

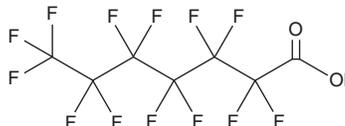
PRODUCT INFORMATION



Perfluoroheptanoic Acid

Item No. 37247

CAS Registry No.: 375-85-9
Formal Name: 2,2,3,3,4,4,5,5,6,6,7,7,7-tridecafluoro-heptanoic acid
Synonym: PFHpA
MF: C₇HF₁₃O₂
FW: 364.1
Purity: ≥90%
Supplied as: A low-melting 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

Perfluoroheptanoic acid (PFHpA) is supplied as a low-melting solid. Aqueous solutions of PFHpA can be prepared by directly dissolving the low-melting solid in aqueous buffers. The solubility of PFHpA in PBS (pH 7.2) is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

PFHpA is a perfluoroalkyl substance (PFAS).¹ Prenatal exposure to PFHpA is positively correlated with preterm birth, primarily in male infants. It is also positively correlated with fat mass, body mass index (BMI), and waist circumference in females but negatively associated with BMI, waist circumference, and weight-to-height ratio in males.²

References

1. Liao, Q., Tang, P., Song, Y., *et al.* Association of single and multiple prefluoroalkyl substances exposure with preterm birth: Results from a Chinese birth cohort study. *Chemosphere* **307(Pt 1)**, 135741 (2022).
2. Zhang, S., Lei, X., Zhang, Y., *et al.* Prenatal exposure to per- and polyfluoroalkyl substances and childhood adiposity at 7 years of age. *Chemosphere* **307(Pt 4)**, 136077 (2022).

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, 12/08/2022

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