

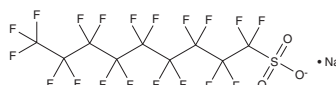
PRODUCT INFORMATION



Perfluorononanesulfonic Acid (sodium salt)

Item No. 38908

CAS Registry No.: 98789-57-2
Formal Name: 1,1,2,2,3,3,4,4,5,5,6,6,7,7,8,8,9,9,9-nonadecafluoro-1-nonanesulfonic acid, monosodium salt
Synonym: PFNS
MF: $C_9F_{19}O_3S \cdot Na$
FW: 572.1
Purity: $\geq 98\%$
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

Perfluorononanesulfonic acid (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the perfluorononanesulfonic acid (sodium salt) in the solvent of choice, which should be purged with an inert gas. Perfluorononanesulfonic acid (sodium salt) is soluble in DMSO and methanol.

Description

Perfluorononanesulfonic acid is a perfluoroalkyl substance (PFAS).^{1,2} It has been found in landfill leachate and marine seafood.

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. Ekperusi, A.O., Bely, N., Pollono, C., *et al.* Prevalence of per- and polyfluoroalkyl substances (PFASs) in marine seafood from the Gulf of Guinea. *Chemosphere* **335**, 139110 (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, 08/28/2023

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