

PRODUCT DATA SHEET

N-Dodecanoyl-NBD-D-*erythro*-dihydrosphingosine

Catalog No: 1625

Common Name: N-C12:0-NBD-Dihydroceramide; N-C12:0-NBD-D-*erythro*-Dihydrosphingosine

Source: synthetic

Solubility: chloroform/methanol (2:1 by vol.)
methanol

CAS No: 474943-05-0

Molecular Formula: C₃₆H₆₃N₅O₆

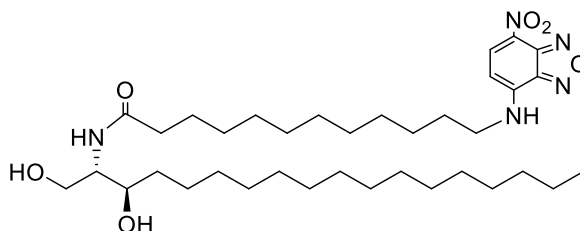
Molecular Weight: 662

Storage: -20°C

Purity: TLC > 98%; identity confirmed by MS

TLC System: chloroform/methanol (90:10 by vol.)

Appearance: solid



Application Notes:

This high purity fluorescent product is ideal for the identification of dihydroceramide in samples and biological systems. 7-nitrobenzofurazan (NBD) has been shown to have only a small influence on lipid adsorption into cells and cellular membranes. This fluorescent analog of natural dihydroceramide is comparable to C12:0-dihydroceramide in many biological functions.^{1,2} Dihydroceramide is a critical intermediate in the synthesis of many complex sphingoid bases. Inhibition of dihydroceramide synthesis by some fungal toxins that have a similar structure causes an increase in dihydroceramide and dihydroceramide-1-phosphate and a decrease in other sphingolipids leading to a number of diseases including oesophageal cancer. Dihydroceramide, synthesized by the acylation of sphinganine, is subsequently converted into ceramide via a desaturase enzyme. N-(4-Hydroxyphenyl) retinamide (4-HPR) has been tested as an anti-cancer agent. It inhibits the dihydroceramide desaturase enzyme in cells resulting in a high concentration of dihydroceramide and dihydro-sphingolipids and this is thought to be the cause of the anti-cancer effects.³ Dihydrosphingosine induces cell death in a number of types of malignant cells.

Selected References:

1. J. Kok et al. "Dihydroceramide Biology STRUCTURE-SPECIFIC METABOLISM AND INTRACELLULAR LOCALIZATION" *Journal of Biological Chemistry*, Vol. 272 pp. 21128-21136, 1997
2. J. Hsu et al. "Enhanced endothelial delivery and biochemical effects of *alpha*-galactosidase by ICAM-1-targeted nanocarriers for Fabry disease" *Journal of Controlled Release*, doi:10.1016/j.jconrel.2010.10.031, 2010
3. W. Zheng "Fenretinide increases dihydroceramide and dihydrosphingolipids due to inhibition of dihydroceramide desaturase" *Georgia Institute of Technology*, 2006

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.