

PRODUCT DATA SHEET

Lyso-Dihydrosphingomyelin

Catalog number: 1913

Common names:

Dihydrosphingosylphosphorylcholine
(mixture of D-erythro and L-threo
isomers)

Source: semisynthetic, bovine buttermilk

Solubility: chloroform/methanol, 2:1

CAS number: N/A

Molecular Formula: C₂₃H₅₁N₂O₅P

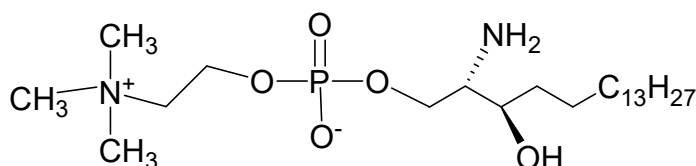
Molecular Weight: 467

Storage: -20°C

Purity: TLC >98%; identity confirmed by MS

TLC System: chloroform/methanol/2.5N
ammonium hydroxide (60:40:9)

Appearance: solid



Application Notes:

This product is the saturated form of the more plentiful sphingomyelin, a lipid that is found in mammalian cell membranes, especially in the membranes of the myelin sheath, and is the most abundant sphingolipid in mammals. Dihydrosphingomyelin has been identified as a minor lipid component in many mammalian tissues but has recently been reported to be present in significant amounts in bovine brain and bovine milk.¹ It is also found in much greater amounts in human lens membranes (half of all the phospholipids) where it has a critical role in ocular function and perhaps in age-related nuclear cataracts.² However, dihydrosphingomyelin has been reported to occur only in small amounts in the lens membranes of other mammals. Dihydrosphingomyelin demonstrates good mixing properties with both sterols and sphingomyelin indicating that it could function as a membrane organizer and this may be the reason it is present in large amounts in human lens membranes where cholesterol is also enriched.³ The enzyme sphingomyelinase is active towards dihydrosphingomyelin and readily converts it to dihydroceramide. Recent evidence has been presented that indicates that dihydrosphingomyelin impairs HIV-1 infection by rigidifying liquid-ordered membrane domains, a finding that could have great potential in providing a therapeutic treatment for this debilitating disease.⁴

Selected References:

1. Wm. Byrdwell and R. Perry "Liquid chromatography with dual parallel mass spectrometry and 31P nuclear magnetic resonance spectroscopy for analysis of sphingomyelin and dihydrosphingomyelin II. Bovine milk sphingolipids" *Journal of Chromatography A*, vol. 1146 pp. 164-185, 2007
2. J. Deeley et al. "Sphingolipid distribution changes with age in the human lens" *Journal of Lipid Research*, vol. 51 pp. 2753-2760, 2010
3. T. Nyholm et al. "A Calorimetric Study of Binary Mixtures of Dihydrosphingomyelin and Sterols, Sphingomyelin, or Phosphatidylcholine" *Biophysical Journal*, vol. 84 pp. 3138-3146, 2003
4. C. Vieira et al. "Dihydrosphingomyelin Impairs HIV-1 Infection by Rigidifying Liquid-Ordered Membrane Domains" *Chemistry and Biology*, vol. 17 pp. 766-775, 2010

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.