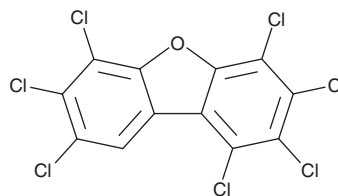


PRODUCT INFORMATION

1,2,3,4,6,7,8-Heptachlorodibenzofuran

Item No. 38617

CAS Registry No.: 67562-39-4
Synonyms: 1,2,3,4,6,7,8-HpCDF, PCDF 131
MF: C₁₂HCl₇O
FW: 409.3
Purity: ≥95%
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

1,2,3,4,6,7,8-Heptachlorodibenzofuran is supplied as a solid. A stock solution may be made by dissolving the 1,2,3,4,6,7,8-heptachlorodibenzofuran in the solvent of choice, which should be purged with an inert gas. 1,2,3,4,6,7,8-Heptachlorodibenzofuran is slightly soluble in chloroform (warmed) and methanol.

Description

1,2,3,4,6,7,8-Heptachlorodibenzofuran is a dioxin-like polychlorinated dibenzofuran (PCDF).¹ It has been found in the soil near municipal waste incinerators, in freshwater fish, and in human breastmilk.¹⁻³

References

1. Harnly, M., Stephens, R., McLaughlin, C., *et al.* Polychlorinated dibenzo-*p*-dioxin and dibenzofuran contamination at metal recovery facilities, open burn sites, and a railroad car incineration facility. *Environ. Sci. Technol.* **29**(3), 667-684 (1995).
2. Zacharewski, T., Safe, L., Safe, S., *et al.* Comparative analysis of polychlorinated dibenzo-*p*-dioxin and dibenzofuran congeners in Great Lakes fish extracts by gas chromatography-mass spectrometry and in-vitro enzyme induction activities. *Environ. Sci. Technol.* **23**(6), 730-735 (1989).
3. Takekuma, M., Saito, K., Ogawa, M., *et al.* Levels of PCDDs, PCDFs and Co-PCBs in human milk in Saitama, Japan, and epidemiological research. *Chemosphere* **54**(1), 127-135 (2004).

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, 04/28/2023

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