PRODUCT INFORMATION



Triticonazole

Item No. 18726

CAS Registry No.: Formal Name:	131983-72-7 5-[(4-chlorophenyl)methylene]- 2,2-dimethyl-1-(1H-1,2,4-triazol- 1-ylmethyl)-cyclopentanol	N N
MF:	C ₁₇ H ₂₀ CIN ₃ O	N
FW:	317.8	∖ он
Purity:	≥95%	CI
UV/Vis.:	λ _{max} : 263 nm	
Supplied as:	A crystalline solid	
Storage:	-20°C	\checkmark \checkmark
Stability:	≥4 years	
1 6 13		

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

Laboratory Procedures

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

Triticonazole is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, triticonazole should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Triticonazole has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Triticonazole is a triazole fungicide used for the control of common soil and seed-borne diseases on cereals and other crops.¹ The antifungal effects of triticonazole are due to its ability to inhibit ergosterol biosynthesis.¹ Formulations containing triticonazole have been used in the control of fungi in agriculture.

Reference

1. Zarn, J.A., Brüschweiler, B.J., and Schlatter, J.R. Azole fungicides affect mammalian steroidogenesis by inhibiting sterol 14 α -demethylase and aromatase. Environ. Health Perspect. 111(3), 255-261 (2003).

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

uyer 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, 07/25/2023

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