

46117 Landing Parkway, Fremont, CA 94538 U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352

www.biotium.com

## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** Biotin-DHPE (N-(biotinoyl)-1,2-dihexadecanoyl-sn-glycero-3-

phosphoethanolamine, triethylammonium salt)

**CATALOG #** 60022

**MOLECULAR**  $C_{53}H_{103}N_4O_{10}PS$  **INFORMATION:** MW: 1019.44

**PROPERTIES:** 

Color & Form White solid

**Soluble** in chloroform

**Absorption/Emission** N/A **Extinction Coefficient** N/A

STORAGE AND HANDLING:

Store at -20 °C. **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:** Biotin-DHPE can be used to couple avidin or streptavidin to cell membranes or

liposomes.

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

**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.