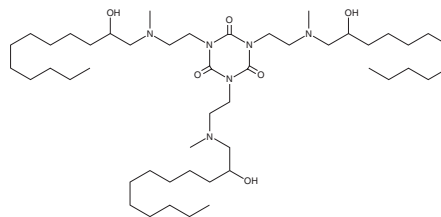


PRODUCT INFORMATION



TNT-b₁₀
Item No. 39294

Formal Name: 1,3,5-tris(2-((2-hydroxydodecyl)(methyl)amino)ethyl)-1,3,5-triazinane-2,4,6-trione
MF: C₄₈H₉₆N₆O₆
FW: 853.3
Purity: ≥95%
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

TNT-b₁₀ is an ionizable cationic lipidoid that has been used in the formation of lipid-like nanoparticles (LLNs).¹ LLNs containing TNT-b₁₀ and encapsulating an mRNA reporter increase luminescence in Hep3B liver cancer cells.

Reference

1. Li, B., Luo, X., Deng, B., *et al.* Effects of local structural transformation of lipid-like compounds on delivery of messenger RNA. *Sci. Rep.* **6**, 22137 (2016).

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, 02/15/2024

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