PRODUCT INFORMATION



ALC-0159

Item No. 34336

CAS Registry No.: 1849616-42-7

Formal Name: α -[2-(ditetradecylamino)-2-oxoethyl]- ω -methoxy-

poly(oxy-1,2-ethanediyl)

Synonyms: N,N-ditetradecylacetamide PEG(2000),

N,N-ditetradecylacetamide Polyethylene Glycol-2000

MF: $(C_2H_4O)_nC_{31}H_{63}NO_2$ Molecular Weight: Average: 2400 Da

≥95% **Purity:** Supplied as: A solid Storage: -20°C Stability: ≥4 years

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

Laboratory Procedures

ALC-0159 is supplied as a solid. A stock solution may be made by dissolving the ALC-0159 in the solvent of choice, which should be purged with an inert gas. ALC-0159 is soluble in organic solvents such as ethanol and dimethyl formamide (DMF). ALC-0159 is slightly soluble in DMSO. The solubility of ALC-0159 in ethanol and DMF is approximately 1 and 10 mg/ml, respectively.

ALC-0159 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ALC-0159 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. ALC-0159 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

ALC-0159 is a lipid excipient. Formulations containing ALC-0159 have been used in the development of lipid nanoparticles (LNPs) for the delivery of mRNA-based vaccines.

Reference

1. Cortés, H., Hernández-Parra, H., Bernal-Chávez, S.A., et al. Non-ionic surfactants for stabilization of polymeric nanoparticles for biomedical uses. Materials (Basel) 14(12), 3197 (2021).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

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

uyer 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, 09/18/2023

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