

# PRODUCT INFORMATION



## 1H,1H,2H,2H-Perfluorodecanesulfonic Acid

Item No. 37259

CAS Registry No.: 39108-34-4

Formal Name: 3,3,4,4,5,5,6,6,7,7,8,8,9,9,10,10,10-heptafluoro-1-decanesulfonic acid

Synonyms: 8:2 Fluorotelomer Sulfonic Acid, 8:2 FtS

MF:  $C_{10}H_5F_{17}O_3S$

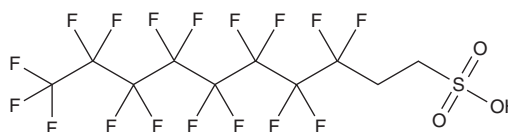
FW: 528.2

Purity:  $\geq 80\%$

Supplied as: A low-melting solid

Storage:  $-20^{\circ}C$

Stability:  $\geq 4$  years



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

### Laboratory Procedures

1H,1H,2H,2H-Perfluorodecanesulfonic acid is supplied as a low-melting solid. A stock solution may be made by dissolving the 1H,1H,2H,2H-perfluorodecanesulfonic acid in the solvent of choice, which should be purged with an inert gas. 1H,1H,2H,2H-Perfluorodecanesulfonic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 1H,1H,2H,2H-perfluorodecanesulfonic acid in these solvents is approximately 10 mg/ml.

### Description

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

### Reference

1. Zhao, Z., Cheng, X., Hua, X., *et al.* Emerging and legacy per- and polyfluoroalkyl substances in water, sediment, and air of the Bohai Sea and its surrounding rivers. *Environ. Pollut.* **263**(Pt. A), 114391 (2020).

#### 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, 03/11/2025

#### 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