

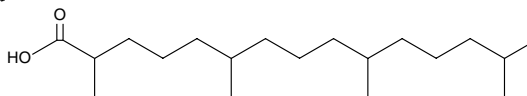
Product Information



Pristanic Acid

Item No. 17264

CAS Registry No.: 1189-37-3
Formal Name: 2,6,10,14-tetramethyl-pentadecanoic acid
MF: C₁₉H₃₈O₂
FW: 298.5
Purity: ≥95%
Stability: ≥1 year at -20°C
Supplied as: A solution in ethanol



Laboratory Procedures

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

Pristanic acid is supplied as a solution in ethanol. To change the solvent, simply evaporate the pristanic acid under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of pristanic acid in these solvents is approximately 20, 10, and 30 mg/ml, respectively.

Pristanic acid is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Pristanic acid is a natural terpenoid acid present at micromolar concentrations in the plasma of healthy individuals. It is obtained from dietary sources or by the α -oxidation of phytanic acid (Item No. 90360), which occurs predominantly in peroxisomes.^{1,2} Pristanic acid is an agonist of PPAR α (EC₅₀ = 40 μ M).³ It is also a substrate of the mammalian peroxisomal ATP-binding cassette (ABC) transporter ABCD3, which facilitates its delivery into peroxisomes for β -oxidation.⁴

References

1. Verhoeven, N.M., Wanders, R.J., Poll-The, B.T., *et al.* The metabolism of phytanic acid and pristanic acid in man: A review. *J. Inherit. Metab. Dis.* **21**, 697-728 (1998).
2. Verhoeven, N.M. and Jakobs, C. Human metabolism of phytanic acid and pristanic acid. *Prog. Lipid Res.* **40**, 453-466 (2001).
3. Zomer, A.W.M., van der Burg, B., Jansen, G.A., *et al.* Pristanic acid and phytanic acid: Naturally occurring ligands for the nuclear receptor peroxisome proliferator-activated receptor α . *J. Lipid Res.* **41**(11), 1801-1807 (2000).
4. Kemp, S., Theodoulou, F.L., and Wanders, R.J.A. Mammalian peroxisomal ABC transporters: From endogenous substrates to pathology and clinical significance. *Br. J. Pharmacol.* **164**(7), 1753-1766 (2011).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/17264

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, 05/11/2015

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