

PRODUCT INFORMATION



Lipid 29 analog 2 (trifluoroacetate salt)

Item No. 43427

Formal Name: ((3-((2-(methylamino)-3,4-dioxocyclobut-1-en-1-yl)amino)propyl)azanediy)bis(hexane-6,1-diyl) bis(2-hexyldecanoate), trifluoroacetate salt

MF: C₅₂H₉₇N₃O₆ • XCF₃COOH

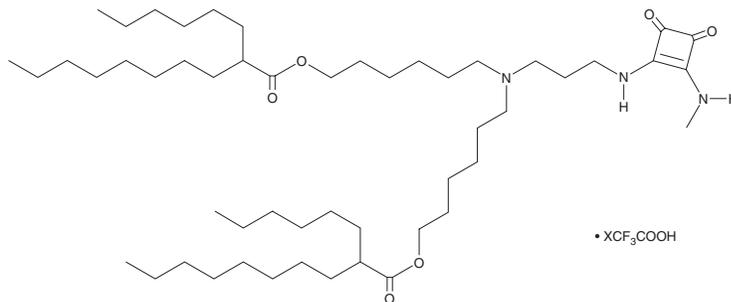
FW: 860.4

Purity: ≥90%

Supplied as: A 10 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 29 analog 2 is an ionizable cationic lipid and a derivative of lipid 29 (Item No. 35337).

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, 12/10/2025

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