

# PRODUCT INFORMATION



## CL15F6

Item No. 40345

**CAS Registry No.:** 2766493-40-5  
**Formal Name:** 1-methyl-4-piperidinecarboxylic acid, 11-[(2-hexyl-1-oxodecyl oxy)-5-[6-[(2-hexyl-1-oxodecyl oxy)hexyl]-5-hydroxyundecyl ester

**MF:** C<sub>56</sub>H<sub>107</sub>NO<sub>7</sub>

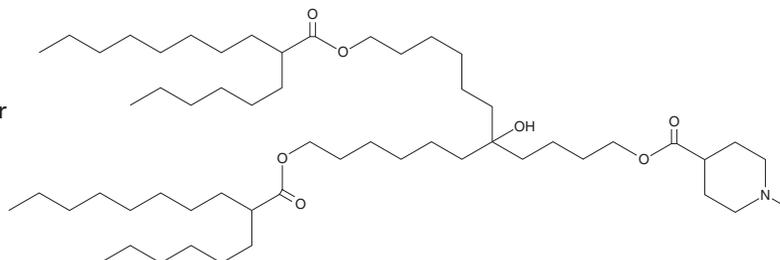
**FW:** 906.5

**Purity:** ≥98%

**Supplied as:** A 10 mg/ml solution in ethanol

**Storage:** -20°C

**Stability:** ≥4 years



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

### Laboratory Procedures

CL15F6 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. Solvents such as chloroform and methanol purged with an inert gas can be used.

### Description

CL15F6 is an ionizable cationic lipid ( $pK_a = 6.75$ ).<sup>1</sup> It has been used in the formation of lipid nanoparticles (LNPs) for the delivery of mRNA and polymer-lipid hybrid nanoparticles for the delivery of plasmid DNA *in vitro*.<sup>1,2</sup>

### References

1. Younis, M.A., Sato, Y., Elewa, Y.H.A., *et al.* Self-homing nanocarriers for mRNA delivery to the activated hepatic stellate cells in liver fibrosis. *J. Control. Release* **353**, 685-698 (2023).
2. Maeki, M., Uno, S., Sugiura, K., *et al.* Development of polymer-lipid hybrid nanoparticles for large-sized plasmid DNA transfection. *ACS Appl. Mater. Interfaces* **16(2)**, 2110-2119 (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, 11/18/2025

#### 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