

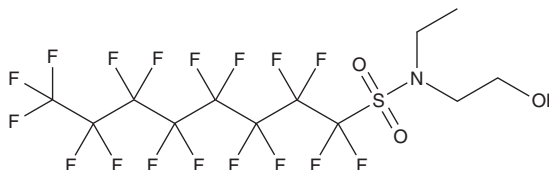
PRODUCT INFORMATION



N-ethyl Perfluorooctanesulfonamidoethanol

Item No. 37916

CAS Registry No.: 1691-99-2
Formal Name: N-ethyl-1,1,2,2,3,3,4,4,5,5,6,6,7,7,8,8,8-heptafluoro-N-(2-hydroxyethyl)-1-octanesulfonamide
Synonym: N-ethyl FOSE
MF: $C_{12}H_{10}F_{17}NO_3S$
FW: 571.2
Purity: ≥85% (mixture of isomers)
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

N-ethyl Perfluorooctanesulfonamidoethanol (N-ethyl FOSE) is supplied as a solid. A stock solution may be made by dissolving the N-ethyl FOSE in the solvent of choice, which should be purged with an inert gas. N-ethyl FOSE is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of N-ethyl FOSE in ethanol is approximately 5 mg/ml and approximately 10 mg/ml in DMSO and DMF.

Description

N-ethyl FOSE is a perfluoroalkyl substance (PFAS).^{1,2} It decreases maternal body weight and food intake in pregnant rats and induces abortions and decreases fetal weight in pregnant rabbits when administered at a dose of 10 mg/kg.³ N-ethyl FOSE has been found in landfill leachate and the atmosphere.^{1,2}

References

1. Capozzi, S.L., Leang, A.L., Rodenburg, L.A., *et al.* PFAS in municipal landfill leachate: Occurrence, transformation, and sources. *Chemosphere* **334**, 138924 (2023).
2. Shoeib, M., Harner, T., and Vlahos, P. Perfluorinated chemicals in the arctic atmosphere. *Environ. Sci. Technol.* **40**(24), 7577-7583 (2006).
3. Case, M.T., York, R.G., and Christian, M.S. Rat and rabbit oral developmental toxicology studies with two perfluorinated compounds. *Int. J. Toxicol.* **20**(2), 101-109 (2001).

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, 02/27/2026

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