

Fluorescein-DHPE

Fluorescein-DHPE is a useful surface pH indicator, and has also been used for measuring lateral diffusion in membranes.



Product attributes

Call us: 800-304-5357

CAS number	87706-98-7
Probe cellular localization	Membrane/cell surface, Membrane/vesicular
Colors	Green
Excitation/Emission	496/519 nm
Conjugation	Fluorescein (FITC)

Product Description

Fluorescein-DHPE

(N-(Fluorescein-5-thiocarbamoyl)-1,2-dihexadecanoyl-sn-glycero-3-phosphoethanolamine, triethylammonium salt) is a useful surface pH indicator (1) and has also been used for measuring lateral diffusion in membranes via fluorescence photobleaching recovery method.

- λεx/λεm (MeOH) = 496/ 519 nm
- Orange solid soluble in chloroform
- Store at -20°C and protect from light
- C₆₄H₁₀₀N₃O₁₃PSMW: 1182.53
- [87706-98-7]

References

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

This datasheet was generated on November 23, 2025 at 03:04:15 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/fluorescein-dhpe-n-fluorescein-5-thiocarbamoyl-12-dihexadecanoyl-sn-glycero-3-phosphoethanolamine-triethylammoniumsalt/