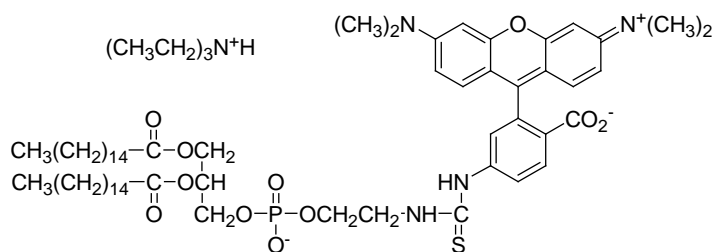


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: **TRITC-DHPE** (*N*-(teramethylrhodamine-6-thiocarbamoyl)-1,2-dihexadecanoyl-*sn*-glycero-3-phosphoethanolamine, triethylammonium salt)

CATALOG # **60028**

MOLECULAR INFORMATION: C₆₈H₁₁₀N₅O₁₁PS
MW: 1236.67

**PROPERTIES:**

Color & Form Dark purple solid
Purity ≥90% by TLC
Solubility Soluble in chloroform
Absorption/Emission 540/566 nm (MeOH)
Extinction Coefficient 93,000

STORAGE AND HANDLING:

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

APPLICATION:

Rhodamine-labeled glycerophosphoethanolamine lipids such as Rhodamine-DHPE (#60026), TRITC-DHPE (#60028) and Texas Red™-DHPE (#60027) have been used as fluorescence energy acceptors in combination with NBD-PE as the donor in membrane fusion assays using the principle of fluorescence energy transfer (FRET).¹⁻³ 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).

RELATED PRODUCTS

Biotin-DHPE(#60022); *Biotin-X-DHPE*(#60023); *Fluorescein-DHPE*(#60024); *NBD-PE* (#60025); *Rhodamine-DHPE*(#60026); *Texas Red-DHPE*(#60027);

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.