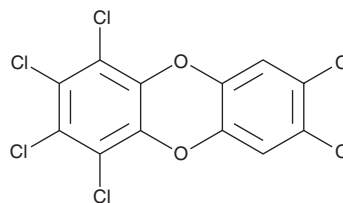


PRODUCT INFORMATION

1,2,3,4,7,8-Hexachlorodibenzo-*p*-dioxin

Item No. 38594

CAS Registry No.: 39227-28-6
Formal Name: 1,2,3,4,7,8-hexachloro-dibenzo[b,e][1,4]dioxin
Synonyms: 1,2,3,4,7,8-HxCDD, PCDD 66
MF: C₁₂H₂Cl₆O₂
FW: 390.9
Purity: ≥98%
Supplied as: A solution in n-nonane:toluene (70:30)
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

1,2,3,4,7,8-Hexachlorodibenzo-*p*-dioxin is a polychlorinated dibenzodioxin (PCDD). It induces the activity of aryl hydrocarbon hydroxylase (AHH) when used at a concentration of 47 pmol/egg in chicken embryo livers.¹ 1,2,3,4,7,8-Hexachlorodibenzo-*p*-dioxin induces toxicity in rats (LD₅₀ = 1.871 mg/kg).² It has been found in freshwater fish and in the air near municipal waste incinerators.^{3,4}

References

- Poland, A. and Glover, E. Studies on the mechanism of toxicity of the chlorinated dibenzo-*p*-dioxins. *Environ. Health Perspect.* **5**, 245-251 (1973).
- Pohjanvirta, R., Unkila, M., and Tuomisto, J. Comparative acute lethality of 2,3,7,8-tetrachlorodibenzo-*p*-dioxin (TCDD), 1,2,3,7,8-pentachlorodibenzo-*p*-dioxin and 1,2,3,4,7,8-hexachlorodibenzo-*p*-dioxin in the most TCDD-susceptible and the most TCDD-resistant rat strain. *Pharmacol. Toxicol.* **73**(1), 52-56 (1993).
- 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).
- 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).

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, 04/23/2026

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