TRITC-DHPE

TRITC-DHPE is a rhodamine-labeled glycerophosphoethanolamine lipid. DHPE conjugates of rhodamine dyes have been used in membrane fusion assays.



Product attributes

Probe cellular localization	Membrane/cell surface, Membrane/vesicular
Colors	Red
Excitation/Emission	540/566 nm
Conjugation	Totromothylrhodomino (TRITC)

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

TRITC-DHPE

(N-(Tetramethylrhodamine-6-thiocarbamoyl)-1,2-dihexadecanoyl-sn-glycero-3-phosphoethanolamine, triethylammonium salt) is a rhodamine-labeled glycerophosphoethanolamine lipid. DHPE conjugates of rhodamine dyes have been used as fluorescence energy acceptors in combination with NBDPE 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).

- $\lambda_{Ex}/\lambda_{Em}(MeOH) = 540/566 \text{ nm}$
- Dark solid soluble in chloroform
- Store at -20°C and protect from light
- C68H110N5O11PS
- MW: 1236.67

This datasheet was generated on November 5, 2025 at 02:16:35 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/tritc-dhpe-n-teramethylrhodamine-6-thiocarbamoyl-12-dihexadecanoyl-sn-glycero-3-phosphoethanolamine-triethylammonium-salt/