

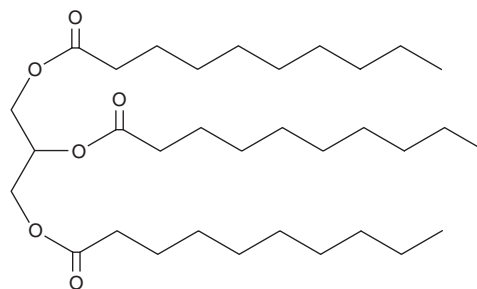
# PRODUCT INFORMATION



## 1,2,3-Tridecanoyl Glycerol

Item No. 26958

**CAS Registry No.:** 621-71-6  
**Formal Name:** decanoic acid, 1,1',1''-(1,2,3-propanetriyl) ester  
**Synonyms:** Glycerol Tricaprate, Glycerol Tridecanoate, TG(10:0/10:0/10:0), Tricaprin, Tridecanoin, Tridecanoylglycerol  
**MF:**  $C_{33}H_{62}O_6$   
**FW:** 554.8  
**Purity:**  $\geq 98\%$   
**Supplied as:** A solid  
**Storage:**  $-20^{\circ}\text{C}$   
**Stability:**  $\geq 4$  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 chloroform at a concentration of approximately 100 mg/ml.

### Description

1,2,3-Tridecanoyl glycerol is a triacylglycerol that contains decanoic acid (Item No. 20838) at the *sn*-1, *sn*-2, and *sn*-3 positions. It has been found in human breast milk.<sup>1</sup> 1,2,3-Tridecanoyl glycerol has been used as a standard for the quantification of triacylglycerols in fish oils by RP-HPLC.<sup>2</sup> Formulations containing 1,2,3-tridecanoyl glycerol have been used in cosmetic products as thickening and skin-conditioning agents.<sup>3</sup>

### References

1. Kim, K.-M., Park, T.-S., and Shim, S.-M. Optimization and validation of HRLC-MS method to identify and quantify triacylglycerol molecular species in human milk. *Anal. Methods* **7**(10), 4362-4370 (2015).
2. Solaesa, Á.G., Bucio, S.L., Sanz, M.T., *et al.* Characterization of triacylglycerol composition of fish oils by using chromatographic techniques. *J. Oleo. Sci.* **63**(5), 449-460 (2014).
3. Johnson, W., Jr. and C.I.R.E. Panel. Final report on the safety assessment of trilaurin, triarachidin, tribehenin, tricaprin, tricapylin, trierucin, triheptanoin, triheptylundecanoin, triisononanoin, triisopalmitin, triisostearin, trilinolein, trimyristin, trioctanoin, triolein, tripalmitin, tripalmitolein, tricinolein, tristearin, triundecanoin, glyceryl triacetyl hydroxystearate, glyceryl triacetyl ricinoleate, and glyceryl stearate diacetate. *Int. J. Toxicol.* **20**(Suppl 4), 61-94 (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, 11/07/2022

#### 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](http://WWW.CAYMANCHEM.COM)