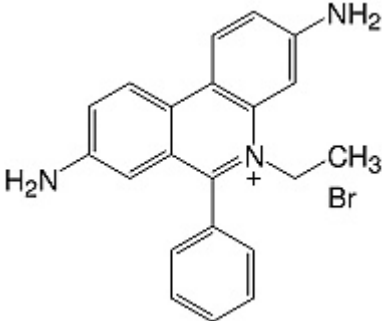


Instruction Manual

Catalog Number	PK-CA707-40042
Description	Because of its toxicity as a potent mutagen, we offer EB in H ₂ O as a safer alternative to the more hazardous powdery form, which is more likely for one to be in contact with. EB is an intercalating DNA-binding dye with little sequence preference. Once bound to nucleic acids, the fluorescence of the dye is enhanced by more than 10 times while the excitation maximum is redshifted by 30-40 nm and emission is blue-shifted by ~15 nm. EB is widely used as a nucleic acid gel stain.
Quantity	10 ml (10 mg/ml solution in H ₂ O)
Excitation / Emission Maxima	$\lambda_{ex} \backslash \lambda_{em} = 518/605$ nm (DNA)
Molecular Structure	
Molecular Weight / Molecular Formula	394.31 g/mol Da; C ₂₁ H ₂₀ BrN ₃
Purity	≥95% (as determined by HPLC)
Appearance / Formulation / Solubility	Red solution. At least 10 mg/ml in H ₂ O.
Storage & Stability	Store desiccated at 4°C. Protect from light!
Applications	Cell stain. Can be used as counterstain for cell nuclei in histological and cytological analysis and in staining of DNA after electrophoresis.
References	<ol style="list-style-type: none"> 1) J.Mol. Biol. 13, 269(1965) 2) Methods Cell Biol. 30, 417(1989) 3) Chromatographia 29, 167(1990).
Caution	Ethidium bromide is a powerful mutagen and moderately toxic. Gloves should be worn when working with solutions that contain this dye. 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, seek medical advice immediately.

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