

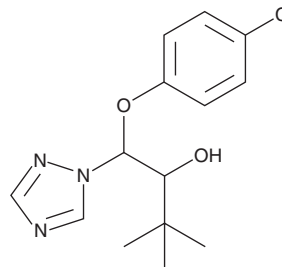
PRODUCT INFORMATION



Triadimenol

Item No. 24216

CAS Registry No.: 55219-65-3
Formal Name: β -(4-chlorophenoxy)- α -(1,1-dimethylethyl)-1H-1,2,4-triazole-1-ethanol
MF: $C_{14}H_{18}ClN_3O_2$
FW: 295.8
Purity: $\geq 95\%$ (mixture of isomers)
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

Triadimenol is supplied as a solid. A stock solution may be made by dissolving the triadimenol in the solvent of choice, which should be purged with an inert gas. Triadimenol is slightly soluble in methanol and DMSO.

Description

Triadimenol is a metabolite of triadimefon (Item No. 18714), a broad-spectrum chiral triazole fungicide, that is formed by reduction of a carbonyl group to the corresponding alcohol.¹ It is teratogenic, inducing cranial nerve and ganglia abnormalities in a rat post-implantation whole embryo culture model when used at concentrations ranging from 12.5 to 125 μM .² *In vivo*, triadimenol induces embryotoxicity in rats and rabbits when administered orally at doses of 100 and 40 mg/kg, respectively. Embryonic exposure to triadimenol (3-3,000 $\mu\text{g/L}$) induces embryonic mortality as well as decreases fertility and increases the number of female offspring in medaka fish (*O. latipes*).³

References

1. Shen, Q., Li, J., Xu, P., et al. Enantioselective metabolism of triadimefon and its chiral metabolite triadimenol in lizards. *Ecotoxicol. Environ. Saf.* **143**, 159-165 (2017).
2. Menegola, E., Broccia, M.L., Di Renzo, F., et al. *In vitro* teratogenic potential of two antifungal triazoles: Triadimefon and triadimenol. *In Vitro Cell. Dev. Biol. Anim.* **36**(2), 88-95 (2000).
3. Chu, S.-H., Liao, P.-H., and Chen, P.-J. Developmental exposures to an azole fungicide triadimenol at environmentally relevant concentrations cause reproductive dysfunction in females of medaka fish. *Chemosphere* **152**, 181-189 (2016).

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, 10/03/2022

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