

# 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:** C<sub>38</sub>H<sub>76</sub>N • Cl

**FW:** 582.5

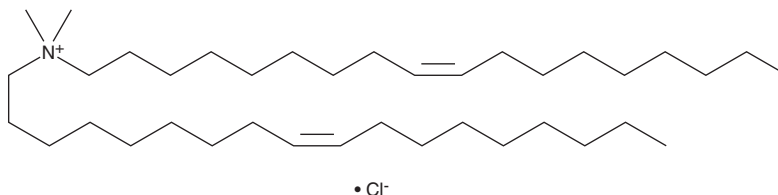
**Purity:** ≥95%

**UV/Vis.:** λ<sub>max</sub>: 220 nm

**Supplied as:** A 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.

### 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.<sup>1</sup> 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, 09/27/2023

#### 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