

PRODUCT DATA SHEET

L-erythro-Dihydrosphingosine

Catalog number: 1846 Common names: L-*erythro*-Sphinganine, C18 chain Source: synthetic

Solubility: chloroform, methanol, ethanol, DMSO CAS number: 6036-76-6 Molecular Formula: C₁₈H₃₉NO₂ Molecular Weight: 301 Storage: -20°C Purity: TLC >98%, GC >98%; identity confirmed by MS TLC System: chloroform/methanol/DI water/ ammonium hydroxide (70:20:1:1) Appearance: solid



Application Notes:

This product is a high purity, well-defined, L-*erythro*-dihydrosphingosine which demonstrates unique properties as compared with the natural D-*erythro* isomer and is therefore ideal for use in studies of dihydrosphingosine. Natural D-*erythro*-dihydrosphingosine is the precursor of dihydroceramide which is then desaturated to form ceramide. It is a critical intermediate in the synthesis of many complex sphingoid bases and ceramide analogs. It has been found that dihydrosphingosine can induce cell death in a number of types of malignant cells and it is being tested for its pharmacological properties.¹ Whereas both D-*threo* and L-*threo*-C2-dihydroceramide induced apoptosis in cells neither D-*erythro*-C2-dihydroceramide showed activity.² One report has concluded that all four of the enantiomers of dihydrosphingosine act as substrates for sphingosine kinase with only the natural D-*erythro*-dihydrosphingosine being metabolized by sphinganine-1-phosphate lyase.³ However, another report concludes that only the *erythro* isomers of dihydrosphingosine act as substrates for this enzyme with both of the *threo* isomers inhibiting its activity.⁴

Selected References:

- 1. W. Zheng "Fenretinide increases dihydroceramide and dihydrosphingolipids due to inhibition of dihydroceramide desaturase" Georgia Institute of Technology, 2006
- 2. A. Bielawska "Selectivity of Ceramide-Mediated Biology Lack of Activity of *erythro*-Dihydroceramide" *Journal of Biological Chemistry*, vol. 268 pp. 26226-26232, *1993*
- 3. W. Stoffel and K. Bister "Stereospecificities in the metabolic reactions of the four isomeric sphinganines (dihydrosphingosines) in rat liver" *Hoppe Seylers Z Physiol Chem*, vol. 354 pp. 169-181, *1973*
- 4. B. Buehrer and R. Bell "Inhibition of Sphingosine Kinase *in Vitro* and in Platelets Implications for Signal Transduction Pathways" *Journal of Biological Chemistry*, vol. 267 pp. 3154-3159, *1992*

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.