

PRODUCT INFORMATION



N-methyl Perfluorooctanesulfonamidoethanol

Item No. 37265

CAS Registry No.: 24448-09-7

Formal Name: 1,1,2,2,3,3,4,4,5,5,6,6,7,7,8,8,8-heptafluoro-N-(2-hydroxyethyl)-N-methyl-1-octanesulfonamide

Synonyms: N-methyl FOSE, N-methyl PFOSE

MF: $C_{11}H_8F_{17}NO_3S$

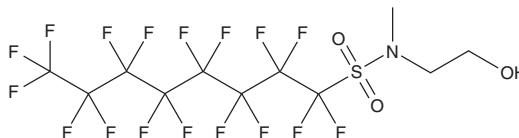
FW: 557.2

Purity: $\geq 90\%$

Supplied as: A low melting 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

N-methyl perfluorooctanesulfonamidoethanol is supplied as a low melting solid. A stock solution may be made by dissolving the N-methyl perfluorooctanesulfonamidoethanol in the solvent of choice, which should be purged with an inert gas. N-methyl perfluorooctanesulfonamidoethanol is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of N-methyl perfluorooctanesulfonamidoethanol in these solvents is approximately 10 mg/ml.

Description

N-methyl Perfluorooctanesulfonamidoethanol is a perfluoroalkyl substance (PFAS).^{1,2} It has been found in landfill leachate and the atmosphere.

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).

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, 11/20/2023

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