

# PRODUCT INFORMATION



## Non-polar Lipid Mixture 2 (plant, ovine)

Item No. 29377

**Supplied as:** A solution in chloroform

**Storage:** -20°C

**Stability:** ≥2 years

**Concentration:** 25 mg/ml

**Special Conditions:** This standard is accurately prepared by gravimetric technique (+/-0.5%)

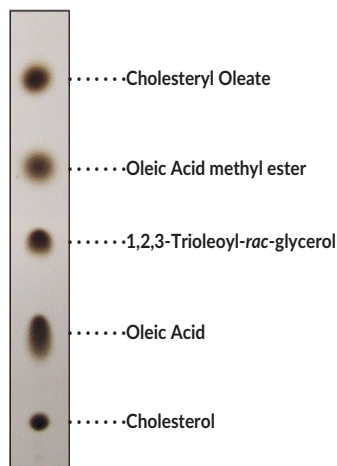
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Description and Data

Non-polar Lipid Mixture 1 contains cholesteryl oleate, oleic acid methyl ester, 1,2,3-trioleoyl-*rac*-glycerol, oleic acid, and cholesterol. It is intended for use as a reference standard for the detection of these non-polar lipids in various sample types by TLC, LC-, or GC-MS.

Included are:

Elution Order*	Item Name
1	Cholesteryl Oleate
2	Oleic Acid methyl ester
3	1,2,3-Trioлеoyl- <i>rac</i> -glycerol
4	Oleic Acid
5	Cholesterol



#### 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, 10/24/2022

#### 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