

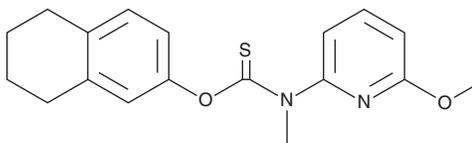
# PRODUCT INFORMATION



## Liranaftate

Item No. 44439

**CAS Registry No.:** 88678-31-3  
**Formal Name:** N-(6-methoxy-2-pyridinyl)-N-methyl-carbamothioic acid, O-(5,6,7,8-tetrahydro-2-naphthalenyl) ester  
**Synonyms:** M-732, Piritetrate  
**MF:** C<sub>18</sub>H<sub>20</sub>N<sub>2</sub>O<sub>2</sub>S  
**FW:** 328.4  
**Purity:** ≥98%  
**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

Liranaftate is supplied as a solid. A stock solution may be made by dissolving the liranaftate in the solvent of choice, which should be purged with an inert gas. Liranaftate is slightly soluble (0.1-1 mg/ml) in DMSO.

### Description

Liranaftate is a thiocarbamate antifungal agent.<sup>1</sup> It is active against *T. rubrum* (MIC = 0.009 g/ml).<sup>1</sup> Liranaftate inhibits lung epithelial cell ferroptosis and malondialdehyde (MDA) production and reduces lung IL-6, IL-1 $\beta$ , and TNF- $\alpha$  levels in a mouse model of acute lung injury induced by paraquat (Item No. 38521).<sup>2</sup>

### References

1. Oku, Y., Sakuma, K., Yokoyama, K., *et al.* Fungicidal activity of liranaftate against *Trichophyton rubrum*. *Nihon Ishinkin Gakkai Zasshi* **43(3)**, 181-187 (2002).
2. Zhu, Y., Zhao, W., Li, W., *et al.* Squalene epoxidase promotes paraquat-induced pulmonary toxicity through endoplasmic reticulum-mediated ferroptosis. *J. Adv. Res.* S2090-1232(25)00385-6 (2025).

#### 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, 01/08/2026

#### 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