

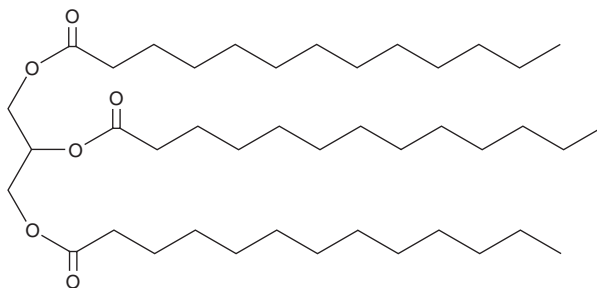
PRODUCT INFORMATION



1,2,3-Tridecanoyl Glycerol

Item No. 26899

CAS Registry No.: 26536-12-9
Formal Name: tridecanoic acid, 1,1',1''-(1,2,3-propanetriyl) ester
Synonyms: Glycerol Tridecanoyl, 13:0 TAG, 13:0/13:0/13:0-TG, TG(13:0/13:0/13:0), Tridecanoin
MF: $C_{42}H_{80}O_6$
FW: 681.1
Purity: $\geq 98\%$
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-Tridecanoyl glycerol is supplied as a solid. A stock solution may be made by dissolving the 1,2,3-tridecanoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,2,3-Tridecanoyl glycerol is soluble in the organic solvent dimethyl formamide at a concentration of approximately 10 mg/ml.

Description

1,2,3-Tridecanoyl glycerol is a synthetic triacylglycerol that contains tridecanoic acid (Item No. 19725) in the *sn*-1, *sn*-2, and *sn*-3 positions.¹ It has been used as a standard for the relative quantification of triacylglycerols in *C. elegans* fat stores. It has also been used as an internal standard for the quantification of triacylglycerols in the serum and liver of adult rat offspring receiving dietary conjugated linoleic acids during and after gestation.²

References

1. Pino, E.C., Webster, C.M., Carr, C.E., *et al.* Biochemical and high throughput microscopic assessment of fat mass in *Caenorhabditis elegans*. *J. Vis. Exp.* **73**, e50180 (2013).
2. Lavandera, J., Gerstner, C.D., Sain, J., *et al.* Maternal conjugated linoleic acid modulates TAG metabolism in adult rat offspring. *Br. J. Nutr.* **118**(11), 906-913 (2017).

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, 10/10/2022

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