

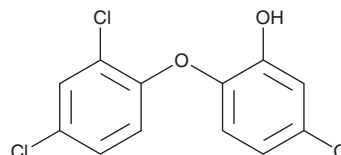
# PRODUCT INFORMATION



## Triclosan

Item No. 20342

**CAS Registry No.:** 3380-34-5  
**Formal Name:** 5-chloro-2-(2,4-dichlorophenoxy)-phenol  
**MF:** C<sub>12</sub>H<sub>7</sub>Cl<sub>3</sub>O<sub>2</sub>  
**FW:** 289.5  
**Purity:** ≥98%  
**Supplied as:** A crystalline solid  
**Storage:** Room temperature  
**Stability:** ≥4 years



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

### Laboratory Procedures

Triclosan is supplied as a crystalline solid. A stock solution may be made by dissolving the triclosan in the solvent of choice, which should be purged with an inert gas. Triclosan is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of triclosan in these solvents is approximately 30 mg/ml.

Triclosan is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, triclosan should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Triclosan 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

Triclosan is an antibiotic and antifungal.<sup>1,2</sup> It is an inhibitor of *E. coli* enoyl-acyl carrier protein reductase (FabI) and FabI containing a glycine-to-valine substitution at position 93 (FabI<sup>G93V</sup>; IC<sub>50</sub>s = 2 and 10 μM, respectively).<sup>1</sup> Triclosan is active against the bacteria *M. epidermidis*, *M. lysodeikticus*, *S. aureus*, *S. lactis*, and *S. haemolyticus* (MICs = 0.03, 0.01, 3, 3, and 1 μg/ml, respectively) and the fungi *A. fumigatus*, *C. albicans*, and *C. tropicalis* (MICs = 10 μg/ml for all).<sup>2</sup> It also decreases reattachment to dentin chips by, reduces the viability of, and increases lactate dehydrogenase (LDH) secretion from, human SG gingival epithelial cells when used at concentrations of 30, 100, and 500 μM, respectively.<sup>3</sup>

### References

1. Heath, R.J., Rubin, J.R., Holland, D.R., *et al.* Mechanism of triclosan inhibition of bacterial fatty acid synthesis. *J. Biol. Chem.* **274**(16), 11110-11114 (1999).
2. Regös, J., Zak, O., Solf, R., *et al.* Antimicrobial spectrum of triclosan, a broad-spectrum antimicrobial agent for topical application. II. Comparison with some other antimicrobial agents. *Dermatologica* **158**(1), 72-79 (1979).
3. Zuckerbraun, H.L., Babich, H., May, R., *et al.* Triclosan: Cytotoxicity, mode of action, and induction of apoptosis in human gingival cells *in vitro*. *Eur. J. Oral Sci.* **106**(2 Pt 1), 628-636 (1998).

#### 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, 07/28/2023

#### 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