



Glowing products for science

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

PMAXx™ Dye, 20 mM in H₂O

Viability PCR dye, a new and improved version of the popular viability dye propidium monoazide (PMA).



Product attributes

Viability dye	PMAXx™
Storage Conditions	Store at -10 to -35 °C
Excitation/Emission	~510/~610 nm (with DNA/RNA, after photolysis), Abs = 464 nm (before photolysis)
Assay type/options	Live/dead discrimination, Viability PCR
Detection method/readout	PCR/qPCR

Product Description

PMAXx™ is a photoreactive DNA-binding dye for viability PCR. It was designed by Biotium scientists to be a superior alternative to the popular viability dye propidium monoazide (PMA).

- Selectively detect viable cells using qPCR
- Better live/dead discrimination of bacteria than PMA
- Dead cell specific dye, binds to DNA
- Covalently attaches to DNA after photoactivation
- 20 mM in water, 100 uL volume
- $\lambda_{Abs} = 464 \text{ nm}$ (before photolysis); $\lambda_{Abs}/\lambda_{Em} = \sim 510/\sim 610 \text{ nm}$ (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](#).

About PMAXx™

PMAXx™ dye is a DNA modifier used for viability PCR, invented by scientists at Biotium. PMAXx™ is a new and improved version of our popular viability dye PMA (propidium monoazide). Like PMA, PMAXx™ is a photo-reactive dye that binds to DNA with high affinity. Upon photolysis with visible light, PMAXx™ dye becomes covalently attached to DNA. This modified DNA cannot be amplified by PCR. PMAXx™ dye is designed to be cell membrane-impermeant. 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 makes PMAXx™ highly useful in selective detection of live bacteria by qPCR.

PMAXx™ was designed by Biotium scientists to be a superior alternative to PMA, and is compatible with all existing PMA protocols, but with superior activity. While PMA is generally effective at differentiating between live and dead bacteria by qPCR, it does not completely eliminate PCR products from dead cell DNA. This could potentially give false positive results. Biotium's new dye PMAXx™ is much more effective at eliminating PCR amplification of dead cell DNA, and therefore provides the best discrimination between live and dead bacteria.

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 June 14, 2026 at 05:17:13 PM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/pmaxx-20-mm-in-h2o/>