

PRODUCT DATA SHEET

Monosialoganglioside GM₂ (Na⁺ salt), bovine

Catalog No: 1554

Common Name: GM₂

Source: semisynthetic, bovine

Solubility: chloroform/methanol/DI water,
(2:1:0.1); forms micellar solution in
water

CAS No: 19600-01-2

Molecular Formula: C₆₇H₁₂₁N₃O₂₆ • Na

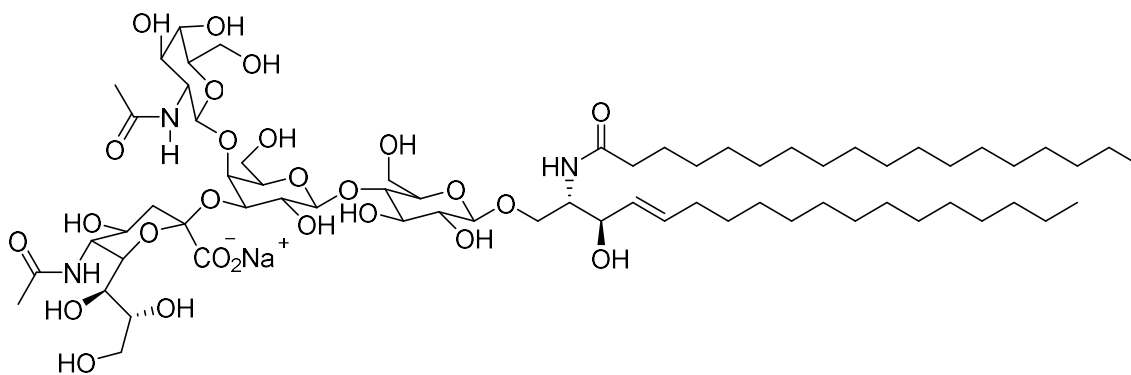
Molecular Weight: 1385+ Na (stearoyl)

Storage: -20°C

Purity: TLC >98%; identity confirmed by MS

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

Appearance: solid



Application notes:

Gangliosides¹ are acidic glycosphingolipids that form lipid rafts in the outer leaflet of the cell plasma membrane, especially in neuronal cells in the central nervous system.² They participate in cellular proliferation, differentiation, adhesion, signal transduction, cell-to-cell interactions, tumorigenesis, and metastasis.³ GM₂ regulates the function of ciliary neurotrophic factor receptors. The accumulation of GM₂ (due to a deficiency in *beta*-hexosaminidase) has characterized Tay-Sachs disease (due to a mutation in the gene *HEXA*) and Sandhoff disease (due to a mutation in the gene *HEXB*). A mutation in the *GM2A* gene results in GM₂ activator deficiency that also leads to accumulation of GM₂.⁴

Selected References:

1. L. Svennerholm, et al. (eds.), *Structure and Function of Gangliosides*, New York, Plenum, 1980
2. T. Kolter, R. Proia, K. Sandhoff "Combinatorial Ganglioside Biosynthesis" *J. Biol. Chem.*, Vol. 277, No. 29, pp. 25859-25862, 2002
3. S. Birkle, G. Zeng, L. Gao, R.K. Yu, and J. Aubry "Role of tumor-associated gangliosides in cancer progression" *Biochimie*, Vol. 85 pp. 455-463, 2003
4. R. Gravel et al., *The Metabolic and Molecular Bases of Inherited Disease* (C. R. Scriver, W. S. Sly, B. Childs, A. L. Beaudet, D. Valle, K. W. Kinzler, and B. Vogelstein, eds) pp. 3827-3876, McGraw-Hill Inc., New York, 2001

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.