

PRODUCT INFORMATION



1,3-Distearoyl-2-Myristoyl Glycerol

Item No. 28860

CAS Registry No.: 57396-99-3

Formal Name: octadecanoic acid,
1,1'-[2-[(1-oxotetradecyl)oxy]-1,3-
propanediyl] ester

Synonyms: 1,2-Stearin-2-Myristin,
TG(18:0/14:0/18:0)

MF: $C_{53}H_{102}O_6$

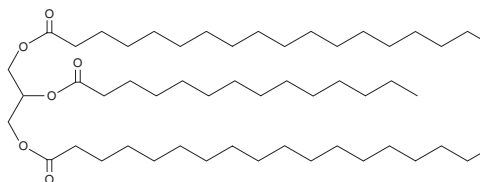
FW: 835.4

Purity: $\geq 95\%$

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

1,3-Distearoyl-2-myristoyl glycerol is supplied as a crystalline solid. A stock solution may be made by dissolving the 1,3-distearoyl-2-myristoyl glycerol in the solvent of choice, which should be purged with an inert gas. 1,3-Distearoyl-2-myristoyl glycerol is soluble in the organic solvent chloroform. The solubility of 1,3-distearoyl-2-myristoyl glycerol in chloroform is approximately 30 mg/ml.

Description

1,3-Distearoyl-2-myristoyl glycerol is a triacylglycerol that contains stearic acid (Item No. 10011298) at the *sn*-1 and *sn*-3 positions and myristic acid (Item No. 13351) at the *sn*-2 position. It has been found in butterfat.¹

Reference

1. Kalo, P., Kemppinen, A., and Ollilainen, V. Determination of triacylglycerols in butterfat by normal-phase HPLC and electrospray-tandem mass spectrometry. *Lipids* **44**(2), 169-195 (2009).

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, 12/12/2022

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