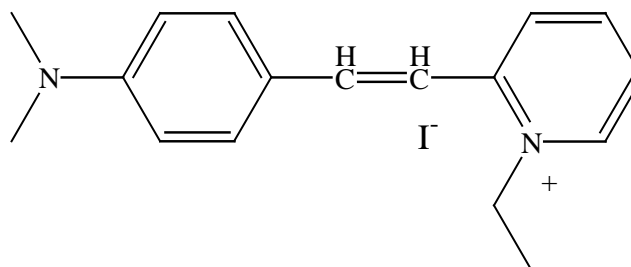


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: **DASPEI** (2-(4-(dimethylamino)styryl)-N-ethylpyridinium iodide)**CATALOG NUMBER:** **70018****MOLECULAR INFORMATION:** C₁₇H₂₁IN₂
Mwt: 380
[3785-01-1]**PROPERTIES:**

Color & Form	Red Solid
Purity	≥ 99% by HPLC
Solubility	Soluble in DMF
$\lambda_{ex}/\lambda_{em}$	461/589 nm (MeOH)
Extinction	39,000
Coefficient (ϵ)	

STORAGE AND HANDLING:

Store at 4 °C and protect from light, especially when in solution.

APPLICATION:

DASPEI is used to stain mitochondria in live cells.

Ref.: 1) *Biochem Biophys Res Comm* **261**, 452 (1999); 2) *J Exp Biol* **202**, 3485 (1999); 3) *J Biol Chem* **272**, 21964 (1997); 4) *Biochim Biophys Acta* **1278**, 205 (1996); 5) *Microscopy Res Technique* **33**, 496(1996).**TOXICITY:**

Unknown

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 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.

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.