

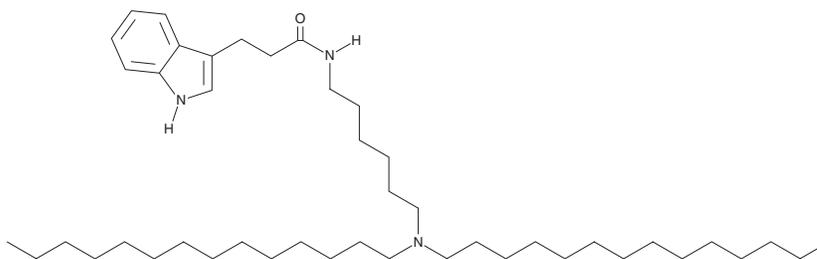
PRODUCT INFORMATION



Lipid TD5

Item No. 44285

CAS Registry No.: 2855180-86-6
Formal Name: N-[6-(ditetradecylamino)hexyl]-1H-indole-3-propanamide
MF: C₄₅H₈₁N₃O
FW: 680.2
Purity: ≥90%
Supplied as: A 10 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥4 years



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

Description

Lipid TD5 is an ionizable cationic lipid ($pK_a = 7.3$) that has been used in the generation of lipid nanoparticles (LNPs) for the delivery of mRNA *in vitro* and *in vivo*.¹ LNPs containing lipid TD5 undergo endocytosis in a serotonin (5-HT) receptor subtype 1A-dependent manner in primary mouse astrocytes. Intrathecal administration of LNPs containing lipid TD5 selectively traffic to the brain over the liver and spleen in mice.

Reference

1. Xue, Y., Wang, C., Li, H., *et al.* Lipid nanoparticles enhance mRNA delivery to the central nervous system upon intrathecal injection. *Adv. Mater.* **37**(27), e2417097 (2025).

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, 09/22/2025

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