

Ethidium Homodimer III (EthD-III)

A red fluorescent dead cell stain for bacteria and mammalian cells that is brighter than Ethidium Homodimer I.



Product Description

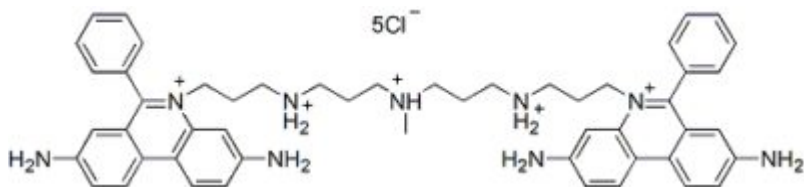
Ethidium Homodimer III, also known as EthD-III, is a red fluorescent dead cell stain for bacteria and mammalian cells. It is a cell membrane-impermeant nucleic acid dye that stains only dead cells with damaged cell membranes. Ethidium Homodimer III was developed by Biotium as a superior alternative to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

- Dead cell-specific DNA stain
- Can be used in mammalian, yeast, and bacteria cells
- 45% brighter than Ethidium Homodimer I
- Provided as a red solid

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, 1 mM solution in DMSO (catalog no. [40051](#)).

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 attributes

| | |
|-----------------------------|--|
| Probe cellular localization | Nucleus & cytoplasm |
| Assay type/options | Real-time imaging |
| Cell permeability | Membrane impermeant |
| Apoptosis/viability marker | Dead cell stain |
| Colors | Red |
| Excitation/Emission | 279, 532/625 nm (with DNA) |
| Storage Conditions | Store at 2 to 8 °C, Protect from light |
| Molecular weight of antigen | ~1000 |

This datasheet was generated on June 21, 2026 at 03:31:19 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/ethidium-homodimer-iii-ethd-iii/>