

# PRODUCT DATA SHEET

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

**Catalog number:** 1828

**Common names:** N-C6:0-L-*threo*-Ceramide

**Source:** synthetic

**Solubility:** chloroform, ethanol, DMSO, DMF  
(up to 5 mg/ml)

**CAS number:** 189894-80-2

**Molecular Formula:** C<sub>24</sub>H<sub>47</sub>NO<sub>3</sub>

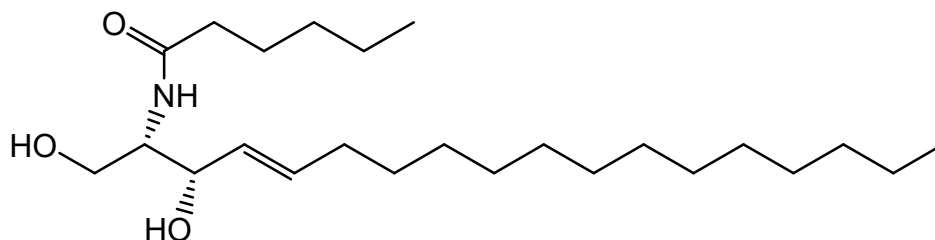
**Molecular Weight:** 398

**Storage:** -20°C

**Purity:** TLC >98%, GC >98%; identity  
confirmed by MS

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

**Appearance:** solid



### Application Notes:

This product is the non-natural L-*threo* stereoisomer of ceramide. Natural D-*erythro* ceramide is a critical compound in cells both as a free ceramide and incorporated into more complex sphingolipids. L-*threo*-ceramides demonstrate a different metabolic functionality from natural ceramides. They have been shown to be metabolized to sphingomyelin but not to glucosylceramide.<sup>1</sup> Another non-natural stereoisomer, L-*erythro* ceramide, is not metabolized to any sphingolipid. In contrast to natural ceramides L-*threo* ceramides are unable to antagonize the disruptive effects of fumonisins B1 on axon growth<sup>2</sup> but it is able to activate intracellular pathways and induces apoptosis.<sup>3</sup> The deacylated form of ceramide, sphingosine, also has many critical cellular functions. L-*threo*-sphingosine, along with other sphingosine isomers, has been found to be an activator of 3-Phosphoinositide-dependent kinase-1<sup>4</sup> and inhibits protein kinase C a little more potently than D-*erythro*-sphingosine.<sup>5</sup>

### Selected References:

1. K. Venkataraman and H. Futerman "Comparison of the metabolism of L-erythro- and L-threo-sphinganine and ceramides in cultured cells and in subcellular fractions" *Biochim Biophys Acta*, vol. 1530 pp. 219-226, 2001
2. A. Schwarz and A. Futerman "Distinct Roles for Ceramide and Glucosylceramide at Different Stages of Neuronal Growth" *The Journal of Neuroscience*, vol. 17 pp. 2929-2938, 1997
3. A. Bielawska et al. "Selectivity of ceramide-mediated biology—lack of activity of erythrodihydroceramide" *J Biol Chem*, vol. 268 pp. 26226–26232, 1993
4. C. King et al. "Sphingosine Is a Novel Activator of 3-Phosphoinositide-dependent Kinase 1" *The Journal of Biological Chemistry*, vol. 275(24) pp. 18108-18113, 2000
5. V. Stevens et al. "Structural requirements for long-chain (sphingoid) base inhibition of protein kinase C in vitro and for the cellular effects of these compounds" *Biochemistry*, vol. 28, 3138-3145, 1989

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.