

PRODUCT INFORMATION



1014

Item No. 38150

Formal Name: 15,19,28-tris(2-hydroxytetradecyl)-
15,19,24,28-tetraazadotetracontane-
13,30-diol

MF: C₈₀H₁₆₆N₄O₅

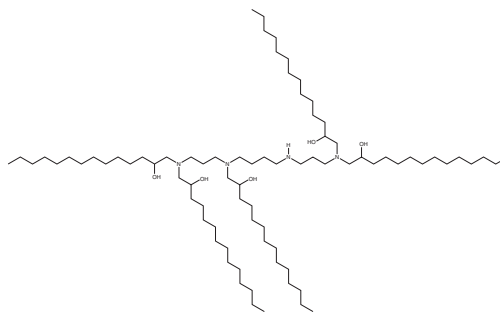
FW: 1,264.2

Purity: ≥95%

Supplied as: A 25 mg/ml 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

1014 is an ionizable cationic lipid that has been used in combination with other lipids in the formation of lipid nanoparticles (LNPs).¹ 1014-containing LNPs have been used for the delivery of IL-1 β siRNA to induce gene silencing and hepatoprotective effects in a mouse model of acute liver injury induced by LPS and galactosamine (GalN).

Reference

1. Zhang, H., Ding, F., Zhu, Z., *et al.* Engineered ionizable lipid nanoparticles mediated efficient siRNA delivery to macrophages for anti-inflammatory treatment of acute liver injury. *Int. J. Pharm.* **631**, 122489 (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, 11/12/2025

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