

PRODUCT INFORMATION



Lipid AX4

Item No. 39070

CAS Registry No.: 2735814-23-8

MF: $C_{83}H_{155}N_3O_{16}$

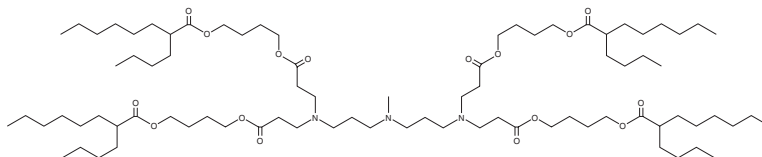
FW: 1,451.2

Purity: $\geq 95\%$

Supplied as: A 10 mg/ml solution in methyl acetate

Storage: $-20^{\circ}C$

Stability: ≥ 3 years



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

Description

Lipid AX4 is an ionizable cationic lipid ($pK_a = 6.89$) that has been used in the formation of lipid nanoparticles (LNPs) for the delivery of mRNA *in vivo*.¹ Administration of LNPs containing lipid AX4 and encapsulating severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) Delta variant mRNA encoding the SARS-CoV-2 spike glycoprotein, also known as surface glycoprotein, receptor-binding domain (RBD) induces the production of antibodies that bind to the SARS-CoV-2 spike RBD in mice.

Reference

1. Huang, K., Li, N., Li, Y., et al. Delivery of circular mRNA via degradable lipid nanoparticles against SARS-CoV-2 Delta variant. *bioRxiv* (2022).

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, 11/13/2025

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