

## **PRODUCT DATA SHEET**

## N-Octanoyl-D-erythro-dihydrosphingosine

Catalog number: 1854

Common Name: N-C8:0-D-*erythro*-Dihydroceramide; N-Octanoyl-D-*erythro*sphinganine

Source: synthetic Solubility: ethanol, DMSO, chloroform CAS number: 145774-33-0 Molecular Formula: C<sub>26</sub>H<sub>53</sub>NO<sub>3</sub> Molecular Weight: 428 Storage: -20°C Purity: TLC >98%, GC >98%; identity confirmed by MS TLC System: chloroform/methanol (95:5) Appearance: solid



## **Application Notes:**

This high purity and well-defined product is ideal as a standard and for biological studies.<sup>1</sup> Dihydroceramide is a critical intermediate in the synthesis of many complex sphingoid bases. Inhibition of dihydroceramide synthesis by some fungal toxins (such as fumonisin B1) 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. Dihydroceramide, synthesized by the acylation of sphinganine, is subsequently converted into ceramide via a desaturase enzyme or into phytosphingosine via the C4-hydrozylase enzyme<sup>2</sup>. 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. Z. Zakeri et al. "Stereospecific Induction of Apoptosis in U937 Cells by N-Octanoyl-Sphingosine Stereoisomers and N-Octyl-Sphingosine" European Journal of Biochemistry, vol. 236 pp. 729-737, 1996

2. Y. Mizutani, A. Kihara, and Y. Igarashi "Identification of the human sphingolipid C4-hydroxylase, hDES2, and its up-regulation during keratinocyte differentiation" *FEBS Letters*, vol. 563 pp. 93-97, 2004

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.