

PRODUCT INFORMATION



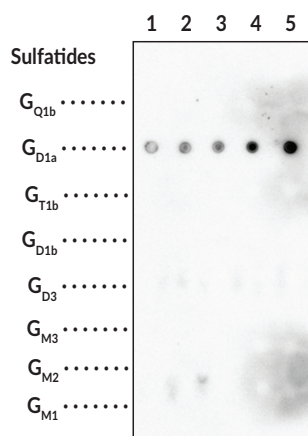
Ganglioside G_{D1a} Monoclonal Antibody (Clone TBG3)

Item No. 38296

Overview and Properties

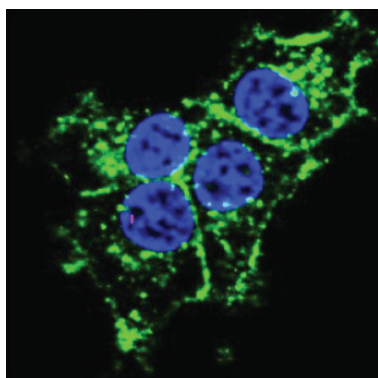
Contents:	This vial contains 100 µg of protein A-purified monoclonal antibody.
Synonyms:	Disialoganglioside G _{D1a} , Ganglioside B ₁ , Ganglioside G ₃
Immunogen:	G _{D1a}
Cross Reactivity:	(+) Ganglioside G _{D1a} ; (-) Ganglioside G _{D1b} , Ganglioside G _{D3} , Ganglioside G _{M1} , Ganglioside G _{M2} , Ganglioside G _{M3} , Ganglioside G _{Q1b} , Sulfatides
Species Reactivity:	Species Independent
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
Clone:	TBG3
Host:	Mouse
Isotype:	IgG3
Applications:	Dot blot, ELISA, Immunofluorescence (IF), and Thin Layer Chromatography (TLC) Immunostaining; the recommended starting dilution for dot blot is 1:500-1:1,000, 1:1,000 for ELISA, 1:10-1:50 for IF, and 1:500 for TLC immunostaining. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Lane 1: 10 ng
Lane 2: 25 ng
Lane 3: 50 ng
Lane 4: 100 ng
Lane 5: 200 ng

Dot blot against gangliosides using Ganglioside G_{D1a} Monoclonal Antibody (Clone TBG3).¹



PC12 cells labeled with Ganglioside G_{D1a} Monoclonal Antibody (Clone TBG3) were incubated at 37°C for 30 min.¹

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

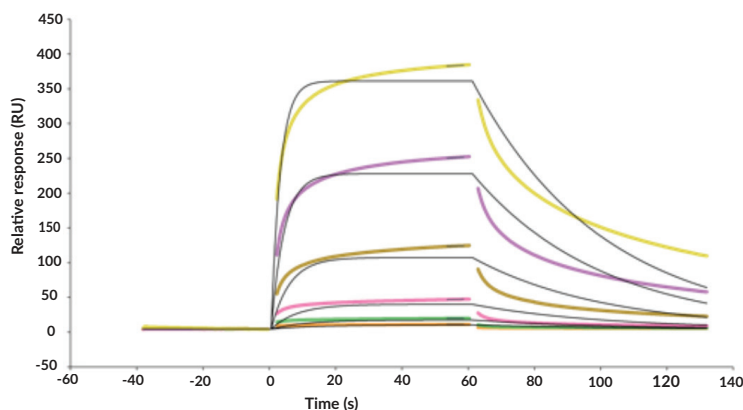
SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your 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, 08/16/2024

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

PRODUCT INFORMATION



Representative image: Ganglioside GD_{1a} Specifically Binds Ganglioside GD_{1a} Monoclonal Antibody (Clone TBG3). Ganglioside GD_{1a} Monoclonal Antibody (Clone TBG3) was immobilized on a Series S Sensor Chip CM5 and SPR analysis was used to determine Ganglioside GD_{1a} (bovine) (ammonium salt) (Item No. 31591) binding affinity on a Biacore 8K, using multi cycle kinetics with six concentrations of Ganglioside GD_{1a} .

Description

Ganglioside G_{D1a} (Item Nos. 31591 | 15585 | 31707) 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.^{2,3} It is formed by the transfer of N-acetylneuraminic acid (Neu5Ac; Item No. 16091) to the precursor ganglioside G_{M1a} by CMP-N-acetylneuraminat- β -galactosamide- α -2,3-sialyltransferase 2 (ST3Gal-II).⁴ 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-B₅) to TLR2.⁵ Ganglioside G_{D1a} also mediates polyomavirus entry into host cells.⁶ Ganglioside G_{D1a} 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, as well as increases VEGF-induced proliferation of human umbilical vein endothelial cells (HUVECs). A monoclonal antibody targeting ganglioside G_{D1a} and ganglioside G_{T1b} (Item Nos. 31712 | 15588 | 31592) inhibits axon regeneration in a mouse model of sciatic nerve crush-induced peripheral nervous system injury.⁷ Ganglioside G_{D1a} accumulates in the CNS of patients with galactosialidosis, a lysosomal storage disorder.⁸ Cayman's Ganglioside G_{D1a} Monoclonal Antibody (Clone TBG3) binds to Ganglioside G_{D1a} (bovine brain) (Item No. 15585) with k_a , k_d , and K_D values of $26,300 \text{ M}^{-1}\text{s}^{-1}$, 0.0342 s^{-1} , and $1.36 \text{ }\mu\text{M}$, respectively, as determined by surface plasmon resonance (SPR). It can be used for dot blot, ELISA, immunofluorescence (IF), and TLC immunostaining applications.

References

1. Fewou, S.N., Rupp, A., Nickolay, L.E., et al. *J. Clin. Invest.* **122**(3), 1037-1051 (2012).
2. Iwamori, M., Shimomura, J., Tsuyuhara, S., et al. *J. Biochem.* **95**(3), 761-770 (1984).
3. Senn, H.J., Orth, M., Fitzke, E., et al. *Eur. J. Biochem.* **181**(3), 657-662 (1989).
4. Cavdarli, S., Groux-Degroote, S., and Delannoy, P. *Biomolecules* **9**(8), 311 (2019).
5. Liang, S., Wang, M., Tapping, R.I., et al. *J. Biol. Chem.* **282**(10), 7532-7542 (2007).
6. Gilbert, J. and Benjamin, T. *J. Virol.* **78**(22), 12259-12267 (2004).
7. Lehmann, H.C., Lopez, P.H.H., Zhang, G., et al. *J. Neurosci.* **27**(1), 27-34 (2007).
8. Walkley, S.U. *Semin. Cell Dev. Biol.* **15**(4), 433-444 (2004).

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