

PRODUCT INFORMATION



CM-DiD

Item No. 45061

Formal Name: 5-((4-(chloromethyl)benzamido)methyl)-2-((1E,3E,5E)-7-((E)-3,3-dimethyl-1-octadecylindolin-2-ylidene)hepta-1,3,5-trien-1-yl)-3,3-dimethyl-1-octadecyl-3H-indol-1-ium, monochloride

MF: $C_{72}H_{109}ClN_3O \bullet Cl$

FW: 1,103.6

Purity: $\geq 90\%$

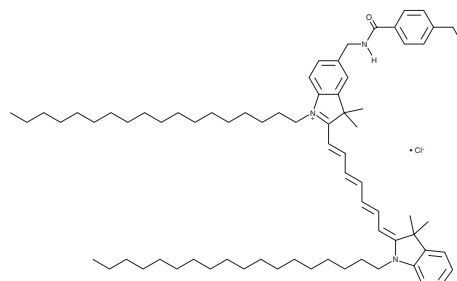
Ex./Em. Max: 643/663 nm

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years

Special Conditions: Light sensitive



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

Laboratory Procedures

CM-DiD is supplied as a solid. A stock solution may be made by dissolving the CM-DiD in the solvent of choice, which should be purged with an inert gas. CM-DiD is soluble in DMSO.

Description

CM-DiD is a hydrophobic, fluorescent probe for cell membranes in culture and in live or fixed tissues. It displays excitation/emission maxima of 643/663 nm, respectively, and is resistant to photobleaching. CM-DiD produces a strong fluorescent signal when incorporated into cell membranes and lipid-containing structures and does not rapidly diffuse within the membrane. It is used to label and trace cell membranes to study cell migration, cell tracking, and axonal projections and is well suited for neuroscience research and fluorescence microscopy.

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, 02/27/2026

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