

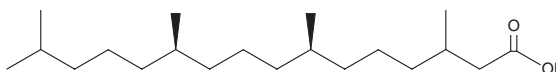
PRODUCT INFORMATION



Phytanic Acid

Item No. 90360

CAS Registry No.: 199484-69-0
Formal Name: (7R,11R)-3,7,11,15-tetramethylhexadecanoic acid
MF: C₂₀H₄₀O₂
FW: 312.5
Purity: ≥96%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥1 year



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

Laboratory Procedures

Phytanic 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 phytanic acid in these solvents is approximately 100 mg/ml.

Phytanic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of phytanic acid should be diluted with the aqueous buffer of choice. Phytanic acid has a solubility of 1 mg/ml in a 0.15 M Tris-HCl (pH 8.5) solution using this method. Therefore, further dilutions of the organic solvent solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that 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.

Description

Phytanic acid is a saturated 20-carbon branched-chain fatty acid which can only be derived from dietary sources. In cases of peroxisomal disorders, elevated levels of phytanic acid have been observed, and this is linked to diseases such as infantile phytanic acid storage disease and Refsum's disease.¹

Reference

1. Wanders, R.J.A., Boltshauser, E., Steinmann, B., *et al.* Infantile phytanic acid storage disease, a disorder of peroxisome biogenesis: A case report. *J. Neurol. Sci.* **98**, 1-11 (1990).

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, 03/24/2022

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