

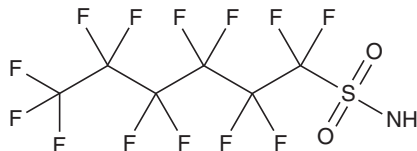
PRODUCT INFORMATION



Perfluorohexanesulfonamide

Item No. 38306

CAS Registry No.: 41997-13-1
Formal Name: 1,1,2,2,3,3,4,4,5,5,6,6,6-tridecafluoro-1-hexanesulfonamide
Synonym: FHxSA
MF: C₆H₂F₁₃NO₂S
FW: 399.1
Purity: ≥75%
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

Perfluorohexanesulfonamide is supplied as a solid. A stock solution may be made by dissolving the perfluorohexanesulfonamide in the solvent of choice, which should be purged with an inert gas. Perfluorohexanesulfonamide is soluble in methanol.

Description

Perfluorohexanesulfonamide is a perfluoroalkyl substance (PFAS).^{1,2} It has been found in surface freshwater foams and in soil from fire-fighting equipment testing sites.

References

1. Schwichtenberg, T., Bogdan, D., Carignan, C.C., *et al.* PFAS and dissolved organic carbon enrichment in surface water foams on a northern U.S. freshwater lake. *Environ. Sci. Technol.* **54**(22), 14455-14464 (2020).
2. Munoz, G., Desrosiers, M., Vetter, L., *et al.* Bioaccumulation of zwitterionic polyfluoroalkyl substances in earthworms exposed to aqueous film-forming foam impacted soils. *Environ. Sci. Technol.* **54**(3), 1687-1697 (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, 11/21/2023

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