

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-12-ANEPPQ

CATALOGUE #: 61015

**MOLECULAR**  $C_{47}H_{77}Br_2N_3$  **INFORMATION:** Mwt: 842.8

 $[CH_{3}(CH_{2})_{11}]_{2}N \xrightarrow{\qquad} CH = CH \xrightarrow{\qquad} N^{+} - CH_{2}CH_{2}CH_{2}N(CH_{3})_{3}$ 

## **PROPERTIES:**

Color & FormDark red solidPurity $\geq$  95% by TLC

**Solubility** Soluble in DMSO or DMF **Absorption/Emission** 522 nm / 720 nm (MeOH)

**Extinction**  $\geq 37,000$ 

Coefficient

STORAGE AND HANDLING:

Store desiccated at -20 °C and protect from light especially in solution.

**APPLICATION:** Di-12-ANEPPQ is a fast-responding membrane potential dye. It is more hydrophobic than

Di-8-ANEPEQ and is useful for potential-sensitive retrograde labeling of neurons.

Ref.: 1) J. Neurosci. Methods. 70, 121(1996); 2) J. Neurosci. Methods 70, 111(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.