

PRODUCT INFORMATION



Dioleoyldimethylammonium (chloride)

Item No. 38797

CAS Registry No.: 7212-69-3
Formal Name: N,N-dimethyl-N-(9Z)-9-octadecen-1-yl-9-octadecen-1-aminium, monochloride
Synonyms: Dimethyldioleoylammonium chloride, DODAC, N,N-Dioleoyl-N,N-dimethylammonium chloride

MF: C₃₈H₇₆N • Cl

FW: 582.5

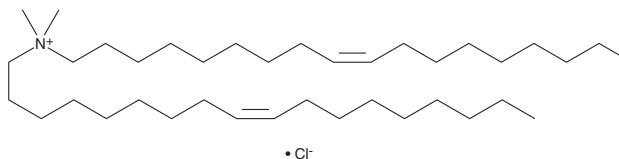
Purity: ≥95%

UV/Vis.: λ_{max}: 220 nm

Supplied as: A 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

Dioleoyldimethylammonium (chloride) (DODAC) is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice, which should be purged with an inert gas. DODAC is soluble in methanol.

Description

DODAC is a cationic lipid that has been used as a transfection reagent.¹ Liposomes containing DODAC and dioleoylphosphatidylethanolamine (DOPE) and encapsulating a plasmid DNA reporter induce reporter gene expression in C2C12 mouse skeletal muscle myoblasts.

Reference

1. Vitiello, L., Bockhold, K., Joshi, P.B., *et al.* Transfection of cultured myoblasts in high serum concentration with DODAC:DOPE liposomes. *Gene Ther.* **5(10)**, 1306-1313 (1998).

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, 04/27/2026

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