

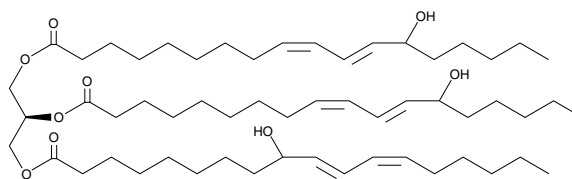
Product Information



Hydroxy Linoleins

Item No. 89420

Purity: ≥98% (A mixture of 132 isomers)
Stability: ≥2 years at -20°C
Supplied as: A solution in ethanol
UV/Vis.: λ_{max} : 234 nm ϵ : 25,000



(only one isomer shown)

Laboratory Procedures

For long term storage, we suggest that hydroxy linoleins be stored as supplied at -20°C. It should be stable for at least two years.

Hydroxy linoleins are supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. Hydroxy linoleins are miscible in these solvents.

Hydroxy linoleins are sparingly soluble in aqueous buffers. An ethanolic solution of hydroxy linoleins at 0.1 mg/ml can be diluted with PBS (pH 7.2) up to 10 times resulting in a colloidal suspension. The suspension should be kept at room temperature and used within two hours or else the compound will separate from the bulk of the liquid.

Hydroxy linoleins are a mixture of 132 possible isomers of mono-, di-, and tri-hydroxy compounds produced by the autoxidation of trilinolein and subsequent reduction of the hydroperoxides. Autoxidation of linoleic acid containing triglycerides (i.e., trilinolein) *in vivo* could result in the formation of hydroperoxides, which could then be reduced by peroxidases to produce hydroxy linoleins. Due to their inherent chemical stability compared to the precursor hydroperoxides, hydroxy linoleins are potential markers of triglyceride lipid peroxidation *in vivo*.

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/89420

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 06/19/2012