

# PRODUCT INFORMATION

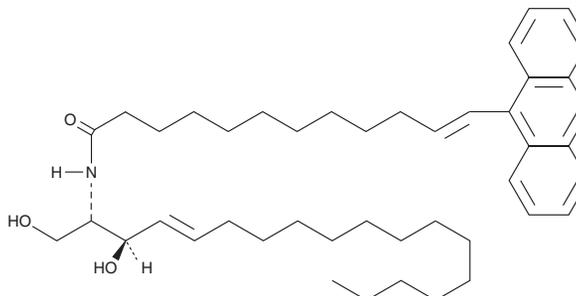


## C10 AV Ceramide (d18:1/10:0)

Item No. 9000753

**CAS Registry No.:** 1263052-40-9  
**Formal Name:** (11E)-12-(9-anthracenyl)-N-[(1S,2R,3E)-2-hydroxy-1-(hydroxymethyl)-3-heptadecen-1-yl]-11-dodecenamide  
**Synonyms:** Anthrylviny Ceramide, AV Ceramide, AV Ceramide (d18:1/10:0), AV Cer(d18:1/10:0)

**MF:** C<sub>44</sub>H<sub>65</sub>NO<sub>3</sub>  
**FW:** 656.0  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 256 nm  
**Supplied as:** A 10 mg/ml 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

C10 AV Ceramide (d18:1/10:0) 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. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of C10 AV ceramide (d18:1/10:0) in these solvents is approximately 10 mg/ml.

### Description

AV-Ceramide is a fluorescently tagged probe consisting of C-10 ceramide with an anthrylviny (AV) group attached to the end of the acyl chain. The AV group, being relatively small and non-polar, readily orients within the central region of a lipid bilayer.<sup>1</sup> The transfer or transport of lipids labeled with AV is commonly evaluated in real time using fluorescence resonance energy transfer, with a second fluorophore.<sup>1-3</sup>

### References

1. Polozov, I., Molotkovsky, J.G., and Bergelson, L.D. Anthrylviny-labeled phospholipids as membrane probes: The phosphatidylcholine-phosphatidylethanolamine system. *Chem. Phys. Lipids* **69(3)**, 209-218 (1994).
2. Mattijus, P., Molotkovsky, J.G., Smaby, J.M., et al. A fluorescence resonance energy transfer approach for monitoring protein-mediated glycolipid transfer between vesicle membranes. *Anal. Biochem.* **268(2)**, 297-304 (1999).
3. Tuuf, J., Kjellberg, M.A., Molotkovsky, J.G., et al. The intermediate ceramide transport catalyzed by CERT is sensitive to the lipid environment. *Biochem. Biophys. Acta* **1808(1)**, 229-235 (2011).

#### 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, 11/24/2025

#### 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