

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: di-8-ANEPPS

CATALOGUE #: 61012

 $^{-}O_{3}S(H_{2}C)_{3} - N$ $CH = CH - (CH_{2})_{7}CH_{3}]_{2}$

PROPERTIES:

Color & Form Purity Solubility Absorption/Emission Extinction Coefficient	Orange red solid. 99% by HPLC Soluble in DMSO or EtOH 498 nm / 713 nm (MeOH) 44,000
STORAGE AND HANDLING:	Store desiccated at \leq 4°C and protect from light especially in solution.
APPLICATION:	ANEPPS dyes are fast responding membrane potential dyes. ¹ Biotium offers both Di-4-ANEPPS and Di-8-ANEPPS. Di-4-ANEPPS has the problem of being internalized in cells over long-term, so it is mainly used for short-term experiments. However, because Di-8-ANEPPS is more lipophilic, the dye is better retained in the outer leaflet of cell plasma membranes and thus more suitable for long-term membrane potential studies. Di-8-ANEPPS is also more photostable and less phototoxic than Di-4-ANEPPS. ^{2,3} Ref: 1) <i>J. Neurosci.</i> 15 , 1392(1995); 2) <i>Neuron</i> 13 , 1187(1994); 3) <i>Neuron</i> , 9 , 393(1992).
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.*