

PRODUCT INFORMATION



Hydroxy Itraconazole

Item No. 22576

CAS Registry No.: 112559-91-8

Formal Name: 4-[4-[4-[[2-(2,4-dichlorophenyl)-2-(1H-1,2,4-triazol-1-ylmethyl)-1,3-dioxolan-4-yl]methoxy]phenyl]-1-piperazinyl]phenyl]-2,4-dihydro-2-(2-hydroxy-1-methylpropyl)-3H-1,2,4-triazol-3-one

Synonyms: OH-ITZ, hydroxy-ITZ

MF: $C_{35}H_{38}Cl_2N_8O_5$

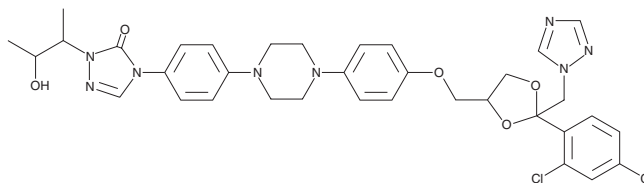
FW: 721.6

Purity: $\geq 95\%$ (mixture of diastereomers)

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.

Description

Hydroxy itraconazole is an active metabolite of the antifungal agent itraconazole (Item No. 13288). Itraconazole is metabolized by the cytochrome P450 isoform 3A4 (CYP3A4) to form hydroxy itraconazole. Hydroxy itraconazole exhibits *in vitro* antifungal activity (MIC = 0.019, 0.078, and 0.078 mg/ml against *C. albicans*, *A. fumigatus*, and *C. neoformans*, respectively).¹ Plasma concentration of hydroxy itraconazole is directly correlated to that of the parent compound and is used as an indicator of prophylaxis for fungal infections.²

References

1. Mikami, Y., Sakamoto, T., Yazawa, K., *et al.* Comparison of *in vitro* antifungal activity of itraconazole and hydroxy-itraconazole by colorimetric MTT assay. *Mycoses* **37**(1-2), 27-33 (1994).
2. Mino, Y., Naito, T., Watanabe, T. *et al.* Hydroxy-itraconazole pharmacokinetics is similar to that of itraconazole in immunocompromised patients receiving oral solution of itraconazole. *Clin. Chim. Acta.* **415**, 128-132 (2013).

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, 11/21/2022

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