

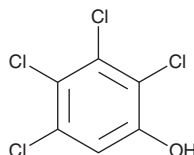
PRODUCT INFORMATION



2,3,4,5-Tetrachlorophenol

Item No. 29743

CAS Registry No.: 4901-51-3
MF: $C_6H_2Cl_4O$
FW: 231.9
Purity: $\geq 95\%$
Supplied as: A crystalline 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

2,3,4,5-Tetrachlorophenol is supplied as a crystalline solid. A stock solution may be made by dissolving the 2,3,4,5-tetrachlorophenol in the solvent of choice, which should be purged with an inert gas. 2,3,4,5-Tetrachlorophenol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 2,3,4,5-tetrachlorophenol in ethanol and DMF is approximately 30 mg/ml and approximately 15 mg/ml in DMSO.

2,3,4,5-Tetrachlorophenol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 2,3,4,5-tetrachlorophenol should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 2,3,4,5-Tetrachlorophenol has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

2,3,4,5-Tetrachlorophenol is a metabolite of the insecticide γ -lindane (Item No. 23952).¹ It is toxic to fathead minnows and rainbow trout ($LC_{50}s = 0.496$ and 0.304 mg/L, respectively, in tank water).²

References

1. Chadwick, R.W. and Freal, J.J. The identification of five unreported lindane metabolites recovered from rat urine. *Bull. Environ. Contam. Toxicol.* **7(2)**, 137-146 (1972).
2. Holcombe, G.W., Phipps, G.L., Knuth, L., et al. The acute toxicity of selected substituted phenols, benzenes and benzoic acid esters to fathead minnows *Pimephales promelas*. *Environ. Pollut.* **35(4)**, 367-381 (1984).

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, 11/18/2022

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