

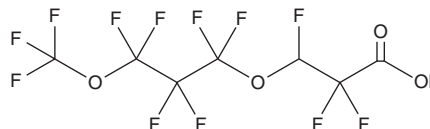
# PRODUCT INFORMATION



## 4,8-Dioxa-3H-perfluorononanoic Acid

Item No. 37267

**CAS Registry No.:** 919005-14-4  
**Formal Name:** 2,2,3-trifluoro-3-[1,1,2,3,3-hexafluoro-3-(trifluoromethoxy)propoxy]propanoic acid  
**MF:** C<sub>7</sub>H<sub>2</sub>F<sub>12</sub>O<sub>4</sub>  
**FW:** 378.1  
**Purity:** ≥90%  
**Supplied as:** A neat oil  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

4,8-Dioxa-3H-perfluorononanoic acid is supplied as a neat oil. A stock solution may be made by dissolving the 4,8-dioxa-3H-perfluorononanoic acid in the solvent of choice, which should be purged with an inert gas. 4,8-Dioxa-3H-perfluorononanoic acid is slightly soluble in chloroform, DMSO, and methanol.

### Description

4,8-Dioxa-3H-perfluorononanoic acid is a perfluoroalkyl substance (PFAS).<sup>1,2</sup> It has been found in bottled water and watersheds.

### References

1. Chow, S.J., Ojeda, N., Jacangelo, J.G., *et al.* Detection of ultrashort-chain and other per- and polyfluoroalkyl substances (PFAS) in U.S. bottled water. *Water Res.* **201**, 117292 (2021).
2. Pétré, M.-A., Salk, K.R., Stapleton, H.M., *et al.* Per- and polyfluoroalkyl substances (PFAS) in river discharge: Modeling loads upstream and downstream of a PFAS manufacturing plant in the Cape Fear watershed, North Carolina. *Sci. Total Environ.* **831**, 154763 (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, 06/26/2023

#### 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