

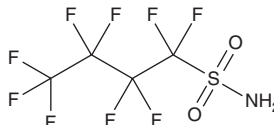
PRODUCT INFORMATION



Perfluorobutylsulfonamide

Item No. 37365

CAS Registry No.: 30334-69-1
Formal Name: 1,1,2,2,3,3,4,4,4-nonafluoro-1-butanesulfonamide
Synonyms: FBSA, Perfluorobutane Sulfonamide
MF: $C_4H_2F_9NO_2S$
FW: 299.1
Purity: $\geq 95\%$
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



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

Laboratory Procedures

Perfluorobutylsulfonamide is supplied as a solid. A stock solution may be made by dissolving the perfluorobutylsulfonamide in the solvent of choice, which should be purged with an inert gas. Perfluorobutylsulfonamide is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of perfluorobutylsulfonamide in these solvents is approximately 20 and 12 mg/ml, respectively. Perfluorobutylsulfonamide is slightly soluble in DMSO.

Description

Perfluorobutylsulfonamide is a perfluoroalkyl substance (PFAS).^{1,2} It has been found as a contaminant in wastewater and consumable Tampa Bay oysters.

References

1. Kim, J., Xin, X., Mamo, B.T., *et al.* Occurrence and fate of ultrashort-chain and other per- and polyfluoroalkyl substances (PFAS) in wastewater treatment plants. *ACS EST Water* **2**(8), 1380–1390 (2022).
2. Lemos, L., Gantiva, L., Kaylor, C., *et al.* American oysters as bioindicators of emerging organic contaminants in Florida, United States. *Sci. Total Environ.* **835**, 155316 (2022).

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, 08/15/2023

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