

PRODUCT DATA SHEET

N-Acetyl-L-*erythro*-sphingosine

Catalog number: 1847

Common names: N-C2:0-L-*erythro*-Ceramide

Source: synthetic

Solubility: chloroform, methanol, ethanol,
DMSO, DMF (up to 5 mg/ml)

CAS number: 150338-90-2

Molecular Formula: C₂₀H₃₉NO₃

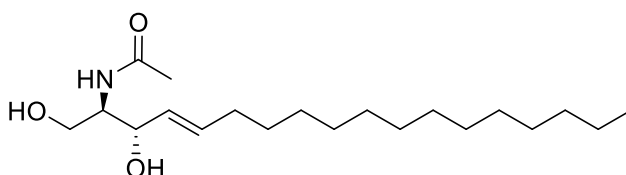
Molecular Weight: 342

Storage: -20°C

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

TLC System: chloroform/methanol (90:10)

Appearance: solid



Application Notes:

This product contains an L-*erythro*-sphingosine containing an acetic amide linkage and is useful for comparison studies against the natural D-*erythro* isomer. This product is also very useful as an internal standard^{1,2} and contains a short-chain fatty acid enabling it to enter easily into cells. L-*erythro*-Sphingosine is an inactive or less active isomer of the naturally occurring D-*erythro*-sphingosine. Natural sphingosine induces dephosphorylation of retinoblastoma gene product and inhibits cell growth while L-*erythro*-sphingosine is 5-8-fold less active. However, the L-*erythro*-sphingosine is taken up by cells to the same extent as the natural sphingosine indicating that cellular uptake was not the factor influencing activity.³ L-*erythro*-sphingosine, along with other sphingosine isomers, has been found to be an activator of 3-Phosphoinositide-dependent kinase 1.

Selected References:

1. T. Cunningham et al. "Product inhibition of secreted phospholipase A2 may explain lysophosphatidylcholines' unexpected therapeutic properties" *Journal of Inflammation*, 5:17 doi:10.1186/1476-9255-5-17, 2008
2. T. Cunningham et al "Uncompetitive Phospholipase A2 Inhibition by CHEC Sequences Including Oral Treatment of Experimental Autoimmune Myeloencephalitis" *The Open Enzyme Inhibition Journal*, vol. 2 pp. 1-7, 2009
3. Y. Hannun et al. "Stereoselectivity of Induction of the Retinoblastoma Gene Product (pRb) Dephosphorylation by D-*erythro*-Sphingosine Supports a Role for pRb in Growth Suppression by Sphingosine" *Biochemistry*, vol. 34 pp. 1885-1892, 1995

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.