

PRODUCT INFORMATION



ω -3 Hydroxy Acid HPLC Mixture

Item No. 34003

Purity: \geq 98% for each compound
Supplied as: A solution in ethanol (5 μ g of each compound)
Storage: -20°C
Stability: \geq 1 year

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

Laboratory Procedures

The ω -3 hydroxy acid HPLC mixture contains the major lipoxygenation products of some ω -3 fatty acids supplied in ethanol. This HPLC mixture contains 5 μ g of the following ω -3 fatty acids: 5(S)-HEPE, 12(S)-HEPE, 15(S)-HEPE, 13(S)-HOTrE, and 15(S)-HETrE.

| Name: | CAS No.: |
|-------------|-------------|
| 5(S)-HEPE | 92008-51-0 |
| 12(S)-HEPE | 116180-17-7 |
| 15(S)-HEPE | 86282-92-0 |
| 13(S)-HOTrE | 87984-82-5 |
| 15(S)-HETrE | 92693-02-2 |

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, 07/25/2022

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