

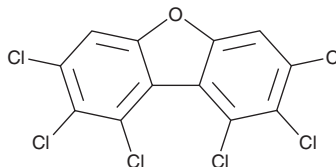
PRODUCT INFORMATION



1,2,3,7,8,9-Hexachlorodibenzofuran

Item No. 38622

CAS Registry No.: 72918-21-9
Synonyms: 1,2,3,7,8,9-HxCDF, PCDF 124
MF: $C_{12}H_2Cl_6O$
FW: 374.9
Purity: $\geq 95\%$
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

1,2,3,7,8,9-Hexachlorodibenzofuran is supplied as a solid. A stock solution may be made by dissolving the 1,2,3,7,8,9-hexachlorodibenzofuran in the solvent of choice, which should be purged with an inert gas. 1,2,3,7,8,9-Hexachlorodibenzofuran is slightly soluble in chloroform and methanol.

Description

1,2,3,7,8,9-Hexachlorodibenzofuran is a dioxin-like polychlorinated dibenzofuran (PCDF).¹ It has been found in the soil near municipal waste incinerators, water leachates from a landfill, and in blood from individuals that lived near or in herbicide-sprayed regions.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. Först, C., Stieglitz, L., and Zwick, G. Isomer-specific determination of PCDD/PCDF in oil extracts from water leachates of a waste landfill. *Chemosphere* **17**(10), 1935-1944 (1988).
3. Vu, T.S., Nishijo, M., Nishino, Y., *et al.* Polychlorinated dibenzo-p-dioxins and polychlorinated dibenzofurans in blood samples from historically herbicide-sprayed areas and rural and industrialized unsprayed areas in Vietnam. *Chemosphere* **326**, 138331 (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, 05/02/2023

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