

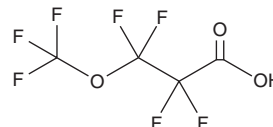
PRODUCT INFORMATION



Perfluoro-3-methoxypropanoic Acid

Item No. 37268

CAS Registry No.: 377-73-1
Formal Name: 2,2,3,3-tetrafluoro-3-(trifluoromethoxy)-propanoic acid
Synonyms: PFMPA, PF4OPeA
MF: $C_4HF_7O_3$
FW: 230.0
Purity: $\geq 90\%$
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-3-methoxypropanoic acid (PF4OPeA) is supplied as a liquid. A stock solution may be made by dissolving the PF4OPeA in the solvent of choice, which should be purged with an inert gas. PF4OPeA is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of PF4OPeA in these solvents is approximately 10 mg/ml.

Description

PF4OPeA is a perfluoroalkyl substance (PFAS).¹ PF4OPeA 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