PRODUCT INFORMATION



C2 Galactosylceramide (d18:1/2:0)

Item No. 24465

CAS Registry No.:	35823-61-1	
Formal Name:	N-[(1S,2R,3E)-1-[(β-D-	H OH
	galactopyranosyloxy)methyl]-2- hydroxy-3-heptadecen-1-yl]-acetamide	
Synonyms:	Galactosylceramide (d18:1/2:0),	
	GalCer(d18:1/2:0), β-Gal-C2-Cer,	
	N-acetyl Psychosine	HO.
MF:	$C_{26}H_{49}NO_8$	o o
FW:	503.7	
Purity:	≥98%	HO
Supplied as:	A solid	
Storage:	-20°C	OH
Stability:	≥4 years	

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

Laboratory Procedures

C2 Galactosylceramide (d18:1/2:0) is supplied as a solid. A stock solution may be made by dissolving the C2 galactosylceramide (d18:1/2:0) in the solvent of choice. C2 Galactosylceramide (d18:1/2:0) is soluble in organic solvents such as chloroform, ethanol, and methanol, which should be purged with an inert gas.

Description

C2 Galactosylceramide is a bioactive sphingolipid.¹⁻³ It inhibits IL-4 production by 23.24% in EL4 T cells when used at a concentration of 10 μ M.¹ C2 Galactosylceramide stimulates NF- κ B activity in C6 glioma cells and activates mast cells in a concentration-dependent manner. It is cytotoxic to U937 cells in vitro $(IC_{50} = 33 \mu M)$. C2 Galactosylceramide also suppresses the allogeneic mixed leukocyte reaction (MLR) response in murine spleen cells in a concentration-dependent manner.³

References

- 1. Park, J., Li, Q., Chang, Y.T., et al. Inhibitory activity of a ceramide library in interleukin-4 production from activated T cells. Bioorg. Med. Chem. 13(7), 2589-2595 (2005).
- 2. Chang, Y.-T., Choi, J., Ding, S., et al. The synthesis and biological characterization of a ceramide library. J. Am. Chem. Soc. 124(9), 1856-1857 (2002).
- 3. Sakai, T., Ueno, H., Natori, T., et al. Effects of α and β -galactosylated C2-ceramides on the immune system. J. Med. Chem. 41(4), 650-652 (1998).

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

SAFFTY 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.

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, 10/14/2022

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