

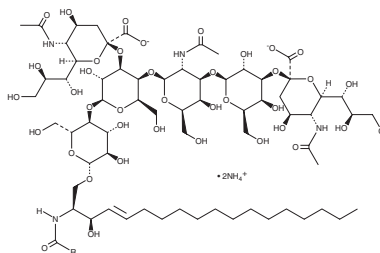
# PRODUCT INFORMATION



## Ganglioside G<sub>D1a</sub> (bovine) (ammonium salt)

Item No. 31591

CAS Registry No.: 12707-58-3  
Synonym: Disialoganglioside G<sub>D1a</sub>  
MF: C<sub>84</sub>H<sub>146</sub>N<sub>4</sub>O<sub>39</sub> • 2NH<sub>4</sub> (for stearoyl)  
FW: 1,872.2  
Purity: ≥98%  
Supplied as: A solid  
Storage: -20°C  
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Ganglioside G<sub>D1a</sub> (bovine) (ammonium salt) is supplied as a solid. A stock solution may be made by dissolving the ganglioside G<sub>D1a</sub> (bovine) (ammonium salt) in the solvent of choice, which should be purged with an inert gas. Ganglioside G<sub>D1a</sub> (bovine) (ammonium salt) is soluble in a 2:1:0.1 solution of chloroform:methanol:DI water.

### Description

Ganglioside G<sub>D1a</sub> is a sialic acid-containing glycosphingolipid that has been found in brain, erythrocytes, bone marrow, testis, spleen, and liver, as well as in serum lipoproteins.<sup>1,2</sup> It is shed into the tumor microenvironment from the surface of tumor cells, where it influences tumor-host cell interactions to promote tumor cell proliferation, invasion, and metastasis.<sup>3</sup> It functions as a toll-like receptor 2 (TLR2) co-receptor in isolated human monocytes, colocalizing with TLR2 and enhancing the binding of type IIb *E. coli* enterotoxin (LT-IIb-B<sub>5</sub>) to TLR2 when used at a concentration of 20 μM.<sup>4</sup> Ganglioside G<sub>D1a</sub> (20 μM) increases VEGF-induced proliferation of human umbilical vein endothelial cells (HUVECs), as well as induces vascularization in a Matrigel™ plug assay in mice when used at a concentration of 20 μM/plug.<sup>3</sup> It accumulates in the CNS of patients with galactosialidosis, a lysosomal storage disorder.<sup>5</sup> This product contains ganglioside G<sub>D1a</sub> molecular species with primarily C18:0 fatty acyl chain lengths. As this product is derived from a natural source, there may be variations in the sphingoid backbone.

### References

1. Iwamori, M., Shimomura, J., Tsuyuhara, S., *et al.* Gangliosides of various rat tissues: Distribution of ganglio-N-tetraose-containing gangliosides and tissue-characteristic composition of gangliosides. *J. Biochem.* **95**(3), 761-770 (1984).
2. Senn, H.J., Orth, M., Fitzke, E., *et al.* Gangliosides in normal human serum. Concentration, pattern and transport by lipoproteins. *Eur. J. Biochem.* **181**(3), 657-662 (1989).
3. Mukherjee, P., Faber, A.C., Shelton, L.M., *et al.* Thematic review series: Sphingolipids. Ganglioside GM3 suppresses the proangiogenic effects of vascular endothelial growth factor and ganglioside GD1a. *J. Lipid Res.* **49**(5), 929-938 (2008).
4. Liang, S., Wang, M., Tapping, R.I., *et al.* Ganglioside GD1a is an essential coreceptor for Toll-like receptor 2 signaling in response to the B subunit of type IIb enterotoxin. *J. Biol. Chem.* **282**(10), 7532-7542 (2007).
5. Walkley, S.U. Secondary accumulation of gangliosides in lysosomal storage disorders. *Semin. Cell Dev. Biol.* **15**(4), 433-444 (2004).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/12/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897  
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM