

PRODUCT INFORMATION



Lipid 20b

Item No. 43121

CAS Registry No.: 2921586-65-2
Formal Name: 4,5,6,7-tetrahydro-2-[[[(9Z,12Z)-1-oxo-9,12-octadecadien-1-yl]amino]-thieno[2,3-c]pyridine-3-carboxylic acid, (9Z,12Z)-9,12-octadecadien-1-yl ester

MF: C₄₄H₇₂N₂O₃S

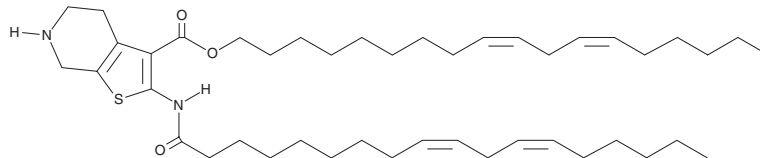
FW: 709.1

Purity: ≥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

Lipid 20b is a thiophene ionizable cationic lipid that has been used in the generation of lipid nanoparticles (LNPs) for the delivery of mRNA *in vivo*.¹ LNPs containing lipid 20b and encapsulating mRNA encoding a luciferase reporter induce bioluminescence in the lungs, spleen, and retina of mice.

Reference

1. Eygeris, Y., Gupta, M., Kim, J., *et al.* Thiophene-based lipids for mRNA delivery to pulmonary and retinal tissues. *Proc. Natl. Acad. Sci. USA* **121**(11), e2307813120 (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