

PRODUCT INFORMATION



FTT5

Item No. 37413

CAS Registry No.: 2803606-43-9
Formal Name: 9,9',9'',9''',9''''',9'''''-[1,3,5-benzenetriyltris (carbonylimino-3,1-propanediyl)nitri]lo]hexakis-nonanoic acid, 1,1',1'',1''',1''''',1'''''-hexakis (1-ethylhexyl) ester

Synonym: Functionalized N¹,N³,N⁵-tris(2-aminoethyl)benzene-1,3,5-Tricarboxamide 5

MF: C₁₂₀H₂₂₂N₆O₁₅

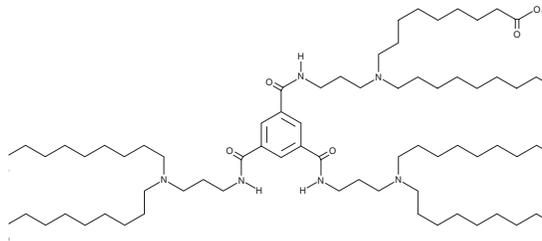
FW: 1,989.1

Purity: ≥98%

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.

Laboratory Procedures

FTT5 is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. FTT5 is soluble in methanol.

Description

FTT5 is a lipid-like compound that has been used in the formation of lipid-like nanoparticles (LLNs) and is a derivative of the ionizable cationic amino lipid TT3 (Item No. 37909).¹ LLNs containing FTT5 and encapsulating an mRNA reporter increase luminescence in Hep3B cells and selectively accumulate in mouse liver. LLNs containing FTT5 and encapsulating mRNA encoding human Factor VIII increase plasma Factor VIII activity in a mouse model of hemophilia A. FTT5-containing LLNs have been used for genome editing in mice.

Reference

1. Zhang, X., Zhao, W., Nguyen, G.N., et al. Functionalized lipid-like nanoparticles for in vivo mRNA delivery and base editing. *Sci. Adv.* **6**(34), eabc2315 (2024).

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, 03/19/2024

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