

# 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)-Trientadecenoyl Glycerol, Tri-10(Z)-Pentadecenoin

**MF:** C<sub>48</sub>H<sub>86</sub>O<sub>6</sub>

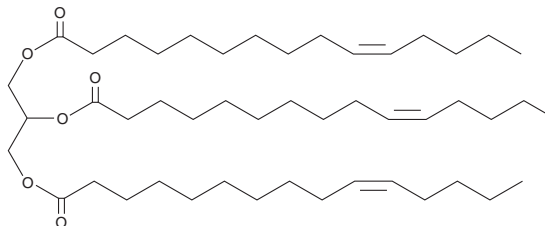
**FW:** 759.2

**Purity:** ≥98%

**Supplied as:** A liquid

**Storage:** -20°C

**Stability:** ≥2 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 liquid. A stock solution may be made by dissolving the 1,2,3-tri-10(Z)-pentadecenoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,2,3-Tri-10(Z)-pentadecenoyl glycerol is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of 1,2,3-tri-10(Z)-pentadecenoyl glycerol in these solvents 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, 06/08/2020

#### 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