

Fluorescein-DHPE

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



Product attributes

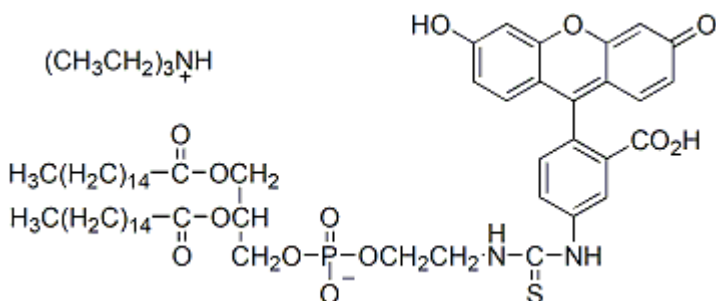
CAS number	87706-98-7
Probe cellular localization	Membrane/cell surface, Membrane/vesicular
Colors	Green
Excitation/Emission	496/519 nm
Conjugation	Fluorescein (FITC)

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.

- $\lambda_{Ex}/\lambda_{Em}$ (MeOH) = 496/ 519 nm
- Orange solid soluble in chloroform
- Store at -20 °C and protect from light
- $C_{64}H_{100}N_3O_{13}PS$
- MW: 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 January 20, 2026 at 03:51:19 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/>