

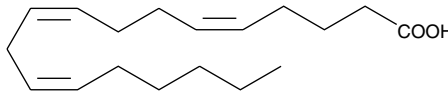
Product Information



Pinolenic Acid

Item No. 10008654

CAS Registry No.: 16833-54-8
Formal Name: 5Z,9Z,12Z-octadecatrienoic acid
MF: C₁₈H₃₀O₂
FW: 278.4
Purity: ≥98%
Stability: ≥1 year at -20°C
Supplied as: A solution in ethanol



Laboratory Procedures

For long term storage, we suggest that pinolenic acid be stored as supplied at -20°C. It should be stable for at least one year.

Pinolenic acid is 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. The solubility of pinolenic acid in these solvents is approximately 30 mg/ml.

Pinolenic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of pinolenic acid should be diluted with the aqueous buffer of choice. Pinolenic acid has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Pinolenic acid is a polyunsaturated fatty acid found in Korean pine (*Pinus orientalis*) and maritime pine (*Pinus pinaster*) seed oils. Both oils have been found to have lipid-lowering properties. A diet containing maritime pine seed oil (MPSO) lowered HDL and ApoA1 levels in transgenic mice expressing human ApoA1. MPSO was found to diminish cholesterol efflux *in vitro*.¹ Korean pine seed oil supplements may help in obesity by reduction of appetite. People taking this oil had an increase in the satiety hormones CCK and GLP-1 and a reduced desire to eat.² The activity of the oil is attributed to pinolenic acid. Pinolenic acid is not converted to arachidonic acid metabolically and can reduce arachidonic acid levels in the phosphatidylinositol fraction of HepG2 cells from 15.9% to 7.0%.³

References

1. Asset, G., Leroy, A., Bauge, E., *et al.* Effects of dietary maritime pine (*Pinus pinaster*)-seed oil on high-density lipoprotein levels and *in vitro* cholesterol efflux in mice expressing human apolipoprotein A-I. *British Journal of Nutrition* **84**, 353-360 (2000).
2. Causey, J.L. Korean pine nut fatty acids induce satiety-producing hormone release in overweight human volunteers. *The 231st ACS National Meeting, Atlanta, GA, March 26-30, 2006.*
3. Tamotsu, T., Tatsunori, T., Morishige, J., *et al.* Non-methylene-interrupted polyunsaturated fatty acids: Effective substitute for arachidonate of phosphatidylinositol. *Biochem. Biophys. Res. Commun.* **264**, 683-688 (1999).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10008654

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

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 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, 05/07/2014

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