

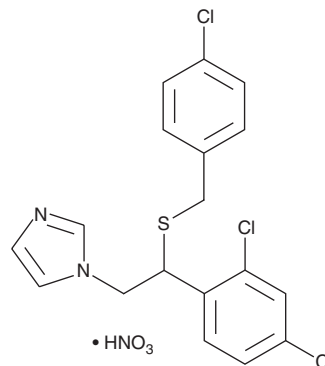
# PRODUCT INFORMATION



## Sulconazole (nitrate)

Item No. 23800

**CAS Registry No.:** 61318-91-0  
**Formal Name:** 1-[2-[[[4-chlorophenyl)methyl]thio]-2-(2,4-dichlorophenyl)ethyl]-1H-imidazole, mononitrate  
**Synonym:** RS 44872  
**MF:**  $C_{18}H_{15}Cl_3N_2S \cdot HNO_3$   
**FW:** 460.8  
**Purity:**  $\geq 95\%$   
**UV/Vis.:**  $\lambda_{max}$ : 219 nm  
**Supplied as:** A crystalline solid  
**Storage:**  $-20^\circ C$   
**Stability:**  $\geq 4$  years



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

### Laboratory Procedures

Sulconazole (nitrate) is supplied as a crystalline solid. A stock solution may be made by dissolving the sulconazole (nitrate) in the solvent of choice, which should be purged with an inert gas. Sulconazole (nitrate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of sulconazole (nitrate) in ethanol is approximately 0.1 mg/ml and approximately 25 mg/ml in DMSO and DMF.

Sulconazole (nitrate) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, sulconazole (nitrate) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Sulconazole (nitrate) has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Sulconazole is an azole antifungal that is active against dermatophyte fungal species, including isolates of various *Candida* species.<sup>1</sup> It acts as a non-specific inhibitor of styrene metabolism by cytochrome P450 (CYP450) enzymes, with  $K_i$  values ranging from 0.008 to 3.5  $\mu M$  against various CYP450 isoforms in human lymphoblast microsomes *in vitro*.<sup>2,3</sup> In rat spleen microsomes, it inhibits the activity of heme oxygenase 1 with an  $IC_{50}$  value of 1.1  $\mu M$ .<sup>4</sup> Sulconazole inhibits DNA damage induced by the metabolism of styrene by bacterial CYP450cam with a  $K_i$  value of 204.2  $\mu M$  in an electrochemical biosensor model.<sup>2</sup> Formulations containing sulconazole have been used for the treatment of fungal infections.

### References

1. Odds, F.C., Webster, C.E., and Abbot, A.B. Antifungal relative inhibition factors: BAY 1-9139, bifonazole, butoconazole, isoconazole, itraconazole (R 51211), oxiconazole, Ro 14-4767/002, sulconazole, terconazole and vibunazole (BAY n-7133) compared *in vitro* with nine established antifungal agents. *J. Antimicrob. Chemother.* **14**(2), 105-114 (1984).
2. Hull, D.O., Bajrami, B., Jansson, I., *et al.* Characterizing metabolic inhibition using electrochemical enzyme/DNA biosensors. *Anal. Chem.* **81**(2), 716-724 (2009).
3. Zhang, W., Ramamoorthy, Y., Kilicarslan, T., *et al.* Inhibition of cytochromes P450 by antifungal imidazole derivatives. *Drug Metab. Dispos.* **30**(3), 314-318 (2002).
4. Kinobe, R.T., Dercho, R.A., Vlahakis, J.Z., *et al.* Inhibition of the enzymatic activity of heme oxygenases by azole-based antifungal drugs. *J. Pharmacol. Exp. Ther.* **319**(1), 277-284 (2006).

#### 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, 10/25/2022

#### 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