

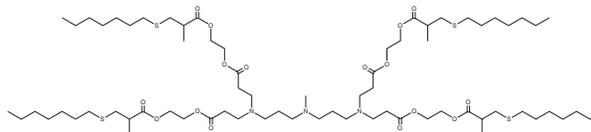
PRODUCT INFORMATION



4A3-SC7

Item No. 43989

CAS Registry No.: 1857340-77-2
MF: $C_{71}H_{131}N_3O_{16}S_4$
FW: 1,411.1
Purity: $\geq 95\%$
Supplied as: A 10 mg/ml solution in methyl acetate
Storage: $-20^{\circ}C$
Stability: ≥ 2 years
Special Conditions: Product degrades in protic solvents



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

Description

4A3-SC7 is an ionizable cationic lipid 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 4A3-SC7 selectively accumulate in mouse liver. LNPs containing 4A3-SC7 have been used to induce gene editing in a mouse model of alpha-1 antitrypsin deficiency.

Reference

1. Kim, M., Song, E.S., Chen, J.C., *et al.* Dual SORT LNPs for multi-organ base editing. *Nat. Biotechnol.* (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, 08/22/2025

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