

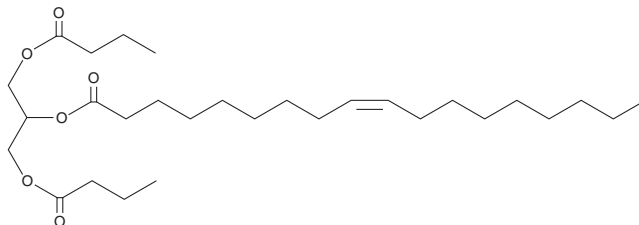
PRODUCT INFORMATION



1,3-Dibutyl-2-Oleoyl Glycerol

Item No. 26985

CAS Registry No.: 175878-34-9
Formal Name: 9Z-octadecenoic acid, 2-(1-oxobutoxy)-1-[(1-oxobutoxy)methyl]ethyl ester
Synonyms: 1,3-Butyryl-2-Olein, TG(4:0/18:1/4:0)
MF: C₂₉H₅₂O₆
FW: 496.7
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

1,3-Dibutyl-2-oleoyl glycerol is supplied as an oil. A stock solution may be made by dissolving the 1,3-dibutyl-2-oleoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,3-Dibutyl-2-oleoyl glycerol is slightly soluble in chloroform.

Description

1,3-Dibutyl-2-oleoyl glycerol is a triacylglycerol that contains butyric acid (Item No. 13121) at the *sn*-1 and *sn*-3 positions and oleic acid (Item Nos. 90260 | 24659) at the *sn*-2 position. It has been found in butterfat.¹

Reference

1. Kalo, P., Kempainen, A., Ollilainen, V., *et al.* Regiospecific determination of short-chain triacylglycerols in butterfat by normal-phase HPLC with on-line electrospray-tandem mass spectrometry. *Lipids* **39**(9), 915-928 (2004).

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, 12/21/2022

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