

PRODUCT INFORMATION



1,2,3-Triheptadecanoyl Glycerol-d₅

Item No. 28918

CAS Registry No.: 944795-74-8

Formal Name: Heptadecanoic acid,
1,1',1''-(1,2,3-propanetriyl-
1,1,2,3,3-d₅) ester

Synonyms: Glycerol Triheptadecanoate-d₅,
TG(17:0/17:0/17:0)-d₅

MF: C₅₄H₉₉D₅O₆

FW: 854.4

Chemical Purity: ≥98% (1,2,3-Triheptadecanoyl
Glycerol)

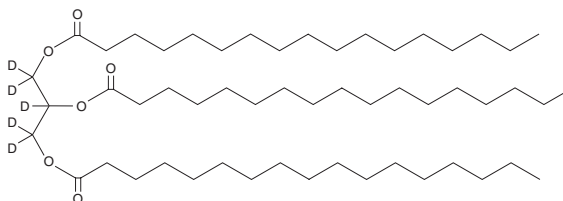
Deuterium

Incorporation: ≥99% deuterated forms (d₁-d₅);
≤1% d₀

Supplied as: A solid

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-Triheptadecanoyl glycerol-d₅ is intended for use as an internal standard for the quantification of 1,2,3-triheptadecanoyl glycerol (Item No. 16489) 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-Triheptadecanoyl glycerol-d₅ is supplied as a solid. A stock solution may be made by dissolving the 1,2,3-triheptadecanoyl glycerol-d₅ in the solvent of choice, which should be purged with an inert gas. 1,2,3-Triheptadecanoyl glycerol-d₅ is soluble in the organic solvent dimethyl formamide at a concentration of approximately 10 mg/ml.

Description

1,2,3-Triheptadecanoyl glycerol is a triacylglycerol that contains heptadecanoic acid (Item No. 19722) at the sn-1, sn-2, and sn-3 positions.

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, 10/27/2022

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