

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: RH 237

CATALOG #: 61018

 $[CH_3(CH_2)_3]_2N$ $(CH=CH)_3$ $N^+(CH_2)_4SO_3$

PROPERTIES:

Color & FormDark red solidSolubilitySoluble in DMSOAbsorption/Emission520/712nm³ (MeOH)Extinction58,000 (MeOH)

Coefficient

STORAGE AND Store desiccated at \leq -20°C and protect from light. **HANDLING:**

APPLICATION: The styryl dye RH237 is a fast-responding potentiometric probe that is primarily used for

functional imaging of neurons.1

Ref.: 1) Neuroscience 31, 613 (1989); 2) The data is for methanol solution of the dye. In cell membranes, spectra of styryl dyes are typically blue-shifted by as much as 20 nm for absorption or excitation and 80 nm for emission.

RELATED PRODUCTS RH414 (#61016); RH421 (#61017); RH795 (#61019)

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.