

PRODUCT INFORMATION

1,2,3-Tri-10(Z)-Pentadecenoyl Glycerol

Item No. 27004

Formal Name: propane-1,2,3-triyl (10Z,10'Z,10''Z)-tris(pentadec-10-enoate)

Synonyms: Glycerol Tri-10(Z)-Pentadecenoate, TG(15:1/15:1/15:1), 1,2,3-10(Z)-Tristadecenoyl Glycerol, Tri-10(Z)-Pentadecenoin

MF: $C_{48}H_{86}O_6$

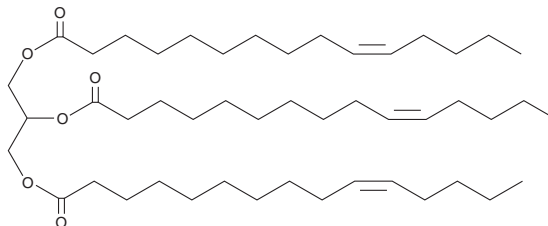
FW: 759.2

Purity: ≥98%

Supplied as: A solution in ethanol

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-Tri-10(Z)-pentadecenoyl glycerol is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as dimethyl formamide purged with an inert gas can be used. The solubility of 1,2,3-tri-10(Z)-pentadecenoyl glycerol in this solvent is approximately 10 mg/ml. 1,2,3-Tri-10(Z)-pentadecenoyl glycerol is also slightly soluble in chloroform.

1,2,3-Tri-10(Z)-pentadecenoyl glycerol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 1,2,3-tri-10(Z)-pentadecenoyl glycerol should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 1,2,3-Tri-10(Z)-pentadecenoyl glycerol has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

1,2,3-Tri-10(Z)-pentadecenoyl glycerol is a triacylglycerol that contains 10(Z)-pentadecenoic acid (Item No. 22362) at the *sn*-1, *sn*-2, and *sn*-3 positions.

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, 01/20/2026

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