

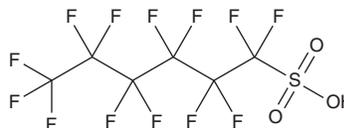
# PRODUCT INFORMATION



## Perfluorohexanesulfonic Acid

Item No. 37248

**CAS Registry No.:** 355-46-4  
**Formal Name:** 1,1,2,2,3,3,4,4,5,5,6,6,6-tridecafluoro-1-hexanesulfonic acid  
**Synonym:** PFHxS  
**MF:** C<sub>6</sub>H<sub>13</sub>O<sub>3</sub>S  
**FW:** 400.1  
**Purity:** ≥95%  
**Supplied as:** A low melting 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

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

### Description

Perfluorohexanesulfonic acid is a perfluoroalkyl substance (PFAS).<sup>1</sup> It activates peroxisome proliferator-activated receptor  $\alpha$  (PPAR $\alpha$ ) in a reporter assay using COS-1 cells expressing the human or mouse receptor when used at concentrations ranging from 5 to 100  $\mu$ M. It has been found in landfill leachate and sea and river water.<sup>2,3</sup>

### References

1. Wolf, C.J., Takacs, M.L., Schmid, J.E., *et al.* Activation of mouse and human peroxisome proliferator-activated receptor alpha by perfluoroalkyl acids of different functional groups and chain lengths. *Toxicol. Sci.* **106**(1), 162-171 (2008).
2. Capozzi, S.L., Leang, A.L., Rodenburg, L.A., *et al.* PFAS in municipal landfill leachate: Occurrence, transformation, and sources. *Chemosphere* **334**, 138924 (2023).
3. Zhao, Z., Cheng, X., Hua, X., *et al.* Emerging and legacy per- and polyfluoroalkyl substances in water, sediment, and air of the Bohai Sea and its surrounding rivers. *Environ. Pollut.* **263**(Pt. A), 114391 (2020).

#### 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, 11/17/2025

#### 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