

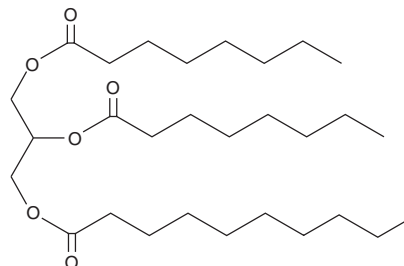
# PRODUCT INFORMATION



## 1,2-Dioctanoyl-3-Decanoyl-*rac*-glycerol

Item No. 28554

**CAS Registry No.:** 69095-45-0  
**Formal Name:** decanoic acid, 2,3-bis[(1-oxooctyl)oxy] propyl ester  
**Synonyms:** 1,2-Octanoin-3-Decanoin, TG(8:0/8:0/10:0)  
**MF:** C<sub>29</sub>H<sub>54</sub>O<sub>6</sub>  
**FW:** 498.7  
**Purity:** ≥95%  
**Supplied as:** A solution in chloroform  
**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-Dioctanoyl-3-decanoyl-*rac*-glycerol is supplied as a solution in chloroform. To change the solvent, simply evaporate the chloroform under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of 1,2-dioctanoyl-3-decanoyl-*rac*-glycerol in these solvents is approximately 30 mg/ml.

### Description

1,2-Dioctanoyl-3-decanoyl-*rac*-glycerol is a triacylglycerol that contains octanoic acid at the *sn*-1 and *sn*-2 positions and decanoic acid (Item No. 20838) at the *sn*-3 position. It has been found in the suberin of wound-healing potatoes.<sup>1</sup>

### Reference

1. Wang, W., Tian, S., and Stark, R.E. Isolation and identification of triglycerides and ester oligomers from partial degradation of potato suberin. *J. Agric. Food Chem.* **58**(2), 1040-1045 (2010).

#### 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, 11/08/2022

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