

Biotin-X-DHPE

Biotin-X-DHPE

(N-((6-(Biotinoyl)amino)hexanoyl)-1,2-dihexadecanoylsn-glycero-3-phosphoethanolamine, triethylammonium salt) can be used to couple avidin or streptavidin to cell membranes or liposomes.



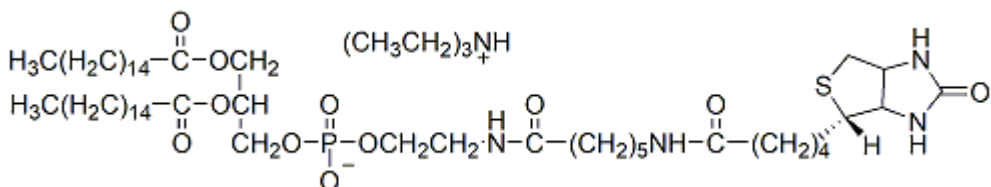
Product attributes

Probe cellular localization	Membrane/cell surface, Membrane/vesicular
Conjugation	Biotin

Product Description

Biotin-X-DHPE (N-((6-(Biotinoyl)amino)hexanoyl)-1,2-dihexadecanoylsn-glycero-3-phosphoethanolamine, triethylammonium salt) can be used to couple avidin or streptavidin to cell membranes or liposomes.

- White solid soluble in chloroform
- Store at -20 °C
- C₅₉H₁₁₄N₅O₁₁PS
- MW: 1132.60



References

1. Immunol Meth 158, 183 (1993).
2. Anal Biochem 207, 341 (1992).
3. Biophys J 59, 387 (1991).

This datasheet was generated on June 22, 2026 at 10:05:14 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/biotin-x-dhpe-n-6-biotinoylaminohexanoyl-12-dihexadecanoylsn-glycero-3-phosphoethanolamine-triethylammonium-salt/>