

# PRODUCT INFORMATION



## AA-T3A-C12

Item No. 38648

**CAS Registry No.:** 2938207-23-7  
**Formal Name:** N-(3-(bis(3-(bis(2-hydroxydodecyl)amino)propyl)amino)propyl)-4-methoxybenzamide

**MF:** C<sub>65</sub>H<sub>126</sub>N<sub>4</sub>O<sub>6</sub>

**FW:** 1,059.7

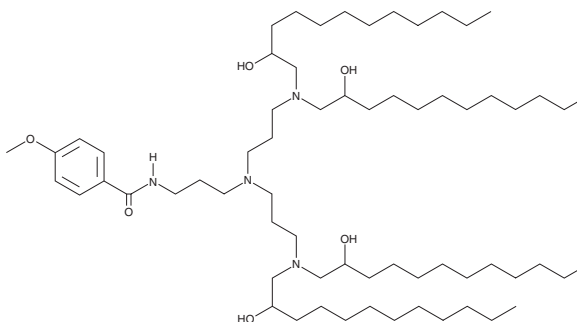
**Purity:** ≥95%

**UV/Vis.:** λ<sub>max</sub>: 254 nm

**Supplied as:** A 10 mg/ml 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

AA-T3A-C12 is an ionizable cationic lipidoid that contains a polyamine core, tethered to the sigma receptor ligand anisamide, and four 12-carbon tails.<sup>1</sup> It has been used in the generation of lipid nanoparticles (LNPs). LNPs containing AA-T3A-C12 and encapsulating siRNA for *Hsp47*, a molecular chaperone whose expression is increased in liver fibrosis, reduce the expression of *Hsp47* in activated NIH3T3 mouse liver fibroblasts and primary mouse hepatic stellate cells, which both express high levels of the sigma receptor. LNPs containing AA-T3A-C12 and encapsulating *Hsp47* siRNA reduce liver fibrosis induced by carbon tetrachloride (CCl<sub>4</sub>) in mice.

### Reference

1. Han, X., Gong, N., Xue, L., et al. Ligand-tethered lipid nanoparticles for targeted RNA delivery to treat liver fibrosis. *Nat. Commun.* **14**(1), 75 (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, 11/11/2025

#### 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