

PRODUCT INFORMATION



NHI-D Mix

Item No. 41447

Supplied as: A neat liquid

Storage: -20°C

Stability: ≥3 years

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

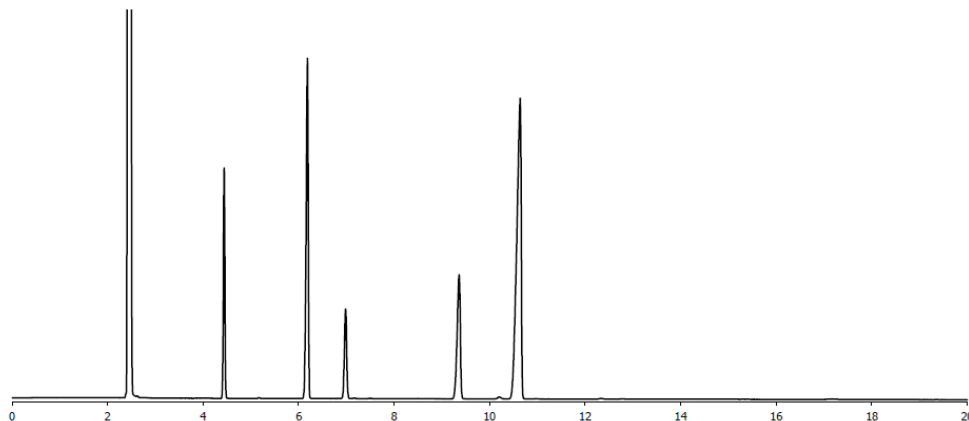
Description and Data

NHI-D Mix contains five fatty acid methyl esters (FAMES): myristic, palmitic, palmitoleic, stearic, and oleic. This product is intended as a reference standard for the detection of these fatty acids in various sample types by GC-FID and GC-MS. This mixture is supplied in an amber ampule in which the headspace has been purged with argon to prevent lipid oxidation.

Included are:

Elution Order*	Item Name	% Concentrations by Weight
1	Methyl tetradecanoate (myristate)	12.0
2	Methyl hexadecanoate (palmitate)	24.0
3	Methyl <i>cis</i> -9-hexadecenoate (palmitoleate)	7.0
4	Methyl octadecanoate (stearate)	13.0
5	Methyl <i>cis</i> -9-octadecenoate (oleate)	44.0

*GC Column Used: SP-2330, 30m x 0.25mm x 0.2 µm



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, 04/29/2026

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