

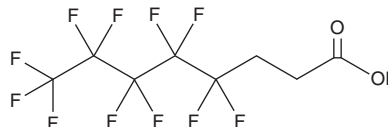
PRODUCT INFORMATION



2H,2H,3H,3H-Perfluorooctanoic Acid

Item No. 37270

CAS Registry No.: 914637-49-3
Formal Name: 4,4,5,5,6,6,7,7,8,8,8-undecafluoro-octanoic acid
Synonyms: 5:3 Fluorotelomer Carboxylic Acid, 5:3 FTCA
MF: C₈H₅F₁₁O₂
FW: 342.1
Purity: ≥90%
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

2H,2H,3H,3H-Perfluorooctanoic acid is supplied as a solid. A stock solution may be made by dissolving the 2H,2H,3H,3H-perfluorooctanoic acid in the solvent of choice, which should be purged with an inert gas. 2H,2H,3H,3H-Perfluorooctanoic acid is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 2H,2H,3H,3H-perfluorooctanoic acid in these solvents is approximately 2 and 5 mg/ml, respectively.

Description

2H,2H,3H,3H-Perfluorooctanoic acid is a polyfluoroalkyl substance (PFAS).¹ It has been found as a contaminant in wastewater.

Reference

1. Kim, J., Xin, X., Mamo, B.T., *et al.* Occurrence and fate of ultrashort-chain and other per- and polyfluoroalkyl substances (PFAS) in wastewater treatment plants. *ACS EST Water* **2**(8), 1380–1390 (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, 10/28/2022

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