

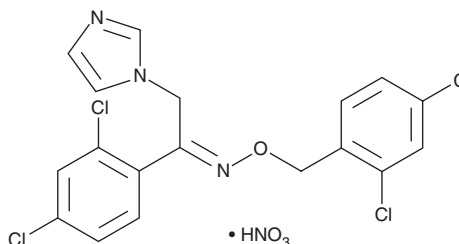
# PRODUCT INFORMATION



## Oxiconazole (nitrate)

Item No. 18724

**CAS Registry No.:** 64211-46-7  
**Formal Name:** (1Z)-O-[(2,4-dichlorophenyl)methyl] oxime 1-(2,4-dichlorophenyl)-2-(1H-imidazol-1-yl)-ethanone, mononitrate  
**Synonym:** Ro 13-8996  
**MF:** C<sub>18</sub>H<sub>13</sub>Cl<sub>4</sub>N<sub>3</sub>O • HNO<sub>3</sub>  
**FW:** 492.1  
**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

Oxiconazole (nitrate) is supplied as a crystalline solid. A stock solution may be made by dissolving the oxiconazole (nitrate) in the solvent of choice, which should be purged with an inert gas. Oxiconazole (nitrate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of oxiconazole (nitrate) in these solvents is approximately 3, 25, and 20 mg/ml, respectively.

Oxiconazole (nitrate) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, oxiconazole (nitrate) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Oxiconazole (nitrate) has a solubility of approximately 0.3 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

Oxiconazole is an antifungal imidazole that inhibits a broad spectrum of pathogenic yeasts and molds, including *A. fumigatus*, *C. neoformans*, *C. albicans*, and *T. mentagrophytes*.<sup>1</sup> Its inhibitory action results from its ability to inhibit the synthesis of ergosterol, the major sterol of the fungal cell membrane.<sup>1</sup>

### Reference

1. Fromtling, R.A. Overview of medically important antifungal azole derivatives. *Clin. Microbiol. Rev.* **1**(2), 187-217 (1988).

#### 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, 10/27/2022

#### 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](http://WWW.CAYMANCHEM.COM)