

# PRODUCT DATA SHEET

## N-Hexanoyl-NBD-L-*threo*-sphingosine

**Catalog number:** 1857; 1857-001

**Synonyms:** N-C6:0-NBD-L-*threo*-Ceramide;  
N-C6:0-NBD-L-*threo*-Sphingosine

**Source:** synthetic

**Solubility:** ethanol, methanol, chloroform

**CAS number:** 114301-95-0

**Molecular Formula:** C<sub>30</sub>H<sub>49</sub>N<sub>5</sub>O<sub>6</sub>

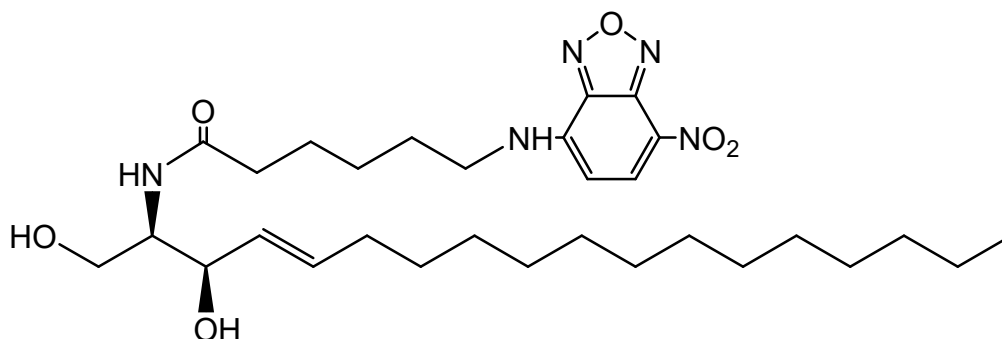
**Molecular Weight:** 576

**Storage:** -20°C

**Purity:** TLC: >98%

**TLC System:** chloroform/methanol (90:10)

**Appearance:** solid



### Application Notes:

This product is a high purity, non-natural L-*threo* ceramide containing a fluorescent NBD label and is ideal as a standard and for biological studies. NBD (7-nitrobenzofurazan) has been shown to have only a small influence on lipid adsorption into cells and cellular membranes especially when the fatty acid is a short chain. The fluorescent analog of natural ceramide is comparable to C6:0-ceramide in many biological functions such as inhibition of VSV-G protein transport<sup>1</sup>, and transport of sphingomyelin and glucocerebroside from the golgi apparatus to the cell surface.<sup>2</sup> Both the natural D-*erythro* and the non-natural L-*erythro* and the D- and L-*threo* ceramides display similar effectiveness in inducing apoptotic damage in cells.<sup>3</sup> The protein phosphatases PP1 and PP2A, which are involved in regulating apoptosis and cell growth, are activated by D-*erythro* ceramide but inhibited by L-*threo*, D-*threo*, and L-*erythro* ceramide.<sup>4</sup> Both D-*erythro* and D-*threo* C2 ceramides have been found to be potent inducers of IL-6 production, while neither the L-*threo* or L-*erythro* stereoisomers of ceramide were effective.<sup>5</sup> D- and L-*erythro* ceramide and D- and L-*threo* ceramide are also comparably effective inhibitors of protein kinase C.<sup>6</sup>

### Selected References:

1. A. Rosenwald and R. Pagano "Inhibition of glycoprotein traffic through the secretory pathway by ceramide" *Journal of Biological Chemistry*, Vol. 268 pp. 4577-4579, 1993
2. N. Lipsky, R. Pagano "Intracellular translocation of fluorescent sphingolipids in cultured fibroblasts: Endogenously synthesized sphingomyelin and glucocerebroside analogues pass through the golgi apparatus en route to the plasma membrane" *Journal of Cell Biology*, Vol. 100 pp. 27-34, 1985
3. W. Jarvis et al. "Induction of Apoptosis and Potentiation of Ceramide-mediated Cytotoxicity by Sphingoid Bases in Human Myeloid Leukemia Cells" *The Journal of Biological Chemistry*, Vol. 271 pp. 8275-8284, 1996
4. C. Chalfant et al. "Long Chain Ceramides Activate Protein Phosphatase-1 and Protein Phosphatase-2A Activation is Stereospecific and Regulated by Phosphatidic Acid" *The Journal of Biological Chemistry*, Vol. 274 pp. 20313-20317, 1999
5. S. Lauderkind et al. "Ceramide Induces Interleukin 6 Gene Expression in Human Fibroblasts" *The Journal of Experimental Medicine*, Vol. 182 pp. 599-604, 1995
6. T. Ariga et al. "Role of sphingolipid-mediated cell death in neurodegenerative diseases" *Journal of Lipid Research*, Vol. 39 pp. 1-16, 1998

This product is to be used for research only. It is not intended for drug or diagnostic use, human consumption or to be used in food or food additives. Matreya assumes no liability for any use of this product by the end user. We believe the information, offered in good faith, is accurate.