

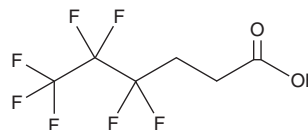
PRODUCT INFORMATION



2H,2H,3H,3H-Perfluorohexanoic Acid

Item No. 37230

CAS Registry No.: 356-02-5
Formal Name: 4,4,5,5,6,6,6-heptafluoro-hexanoic acid
Synonyms: 3:3 Fluorotelomer Carboxylic Acid, 3:3 FTCA, 3-Perfluoropropyl Propanoic Acid, FPrPA
MF: C₆H₅F₇O₂
FW: 242.1
Purity: ≥95%
Supplied as: A solution in ethanol
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-Perfluorohexanoic acid is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. 2H,2H,3H,3H-Perfluorohexanoic acid is soluble in chloroform and methanol.

Description

2H,2H,3H,3H-Perfluorohexanoic acid is a polyfluoroalkyl substance (PFAS).¹ It has been found as a contaminant in sea and river water.

Reference

1. Zhao, Z., Cheng, X., Hua, X., *et al.* Emerging and legacy per- and polyfluoroalkyl substances in water, sediment, and air of the Bohai Sea and its surrounding rivers. *Environ. Pollut.* **263(Pt. A)**, 114391 (2020).

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, 03/25/2026

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