

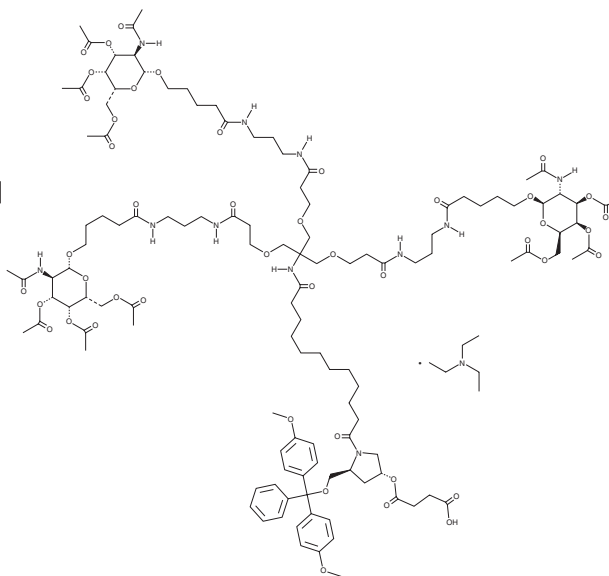
PRODUCT INFORMATION



GalNac-L96 (triethylamine salt)

Item No. 40456

CAS Registry No.: 1159408-62-4
Formal Name: butanedioic acid, 1-[[[3R,5S)-5-[[[bis(4-methoxyphenyl)phenylmethoxy]methyl]-1-[1,12,19,25-tetraoxo-14,14-bis[[[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]-29-[[[3,4,6-tri-O-acetyl-2-(acetylamino)-2-deoxy-β-D-galactopyranosyl]oxy]-16-oxa-13,20,24-triazanonacos-1-yl]-3-pyrrolidinyl] ester, compd. with N,N-diethylethanamine
Synonyms: N-Acetylgalactosamine-L96, N-acetyl-D-Galactosamine-L96
MF: C₁₂₁H₁₇₉N₁₁O₄₅ • C₆H₁₅N
FW: 2,609.0
Purity: ≥95%
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

GalNac-L96 (triethylamine salt) is supplied as a solid. A stock solution may be made by dissolving the GalNac-L96 (triethylamine salt) in the solvent of choice, which should be purged with an inert gas. GalNac-L96 (triethylamine salt) is soluble in acetonitrile.

Description

GalNac-L96 is a triantennary N-acetylgalactosamine (GalNac) asialoglycoprotein receptor (ASGPR) ligand.¹ It has been used in the synthesis of siRNA-GalNac conjugates for RNA interference-mediated gene silencing *in vitro* and in mice.

Reference

1. Nair, J.K., Willoughby, J.L., Chan, A., *et al.* Multivalent N-acetylgalactosamine-conjugated siRNA localizes in hepatocytes and elicits robust RNAi-mediated gene silencing. *J. Am. Chem. Soc.* **136**(49), 16958-16961 (2014).

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, 06/19/2024

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