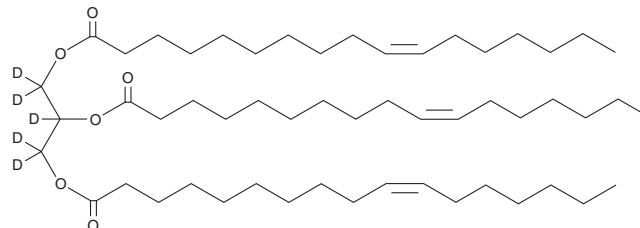


PRODUCT INFORMATION



1,2,3-Tri-10(Z)-Heptadecenoyl Glycerol-d₅ Item No. 28920

Formal Name: propane-1,2,3-triyl-d₅ (10Z,10'Z,10''Z)-tris(heptadec-10-enoate)
Synonyms: Glycerol Tri-10(Z)-Heptadecenoate-d₅, TG(17:1/17:1/17:1)-d₅, Tri-10(Z)-Heptadecenoin-d₅
MF: C₅₄H₉₃D₅O₆
FW: 848.4
Chemical Purity: ≥98% (1,2,3-Tri-10(Z)-Heptadecenoyl Glycerol)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀
Supplied as: An oil
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

1,2,3-Tri-10(Z)-heptadecenoyl glycerol-d₅ is intended for use as an internal standard for the quantification of 1,2,3-tri-10(Z)-heptadecenoyl glycerol by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

1,2,3-Tri-10(Z)-heptadecenoyl glycerol-d₅ is supplied as an oil. A stock solution may be made by dissolving the 1,2,3-tri-10(Z)-heptadecenoyl glycerol-d₅ in the solvent of choice, which should be purged with an inert gas. 1,2,3-Tri-10(Z)-heptadecenoyl glycerol-d₅ is miscible in organic solvents such as ethanol, DMSO, and dimethyl formamide.

Description

1,2,3-Tri-10(Z)-heptadecenoyl glycerol is a triacylglycerol that contains 10(Z)-heptadecenoic acid (Item No. 19748) at the *sn*-1, *sn*-2, and *sn*-3 positions. It has been used as an internal standard for the quantification of triglycerides in green microalgae and phospholipid bilayers.^{1,2}

References

1. Allen, J.W., DiRusso, C.C., and Black, P.N. Triglyceride quantification by catalytic saturation and LC-MS/MS reveals an evolutionary divergence in regioisometry among green microalgae. *Algal Res.* **5**(1), 23-31 (2014).
2. Duelund, L., Jensen, G.V., Hannibal-Bach, H.K., *et al.* Composition, structure and properties of POPC-triolein mixtures. Evidence of triglyceride domains in phospholipid bilayers. *Biochim. Biophys. Acta* **1828**(8), 1909-1917 (2013).

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, 11/08/2022

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