

PRODUCT INFORMATION



98N12-5

Item No. 37651

CAS Registry No.: 917572-74-8
Formal Name: N1,N16-didodecyl-4,7,13-tris[3-(dodecylamino)-3-oxopropyl]-4,7,10,13-tetraazahexadecanediamide

Synonym: ND98

MF: C₈₁H₁₆₃N₉O₅

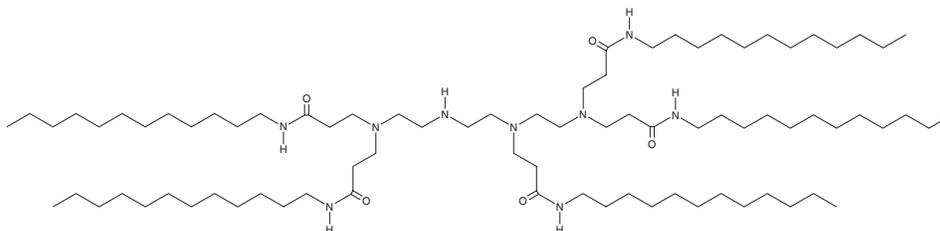
FW: 1,343.2

Purity: ≥95%

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥2 years



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

Laboratory Procedures

98N12-5 is supplied as a crystalline solid. A stock solution may be made by dissolving the 98N12-5 in the solvent of choice, which should be purged with an inert gas. 98N12-5 is soluble in the organic solvent chloroform at a concentration of approximately 30 mg/ml.

Description

98N12-5 is an ionizable cationic lipid.¹ It has been used in combination with other lipids in the generation of lipid nanoparticles (LNPs). LNPs containing 98N12-5 and encapsulating proprotein convertase subtilisin kexin type 9 (PCSK9) siRNA selectively accumulate in the liver and reduce total serum cholesterol levels in mice and rats and serum LDL levels in cynomolgus monkeys.²

References

1. Akinc, A., Goldberg, M., Qin, J., *et al.* Development of lipidoid-siRNA formulations for systemic delivery to the liver. *Mol. Ther.* **17**(5), 872-879 (2009).
2. Frank-Kamenetsky, M., Grefhorst, A., Anderson, N.N., *et al.* Therapeutic RNAi targeting PCSK9 acutely lowers plasma cholesterol in rodents and LDL cholesterol in nonhuman primates. *Proc. Natl. Acad. Sci. USA* **105**(33), 11915-11920 (2008).

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, 02/15/2023

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