

EthD-2 [Ethidium Homodimer-2]

Catalog number: 17782

Unit size: 1 mg

Product Details

Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Solid
Molecular Weight	1292.71
Soluble In	DMSO

Applications

Ethidium Homodimer-2 (EthD-2) is usually used as a cell-impermeant viability indicator. It is a high-affinity nucleic acid stain that is weakly fluorescent until bound to DNA and emits red fluorescence (excitation/emission maxima ~535/624). It is often used together with Calcein AM. The viability of cells can be readily assessed by staining with calcein acetoxymethyl ester (Calcein AM) and EthD-2. The live cells (green, stained with Calcein AM) and dead cells (red, stained with EthD-2) can be easily imaged and even quantified from the images using the ImageJ program. It might also be combined with our other cell viability stains such as the superior Calcein Ultra Blue, Calcein Ultra Green, Cytocalcein, Cytotrace dyes and CytoTell dyes. AAT Bioquest offers the largest collection of fluorescent stains for monitoring cell viabilities.