

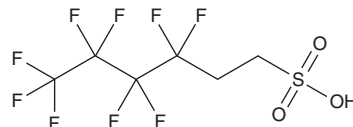
# PRODUCT INFORMATION



## 1H,1H,2H,2H-Perfluorohexanesulfonic Acid

Item No. 37261

**CAS Registry No.:** 757124-72-4  
**Formal Name:** 3,3,4,4,5,5,6,6,6-nonafluoro-1-hexanesulfonic acid  
**Synonyms:** 4:2 Fluorotelomer Sulfonic Acid, 4:2 FtS  
**MF:** C<sub>6</sub>H<sub>5</sub>F<sub>9</sub>O<sub>3</sub>S  
**FW:** 328.2  
**Purity:** ≥85%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

1H,1H,2H,2H-Perfluorohexanesulfonic acid is supplied as a solid. A stock solution may be made by dissolving the 1H,1H,2H,2H-perfluorohexanesulfonic acid in the solvent of choice, which should be purged with an inert gas. 1H,1H,2H,2H-Perfluorohexanesulfonic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 1H,1H,2H,2H-perfluorohexanesulfonic acid 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 1H,1H,2H,2H-perfluorohexanesulfonic acid can be prepared by directly dissolving the solid in aqueous buffers. The solubility of 1H,1H,2H,2H-perfluorohexanesulfonic 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

1H,1H,2H,2H-Perfluorohexanesulfonic acid is a polyfluoroalkyl substance (PFAS).<sup>1</sup> It has been found as a contaminant in drinking water.

### Reference

1. Hall, S.M., Zhang, S., Tait, G.H., *et al.* PFAS levels in paired drinking water and serum samples collected from an exposed community in Central North Carolina. *Sci. Total Environ.* **895**, 165091 (2023).

#### 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, 08/16/2023

#### 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