Gram-Negative Bacteria, a reagent that improves live/dead discrimination for gram-negative bacteria (note: Enhancer is not recommended for use with gram-positive bacteria or samples with mixed strains).

To learn more about the advantages of determining microbial or cell viability using viability PCR, visit the <u>Viability PCR Technology Page</u> and see the <u>PMA/PMAxx™ Reference List</u> and <u>Validated Bacterial</u> <u>Strains List</u>.

PMAxx technology is covered by granted and/or pending US and international patents.

References

Find a list of PMA and PMAxx™ references and a list of validated bacterial strains under Supporting Documents.

This datasheet was generated on November 23, 2025 at 10:51:34 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/viability-pcr-starter-kits/