

PRODUCT INFORMATION



FO-35

Item No. 42923

Formal Name: bis((4-(prop-1-en-2-yl)cyclohex-1-en-1-yl)methyl) 5-(butylcarbamoyl)-5-(N-(4-(diethylamino)butyl)-6-(heptadecan-9-yloxy)-6-oxohexanamido) nonanedioate

MF: C₆₅H₁₁₃N₃O₈

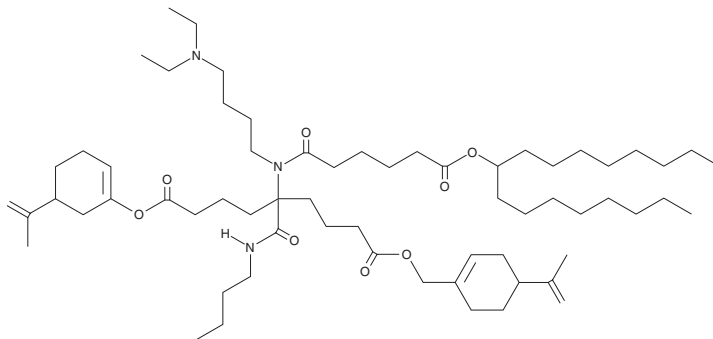
FW: 1,064.6

Purity: ≥98%

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

Storage: -20°C

Stability: ≥2 years



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

Description

FO-35 is an ionizable cationic lipid ($pK_a = 7.67$) that has been used in the generation of lipid nanoparticles (LNPs) for the delivery of mRNA *in vivo*.¹ Intranasal administration of LNPs containing FO-35 and encapsulating reporter mRNA accumulate in nasal epithelial tissue in mice. Nebulized administration of FO-35-containing LNPs encapsulating Cre mRNA induces the expression of tdTomato in the lungs of ferrets engineered to express tdTomato upon Cre-mediated recombination.

Reference

1. Witten, J., Raji, I., Manan, R.S., et al. Artificial intelligence-guided design of lipid nanoparticles for pulmonary gene therapy. *Nat. Biotechnol.* (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, 11/24/2025

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