

PRODUCT DATA SHEET

N-Acetyl-D-erythro-dihydrosphingosine

Catalog number: 1834 Common names: N-Acetyl-D-*erythro*sphinganine; N-C2:0-D*erythro*-Dihydroceramide

Source: synthetic Solubility: chloroform, ethanol, methanol CAS number: 13031-64-6 Molecular Formula: C₂₀H₄₁NO₃ Molecular Weight: 344 Storage: -20°C Purity: TLC: 98%, GC >98%; identity confirmed by MS TLC System: chloroform/methanol (90:10) Appearance: solid



Application Notes:

N-Acetyl-D-*erythro*-dihydrosphingosine is a non-natural analog of the ceramide precursor dihydroceramide. This acetyl dihydroceramide has properties and functions that are different from natural dihydroceramide and is therefore a useful tool for studying dihydroceramides and ceramides. Whereas short chain ceramides potently induce apoptosis in cells neither short nor long-chain dihydroceramides are able to do so.¹ N-Acetyl-dihydroceramide has been shown to be incorporated into mitochondria membranes at the same rate as N-acetyl-ceramide and at a greater rate than N-palmitoyl-ceramide. This demonstrates that both dihydroceramide and ceramide insert into the mitochondrial membrane at the same rate indicating that the inability of dihydroceramide to induce apoptosis is not due to a lack of its insertion into the membrane.² The presence of the 4–5 *trans* double bond is essential for ceramide-channel formation. Dihydroceramides are unable to form these ceramide-channels due to their lack of a 4-5 double bond and it is suggested that this is the reason for its lack of apoptotic activity.³ Natural 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 dihydrosphingosine and dihydrosphingosine-1-phosphate and a decrease in other sphingolipids leading to a number of diseases including oesophageal cancer. 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.⁴

Selected References:

1. F. Bi et al. "A Conserved Cysteine Motif Is Critical for Rice Ceramide Kinase Activity and Function" PLoS ONE 6(3): e18079.

- doi:10.1371/journal.pone.0018079, 2011
- 2. L. Siskind et al. "Ceramide forms channels in mitochondrial outer membranes at physiologically relevant concentrations" *Mitochondrion*, vol. 6 pp. 118-125, 2006
- 3. T. Goldkorn et al. "H2O2 acts on cellular membranes to generate ceramide signaling and initiate apoptosis in tracheobronchial epithelial cells" *Journal of Cell Science*, vol. 111 pp. 3209-3220, *1998*
- 4. 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.