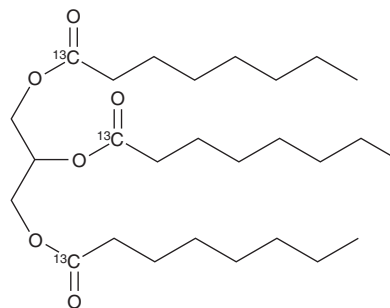


PRODUCT INFORMATION



1,2,3-Trioctanoyl Glycerol-¹³C₃ Item No. 29464

CAS Registry No.: 65402-55-3
Formal Name: octanoic-1-¹³C acid
1,1',1''-(1,2,3-propanetriyl) ester
Synonyms: Glycerol Trioctanoate-¹³C₃,
TG(8:0/8:0/8:0)-¹³C₃, Trioctanoin-¹³C₃
MF: C₂₄[¹³C₃]H₅₀O₆
FW: 473.7
Purity: ≥95%
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-Trioctanoyl glycerol-¹³C₃ is supplied as an oil. A stock solution may be made by dissolving the 1,2,3-trioctanoyl glycerol-¹³C₃ in the solvent of choice, which should be purged with an inert gas. 1,2,3-Trioctanoyl glycerol-¹³C₃ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 1,2,3-trioctanoyl glycerol-¹³C₃ in ethanol is approximately 20 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Description

1,2,3-Trioctanoyl glycerol-¹³C₃ is intended for use as an internal standard for the quantification of 1,2,3-trioctanoyl glycerol by GC- or LC-MS. 1,2,3-Trioctanoyl glycerol is a triacylglycerol that contains octanoic acid at the *sn*-1, *sn*-2, and *sn*-3 positions. Dietary administration of 1,2,3-trioctanoyl glycerol increases hippocampal levels of the glycolytic metabolites glucose 6-phosphate, fructose 6-phosphate, and β-hydroxybutyrate and the seizure threshold in the 6 Hz psychomotor seizure test in mice.¹ Formulations containing 1,2,3-trioctanoyl glycerol have been used in cosmetic products as thickening and skin-conditioning agents.

Reference

1. McDonald, T.S., Tan, K.N., Hodson, M.P., *et al.* Alterations of hippocampal glucose metabolism by even versus uneven medium chain triglycerides. *J. Cereb. Blood Flow Metab.* **34**(1), 153-160 (2014).

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, 11/22/2022

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