

# PRODUCT DATA SHEET

## N-Hexanoyl-NBD-D-erythro-dihydrosphingosine

**Catalog number:** 1626

**Common Name:** N-C6:0-NBD-Dihydroceramide; N-C6:0-NBD-D-erythro-Dihydrosphingosine

**Source:** synthetic

**Solubility:** chloroform/methanol (2:1), methanol

**CAS number:** 114301-97-2

**Molecular Formula:** C<sub>30</sub>H<sub>51</sub>N<sub>5</sub>O<sub>6</sub>

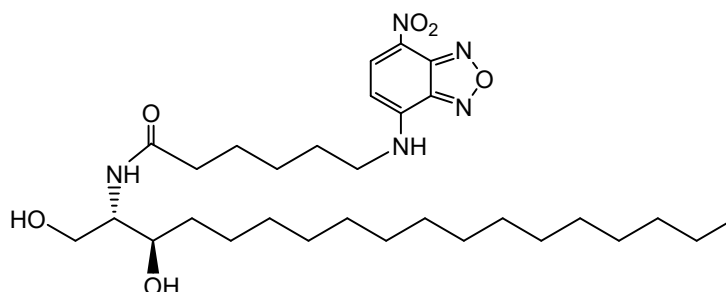
**Molecular Weight:** 578

**Storage:** -20°C

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

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

**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 C6:0-dihydroceramide in many biological functions.<sup>1,2</sup> 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.<sup>3</sup> Dihydrosphingosine induces cell death in a number of types of malignant cells.

### Selected References:

1. G van Meer et al. "Epithelial sphingolipid sorting allows for extensive variation of the fatty acyl chain and the sphingosine backbone" *Journal of Biochemistry*, vol. 283 pp. 913-917, 1992
2. A. Merrill, Jr. et al. "Dihydroceramide Biology STRUCTURE-SPECIFIC METABOLISM AND INTRACELLULAR LOCALIZATION" *Journal of Biological Chemistry*, vol. 272 pp. 21128-21136, 1997
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.