

PRODUCT INFORMATION



1,2,3-Trielaidoyl Glycerol

Item No. 26950

CAS Registry No.: 537-39-3

Formal Name: (9E,9'E,9''E)-9-octadecenoic acid, 1,1,1''-(1,2,3-propanetriyl) ester

Synonyms: Glycerol Trielaidate, TG(18:1(9E)/18:1(9E)/18:1(9E)), Trielaidin

MF: $C_{57}H_{104}O_6$

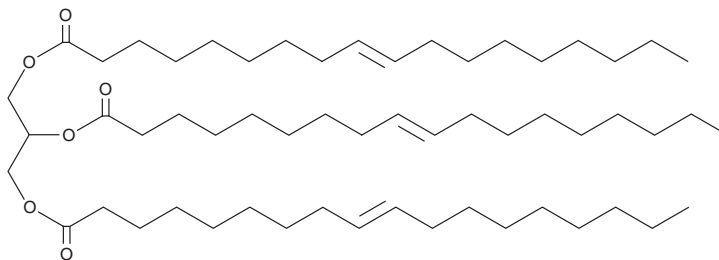
FW: 885.4

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

1,2,3-Trielaidoyl glycerol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 1,2,3-trielaidoyl glycerol should first be dissolved in chloroform and then diluted with the aqueous buffer of choice. 1,2,3-Trielaidoyl glycerol has a solubility of approximately 0.25 mg/ml in a 1:3 solution of chloroform:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

1,2,3-Trielaidoyl glycerol is a triacylglycerol that contains elaidic acid (Item No. 90250) at the *sn*-1, *sn*-2, and *sn*-3 positions. Dietary administration of 1,2,3-trielaidoyl glycerol reduces serum cholesterol levels in rats compared to corn oil-fed control animals.¹ 1,2,3-Trielaidoyl glycerol has been used as a substrate to determine the substrate specificity of *S. rimosus* extracellular lipase.²

References

1. Clifford, A.J., Smith, L.M., Creveling, R.K., et al. Effects of dietary triglycerides on serum and liver lipids and sterol excretion of rats. *J. Nutr.* **116**(6), 944-956 (1986).
2. Lešćić, I., Vukelić, B., Majerić-Elenkov, M., et al. Substrate specificity and effects of water-miscible solvents on the activity and stability of extracellular lipase from *Streptomyces rimosus*. *Enzyme Microb. Technol.* **29**(8-9), 548-553 (2001).

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, 12/05/2022

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