

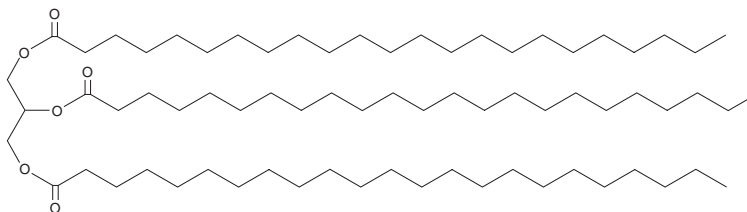
PRODUCT INFORMATION



1,2,3-Trtricosanoyl Glycerol

Item No. 26929

CAS Registry No.: 86850-72-8
Formal Name: tricosanoic acid,
1,2,3-propanetriyl ester
Synonyms: Glycerol Trtricosanoate,
TG(23:0/23:0/23:0),
Trtricosanoin
MF: C₇₂H₁₄₀O₆
FW: 1,101.9
Purity: ≥98%
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-Trtricosanoyl glycerol is supplied as a solid. A stock solution may be made by dissolving the 1,2,3-trtricosanoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,2,3-Trtricosanoyl glycerol is soluble in the organic solvent dimethyl formamide at a concentration of approximately 10 mg/ml.

Description

1,2,3-Trtricosanoyl glycerol is a triacylglycerol that contains tricosanoic acid (Item No. 26867) at the *sn*-1, *sn*-2, and *sn*-3 positions. It has been used as an internal standard for the quantification of fatty acids in the triglyceride component of human aortic endothelial cells (HAECs) grown in media supplemented with stearic and/or oleic acid.¹

Reference

1. Harvey, K.A., Walker, C.L., Xu, Z., *et al.* Oleic acid inhibits stearic acid-induced inhibition of cell growth and pro-inflammatory responses in human aortic endothelial cells. *J. Lipid Res.* **51(12)**, 3470-3480 (2010).

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, 10/25/2022

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