

# PRODUCT INFORMATION



**TT3**

Item No. 37909

**CAS Registry No.:** 1821214-50-9  
**Formal Name:** N<sup>1</sup>,N<sup>3</sup>,N<sup>5</sup>-tris[3-(didodecylamino)propyl]-1,3,5-benzenetricarboxamide

**MF:** C<sub>90</sub>H<sub>174</sub>N<sub>6</sub>O<sub>3</sub>

**FW:** 1,388.4

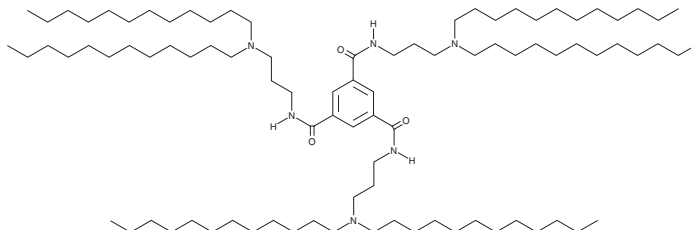
**Purity:** ≥95%

**UV/Vis.:** λ<sub>max</sub>: 210 nm

**Supplied as:** A solution in ethanol

**Storage:** -20°C

**Stability:** ≥2 years



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

## Description

TT3 is an ionizable cationic amino lipid that has been used in combination with other lipids in the formation of lipid-like nanoparticles (LLNs).<sup>1</sup> Administration of LLNs containing TT3 and encapsulating mRNA encoding human coagulation Factor IX induces human coagulation Factor IX expression in the plasma of mice.

## Reference

1. Li, B., Luo, X., Deng, B., *et al.* An orthogonal array optimization of lipid-like nanoparticles for mRNA delivery in vivo. *Nano Lett.* **15**(12), 8099-8107 (2015).

### 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, 02/14/2023

## 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](http://WWW.CAYMANCHEM.COM)