

PRODUCT INFORMATION



Tri-GalNAc COOH

Item No. 40934

CAS Registry No.: 1953146-81-0

Formal Name: N-[3-[[[5-[[2-(acetylamino)-2-deoxy-β-D-galactopyranosyl]oxy]-1-oxopentyl]amino]propyl]-6,6-bis[[[3-[[[5-[[2-(acetylamino)-2-deoxy-β-D-galactopyranosyl]oxy]-1-oxopentyl]amino]propyl]amino]-3-oxopropoxy]methyl]-25-hydroxy-8-oxo-4,11,14,17,20,23-hexaoxa-7-azapentacosanamide

Synonym: SGN-46810

MF: C₇₅H₁₃₄N₁₀O₃₅

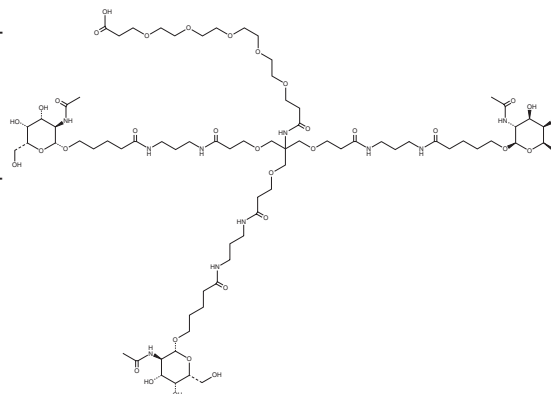
FW: 1,735.9

Purity: ≥95%

Supplied as: A solution in ethanol

Storage: -80°C

Stability: ≥4 years



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

Laboratory Procedures

Tri-GalNAc COOH is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Tri-GalNAc COOH is sparingly soluble (1-10 mg/ml) in methanol and chloroform purged with an inert gas can be used.

Description

Tri-GalNAc COOH is an intermediate in the synthesis of lysosome-targeting chimeras (LYTACs) containing a triad of the galactose amino sugar derivative GalNAc (Item No. 31728) connected by acyl-amino chains to an ether linker with a terminal carboxylic acid that can be functionalized with antibodies or biotin.¹ It has been used in the synthesis of LYTACs functionalized with anti-EGFR antibodies that reduce EGFR levels without affecting total protein levels in HepG2 hepatocellular carcinoma cells.

Reference

1. Zhou, Y., Teng, P., Montgomery, N.T., *et al.* Development of triantennary N-acetylgalactosamine conjugates as degraders for extracellular proteins. *ACS Cent. Sci.* **7**(3), 499-506 (2021).

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, 05/11/2026

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