

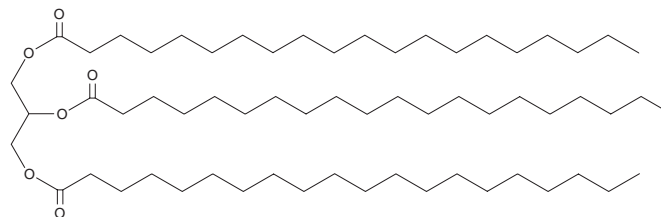
# PRODUCT INFORMATION



## 1,2,3-Triarachidoyl Glycerol

Item No. 23333

**CAS Registry No.:** 620-64-4  
**Formal Name:** eicosanoic acid, 1,2,3-propanetriyl ester  
**Synonym:** Dynasan 120, TG(20:0/20:0/20:0),  
Trieicosanoate  
**MF:**  $C_{63}H_{122}O_6$   
**FW:** 975.6  
**Purity:**  $\geq 98\%$   
**Supplied as:** A crystalline 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-Triarachidoyl glycerol is supplied as a crystalline solid. A stock solution may be made by dissolving the 1,2,3-triarachidoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,2,3-Triarachidoyl glycerol is soluble in the organic solvent chloroform at a concentration of approximately 10 mg/ml.

### Description

1,2,3-Triarachidoyl glycerol is a triacylglycerol containing three arachidic acid (Item No. 9000339 | 21906) groups. It has been found as a component of tomato seed oil.<sup>1</sup>

### Reference

1. Ma, Y., Ma, J., Yang, T., *et al.* Components, antioxidant and antibacterial activity of tomato seed oil. *Food Sci. Technol. Res.* **20(1)**, (2014).

#### 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, 12/02/2022

#### 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)