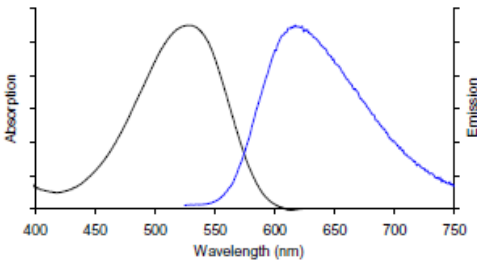


Instruction Manual

Catalog Number	PK-CA707-40051
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 40% 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). It is also included in PromoKine's Live/Dead Cell Staining Kit II PK-CA707-30002).
Quantity	200 μ l
Excitation / Emission Maxima	$\lambda_{ex} \backslash \lambda_{em}$ (DNA-bound) = ~530/~620 nm; $\lambda_{ex} \backslash \lambda_{em}$ (no DNA) = 493/- nm Extinction coefficient: $\geq 8,500$
	 <p>Absorption and emission of EthD-III bound to DNA.</p>
Molecular Structure	proprietary
Molecular Weight / Molecular Formula	~1000 Da; Structure: proprietary
Purity	>90% (as determined by TLC)
Appearance / Formulation / Solubility	1 mM red solution in DMSO, ready-to-use.
Storage & Stability	Stored at 4°C. Protect from light.
Applications	Nucleus and dead cell stain (nucleic acid stain)
References	NA
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.