

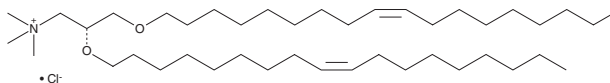
PRODUCT INFORMATION



(R)-DOTMA (chloride)

Item No. 45535

CAS Registry No.: 2101981-60-4
Formal Name: (2R)-N,N,N-trimethyl-2,3-bis[(9Z)-9-octadecen-1-yloxy]-1-propanaminium, monochloride
MF: $C_{42}H_{84}NO_2 \cdot Cl$
FW: 670.6
Purity: $\geq 98\%$
Supplied as: A 50 mg/ml solution in ethanol
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Description

(R)-DOTMA is a cationic lipid that has been used in the formation of liposomes with a net cationic charge.¹ (R)-DOTMA-containing liposomes have been complexed with mRNA to form RNA-lipoplexes (RNA-LPXs).

Reference

1. Meibom, D., Albrecht-Küpper, B., Diedrichs, N., *et al.* Neladenoson bialanate hydrochloride: A prodrug of a partial adenosine A_1 receptor agonist for the chronic treatment of heart diseases. *ChemMedChem* **12**(10), 728-737 (2017).

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, 04/17/2026

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