

PRODUCT INFORMATION



Disialoganglioside Mixture

Item No. 29360

Supplied as: A solution in chloroform:methanol:water (2:1:0.1)

Storage: -20°C

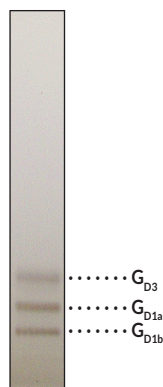
Stability: ≥2 years

Concentration: 500 µg/ml

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

Description and Data

Disialogangliosides are membrane sphingolipids that contain two sialic acid residues linked to a galactose moiety.¹ They have roles in cell proliferation, immune regulation, tumor cell invasion and metastasis, intra- and intercellular signaling, and neurogenesis.¹⁻⁵ Disialoganglioside Mixture contains the disialogangliosides G_{D1a} (Item No. 31591), G_{D1b} (Item No. 31708), and G_{D3} (Item No. 31709) and is intended for use as a reference standard for the detection of these disialogangliosides in various sample types by TLC or LC-MS. This mixture is supplied in an amber ampule in which the headspace has been purged with argon to prevent lipid oxidation.



References

1. 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).
2. 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).
3. Kolter, T. Ganglioside biochemistry. *ISRN Biochem.* (2012).
4. De Maria, R., Lenti, L., Malisan, F., *et al.* Requirement for G_{D3} ganglioside in CD95- and ceramide-induced apoptosis. *Science* **277**(5332), 1652-1655 (1997).
5. Wang, J., Cheng, A., Wakade, C., *et al.* Ganglioside GD3 is required for neurogenesis and long-term maintenance of neural stem cells in the postnatal mouse brain. *J. Neurosci.* **34**(41), 13790-13800 (2014).

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, 08/07/2026

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