

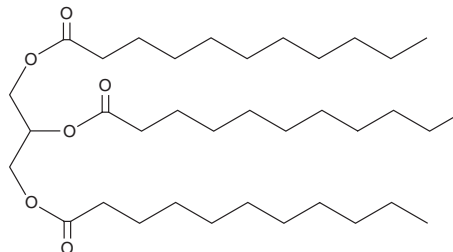
# PRODUCT INFORMATION



## 1,2,3-Triundecanoyl Glycerol

Item No. 27008

**CAS Registry No.:** 13552-80-2  
**Formal Name:** undecanoic acid, 1,1',1''-(1,2,3-propanetriyl) ester  
**Synonyms:** Glycerol Triundecanoate, TG(11:0/11:0/11:0), Triundecanoin, Triundecanoyl Glycerol  
**MF:** C<sub>36</sub>H<sub>68</sub>O<sub>6</sub>  
**FW:** 596.9  
**Purity:** ≥95%  
**Supplied as:** A crystalline 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-Triundecanoyl glycerol is supplied as a crystalline solid. A stock solution may be made by dissolving the 1,2,3-triundecanoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,2,3-Triundecanoyl glycerol is soluble in ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 1,2,3-triundecanoyl glycerol in these solvents is approximately 25 mg/ml in ethanol and 20 mg/ml in DMSO and DMF.

### Description

1,2,3-Triundecanoyl glycerol is a triacylglycerol that contains undecanoic acid (Item No. 19721) at the *sn*-1, *sn*-2, and *sn*-3 positions. Dietary administration of a 7:3 (w/w) mixture of 1,2,3-triundecanoyl glycerol and corn oil prevents fasting-induced decreases in plasma glucose, alanine, and insulin levels, as well as the ratio of insulin to glucagon, in rats compared with dietary administration of corn oil alone.<sup>1</sup>

### Reference

1. Pi-Sunyer, F.X., Conway, J.M., Lavau, M., *et al.* Glucagon, insulin, and gluconeogenesis in fasted odd carbon fatty acid-enriched rats. *Am. J. Physiol.* **231(2)**, 366-369 (1976).

#### 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, 12/08/2022

#### 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