

PRODUCT INFORMATION



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

Item No. 24623

Formal Name: 2-((3S,5S,7S)-adamantan-1-yl)-N-((2S,3R,4E,8E)-3-hydroxy-1-(((2R,3R,4S,5S,6R)-3,4,5-trihydroxy-6-(hydroxymethyl)tetrahydro-2H-pyran-2-yl)oxy)octadeca-4,8-dien-2-yl)acetamide

Synonyms: N-(1-Adamantaneacetyl) Glucosylceramide, Adamantanyl GluCer(d18:1/2:0), Adamantanyl Glucosylceramide (d18:1/2:0), AdaGluCer(d18:1/2:0)

MF: C₃₆H₆₁NO₈

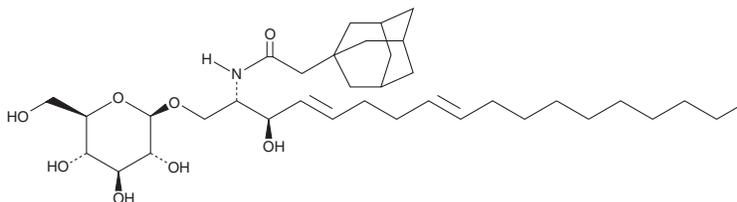
FW: 635.9

Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

C2 Adamantanyl glucosylceramide (d18:1/2:0) (AdaGluCer) is supplied as a solid. A stock solution may be made by dissolving the AdaGluCer in the solvent of choice, which should be purged with an inert gas. AdaGluCer is soluble in chloroform and methanol and a 9:1 solution of chloroform:methanol.

Description

AdaGluCer is a bioactive sphingolipid.¹ It increases levels of glycosphingolipids in Vero cells at low concentrations (≤20 μM) but decreases them at high concentrations (≥40 μM). Pretreatment of microsomes with AdaGluCer increases production of adamantanyl lactosylceramide (AdaLacCer) and reduces synthesis of LacCer from exogenous GlcCer. AdaGluCer inhibits glucocerebrosidase activity *in vitro* in a pH-dependent manner.

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, 12/17/2022

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