

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 421 (*N*-(4-sulfobutyl)-4-(4-(dipentylamino)phenyl)butadienyl) pyridinium, inner salt)

CATALOGUE #: 61017

MOLECULAR $C_{29}H_{42}N_2O_3S$. **INFORMATION:** Mwt: 498.72.

[107610-19-5]

N—(CH=CH)—N

PROPERTIES:

Color & Form Red solid.

Soluble in DMSO, EtOH and DMF

Absorption/Emission 515nm / 704nm (MeOH)

Extinction 50,000 (MeOH)

Coefficient

STORAGE AND

HANDLING:

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

APPLICATION: RH421 is a fast responding membrane potential dye, exhibiting >20% fluorescence change

per 100 mV on neuroblastoma cells.

Ref.: Biophys. J. 42, 195(1983)

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.