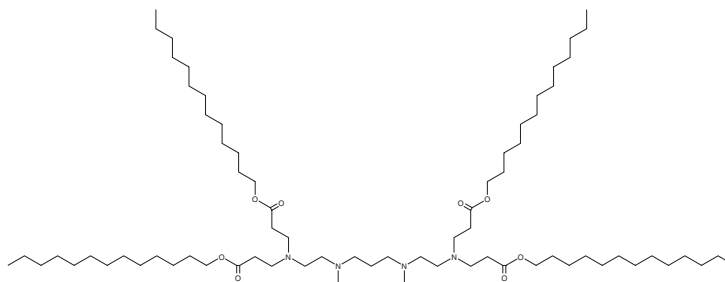


# PRODUCT INFORMATION



**503O<sub>13</sub>**  
Item No. 38628

**CAS Registry No.:** 1566560-20-0  
**Formal Name:** 7,11-dimethyl-4,14-bis[3-oxo-3-(tridecyloxy)propyl]-4,7,11,14-tetraazaheptadecanedioic acid, 1,17-ditridecyl ester  
**MF:** C<sub>73</sub>H<sub>144</sub>N<sub>4</sub>O<sub>8</sub>  
**FW:** 1,206.0  
**Purity:** 90%  
**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.

## Laboratory Procedures

503O<sub>13</sub> 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. 503O<sub>13</sub> is soluble in methanol.

## Description

503O<sub>13</sub> is an ionizable cationic lipidoid that has been used in the generation of lipid nanoparticles (LNPs) for the delivery of siRNA *in vivo*.<sup>1</sup> LNPs containing 503O<sub>13</sub> and encapsulating siRNA targeting mRNA encoding Factor VII decrease serum Factor VII levels in mice.

## Reference

1. Whitehead, K.A., Dorkin, J.R., Vegas, A.J., *et al.* Degradable lipid nanoparticles with predictable *in vivo* siRNA delivery activity. *Nat. Commun.* **5**, 4277 (2014).

### 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, 09/03/2025

## 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](http://WWW.CAYMANCHEM.COM)