



Glowing products for science

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

Viability PCR Starter Kits

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



Product attributes

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 PMAxx™, and perform 200 PCR reactions. There is an option to choose either PMA or PMAxx™ 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. PMAxx™ 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 Enhancer 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 [Viability PCR Technology Page](#) and see the [PMA/PMAxx™ Reference List](#) and [Validated Bacterial Strains List](#).

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 August 27, 2026 at 03:55:20 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/viability-pcr-starter-kits/>