

# PRODUCT INFORMATION



## Lipid 29

Item No. 35337

**CAS Registry No.:** 2244716-55-8  
**Formal Name:** 8-[[8-[(1-ethylonyl)oxy]-8-oxooctyl]  
[3-[[2-(methylamino)-3,4-dioxo-1-  
cyclobuten-1-yl]amino]propyl]amino]-  
octanoic acid, 1-octylonyl ester

**MF:** C<sub>52</sub>H<sub>97</sub>N<sub>3</sub>O<sub>6</sub>

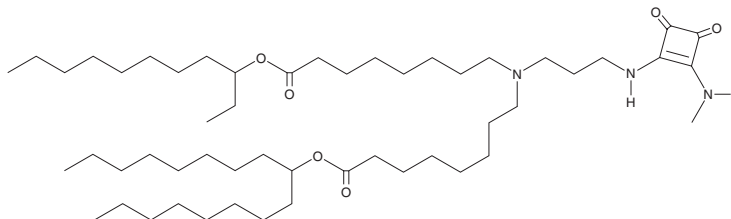
**FW:** 860.4

**Purity:** ≥98%

**Supplied as:** A 25 mg/ml 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

Lipid 29 is an ionizable cationic amino lipid ( $pK_a = 6.91$ ) that has been used in combination with other lipids in the formation of lipid nanoparticles (LNPs).<sup>1</sup> Administration of human erythropoietin (EPO) mRNA in lipid 29-containing LNPs increases serum EPO levels in mice.

### Reference

1. Cornebise, M., Narayanan, E., Xia, Y., *et al.* Discovery of a novel amino lipid that improves lipid nanoparticle performance through specific interactions with mRNA. *Adv. Funct. Mater.* **32(8)**, 2106727(1-12) (2021).

#### 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, 07/29/2025

#### 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