

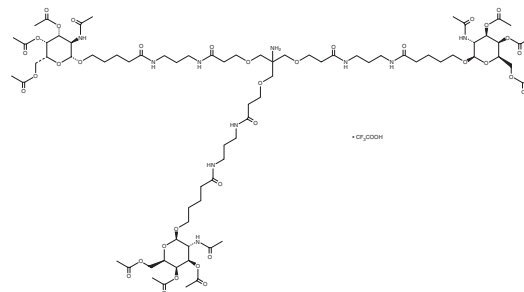
PRODUCT INFORMATION



Tri-GalNAc(OAc)3 (trifluoroacetate salt)

Item No. 40846

CAS Registry No.: 1159408-65-7
Formal Name: 6-amino-11,17-dioxo-6-[[[3-oxo-3-[[[3-[[[1-oxo-5-[[[3,4,6-tri-O-acetyl-2-(acetylamino)-2-deoxy-β-D-galactopyranosyl]oxy]pentyl]amino]propyl]amino]propoxy]methyl]-N-[3-[[[1-oxo-5-[[[3,4,6-tri-O-acetyl-2-(acetylamino)-2-deoxy-β-D-galactopyranosyl]oxy]pentyl]amino]propyl]-21-[[[3,4,6-tri-O-acetyl-2-(acetylamino)-2-deoxy-β-D-galactopyranosyl]oxy]-4,8-dioxo-12,16-diazaheneicosanamide, trifluoroacetate salt
MF: C₇₉H₁₂₈N₁₀O₃₆ • CF₃COOH
FW: 1,907.9
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

Tri-GalNAc(OAc)3 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the tri-GalNAc(OAc)3 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Tri-GalNAc(OAc)3 (trifluoroacetate salt) is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of tri-GalNAc(OAc)3 (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. Tri-GalNAc(OAc)3 (trifluoroacetate salt) is soluble (≥10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Tri-GalNAc(OAc)3 is an intermediate in the synthesis of lysosome-targeting chimeras (LYTACs) with anticancer activity.¹

Reference

1. Ahn, G., Banik, S.M., Miller, C.L., *et al.* LYTACs that engage the asialoglycoprotein receptor for targeted protein degradation. *Nat. Chem. Biol.* **17**(9), 937-946 (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, 06/03/2024

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