

# PRODUCT DATA SHEET

## N-Pentadecanoyl-D-erythro-sphingosine

**Catalog number:** 2037

**Common Name:** N-C15:0-D-erythro-Ceramide

**Source:** synthetic

**Solubility:** chloroform, warm ethanol, warm methanol

**CAS number:** 67492-15-3

**Molecular Formula:** C<sub>33</sub>H<sub>65</sub>NO<sub>3</sub>

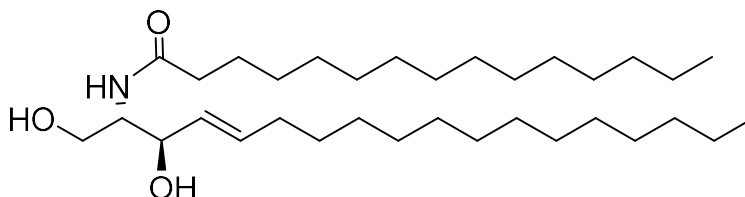
**Molecular Weight:** 524

**Storage:** -20°C

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

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

**Appearance:** solid



### Application Notes:

This product is a high purity ceramide containing an uncommon C15:0 fatty acid acylated to sphingosine making it ideal as an internal standard and for biological studies.<sup>1</sup> Ceramide is a fatty acid amide of sphingosine. 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-1-phosphate and glucosylceramide, produce cell proliferation and other cellular functions.<sup>2</sup> Ceramide exerts numerous biological effects, including induction of cell maturation, cell cycle arrest, terminal cell differentiation, cell senescence, and cell death.<sup>3</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>4</sup> Other effects include producing reactive oxygen in mitochondria (followed by apoptosis) and stimulating phosphorylation of certain proteins (especially mitogen activated protein). It also stimulates some protein phosphatases (especially protein phosphatase 2A) making it an important controller of protein activity.

### Selected References:

1. C. Vieu et al. "Coupled assay of sphingomyelin and ceramide molecular species by gas liquid chromatography" *Journal of Lipid Research*, vol. 43 pp. 510-522, 2002
2. J. M. Hauser, B. M. Buehrer, and R. M. Bell "Role of ceramide in mitogenesis induced by exogenous sphingoid bases." *Journal of Biological Chemistry* Vol. 269 pp. 6803, 1994
3. N. S. Radin, "Killing tumours by ceramide-induced apoptosis: a critique of available drugs" *Biochemical Journal*, Vol. 371 pp. 243-256, 2003
4. N. S. Radin, "Designing anticancer drugs via the achilles heel: ceramide, allylic ketones, and mitochondria" *Bioorganic and Medicinal Chemistry*, Vol. 11(10) pp. 2123-2142, 2003

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.