

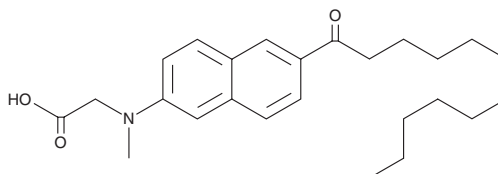
# PRODUCT INFORMATION



## C-Laurdan

Item No. 43531

**CAS Registry No.:** 959839-06-6  
**Formal Name:** N-methyl-N-[6-(1-oxododecyl)-2-naphthalenyl]-glycine  
**MF:** C<sub>25</sub>H<sub>35</sub>NO<sub>3</sub>  
**FW:** 397.6  
**Purity:** ≥98%  
**EX. Max:** 820 nm in water  
**Abs./Em. Max:** 384/522 nm in water  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥3 years  
**Special Conditions:** Light sensitive



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

### Laboratory Procedures

C-Laurdan is supplied as a solid. A stock solution may be made by dissolving the C-laurdan in the solvent of choice, which should be purged with an inert gas. C-Laurdan is soluble (≥10 mg/ml) in DMSO.

### Description

C-Laurdan is a polarity-sensitive fluorescent probe for lipid rafts and a derivative of the fluorescent membrane probe laurdan (Item No. 19706).<sup>1</sup> It displays a two-photon excitation maximum of 820 nm and one-photon absorption/emission maxima of 384/522 nm, respectively, in water. C-Laurdan fluorescence exhibits polarity-dependent spectral shifts, with emission at longer wavelengths in polar environments and shorter wavelengths in non-polar environments. It has been used in the study of lipid rafts and membrane lateral organization.<sup>1,2</sup>

### References

1. Kim, H.M., Choo, H.-J., Jung, S.-Y., *et al.* A two-photon fluorescent probe for lipid raft imaging: C-laurdan. *Chembiochem* **8**(5), 553-559 (2007).
2. Dodes Traian, M.M., Flecha, F.L.G., and Levi, V. Imaging lipid lateral organization in membranes with C-laurdan in a confocal microscope. *J. Lipid Res.* **53**(3), 609-616 (2012).

#### 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, 12/18/2025

#### 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](http://WWW.CAYMANCHEM.COM)