

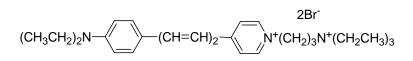
46117 Landing Parkway, Fremont, CA 94538 U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352

www.biotium.com

## **PRODUCT AND SAFETY DATA SHEET**

<b>PRODUCT NAME:</b>	RH 414
	TTT TTT

CATALOG #: 61016



## **PROPERTIES:**

Color & Form Solubility Absorption/Emission Extinction Coefficient	Orange red solid. Soluble in H <sub>2</sub> O or DMSO 532 / 716nm (MeOH) 42,000 (MeOH)
STORAGE AND HANDLING:	Store desiccated at -20 °C and protect from light especially in solution.
APPLICATION:	The styryl dye RH414 is a fast-responding potentiometric probe that is primarily used for functional imaging of neurons. 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. Ref.: <i>Neuroscience</i> <b>31</b> , 613 (1989)
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.*