

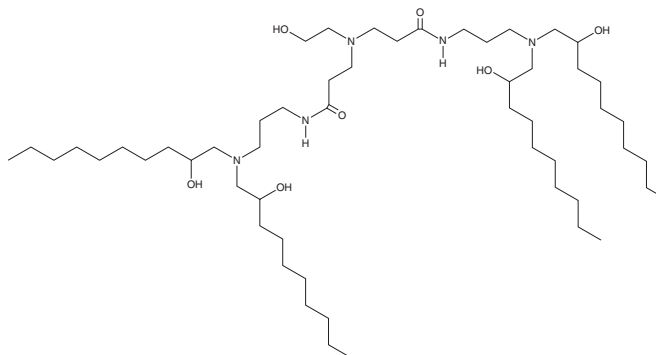
PRODUCT INFORMATION



OC2-K3-E10

Item No. 38243

CAS Registry No.: 2933216-12-5
Formal Name: 3,3'-((2-hydroxyethyl)azanediyl)bis(N-(3-(bis(2-hydroxydecyl)amino)propyl)propanamide)
Synonym: I-28
MF: C₅₄H₁₁₁N₅O₇
FW: 942.5
Purity: ≥95%
Supplied as: A 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

OC2-K3-E10 is an ionizable cationic lipid ($pK_a = 6.8$).¹ It has been used in the formation of lipid nanoparticles (LNPs) for the delivery of CRISPR complementary single-guide RNA (sgRNA) and Cas9 mRNA for genome editing in transgenic mice.

Reference

1. Khan, O.F., Tilstra, G., Manning, A.M., *et al.* Multi-motif dendrons and their supramolecular structures and uses thereof *The Governing Council of the University of Toronto WO 2023/092242 AI* (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, 12/11/2023

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