

PRODUCT DATA SHEET

N-Hexanoyl-biotin-disialoganglioside GD₃

Catalog number: 2055

Synonyms: Biotin-C6:0-GD₃

Source: semisynthetic, bovine buttermilk

Solubility: chloroform/methanol/DI water,
2:1:0.1

CAS number: N/A

Molecular Formula: C₆₈H₁₁₆N₆O₃₁S • 2NH₃

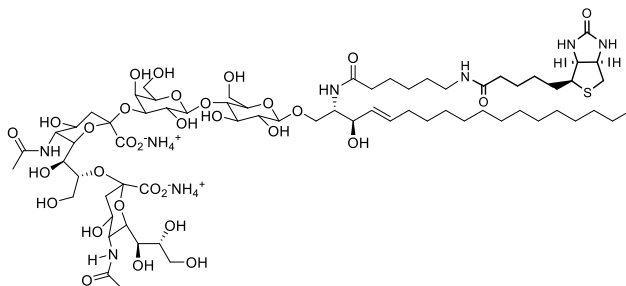
Molecular Weight: 1546 + 2NH₃

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 ganglioside GD₃ analogue contains a biotin unit attached to the amine of the sphingosine moiety via a hexanoic acid linker and is ideal for use in ganglioside studies. The biotin structure allows for attachment of the ganglioside to streptavidin and avidin substrates making it extremely useful for binding to substrates and for toxin detection.¹

Gangliosides are acidic glycosphingolipids containing one or more sialic acids that generally form lipid rafts in the outer leaflet of the cell plasma membrane, especially in neuronal cells in the central nervous system.^{2,3} They participate in many cellular activities including proliferation, differentiation, adhesion, signal transduction, cell-to-cell interactions, tumorigenesis, and metastasis.⁴ The accumulation of gangliosides has been linked to several diseases including Tay-Sachs and Sandhoff disease while an autoimmune response against gangliosides can lead to Guillain-Barre syndrome. Gangliosides act as receptors for various toxins and bacteria, accumulate in various tumors, and aid in many neuronal functions.

Disialoganglioside GD₃ is predominantly expressed during neuronal development and its expression becomes very limited in adult tissues. GD₃ expression is unusually high in basal cell carcinomas and malignant melanomas and is thought to be a human melanoma-specific antigen. Although GD₃ is not immunogenic it has been investigated as a tool for immunotargeting human melanoma cells.⁴ Over expression of GD₃ has led to apoptosis by recruiting mitochondria to apoptotic pathways and suppressing NF-κB activation and subsequent κB-dependent gene induction.⁵ Increased levels of GD₃ have also been found to be associated with proliferative diseases, such as atherosclerosis.

Selected References:

1. A. Pukin et al. Chemoenzymatic synthesis of biotin-appended analogues of gangliosides GM₂, GM₁, GD_{1a} and GalNAc-GD_{1a} for solid-phase applications and improved ELISA tests. *Org. Biomol. Chem.*, 9(16):5809-5815, 2011
2. L. Svennerholm, et al. (eds.), *Structure and Function of Gangliosides*, New York, Plenum, 1980
3. T. Kolter, R. Proia, K. Sandhoff, Combinatorial Ganglioside Biosynthesis. *J. Biol. Chem.*, July Vol. 277, No. 29, pp. 25859-25862, 2002
4. H. Jennings et al. "Bioengineering of Surface GD₃ Ganglioside for Immunotargeting Human Melanoma Cells" *Journal of Biological Chemistry*, Vol. 279:24 pp. 25390, 2004
5. J. Fernández-Checa et al. "Ganglioside GD₃ Sensitizes Human Hepatoma Cells to Cancer Therapy" *Journal of Biological Chemistry*, Vol. 277:51 pp. 49870, 2002

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.