

## PMA (Propidium Monoazide)

A photo-reactive DNA-binding dye typically used in viability PCR (v-PCR) of microorganisms like bacteria, viruses, and fungi.



## Product Description

PMA (propidium monoazide) is a photo-reactive DNA-binding dye used in viability PCR (v-PCR) of microorganisms like bacteria, viruses, and fungi. PMA is available as a 1 mg lyophilized solid or 20 mM solution in water. Also try PMAxx™ ([40069](#)), a superior alternative to PMA.

- Selectively detect viable cells using qPCR
- Validated in hundreds of publications
- Dead cell specific dye, binds to DNA
- Covalently attaches to DNA after photoactivation
- Available as 1 mg lyophilized solid or 20 mM solution in water
- $\lambda_{\text{Abs}} = 464 \text{ nm}$  (before photolysis);  $\lambda_{\text{Abs}}/\lambda_{\text{Em}} = \sim 510/\sim 610 \text{ nm}$  (with DNA/RNA, after photolysis)

To learn more about the advantages of determining microbial or cell viability using viability PCR, visit the [Viability PCR Technology Page](#).

| Product                            | Catalog Number | Unit Size               | Format            |
|------------------------------------|----------------|-------------------------|-------------------|
| PMA Dye                            | 40013          | 1 mg                    | Lyophilized solid |
| PMA Dye, 20 mM in H <sub>2</sub> O | 40019          | 100 uL (20 mM in water) | Solution          |

## About PMA

PMA dye is a DNA modifier invented by scientists at Biotium. It is a photo-reactive dye that binds to DNA with high affinity. Upon photolysis with visible light, PMA dye becomes covalently attached to DNA. This modified DNA cannot be amplified by PCR. The dye is designed to be cell membrane-impermeable. Thus, in a population of live and dead cells, only dead cells are susceptible to DNA modification due to compromised cell membranes. This unique feature of PMA dye makes it highly useful in selective detection of live bacteria by qPCR.

Since Biotium first developed PMA dye, there have been hundreds of publications on the use of the dye in many sample types including dozens of bacterial strains, biofilms, yeast, fungi, viruses, and eukaryotic cells. It has been used in such applications as food and water safety and environmental testing, and has been used in conjunction with qPCR, NextGen Sequencing (NGS), Sanger sequencing, and Loop-mediated Isothermal Amplification (LAMP).

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

## References

Download list of curated [PMA and PMAxx™ References](#) and a list of [PMA and PMAxx™ Validated Bacterial Strains](#).

This datasheet was generated on December 29, 2025 at 12:15:33 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/pma-propidium-monoazide/>

## Product attributes

|                          |   |
|--------------------------|---|
| Viability dye            | PMA   |
| Molecular weight         | 511   |
| Excitation/Emission      | Abs = 464 nm (before photolysis), $\sim 510/\sim 610 \text{ nm}$ (with DNA/RNA, after photolysis) |
| Assay type/options       | Live/dead discrimination, Viability PCR   |
| Storage Conditions       | 1 mg solid, store at 2 to 8 °C, 100 uL in water, store at -10 to -35 °C                           |
| Detection method/readout | PCR/qPCR  |