

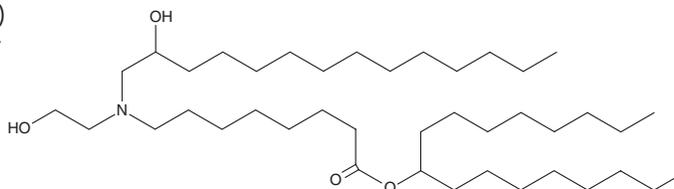
PRODUCT INFORMATION



Lipid HTO12

Item No. 41202

CAS Registry No.: 3047222-45-4
Formal Name: 8-[(2-hydroxyethyl)(2-hydroxytetradecyl) amino]-octanoic acid, 1-octylnonyl ester
MF: C₄₁H₈₃NO₄
FW: 654.1
Purity: ≥98%
Supplied as: A 50 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 HTO12 is an ionizable cationic lipid ($pK_a = 6.432$) that has been used in combination with other lipids in the formation of lipid nanoparticles (LNPs) for the delivery of mRNA *in vitro* and *in vivo*.¹ LNPs containing lipid HTO12 and encapsulating adenine base editor mRNA targeting *Pcsk9* reduce hepatic *Pcsk9* levels in mice at a reduced ionizable lipid/total lipids-to-mRNA ratio than LNPs containing SM-102 (Item No. 33474).

Reference

1. Chen, Q., Wang, X., Zhang, Y., *et al.* Minimizing the ratio of ionizable lipid in lipid nanoparticles for *in vivo* base editing. *Natl. Sci. Rev.* **11(6)**, nwae135 (2024).

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, 11/21/2025

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