

PRODUCT INFORMATION



2-(N-Ethylperfluorooctanesulfonamido)acetic Acid

Item No. 37363

CAS Registry No.: 2991-50-6

Formal Name: N-ethyl-N-[(1,1,2,2,3,3,4,4,5,5,6,6,7,7,8,8,8-heptafluorooctyl)sulfonyl]-glycine

Synonyms: EtFOSAA,
N-Ethylperfluorooctanesulfonamidoacetate,
Et-PFOSA-AcOH

MF: $C_{12}H_8F_{17}NO_4S$

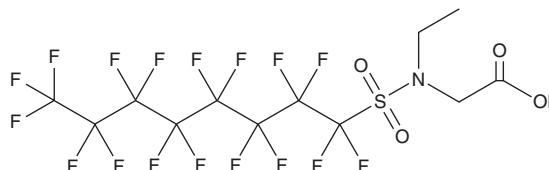
FW: 585.2

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

2-(N-Ethylperfluorooctanesulfonamido)acetic acid (Et-PFOSA-AcOH) is supplied as a solid. A stock solution may be made by dissolving the Et-PFOSA-AcOH in the solvent of choice, which should be purged with an inert gas. Et-PFOSA-AcOH is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of Et-PFOSA-AcOH in these solvents is approximately 2 mg/ml.

Description

Et-PFOSA-AcOH is a perfluoroalkyl substance (PFAS).¹ It has been found as a contaminant in wastewater samples. Et-PFOSA-AcOH is associated with decreased plasma levels of maternal and fetal free thyroxine (T_4).²

References

1. Worley, R.R., Moore, S.M., Tierney, B.C., *et al.* Per- and polyfluoroalkyl substances in human serum and urine samples from a residentially exposed community. *Environ. Int.* **106**, 135-143 (2017).
2. Rickard, B.P., Rizvi, I., and Fenton, S.E. Per- and poly-fluoroalkyl substances (PFAS) and female reproductive outcomes: PFAS elimination, endocrine-mediated effects, and disease. *Toxicology* **465:153031**, (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, 02/14/2023

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