

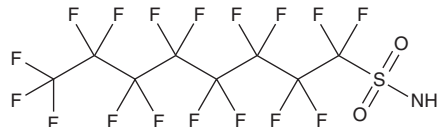
PRODUCT INFORMATION



Perfluorooctanesulfonamide

Item No. 37258

CAS Registry No.: 754-91-6
Formal Name: 1,1,2,2,3,3,4,4,5,5,6,6,7,7,8,8,8-heptadecafluoro-1-octanesulfonamide
Synonym: PFOSA
MF: C₈H₂F₁₇NO₂S
FW: 499.1
Purity: ≥85%
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

Perfluorooctanesulfonamide (PFOSA) is supplied as a solid. A stock solution may be made by dissolving the PFOSA in the solvent of choice, which should be purged with an inert gas. PFOSA is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of PFOSA in these solvents is approximately 10 mg/ml.

Description

PFOSA is a perfluoroalkyl substance (PFAS).¹ It induces hepatic macrophage and neutrophil accumulation and fibrosis in zebrafish larvae when administered in tank water at concentrations ranging from 0.1-100 µg/L. PFOSA has been found in municipal wastewater networks.²

References

1. Xuan, R., Qiu, X., Wang, J., *et al.* Hepatotoxic response of perfluorooctane sulfonamide (PFOSA) in early life stage zebrafish (*Danio rerio*) is greater than perfluorooctane sulfonate (PFOS). *J. Hazard Mater.* **461**, 132552 (2024).
2. Gobelius, L., Glimstedt, L., Olsson, J., *et al.* Mass flow of per- and polyfluoroalkyl substances (PFAS) in a Swedish municipal wastewater network and wastewater treatment plant. *Chemosphere* **336**, 139182 (2023).

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, 11/28/2023

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