

www.biotium.com

Revised: August 5, 2020

Product Information

Ethidium Homodimer III (EthD-III)

Catalog Number

40050: 1 mg solid 40051: 1 mM in DMSO

Storage and Handling

The solid is soluble in DMSO, methanol, or water. Store the solid and the solution at 4°C, protected from light. Product is stable for at least 5 years from date of receipt when stored as recommended. Ethidium Homodimer III dye binds to nucleic acids. The mutagenicity or toxicity of EthD-III is currently unknown. It should be handled using universal laboratory safety precautions.

Spectral Properties

Ex/Em: 532/625 nm* (with DNA) *Ethidium Homodimer III also has a strong UV absorbance peak at 279 nm

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 only stains 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 useful for detecting dead bacteria or mammalian cells by fluorescence microscopy or flow cytometry. It can be used in combination with cell permeable stains (such as: Hoechst, DMAO, or BactoView[™] Live Green for bacteria; and Hoechst, DAPI, Calcein AM, or ViaFluor® SE for mammalian cells) to simultaneously label live and dead cells.

Ethidium Homodimer III is not fixable. If a mixture of live and stained dead cells is fixed, dye will leak from the dead cells into the live cells. If you are looking for a fixable, nuclear dead cell stain, try Live-or-Dye NucFix™ Red (32010).

We also offer Ethidium Homodimer III in several combination kits for added convenience, including the Viability/Cytotoxicity Assay Kit for Bacteria (30027) and the Viability/Cytotoxicity Assay Kit for Animal Cells (30002). For more information visit our website.



Figure 1. Normalized absorption and emission of Ethidium Homodimer III

General Staining Protocol

This protocol is a general guideline for staining mammalian tissue culture cells and laboratory bacteria strains. Optimization may be needed for other sample types.

Note: EthD-III selectively stains dead cells based on membrane integrity, and must be used on unfixed cells for live/dead discrimination. If used on cells after fixation, it will stain nuclei and cytoplasm of all cells.

- 1. Grow cells in the appropriate growth medium and growth conditions.
- 2. Optional positive control (dead) cells preparation: For bacteria, incubate cells at 90°C for 5 minutes and allow to cool to room temperature. For adherent mammalian cells, to kill a subset of cells incubate with 15% ethanol in buffer or medium for 10 minutes at room temperature then wash. To prepare a uniformly killed mammalian cell population, incubate at 56°C for 45 minutes and then cool.
- If desired, collect the cells by centrifugation (or wash adherent cells) and resuspend in a buffer for staining. EthD-III can stain cells in growth medium, as well as PBS, HBSS, and 150 mM NaCl.
- Add dye to the cells at a final concentration of 2.5-5 uM (for mammalian cells) or 5 uM (bacterial cells) and mix well. Dye concentration may need to be optimized for different cell or sample types. Other stains may be added simultaneously.
- 5. Incubate at room temperature or 37°C for 15-30 minutes, in the dark.
- 6. Optional wash: Collect cells by centrifugation (or wash adherent cells) and resuspend in fresh buffer of your choice.

Note: Ethidium Homodimer III is not fixable. If a mixture of live and dead cells are fixed after staining, dye will leak from the dead cells and into the live cells.

 For fluorescence microscopy, image cells with Cy®3 or Texas Red® bandpass filter sets. For flow cytometry, detect cells in the PE channel.

Related Products

Catalog number	Product
40101	BactoView™ Live Red
40102	BactoView™ Live Green
30002	Viability/Cytotoxicity Assay Kit for Animal Live & Dead Cells
30066	Apoptotic, Necrotic & Healthy Cells Quantitation Kit Plus
30027	Viability/Cytotoxicity Assay Kit for Bacteria Live and Dead Cells
32001	Bacterial Viability and Gram Stain Kit
32010	Live-or-Dye NucFix™ Red
32002- 32009	Live-or-Dye™ Fixable Viability Staining Kits
40069	PMAxx [™] Dye for Viability PCR, 20 mM in Water
30068	ViaFluor® 405 SE Cell Proliferation Kit
30086	ViaFluor® 488 SE Cell Proliferation Kit

Please visit our website at www.biotium.com for information on our life science research products, including kits for cell biology and microbiology research.

Texas Red is a registered trademark of Thermo Fisher. Cy is a registered trademark of GE Healthcare.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.