

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: Rhodamine-DHPE (N-(Lissamine Rhodamine B sulfonyl)-1,2-dihexadecanoyl-

sn-glycero-3-phosphoethanolamine, triethylammonium salt)

CATALOG# 60026

MOLECULAR $C_{70}H_{117}N_4O_{14}PS_2$ MW: 1333.80 **INFORMATION:**

(CH₃CH₂)₂N $N^+(CH_2CH_3)_2$ (CH₃CH₂)₃N⁺H

PROPERTIES:

Color & Form Dark red solid **Purity** ≥90% by TLC

Solubility Soluble in chloroform Absorption/Emission 560/581 nm (MeOH)

Extinction Coefficient >85,000

STORAGE AND Store at -20°C and protect from light. Please note: Chloroform may degrade the O-HANDLING:

ring in the cap on the product vial. For long term storage after reconstitution in

chloroform, transfer the solution to a vial without an O-ring.

Rhodamine-labeled glycerophosphoethanolamine lipids such as Rhodamine-DHPE APPLICATION:

(#60026), TRITC-DHPE (#60028) and Texas Red TM -DHPE (#60027) have been used as fluorescence energy acceptors in combination with NBD-PE as the donor in membrane fussion assays using the principle of fluorescence energy transfer (FRET). These probes have also been used for following membrane trafficking

during endocytosis. 4,5

Ref.: 1) Biochemistry 24, 6390 (1985); 2) Biochemistry 21, 1720(1982); 3) Biochemistry 20, 4093(1981); 4) Biochim.

Biophys. Acta 1103, 185(1992); 5) Eur. J. Cell Biol. 53, 173(1990).

Biotin-DHPE(#60022); Biotin-X-DHPE(#60023); Fluorescein-DHPE(#60024); RELATED PRODUCTS

NBD-PE (#60025); Texas Red-DHPE(#60027); TRITC-DHPE (#60028)

TOXICITY: Not established. Not listed by NTP, IARC or OSHA.

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.