

PRODUCT DATA SHEET

L-threo-Sphingosine

Catalog number: 1806

Common Name: L-threo-Sphingosine, C18 chain

Source: synthetic

Solubility: ethanol, methanol, chloroform, DMSO

CAS number: 25695-95-8

Molecular Formula: C₁₈H₃₇NO₂

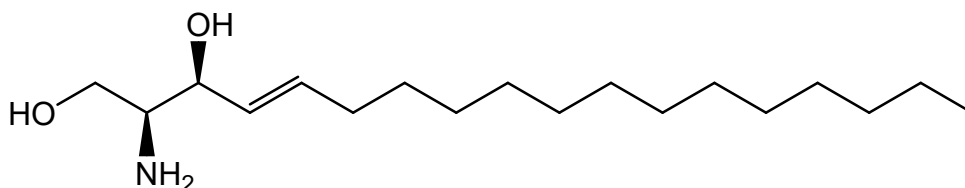
Molecular Weight: 300

Storage: -20°C

Purity: TLC >98%, GC >98%

TLC System: chloroform/methanol/DI water/ammonium hydroxide (70:20:1:1 by vol.)

Appearance: solid



Application Notes:

L-threo-Sphingosine is an inactive or less active isomer of the naturally occurring D-erythro-sphingosine. Natural sphingosine induces dephosphorylation of retinoblastoma gene products and inhibits cell growth while L-threo-sphingosine is less active. However, the L-threo-sphingosine is taken up by cells to the same extent as the natural sphingosine indicating that cellular uptake was not the factor influencing activity.¹ L-threo-sphingosine, along with other sphingosine isomers, has been found to be an activator of 3-Phosphoinositide-dependent kinase-1.² Natural D-erythro-sphingosine is a positive regulator of cell growth in fibroblasts whereas L-threo-sphingosine has no regulatory effect.³ However, non-natural stereoisomers of sphingosine are not always inactive; L-threo-sphingosine has been shown to inhibit protein kinase C a little more potently than D-erythro-sphingosine.⁴

Sphingosine is a characteristic structural unit of many sphingolipids such as ceramides, gangliosides, globosides, sulfatides, sphingomyelin, and others. It is most abundant in nervous tissue and cell membranes. Sphingosine with an 18-carbon chain and a double bond at carbon 4 is the most abundant sphingosine in animal tissues. Lysosphingolipids inhibit protein kinase C activity resulting in the pathogenesis of sphingolipidoses such as Krabbe's disease and Gaucher's disease. Sphingosine can be phosphorylated via two kinases to form sphingosine-1-phosphate, which has important signaling functions. While sphingosines and ceramides can induce apoptosis, sphingosine-1-phosphate can promote cell survival or proliferation. Sphingosine has been shown to cause an increase in the cytoplasmic calcium level of cells.

Selected References:

1. Y. Hannun et al. "Stereo-selectivity of Induction of the Retinoblastoma Gene Product (pRb) Dephosphorylation by D-erythro-Sphingosine Supports a Role for pRb in Growth Suppression by Sphingosine" *Biochemistry*, vol. 34 pp. 1885-1892, 1995
2. 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
3. S. Spiegel et al. "Stereo-specificity of Sphingosine-induced Intracellular Calcium Mobilization and Cellular Proliferation" *Journal of Biological Chemistry*, vol. 269 pp. 17924-17930, 1994
4. 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.