

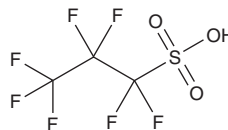
PRODUCT INFORMATION



Perfluoropropanesulfonic Acid

Item No. 38156

CAS Registry No.: 423-41-6
Formal Name: 1,1,2,2,3,3,3-heptafluoro-1-propanesulfonic acid
Synonyms: Heptafluoropropanesulfonic Acid, PFPrS
MF: C₃HF₇O₃S
FW: 250.1
Purity: ≥90%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Perfluoropropanesulfonic acid is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Perfluoropropanesulfonic acid is soluble in methanol.

Description

Perfluoropropanesulfonic acid is a perfluoroalkyl substance (PFAS).¹ It has been found as a contaminant in groundwater in areas treated with aqueous film-forming foam fire extinguishing agents.

Reference

1. Barzen-Hanson, K.A. and Field, J.A. Discovery and implications of C₂ and C₃ perfluoroalkyl sulfonates in aqueous film-forming foams and groundwater. *Environ. Sci. Technol.* **2(4)**, 95-99 (2015).

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, 02/05/2024

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