

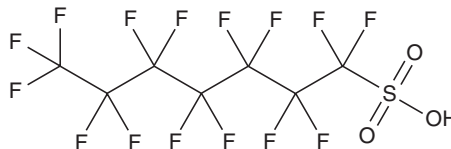
PRODUCT INFORMATION



Perfluoroheptanesulfonic Acid

Item No. 37246

CAS Registry No.: 375-92-8
Formal Name: 1,1,2,2,3,3,4,4,5,5,6,6,7,7,7-pentadecafluoro-1-heptanesulfonic acid
Synonym: PFHpS
MF: C₇HF₁₅O₃S
FW: 450.1
Purity: ≥95%
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

Perfluoroheptanesulfonic acid (PFHpS) is supplied as a solid. A stock solution may be made by dissolving the PFHpS in the solvent of choice, which should be purged with an inert gas. PFHpS is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of PFHpS in ethanol is approximately 5 mg/ml and approximately 10 mg/ml in DMSO and DMF.

Description

PFHpS is a perfluoroalkyl substance (PFAS).¹ PFHpS induces malformations in zebrafish larvae (EC₅₀ = 168.1 μM). It has been found in landfill leachate, and prenatal exposure to PFHpS is associated with decreased birth weight.^{2,3}

References

1. Gaballah, S., Swank, A., Sobus, J.R., *et al.* Evaluation of developmental toxicity, developmental neurotoxicity, and tissue dose in zebrafish exposed to GenX and other PFAS. *Environ. Health Perspect.* **128**(4), 47005 (2020).
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. Gui, S.Y., Chen, Y.-N., Wu, K.-J., *et al.* Association between exposure to per- and polyfluoroalkyl substances and birth outcomes: A systematic review and meta-analysis. *Front. Public Health* **10**, 855348 (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, 11/14/2023

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