

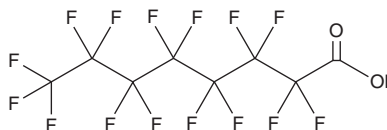
PRODUCT INFORMATION



Pentadecafluorooctanoic Acid

Item No. 37232

CAS Registry No.: 335-67-1
Formal Name: 2,2,3,3,4,4,5,5,6,6,7,7,8,8,8-pentadecafluorooctanoic acid
Synonyms: A-5717, NSC 95114, Perfluorooctanoic Acid, PFOA
MF: C₈HF₁₅O₂
FW: 414.1
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

Pentadecafluorooctanoic acid is supplied as a solid. A stock solution may be made by dissolving the pentadecafluorooctanoic acid in the solvent of choice, which should be purged with an inert gas. Pentadecafluorooctanoic acid is slightly soluble in chloroform and methanol.

Description

Pentadecafluorooctanoic acid is a perfluoroalkyl substance (PFAS).¹ Chronic administration of pentadecafluorooctanoic acid (5 and 20 mg/kg per day) leads to accumulation in the kidney, liver, lung, heart, testes, and brain in rats.² Pentadecafluorooctanoic acid (20 mg/kg per day) is toxic to rats, inducing weight loss, sensitivity to external stimuli, lethargy, and hepatic inflammation. It has been found in marine life, including lake trout, salmon, and whitefish, and as an environmental contaminant.¹ Formulations containing pentadecafluorooctanoic acid have been used as emulsifiers in fluoropolymer polymerization.

References

1. Sinclair, E., Taniyasu, S., Yamashita, N., *et al.* Perfluorooctanoic acid and perfluorooctane sulfonate in Michigan and New York waters. *Organohalo. Comp.* **66**, 4019-4023 (2004).
2. Cui, L., Zhou, Q.-F., Liao, C.-Y., *et al.* Studies on the toxicological effects of PFOA and PFOS on rats using histological observation and chemical analysis. *Arch. Environ. Contam. Toxicol.* **56(2)**, 338-349 (2009).

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, 11/01/2022

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