

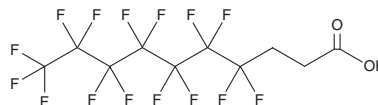
PRODUCT INFORMATION



2H,2H,3H,3H-Perfluorodecanoic Acid

Item No. 37271

CAS Registry No.: 812-70-4
Formal Name: 4,4,5,5,6,6,7,7,8,8,9,9,10,10,10-pentadecafluorodecanoic acid
Synonyms: 7:3 FTCA,
2H,2H,3H,3H-Pentadecafluorodecanoic Acid,
2H,2H,3H,3H-PFDA
MF: $C_{10}H_5F_{15}O_2$
FW: 442.1
Purity: $\geq 80\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

2H,2H,3H,3H-Perfluorodecanoic acid is supplied as a solid. A stock solution may be made by dissolving the 2H,2H,3H,3H-perfluorodecanoic acid in the solvent of choice, which should be purged with an inert gas. 2H,2H,3H,3H-Perfluorodecanoic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 2H,2H,3H,3H-perfluorodecanoic acid in these solvents is approximately 10 mg/ml.

Description

2H,2H,3H,3H-Perfluorodecanoic acid is a polyfluoroalkyl substance (PFAS).^{1,2} It has been found as a contaminant in freshwater reserves and marine seafood.

References

1. Riaz, R., Junaid, M., Rehman, M.Y.A., *et al.* Spatial distribution, compositional profile, sources, ecological and human health risks of legacy and emerging *per*- and polyfluoroalkyl substances (PFASs) in freshwater reservoirs of Punjab, Pakistan. *Sci. Total Environ.* **856(Pt 2)**, 159144 (2023).
2. Ekperusi, A.O., Bely, N., Pollono, C., *et al.* Prevalence of *per*- and polyfluoroalkyl substances (PFASs) in marine seafood from the Gulf of Guinea. *Chemosphere* **335**, 139110 (2023).

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, 08/17/2023

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