

PRODUCT INFORMATION



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

Item No. 26935

CAS Registry No.: 93824-29-4

Formal Name: 10-undecenoic acid, 1,2,3-propanetriyl ester

Synonyms: Glycerol Tri-10(Z)-Undecenoate, TG(11:1/11:1/11:1), Tri-10(Z)-Undecenoyl Glycerol, Triundecylenoyl Glycerol, 10(Z)-Undecenoin

MF: $C_{36}H_{62}O_6$

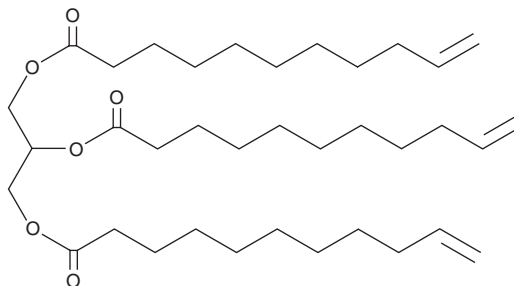
FW: 590.9

Purity: $\geq 98\%$

Supplied as: A liquid

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)-undecenoyl glycerol is supplied as a liquid. A stock solution may be made by dissolving the 1,2,3-tri-10(Z)-undecenoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,2,3-Tri-10(Z)-undecenoyl glycerol is slightly soluble in chloroform.

Description

1,2,3-Tri-10(Z)-undecenoyl glycerol is a triacylglycerol that contains 10(Z)-undecenoic acid 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, 01/19/2026

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