

# PRODUCT INFORMATION



## Polyunsaturated Fatty Acid LC-MS Mixture

Item No. 17941

**Purity:** ≥98% for each compound  
**Supplied as:** A solution in ethanol (10 µg/ml of each compound)  
**Storage:** -20°C  
**Stability:** ≥5 years

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

## Description

This mixture contains various polyunsaturated fatty acids, including many important essential fatty acids. The mixture is supplied in an amber ampule in which the headspace has been purged with argon to prevent lipid oxidation. This product has been designed for direct use in LC-MS applications. The solution may be serially diluted for preparation of calibrators and QC standards and/or used directly as a system suitability standard or tuning standard. After opening, we recommend that the mixture be transferred immediately to a 1 ml glass screw cap vial, to prevent solvent evaporation, and stored at -20°C. The mixture should be discarded after multiple freeze/thaw cycles.

The polyunsaturated fatty acids represented in this mixture are composed of 18-22 carbons and include many of the key  $\omega$ -3 and  $\omega$ -6 fatty acids. These compounds are either produced endogenously or found naturally in plants and marine organisms and most have significant biological roles as eicosanoid precursors. Some compounds have also demonstrated protective effects against inflammatory diseases, coronary heart disease, thrombosis, or ischemic brain injury.

## Contents

Item Number	Item Name
90300	Adrenic Acid
90010	Arachidonic Acid
90230	Dihomo- $\gamma$ -Linolenic Acid
90310	Docosahexaenoic Acid
90165	Docosapentaenoic Acid
90110	Eicosapentaenoic Acid
90150	Linoleic Acid
90210	$\alpha$ -Linolenic Acid
90220	$\gamma$ -Linolenic Acid
90320	Stearidonic Acid

### 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, 03/24/2022

### 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