

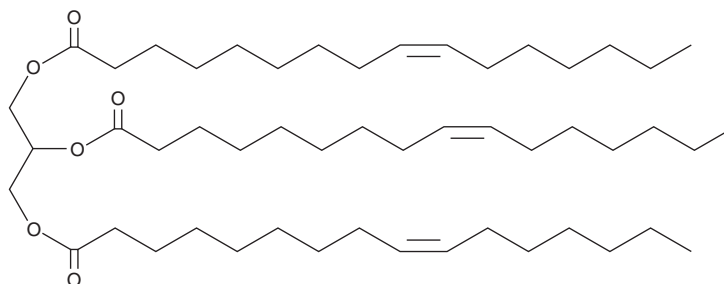
PRODUCT INFORMATION



1,2,3-Tripalmitoleoyl Glycerol

Item No. 26897

CAS Registry No.: 20246-55-3
Formal Name: (9Z,9'Z,9''Z)-9-hexadecenoic acid, 1,1',1''-(1,2,3-propanetriyl) ester
Synonyms: Glycerol Tripalmitoleate, TG(16:1/16:1/16:1), Tripalmitolein
MF: C₅₁H₉₂O₆
FW: 801.3
Purity: ≥98%
Supplied as: A liquid
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-Tripalmitoleoyl glycerol is supplied as a liquid. A stock solution may be made by dissolving the 1,2,3-tripalmitoleoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,2,3-Tripalmitoleoyl glycerol is soluble in the organic solvent chloroform.

Description

1,2,3-Tripalmitoleoyl glycerol is a triacylglycerol that contains palmitoleic acid (Item Nos. 10009871 | 21911) at the *sn*-1, *sn*-2, and *sn*-3 positions. It reduces red blood cell deformability in a concentration-dependent manner in a Reid's filtration assay.¹ Hepatic levels of 1,2,3-tripalmitoleoyl glycerol are increased in the JAK2L mouse model of hepatic steatosis.² 1,2,3-Tripalmitoleoyl glycerol plasma levels are decreased in renal patients before dialysis.³

References

1. Hong, C.Y., Lai, L.J., Shiao, M.S., *et al.* Effect of triacylglycerols on erythrocyte deformability in vitro. *Prostaglandins Leukot. Essent. Fatty Acids* **48(5)**, 351-353 (1993).
2. Wilson, C.G., Tran, J.L., Erion, D.M., *et al.* Hepatocyte-specific disruption of CD36 attenuates fatty liver and improves insulin sensitivity in HFD-fed mice. *Endocrinology* **157(2)**, 570-585 (2016).
3. Lee, D.M., Jackson, K.W., Knowlton, N., *et al.* Oxidative stress and inflammation in renal patients and healthy subjects. *PLoS One* **6(7)**, e22360 (2011).

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, 01/16/2026

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