

# PRODUCT INFORMATION



## A12-Iso5-2DC18

Item No. 38586

**CAS Registry No.:** 2412492-06-7  
**Formal Name:** 1-[3-(dimethylamino)propyl]-5,5-di-(8Z)-8-heptadecen-1-yl-2,5-dihydro-1H-imidazole-2-carboxylic acid, ethyl ester

**MF:** C<sub>45</sub>H<sub>85</sub>N<sub>3</sub>O<sub>2</sub>

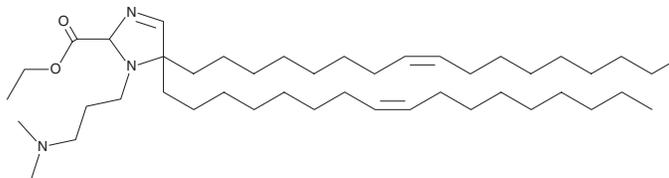
**FW:** 700.2

**Purity:** ≥95%

**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.

### Description

A12-Iso5-2DC18 is an ionizable cationic lipidoid that has been used in the formation of lipid nanoparticles (LNPs).<sup>1</sup> LNPs containing A12-Iso5-2DC18 and encapsulating mRNA encoding luciferase have been used to induce luciferase expression in HeLa cells.

### Reference

1. Miao, L., Li, L., Huang, Y., *et al.* Delivery of mRNA vaccines with heterocyclic lipids increases anti-tumor efficacy by STING-mediated immune cell activation. *Nat. Biotechnol.* **37**(10), 1174-1185 (2019).

#### 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, 05/25/2023

#### 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