

PRODUCT INFORMATION



ATX-001

Item No. 39037

CAS Registry No.: 1777792-33-2
Formal Name: 8,8'-[[2-[[2-(dimethylamino)ethyl]thio]acetyl]imino]bis-octanoic acid, 1,1'-di-(2Z)-2-nonen-1-yl ester

MF: C₄₀H₇₄N₂O₅S

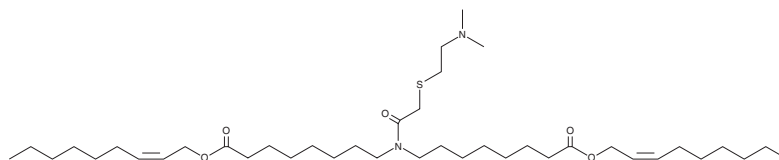
FW: 695.1

Purity: ≥98%

Supplied as: A 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.

Laboratory Procedures

ATX-001 is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. ATX-001 is soluble in the organic solvent DMSO.

Description

ATX-001 is an ionizable cationic lipid.¹ It has been used in combination with other lipids in the formation of lipid nanoparticles (LNPs) for the delivery of mRNA *in vivo*.

Reference

1. Karmali, P., Chivukula, P., Payne, J.E., *et al.* Method of making lipid-encapsulated RNA nanoparticles. *Arcturus Therapeutics*. **WO 2020/191103 A1** (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, 06/15/2023

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