

PRODUCT INFORMATION



Cannabis Lipid Residue Mixture

Item No. 41690

Supplied as: A solution in methanol (1 mg/ml of each compound)

Storage: -20°C

Stability: ≥3 years

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

Description

Cannabis Lipid Residue Mixture (Item No. 41690) is a mixture of analytical reference standards containing compounds that can be isolated from plants in the genus *Cannabis*. The mixture contains the medium-chain triglycerides (MCTs) 1,2,3-trihexanoyl glycerol, 1,2,3-trioctanoyl glycerol, 1,2,3-tridecanoyl glycerol, and 1,2,3-trilauroyl glycerol, as well as vitamin E acetate, also known as (+)-α-tocopherol acetate. The mixture is supplied in an amber ampule in which the headspace has been purged with argon. It is designed for use in method development and as a reference standard for these substances during routine analytical checks using HPLC and GC-MS instrumentation. This product is intended for research and forensic applications.

Contents

Item No. 41690	Cannabis Lipid Residue Mixture	
Item Name	CAS Number	Concentration (mg/ml)
1,2,3-Trihexanoyl Glycerol	621-70-5	1.0
1,2,3-Trioctanoyl Glycerol	538-23-8	1.0
1,2,3-Tridecanoyl Glycerol	621-71-6	1.0
1,2,3-Trilauroyl Glycerol	538-24-9	1.0
(+)-α-Tocopherol Acetate	58-95-7	1.0

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, 08/29/2024

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