

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: NBD C6-Sphingomyelin (6-((N-(7-nitrobenz-2-oxa-1,3-diazol-4-yl)

amino)hexanoyl)sphingosyl phosphocholine)

CATALOG #: 60031

 $CH_{3}(CH_{2})_{12}CH=CH-CHOH$ $NH(CH_{2})_{5}-C-NHCH$ O $CH_{2}O-P-OCH_{2}CH_{2}N^{+}(CH_{3})_{5}$ O O

PROPERTIES:

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

Soluble in chloroform or DMSO

 $\lambda_{ex} \setminus \lambda_{em}$ 466/ 536nm (MeOH)

STORAGE AND HANDLING:

Store at -20 °C and protect from light.

APPLICATION: NBD C6-Sphingomyelin has been used for following sphingolipid metabolism in cells.

Ref.:1) Methods Cell Biol. 38, 221(1993); 2) Biochim Biophys Acta 1113, 277 (1992); 3) Adv. Cell Mol. Biol. Membranes 1, 199(1993); 4) Biochim Biophys Acta 1082, 113 (1991).

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.