

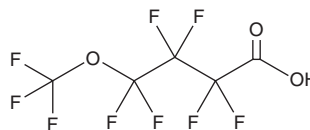
PRODUCT INFORMATION



Perfluoro-4-methoxybutanoic Acid

Item No. 37269

CAS Registry No.: 863090-89-5
Formal Name: 2,2,3,3,4,4-hexafluoro-4-(trifluoromethoxy)-butanoic acid
Synonyms: PF5OHxA, PFMOMA
MF: C₅HF₉O₃
FW: 280.0
Purity: ≥90%
Supplied as: A liquid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Perfluoro-4-methoxybutanoic acid (PF5OHxA) is supplied as a liquid. A stock solution may be made by dissolving the PF5OHxA in the solvent of choice, which should be purged with an inert gas. PF5OHxA is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of PF5OHxA in these solvents is approximately 10 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of PF5OHxA can be prepared by directly dissolving the liquid in aqueous buffers. The solubility of PF5OHxA in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

PF5OHxA is a perfluoroalkyl substance (PFAS).¹ PF5OHxA 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, 01/25/2024

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