

PRODUCT INFORMATION



1-Palmitoyl-2-11(Z)-Octadecenoyl-3-Butyroyl-*rac*-glycerol

Item No. 28558

CAS Registry No.: 2692623-15-5
Formal Name: 11Z-octadecenoic acid, 1-[(1-oxobutoxy)methyl]-2-[(1-oxohexadecyl)oxy]ethyl ester

Synonyms: 1-Palmitin-2-Vaccenate-3-Butyrin,
1-Palmitoyl-2-*cis*-Vaccenoyl-3-Butyroyl-*rac*-glycerol,
TG(16:0/18:1(11Z)/4:0), TG(16:0/18:1/4:0)

MF: C₄₁H₇₆O₆

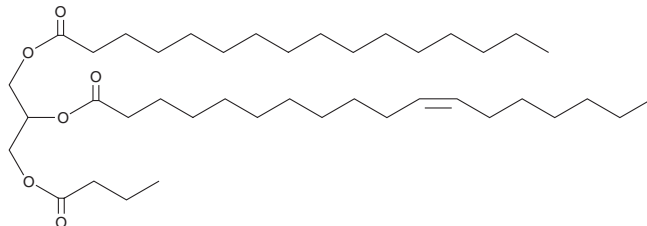
FW: 665.1

Purity: ≥98%

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-Palmitoyl-2-11(Z)-octadecenoyl-3-butyroyl-*rac*-glycerol is supplied as an oil. A stock solution may be made by dissolving the 1-palmitoyl-2-11(Z)-octadecenoyl-3-butyroyl-*rac*-glycerol in the solvent of choice, which should be purged with an inert gas. 1-Palmitoyl-2-11(Z)-octadecenoyl-3-butyroyl-*rac*-glycerol is soluble in chloroform and methanol.

Description

1-Palmitoyl-2-11(Z)-octadecenoyl-3-butyroyl-*rac*-glycerol is a triacylglycerol that contains palmitic acid (Item No. 10006627) at the *sn*-1 position, *cis*-vaccenic acid (Item No. 20023) at the *sn*-2 position, and butyric acid (Item No. 13121) at the *sn*-3 position.

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, 03/07/2023

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