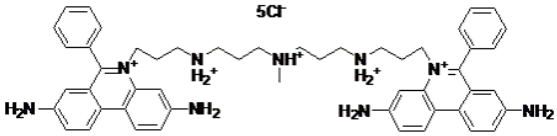


Dead cell stain (also known as EthD-1, or EtDi)

Instruction Manual

Catalog Number	PK-CA707-40050
Description	Ethidium Homodimer III was developed as an alternative to Ethidium Homodimer I. It has absorption and emission spectra similar to those of Ethidium Homodimer I. However, the dye stains DNA 70% brighter than Ethidium Homodimer I. The dye is also a component of PromoKine's Apoptotic/Necrotic Cell Detection Kit and Apoptotic/Necrotic/Healthy Cell Detection Kit for detecting apoptotic, necrotic (and healthy) cells in the same population (PK-CA707-30017 & PK-CA707-30018).
Quantity	1 mg
Excitation / Emission Maxima	$\lambda_{ex}\backslash\lambda_{em}$ (DNA-bound) = ~530/~620 nm; $\lambda_{ex}\backslash\lambda_{em}$ (no DNA) = 493/- nm Extinction coefficient: $\geq 8,500$
Molecular Structure	
Molecular Weight / Molecular Formula	~1000 Da; NA
Purity	>90% (as determined by HPLC)
Appearance / Formulation / Solubility	Red solid; soluble in DMSO, MeOH, or water.
Storage & Stability	Stored desiccated at 4°C. Protect from light, especially when in solution.
Applications	Nucleus stain and dead cell stain (nucleic acid stain)
References	1. Mol Cancer Res 6, 965 (2008). 2. J Pharmacol Exp Therap 332, 738 (2010).
Caution	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.