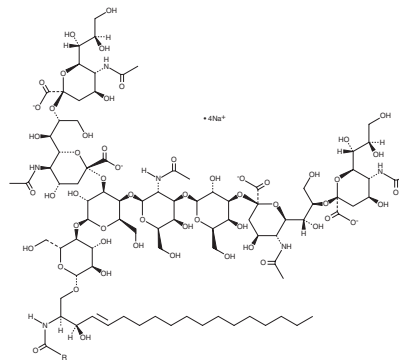


PRODUCT INFORMATION

Ganglioside G_{Q1b} Mixture (sodium salt) Item No. 15589

CAS Registry No.: 68652-37-9
MF: C₁₀₆H₁₇₈N₆O₅₅ • 4Na
 (for stearoyl)
FW: 2,508.5
Purity: ≥98%
Supplied as: A lyophilized solid
Storage: -20°C
Stability: ≥4 years
Special Conditions: Hygroscopic. Protect from moisture.



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

Laboratory Procedures

Ganglioside G_{Q1b} mixture (sodium salt) is supplied as a lyophilized solid. A stock solution may be made by dissolving the ganglioside G_{Q1b} mixture (sodium salt) in the solvent of choice. Ganglioside G_{Q1b} mixture (sodium salt) is soluble in 2:1 chloroform:methanol, which should be purged with an inert gas. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Ganglioside G_{Q1b} is a tetrasialoganglioside that contains two sialic acid residues linked to an inner galactose unit. It stimulates phosphorylation of several ecto-type protein kinase substrates on the surface of GOTO human neuroblastoma cells when used at a concentration of 5 nM.¹ Ganglioside G_{Q1b} promotes differentiation of murine embryonic stem cells (mESCs) to neuronal precursor and glial cells *via* activation of the ERK1/2 pathway.² It also induces differentiation of murine keratinocytes through phosphoinositide turnover.³ Ganglioside G_{Q1b} mixture contains ganglioside G_{Q1b} molecular species with C18:1 and C20:1 sphingoid backbones.

References

1. Tsuji, S., Yamashita, T., and Nagai, Y. A novel, carbohydrate signal-mediated cell surface protein phosphorylation: Ganglioside GQ1b stimulates ecto-protein kinase activity on the cell surface of a human neuroblastoma cell line, GOTO. *J. Biochem.* **104(4)**, 498-503 (1988).
2. Kwak, D.H., Jin, J.W., Ryu, J.S., *et al.* Regulatory roles of ganglioside GQ1b in neuronal cell differentiation of mouse embryonic stem cells. *BMB Rep.* **44(12)**, 799-804 (2011).
3. Yada, Y., Okano, Y., and Nozawa, Y. Ganglioside GQ1b-induced terminal differentiation in cultured mouse keratinocytes. Phosphoinositide turnover forms the onset signal. *Biochem. J.* **279(Pt 3)**, 665-670 (1991).

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, 11/11/2022

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