

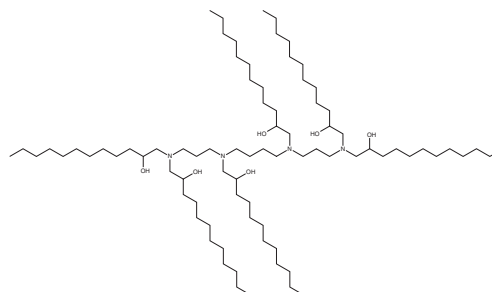
# PRODUCT INFORMATION



## C12-SPM

Item No. 38784

**CAS Registry No.:** 2055647-81-7  
**Formal Name:** 13,17,22,26-tetrakis(2-hydroxydodecyl)-13,17,22,26-tetraazaoctatriacontane-11,28-diol  
**Synonym:** C12-Spermine  
**MF:**  $C_{82}H_{170}N_4O_6$   
**FW:** 1,308.3  
**Purity:**  $\geq 95\%$  (mixture of isomers)  
**Supplied as:** A 25 mg/ml solution in methanol  
**Storage:**  $-20^{\circ}C$   
**Stability:**  $\geq 3$  years



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

### Description

C12-SPM is a polyamine branched-chain lipidoid.<sup>1</sup> It has been used in combination with other lipids in the formation of lipid nanoparticles (LNPs) for the delivery of siRNA targeting the host protein kinase protein kinase C-related kinase 2 (PRK2) to inhibit hepatitis C virus (HCV) replication in Huh7 cells and a mouse model of HCV infection.

### Reference

1. Park, H.-J., Jeon, E.J., Lee, J.S., *et al.* Galactosylated lipidoid nanoparticles for delivery of small interfering RNA to inhibit hepatitis C viral replication in vivo. *Adv. Healthc. Mater.* **5(22)**, 2931-2941 (2016).

#### 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, 01/26/2026

#### 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)