

PRODUCT INFORMATION

Lipid OEs-K

Item No. 43862

Formal Name: (1,4,7,10-tetraazacyclododecane-1,4,7,10-tetrayl)tetrakis(2-hydroxypropane-3,1-diyl)tetrakis(2-hexyldecanoate)

Synonyms: Lipidoid OEs-K, OEs-K

MF: $C_{84}H_{164}N_4O_{12}$

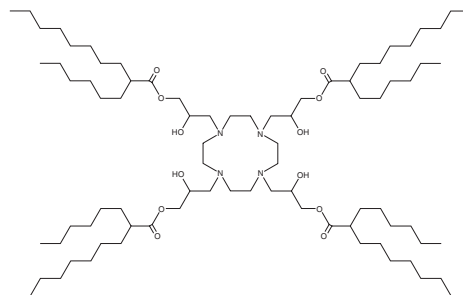
FW: 1,422.2

Purity: $\geq 95\%$ (mixture of diastereomers)

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.

Laboratory Procedures

Lipid OEs-K is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice, which should be purged with an inert gas. Lipid OEs-K is sparingly soluble (1-10 mg/ml) in 2:1 solution of chloroform:methanol.

Description

Lipid OEs-K is a cyclen-based ionizable cationic lipidoid that has been used in the formation of lipid nanoparticles (LNPs) for the delivery of mRNA *in vivo*.¹ LNPs containing lipid OEs-K accumulate in the lymph nodes of mice. LNPs containing lipid OEs-K and encapsulating ovalbumin (OVA) mRNA increase the uptake of dendritic cells by the lymph nodes and inhibit tumor growth in a murine E.G7-OVA model of lymphoma.

Reference

1. Zhang, L., Chen, P., Tian, X.-L., *et al.* Cyclen-based lipidoids for mRNA delivery and immunotherapy. *Biomater. Sci.* **13**(10), 2783-2793 (2025).

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, 04/07/2026

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