

PRODUCT DATA SHEET

D-erythro-C20-Dihydrosphingosine

Catalog number: 1845

Common Name: D-erythro-Sphinganine, C20 chain

Source: synthetic

Solubility: chloroform/methanol (5:1 by vol.)
warm ethanol

CAS number: 24006-62-0

Molecular Formula: C₂₀H₄₃NO₂

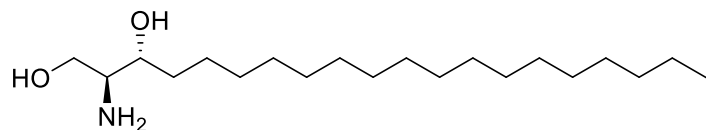
Molecular Weight: 330

Storage: -20°C

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

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

Appearance: solid



Application Notes:

This product is the natural D-erythro isomer of dihydrosphingosine. Sphinganine (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 sphinganine can induce cell death in a number of types of malignant cells and is being tested for its pharmacological properties.¹ Inhibition of dihydroceramide synthesis by some fungal toxins that have a similar structure causes an increase in sphinganine and sphinganine-1-phosphate and a decrease in other sphingolipids leading to a number of diseases including oesophageal cancer.² Sphinganine has been found to mediate fumonisin (a toxic sphinganine analog) induced hypotension.³ In yeast the amount of C20-dihydrosphingosine increases as a response to heat stress along with other sphingolipids, indicating that it is involved in heat stress adaptation.

Selected References:

1. W. Zheng "Fenretinide increases dihydroceramide and dihydrosphingolipids due to inhibition of dihydroceramide desaturase" Georgia Institute of Technology, 2006
2. L. van der Westhuizen et al. "Sphingoid base levels in humans consuming fumonisin-contaminated maize in rural areas of the former Transkei, South Africa: a cross-sectional study" *Food Additives and Contaminants*, Vol. 25(11), pages 1385 – 1391, 2008
3. Shih-Hsuan Hsiao et al. "Effects of Exogenous Sphinganine, Sphingosine, and Sphingosine-1-Phosphate on Relaxation and Contraction of Porcine Thoracic Aortic and Pulmonary Arterial Rings" *Toxicological Sciences*, Vol. 86(1) Pp. 194-199, 2005

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.