

PRODUCT INFORMATION



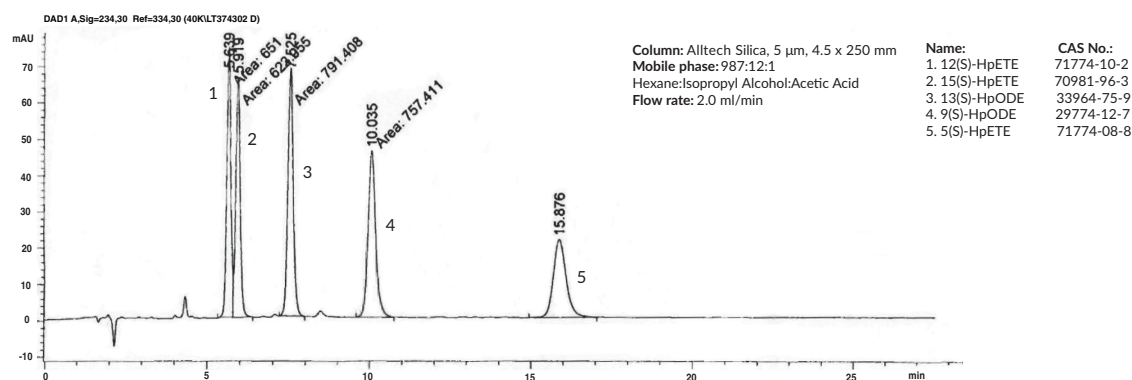
Hydroperoxy HPLC Mixture Item No. 44001

Purity: ≥98% for each compound
Supplied as: A solution in ethanol
Storage: -80°C
Stability: ≥6 months

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

Description and Data

The hydroperoxy HPLC mixture contains several of the principle lipoxygenase-derived hydroperoxides of arachidonic acid and linoleic acid supplied in ethanol. This HPLC mixture contains 5 µg of the following fatty acid hydroperoxides: 5(S)-HpETE, 12(S)-HpETE, 15(S)-HpETE, 9(S)-HpODE, and 13(S)-HpODE.



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, 07/29/2022

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