

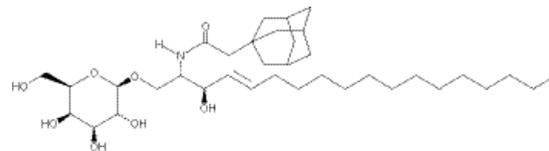
PRODUCT INFORMATION



C2 Adamantanyl Galactosylceramide (d18:1/2:0)

Item No. 24472

CAS Registry No.: 574738-16-2
Formal Name: N-[(1S,2R,3E)-1-[(β-D-galactopyranosyloxy)methyl]-2-hydroxy-3-heptadecen-1-yl]-tricyclo[3.3.1.1^{3,7}]decane-1-acetamide
Synonyms: AdaGalCer(d18:1/2:0), Admantanyl Galactosylceramide (d18:1/2:0), Admantanyl GalCer(d18:1/2:0), N-(1-adamantaneacetyl) Galactosylceramide
MF: C₃₆H₆₃NO₈
FW: 637.9
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

C2 Adamantanyl galactosylceramide (d18:1/2:0) (AdaGalCer(d18:1/2:0)) is supplied as a solid. A stock solution may be made by dissolving the AdaGalCer(d18:1/2:0) in the solvent of choice. AdaGalCer(d18:1/2:0) is soluble in chloroform, methanol, and a 9:1 solution of chloroform:methanol.

Description

AdaGalCer is a bioactive sphingolipid.¹ It reduces globotriaosylceramide 3 synthesis from exogenous lactosylceramide in microsomes. AdaGalCer stimulates recombinant glucocerebrosidase activity in a pH-dependent manner. It activates glucocerebrosidase to decrease glucosylceramide accumulation in fibroblasts and lymphoblasts isolated from patients with Gaucher and Fabry disease, respectively.

Reference

1. Kamani, M., Mylvaganam, M., Tian, R., *et al.* Adamantyl glycosphingolipids provide a new approach to the selective regulation of cellular glycosphingolipid metabolism. *J. Biol. Chem.* **286**(24), 21413-21426 (2011).

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, 08/11/2025

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