

PRODUCT INFORMATION



Fatty Acid ethyl ester Standard Pack

Item No. 10008188

Purity:	≥98%
Supplied as:	A solution in ethanol
Storage:	-20°C
Stability:	≥1 year

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

Description

The fatty acid ethyl ester standard pack contains nine vials each containing 100 mg of an individual fatty acid ethyl ester.

This pack contains stearic acid ethyl ester (Item No. 10008196), myristic acid ethyl ester (Item No. 10008197), linoleic acid ethyl ester (Item No. 10008198), α -linolenic acid ethyl ester (Item No. 10008199), arachidonic acid ethyl ester (Item No. 10008200), oleic acid ethyl ester (Item No. 10008201), palmitic acid ethyl ester (Item No. 10008202), lauric acid ethyl ester (Item No. 10008203), and palmitoleic acid ethyl ester (Item No. 10008204).

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, 07/16/2019

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