

# Ethidium Homodimer III (EthD-III), 1 mM in DMSO

A red fluorescent dead cell stain for bacteria and mammalian cells



## Product Description

Ethidium Homodimer III, also known as EthD-III, is a red fluorescent dead cell stain for bacteria and mammalian cells. It is a cell membrane-impermeant nucleic acid dye that stains only dead cells with damaged cell membranes. Ethidium Homodimer III was developed by Biotium as a superior alternative to Ethidium Homodimer I. The absorption and emission spectra are similar, but EthD-III is 45% brighter.

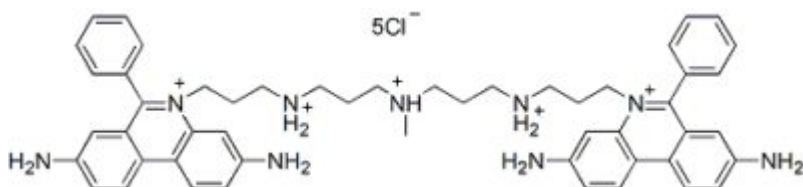
Ethidium Homodimer III is membrane impermeant and useful for selectively staining dead cells with damaged plasma membranes. It also is a component of several combination viability assay kits for detecting both live and dead cells in the same population, such as our Viability/Cytotoxicity Assay Kit for Animal Cells (Cat. No. [30002](#)) and our Viability/Cytotoxicity Assay Kit for Bacteria (Cat. No. [30027](#)).

We also offer Ethidium Homodimer III in solid form (catalog no. [40050](#)).

EthD-III is dead cell specific in all cell types, including mammalian cells, bacteria and yeast. See our [Cellular Stains Table](#) for more information on how our dyes stain various organisms.

### Product Features:

- $\lambda_{Ex}/\lambda_{Em}$  (with DNA) = 522/593 nm\* (EthD-III also has a large absorption peak at 279 nm)
- 1 mM solution in DMSO
- Store at 4 °C and protect from light
- MW: ~1000



This datasheet was generated on January 7, 2026 at 12:21:26 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/ethidium-homodimer-iii-1mm-in-dmsoethd-iii-in-dmso/>

## Product attributes

Probe cellular localization	Nucleus & cytoplasm
For live or fixed cells	For live/intact cells
Assay type/options	Real-time imaging
Cell permeability	Membrane impermeant
Apoptosis/viability marker	Dead cell stain
Colors	Red
Excitation/Emission	279, 522/593 nm (with DNA)