

PRODUCT INFORMATION



Perfluorotetradecanoic Acid

Item No. 37255

CAS Registry No.: 376-06-7
Formal Name: 2,2,3,3,4,4,5,5,6,6,7,7,8,8,9,9,10,10,11,11,12,12,13,13,14,14,14-heptacosafuoro-tetradecanoic acid

Synonyms: Perfluoromyristic Acid, PFMA, PFTeDA

MF: C₁₄HF₂₇O₂

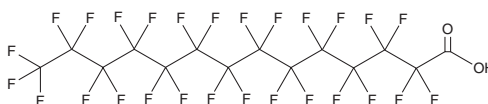
FW: 714.1

Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

Perfluorotetradecanoic acid (PFTeDA) is supplied as a solid. A stock solution may be made by dissolving the PFTeDA in the solvent of choice, which should be purged with an inert gas. PFTeDA is soluble in the organic solvent ethanol at a concentration of approximately 5 mg/ml.

Description

PFTeDA is a perfluoroalkyl substance (PFAS).¹ PFTeDA (20 mg/kg) decreases serum testosterone levels and epididymal sperm count at 56 days post-Leydig cell ablation in a rat model of Leydig cell regeneration induced by ethane dimethyl sulfonate (EDS). It has been found in landfill leachate and shellfish.^{2,3}

References

1. Zhang, S., Wen, Z., Li, X., *et al.* Short-term exposure to perfluorotetradecanoic acid affects the late-stage regeneration of Leydig cells in adult male rats. *Toxicol. Appl. Pharmacol.* **433**, 115777 (2021).
2. Capozzi, S.L., Leang, A.L., Rodenburg, L.A., *et al.* PFAS in municipal landfill leachate: Occurrence, transformation, and sources. *Chemosphere* **334**, 138924 (2023).
3. Catherine, M., Nadège, B., Charles, P., *et al.* Perfluoroalkyl substances (PFASs) in the marine environment: Spatial distribution and temporal profile shifts in shellfish from French coasts. *Chemosphere* **228**, (2019).

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, 09/29/2023

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