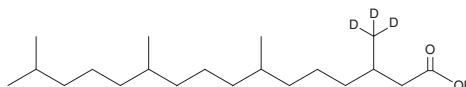


PRODUCT INFORMATION



Phytanic Acid-d₃ Item No. 29068

CAS Registry No.: 1383920-40-8
Formal Name: 7,11,15-trimethyl-3-(methyl-d₃)
hexadecanoic acid
MF: C₂₀H₃₇D₃O₂
FW: 315.6
Chemical Purity: ≥95% (Phytanic Acid)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Phytanic acid-d₃ is intended for use as an internal standard for the quantification of phytanic acid (Item No. 90360) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Phytanic acid-d₃ 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 chloroform, and methanol purged with an inert gas can be used.

Description

Phytanic acid is a saturated 20-carbon branched-chain fatty acid which can only be derived from dietary sources.¹ 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.

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)**, 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, 11/21/2022

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