

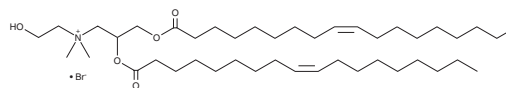
PRODUCT INFORMATION



DORI

Item No. 38981

CAS Registry No.: 153312-59-5
Formal Name: N-(2-hydroxyethyl)-N,N-dimethyl-2,3-bis[[[(9Z)-1-oxo-9-octadecen-1-yl]oxy]-1-propanaminium, monobromide
Synonym: 1,2-Dioleoyl-3-dimethyl-hydroxyethylammonium bromide
MF: C₄₃H₈₂NO₅ • Br
FW: 773.0
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

DORI is supplied as a crystalline solid. A stock solution may be made by dissolving the DORI in the solvent of choice, which should be purged with an inert gas. DORI is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of DORI in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMF. DORI is slightly soluble in DMSO.

Description

DORI is a cationic lipid that has been used in the formation of liposomes for the delivery of plasmid DNA *in vitro*.¹

Reference

1. Felgner, J.H., Kumar, R., Sridhar, C.N., *et al.* Enhanced gene delivery and mechanism studies with a novel series of cationic lipid formulations. *J. Biol. Chem.* **269**(4), 2550-2561 (1994).

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, 10/09/2025

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