

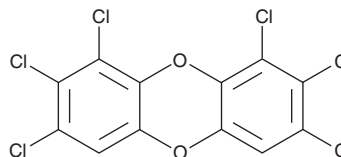
PRODUCT INFORMATION



1,2,3,7,8,9-Hexachlorodibenzo-*p*-dioxin

Item No. 38596

CAS Registry No.: 19408-74-3
Formal Name: 1,2,3,7,8,9-hexachloro-dibenzo[b,e][1,4]dioxin
Synonyms: 1,2,3,7,8,9-HCDD,
1,2,3,7,8,9-HxCDD
MF: C₁₂H₂Cl₆O₂
FW: 390.9
Purity: ≥90%
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

1,2,3,7,8,9-Hexachlorodibenzo-*p*-dioxin is supplied as a solid. A stock solution may be made by dissolving the 1,2,3,7,8,9-hexachlorodibenzo-*p*-dioxin in the solvent of choice, which should be purged with an inert gas. 1,2,3,7,8,9-Hexachlorodibenzo-*p*-dioxin is slightly soluble in chloroform and ethyl acetate (heated).

Description

1,2,3,7,8,9-Hexachlorodibenzo-*p*-dioxin is a polychlorinated dibenzodioxin (PCDD).¹ It induces the formation of liver adenomas or carcinomas in rats when administered at a dose of 2.5 µg/kg once per week in combination with 1,2,3,6,7,8-hexachlorodibenzo-*p*-dioxin. 1,2,3,7,8,9-Hexachlorodibenzo-*p*-dioxin has been found in the air near municipal waste incinerators and in blood from individuals that lived near or in herbicide-sprayed regions.^{2,3}

References

1. Huff, J.E., Salmon, A.G., Hooper, N.K., *et al.* Long-term carcinogenesis studies on 2,3,7,8-tetrachlorodibenzo-*p*-dioxin and hexachlorodibenzo-*p*-dioxins. *Cell Biol. Toxicol.* **7**(1), 67-94 (1991).
2. Hu, S.-W., ChangChien, G.-P., and Chan, C.-C. PCDD/Fs levels in indoor environments and blood of workers of three municipal waste incinerators in Taiwan. *Chemosphere* **55**(4), 611-620 (2004).
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, 05/01/2023

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