

PRODUCT INFORMATION



9A1P9

Item No. 37276

CAS Registry No.: 2760467-57-8
Formal Name: 2-(dioctylamino)ethyl nonyl hydrogen phosphate

MF: C₂₇H₅₈NO₄P

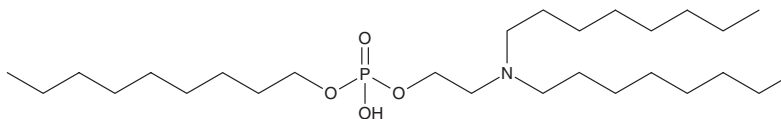
FW: 491.7

Purity: ≥95%

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

9A1P9 is supplied as a solution in ethanol. To change the solvent, simply evaporate the 9A1P9 under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of 9A1P9 in DMF is approximately 2 mg/ml.

Description

9A1P9 is a multi-tail ionizable cationic phospholipid.¹ It induces membrane destabilization and has been used for organ-selective mRNA delivery to the spleen and liver and CRISPR-Cas9 gene editing in mice.

Reference

1. Liu, S., Cheng, Q., Wei, T., *et al.* Membrane-destabilizing ionizable phospholipids for organ-selective mRNA delivery and CRISPR-Cas gene editing. *Nat. Mater.* **20(5)**, 701-710 (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, 11/06/2025

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