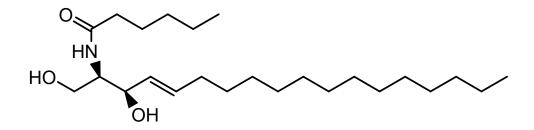


## **PRODUCT DATA SHEET**

## N-Hexanoyl-D-threo-sphingosine

Catalog number: 1809 Common Name: N-C6:0-D-threo-Ceramide Source: synthetic Solubility: chloroform, methanol, DMSO (up to 5mg/ml) CAS number: 189894-79-9 Molecular Formula: C<sub>24</sub>H<sub>47</sub>NO<sub>3</sub> Molecular Weight: 398 Storage: -20°C Purity: TLC > 98%, GC > 98%; identity confirmed by MS TLC System: chloroform/methanol (95:5 by vol.) Appearance: solid



## **Application Notes:**

Ceramide is a fatty acid amide of sphingosine. This product is a well-defined ceramide with a hexanoyl acyl group. Ceramide functions as a precursor in the synthesis of sphingomyelin, glycosphingolipids, and of free sphingosine and fatty acids. The sphingosine can be phosphorylated to form sphingosine-1-phosphate. Two of ceramide's metabolites, sphingosine-1phosphate and glucosylceramide, produce cell proliferation and other cellular functions.<sup>1</sup> Ceramide exerts numerous biological effects, including induction of cell maturation, cell cycle arrest, terminal cell differentiation, cell senescence, and cell death.<sup>2</sup> Because of these effects ceramide has been investigated for its use in cancer treatment and many potential approaches to cancer therapy have been presented.<sup>3</sup> Other effects include producing reactive oxygen in mitochondria (followed by apoptosis) and stimulating phosphorylation of certain proteins (especially mitogen activated protein).<sup>4</sup> It also stimulates some protein phosphatases (especially protein phosphatase 2A) making it an important controller of protein activity. Ceramides with short side chains have been shown to enter easily into cells where they are biologically active. Ceramides with longer side chains also enter cells if dissolved in dodecane-isopropanol first.

## **Selected References:**

- 1. J. Hauser, B. Buehrer, and R. Bell "Role of ceramide in mitogenesis induced by exogenous sphingoid bases." *Journal of Biological Chemistry* Vol. 269 pp. 6803, *1994*
- 2. N. Radin, "Killing tumours by ceramide-induced apoptosis: a critique of available drugs" Biochemical Journal, Vol. 371 pp. 243-256, 2003
- 3. N. Radin, "Designing anticancer drugs via the achilles heel: ceramide, allylic ketones, and mitochondria" *Bioorganic and Medicinal Chemistry*, Vol. 11(10) pp. 2123-2142, *2003*
- 4. M. Yano, E. Kishida, Y. Muneyuki, and Y. Masuzawa "Quantitative analysis of ceramide molecular species by high performance liquid chromatography" *Journal of Lipid Research*, Vol. 39 pp. 2091-2098, *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.