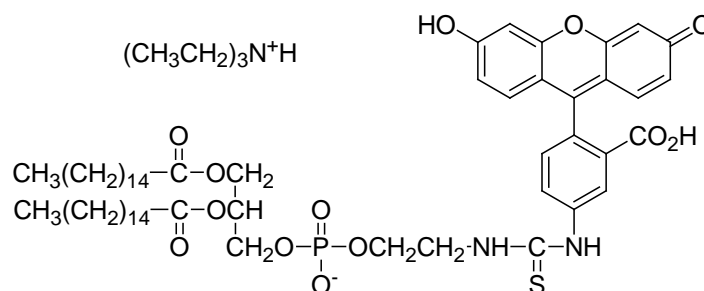


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: **Fluorescein-DHPE** (*N*-(fluorescein-5-thiocarbamoyl)-1,2-dihexadecanoyl-*sn*-glycero-3-phosphoethanolamine, triethylammonium salt)

CATALOG # **60024**

MOLECULAR INFORMATION: C₆₄H₁₀₀N₃O₁₃PS
MW: 1182.53

**PROPERTIES:**

Color & Form Orange solid
Purity ≥95% by TLC
Solubility Soluble in chloroform
Absorption/Emission 496/519 nm (MeOH)
Extinction Coefficient ≥80,000

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:

Fluorescein-DHPE is a useful surface pH indicator¹ and has also been used for measuring lateral diffusion in membranes via fluorescence photobleaching recovery method

Ref.: 1) *Biochim. Biophys. Acta* **939**, 289(1988); 2) *Biophys. J.* **66**, 25(1994); 3) *J. Cell Biol.* **103**, 807(1986).

RELATED PRODUCTS

Biotin-DHPE(#60022); *Biotin-X-DHPE*(#60023); *NBD-DHPE*(#60025); *Rhodamine-DHPE*(#60026); *Texas Red-DHPE*(#60027); *TRIC 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.