

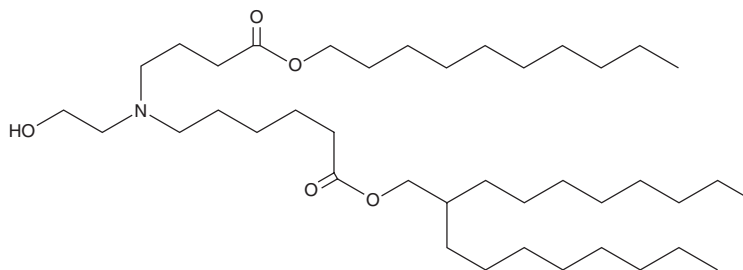
PRODUCT INFORMATION



YK-009

Item No. 38282

CAS Registry No.: 2761458-86-8
Formal Name: 6-[[4-(decyloxy)-4-oxobutyl] (2-hydroxyethyl)amino]-hexanoic acid, 2-octyldecyl ester
MF: C₄₀H₇₉NO₅
FW: 654.1
Purity: ≥95%
Supplied as: A 10 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥3 years



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

Description

YK-009 is an ionizable cationic lipid (pKa = 6.512) that has been used in combination with other lipids in the formation of lipid nanoparticles (LNPs).¹ Administration of severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) Omicron variant mRNA in YK-009-containing LNPs induces the production of IgG that binds to the SARS-CoV-2 receptor-binding domain (RBD) in mice, with a boost in IgG titers 10 days after a second dose.

Reference

1. Long, J., Yu, C., Zhang, H., *et al.* Novel ionizable lipid nanoparticles for SARS-CoV-2 omicron mRNA delivery. *Adv. Healthc. Mater.* **12**(13), 2202590 (2023).

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, 01/26/2026

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