

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-12 SE (4-((6-((biotinoyl)amino)hexanoyl)amino)butyric acid succinimidyl

ester)

CATALOG #: 90083

MOLECULAR $C_{24}H_{37}N_5O_7S$ **INFORMATION:** MW: 539.64

HN NH
H O O O
S
$$(CH_2)_4$$
 $(CH_2)_5$
 $(CH_2)_5$

PROPERTIES:

Color & FormWhite solidPurity $\geq 95\%$ by TLC

Soluble in DMSO or DMF

STORAGE AND HANDLING:

Store desiccated at -20 °C.

APPLICATION: Biotin-12 SE is an amine-reactive biotin derivative with a long spacer group. This

product can be conjugated to AA-dUTP or AA-UTP to make biotin-16-dUTP or

biotin-16-UTP, two important products used in genomic studies.

TOXICITY: Unknown.

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.