

# PRODUCT INFORMATION



## 1,2,3-Triarachidonoyl Glycerol

Item No. 10009826

CAS Registry No.: 23314-57-0

Formal Name: (5Z,5'Z,5''Z,8Z,8'Z,8''Z,11Z,11'Z,11''Z,14Z,14'Z,14''Z)-5,8,11,14-eicosatetraenoic acid, 1,1,1''-(1,2,3-propanetriyl) ester

Synonyms: Glycerol Triarachidonoyl, Glycerol Trieicosatetraenoate, Optimar, TG(20:4/20:4/20:4), Triarachidonin

MF:  $C_{63}H_{98}O_6$

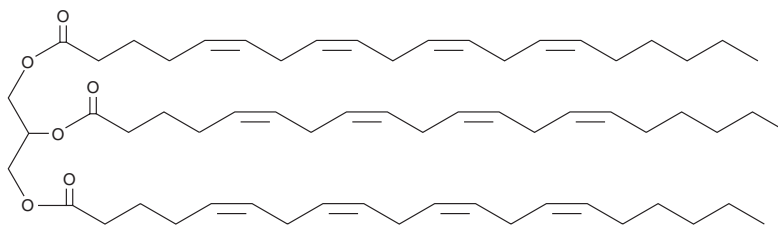
FW: 951.5

Purity:  $\geq 98\%$

Supplied as: A solution in ethanol

Storage:  $-20^{\circ}C$

Stability:  $\geq 2$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

1,2,3-Triarachidonoyl glycerol is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. The solvent dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of 1,2,3-triarachidonoyl glycerol in DMF is approximately 10 mg/ml.

1,2,3-Triarachidonoyl glycerol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of 1,2,3-triarachidonoyl glycerol should be diluted with the aqueous buffer of choice. 1,2,3-Triarachidonoyl glycerol has a solubility of approximately 500  $\mu\text{g/ml}$  in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

1,2,3-Triarachidonoyl glycerol is a polyunsaturated triacylglycerol with arachidonoyl (20:4) side-chains attached at the C-1, C-2, and C-3 positions that is useful in lipid research.

#### 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