

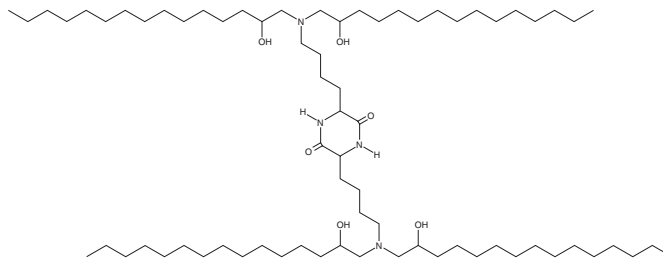
PRODUCT INFORMATION



cKK-E15

Item No. 38211

CAS Registry No.: 1432494-71-7
Formal Name: 3,6-bis[4-[bis(2-hydroxypentadecyl) amino]butyl]-2,5-piperazinedione
MF: C₇₂H₁₄₄N₄O₆
FW: 1,161.9
Purity: ≥95%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥1 year



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

Description

cKK-E15 is an ionizable cationic lipid and a derivative of cKK-E12 (Item No. 36700) that has been used in the generation of lipid nanoparticles (LNPs).¹ LNPs containing cKK-E15 and encapsulating Cre mRNA induce the expression of Cre in Kupffer cells, endothelial cells, and hepatocytes in Ai14 mice engineered to express the fluorescent protein tdTomato upon translation of Cre. Translation of mRNA delivered to the liver by LNPs containing cKK-E15 is inhibited by the toll-like receptor 4 (TLR4) agonist LPS in Ai14 mice but not by low-grade chronic inflammation in a B16/F10 murine model of cancer-induced inflammation. Unlike cKK-E12, the delivery of LNPs containing cKK-E15 is not dependent upon *ApoE* or *Ldlr* expression in mice.²

References

1. Lokugamage, M.P., Gan, Z., Zurla, C., *et al.* Mild innate immune activation overrides efficient nanoparticle-mediated RNA delivery. *Adv. Mater.* **32(1)**, e1904905 (2020).
2. Paunovska, K., Da Silva Sanchez, A.J., Lokugamage, M.P., *et al.* The extent to which lipid nanoparticles require apolipoprotein E and low-density lipoprotein receptor for delivery changes with ionizable lipid structure. *Nano. Lett.* **22(24)**, 10025-10033 (2022).

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, 10/13/2023

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