

Biotin-DHPE

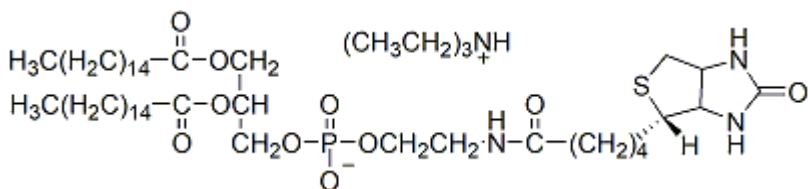


The image shows the Biotium logo, which consists of the word "Biotium" in a sans-serif font next to a stylized graphic of blue and white circles. Below the logo is a small vial of Biotium product, which is a clear glass vial with a black cap and a white label. The vial is placed on a white surface, and its reflection is visible below it.

Product Description

Biotin-DHPE (N-(Biotinoyl)-1,2-dihexadecanoyl-sn-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
- C₅₃H₁₀₃N₄O₁₀PS
- MW: 1019.44
- [136235-58-0]



References

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

This datasheet was generated on January 6, 2026 at 08:50:47 PM. Visit [product page](#) to check for updated information before use.

Product link: <https://biotium.com/product/biotin-dhpe-n-biotinoyl-12-dihexadecanoyl-sn-glycero-3-phosphoethanolamine-triethylammonium-salt/>