

PRODUCT INFORMATION



1,2,3-Trilignoceroyl Glycerol

Item No. 29055

CAS Registry No.: 56983-26-7

Formal Name: tetracosanoic acid,
1,1',1''-(1,2,3-propanetriyl) ester

Synonyms: Glycerol Trilignoceranoate,
TG(24:0/24:0/24:0),
1,2,3-Trilignocerin

MF: $C_{75}H_{146}O_6$

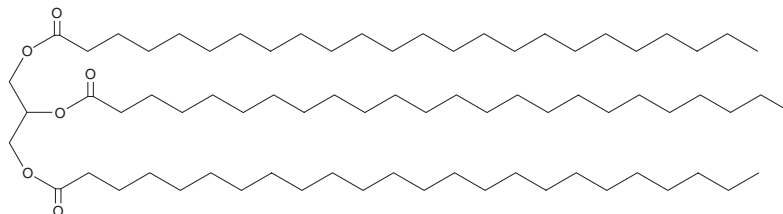
FW: 1,144.0

Purity: $\geq 95\%$

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-Trilignoceroyl glycerol is supplied as a solid. A stock solution may be made by dissolving the 1,2,3-trilignoceroyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,2,3-Trilignoceroyl glycerol is soluble in the organic solvent dimethyl formamide at a concentration of approximately 10 mg/ml.

Description

1,2,3-Trilignoceroyl glycerol is a triacylglycerol that contains lignoceric acid (Item No. 13353) at the *sn*-1, *sn*-2, and *sn*-3 positions. It has been found in goat milk fat.¹

Reference

1. Fontecha, J., Díaz, V., Fraga, M.J., *et al.* Triglyceride analysis by gas chromatography in assessment of authenticity of goat milk fat. *J. Am Chem. Soc.* **75**(12), 1893-1896 (1998).

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, 11/08/2022

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