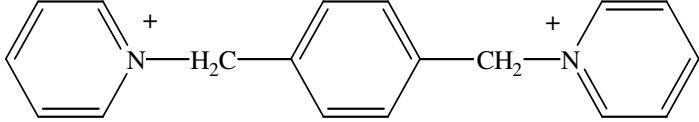


p-xylene-bis-pyridinium bromide

Instruction Manual

Catalog Number	PK-CA707-80012
Description	A positively charged quencher that is often used with ANTS to study membrane fusion or permeability. See information for ANTS (Cat.No. PK-CA707-90010).
Quantity	500 mg
Excitation / Emission Maxima	Absorption at 259 nm; Extinction coefficient = 8,800
Molecular Structure	<p style="text-align: center;">2Br^-</p> 
Molecular Weight / Molecular Formula	422 Da; $\text{C}_{18}\text{H}_{18}\text{Br}_2\text{N}_2$
Purity	>98% (as determined by TLC)
Appearance / Formulation / Solubility	Off-white solid; soluble in water.
Storage & Stability	Store desiccated at $\leq 4^\circ\text{C}$. Protect from light, especially when in solution.
Applications	See Description
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.