

46117 Landing Parkway, Fremont, CA 94538 U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352

www.biotium.com

## **PRODUCT AND SAFETY DATA SHEET**

**PRODUCT NAME:** 

**DPX** (p-xylene-bis-pyridinium bromide)

CATALOG #:

80012

 $2Br^{-}$ 

<b>PROPERTIES:</b>	
Color & Form	Off-white solid
Purity	$\geq$ 98% by TLC
Solubility	Soluble in water
Absorption	259 nm
Extinction	
Coefficient	8,800
STORAGE AND	Store desiccated at 4 °C.
HANDLING:	
SHELF LIFE	Minimum 5 years at -20 °C
APPLICATION:	A monitorial allowed quanches that is often used with ANTE to study membrane fusion
APPLICATION:	A positively charged quencher that is often used with ANTS to study membrane fusion or permeability. See information on <b>ANTS</b> (90010).
	or permeability. See information on AINIS (90010).
TOXICITY:	Unknown.
IOMCIII.	Clikiowi.
FIRST AID:	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

**Disclaimer:** Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.

move individual to fresh air and seek medical advice immediately.

plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing,