

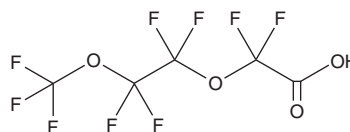
PRODUCT INFORMATION



Nonafluoro-3,6-dioxaheptanoic Acid

Item No. 37917

CAS Registry No.: 151772-58-6
Formal Name: 2,2-difluoro-2-[1,1,2,2-tetrafluoro-2-(trifluoromethoxy)ethoxy]-acetic acid
Synonyms: NFDHA, Perfluoro-3,6-dioxaheptanoic Acid, PFO2HpA
MF: $C_5HF_9O_4$
FW: 296.0
Purity: $\geq 95\%$
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

Nonafluoro-3,6-dioxaheptanoic acid (NFDHA) is supplied as a liquid. A stock solution may be made by dissolving the NFDHA in the solvent of choice, which should be purged with an inert gas. NFDHA is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of NFDHA in ethanol is approximately 5 mg/ml and approximately 10 mg/ml in DMSO and DMF.

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 NFDHA can be prepared by directly dissolving the liquid in aqueous buffers. The solubility of nonafluoro-3,6-dioxaheptanoic acid in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

NFDHA is a perfluoroalkyl ether carboxylic acid (PFECA).¹ It has been found as an environmental contaminant in watersheds.

Reference

1. Liu, Z., Zhou, J., Xu, Y., *et al.* Distributions and sources of traditional and emerging per- and polyfluoroalkyl substances among multiple environmental media in the Qiantang River watershed, China. *RSC Adv.* **12**(33), 21247-2125 (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, 12/16/2024

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