

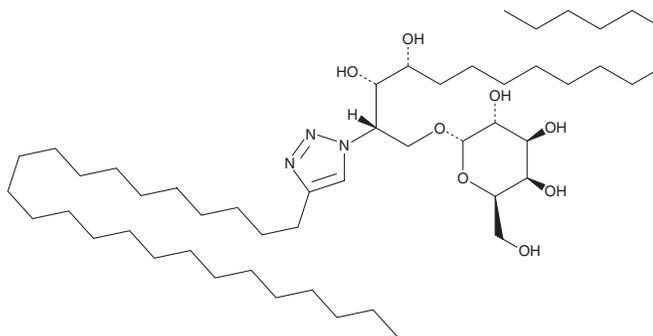
PRODUCT INFORMATION



α -GalCer analog 8

Item No. 26660

CAS Registry No.: 922727-14-8
Formal Name: (2S,3S,4R)-3,4-dihydroxy-2-(4-tetracosyl-1H-1,2,3-triazol-1-yl) octadecyl, α -D-galactopyranoside
Synonym: α -Galactosylceramide analog 8
MF: C₅₀H₉₇N₃O₈
FW: 868.3
Purity: \geq 95%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

α -Galactosylceramide analog 8 (α -GalCer analog 8) is supplied as a solid. A stock solution may be made by dissolving the α -GalCer analog 8 in the solvent of choice. α -GalCer analog 8 is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide, which should be purged with an inert gas.

Description

α -GalCer analog 8 is a triazole derivative of α -galactosylceramide.¹ It increases IL-2 secretion by DN32.D3 NKT hybridoma cells when co-cultured with CD1d-transfected RBL cells pre-loaded with α -GalCer analog 8 at a concentration of 32 ng/ml. α -GalCer analog 8 (32 ng/ml) induces IL-4 secretion to a greater extent than the synthetic α -GalCer KRN 7000 (Item No. 11208) in mouse splenocytes *in vitro* and in mouse serum following administration of a 1 μ g per animal dose, indicating a Th2 response.

Reference

1. Lee, T., Cho, M., Ko, S.-Y., *et al.* Synthesis and evaluation of 1,2,3-triazole containing analogues of the immunostimulant α -GalCer. *J. Med. Chem.* **50**(3), 585-589 (2007).

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, 12/20/2022

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