

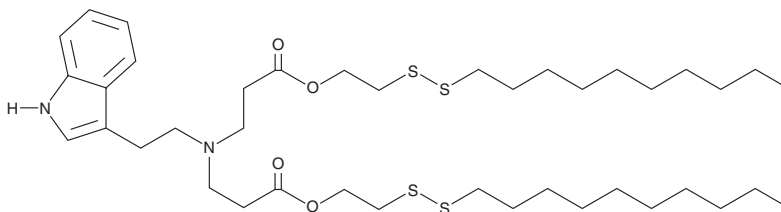
PRODUCT INFORMATION



NT1-O14B

Item No. 37095

CAS Registry No.: 2739805-64-0
MF: $C_{40}H_{68}N_2O_4S_4$
FW: 769.2
Purity: $\geq 98\%$
Supplied as: A 50 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥ 2 years



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

Description

NT1-O14B is a tryptamine-containing ionizable cationic lipidoid.¹ It has been used in combination with other lipids in the formation of lipid nanoparticles (LNPs). Intravenous administration of LNPs containing NT1-O14B and the cationic lipidoid 306-O12B-3 (Item No. 37096) encapsulating antisense oligonucleotides (ASOs) against tau decreases tau brain levels in mice.

Reference

1. Ma, F., Yang, L., Sun, Z., *et al.* Neurotransmitter-derived lipidoids (NT-lipidoids) for enhanced brain delivery through intravenous injection. *Sci. Adv.* **6**(30), eabb4429 (2020).

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, 11/06/2025

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