

PRODUCT INFORMATION



ZA3-Ep10

Item No. 40764

CAS Registry No.: 2090299-48-0
Formal Name: 10-[2-[bis(2-hydroxydecyl)amino]ethyl]-15-hydroxy-7,13-bis(2-hydroxydecyl)-N,N-dimethyl-4-oxo-N-(3-sulfopropyl)-3,7,10,13-tetraazatricosan-1-aminium, inner salt

MF: C₆₆H₁₃₈N₆O₉S

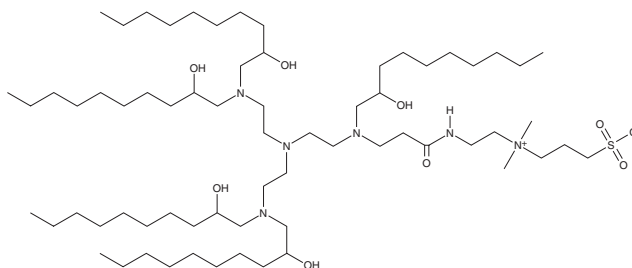
FW: 1,191.9

Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

ZA3-Ep10 is supplied as a solid. A stock solution may be made by dissolving the ZA3-Ep10 in the solvent of choice, which should be purged with an inert gas. ZA3-Ep10 is slightly soluble (0.1-1 mg/ml) in chloroform and methanol.

Description

ZA3-Ep10 is a zwitterionic aminolipid that has been used in the generation of lipid nanoparticles (LNPs) for the delivery of mRNA and single-guide RNA (sgRNA) *in vitro* and *in vivo*.¹ LNPs containing ZA3-Ep10 selectively accumulate in the liver, lung, and kidneys over the heart and spleen in mice. ZA3-Ep10-containing LNPs encapsulating sgRNA targeting DNA encoding luciferase eliminate luciferase activity in HeLa-Luc-Cas9 cells.

Reference

1. Miller, J.B., Zhang, S., Kos, P., *et al.* Non-viral CRISPR/Cas gene editing *in vitro* and *in vivo* enabled by synthetic nanoparticle co-delivery of Cas9 mRNA and sgRNA. *Angew. Chem. Int. Ed. Engl.* **56**(4), 1059-1063 (2017).

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, 01/23/2025

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