

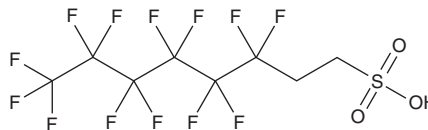
PRODUCT INFORMATION



1H,1H,2H,2H-Perfluorooctanesulfonic Acid

Item No. 37260

CAS Registry No.: 27619-97-2
Formal Name: 3,3,4,4,5,5,6,6,7,7,8,8,8-tridecafluoro-1-octanesulfonic acid
Synonyms: 6:2 Fluorotelomer Sulfonate, 6:2 Fluorotelomer Sulfonic Acid, 6:2 FTS, 6:2 FTSA
MF: C₈H₅F₁₃O₃S
FW: 428.2
Purity: ≥90%
Supplied as: A low-melting 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-Perfluorooctanesulfonic acid is supplied as a low-melting solid. A stock solution may be made by dissolving the 1H,1H,2H,2H-perfluorooctanesulfonic acid in the solvent of choice, which should be purged with an inert gas. 1H,1H,2H,2H-Perfluorooctanesulfonic acid is slightly soluble in acetone, chloroform, DMSO and methanol.

Description

1H,1H,2H,2H-Perfluorooctanesulfonic acid is a polyfluoroalkyl substance (PFAS).¹ 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, 12/09/2022

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