

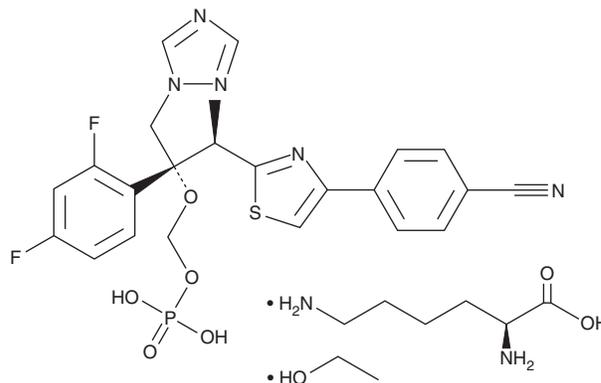
PRODUCT INFORMATION



Fosravuconazole (L-lysine ethanolate)

Item No. 42133

CAS Registry No.: 914361-45-8
Formal Name: L-Lysine, compd. with 4-[2-[(1R,2R)-2-(2,4-difluorophenyl)-1-methyl-2-[(phosphonoxy)methoxy]-3-(1H-1,2,4-triazol-1-yl)propyl]-4-thiazolyl]benzotrile and ethanol (1:1:1)
Synonyms: BMS 379224, E1224, F-RVCZ
MF: C₂₃H₂₀F₂N₅O₅PS • C₆H₁₄N₂O₂ • C₂H₆O
FW: 739.7
Purity: ≥95%
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

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

Fosravuconazole (L-lysine ethanolate) is slightly soluble (0.1-1 mg/ml) in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Fosravuconazole is a prodrug form of the triazole antifungal agent ravuconazole (Item No. 18750).¹ It increases survival in a mouse model of systemic *C. albicans* infection. Formulations containing fosravuconazole have been used in the treatment of onychomycosis.

Reference

1. Ueda, Y., Matiskella, J.D., Golik, J., *et al.* Phosphonooxymethyl prodrugs of the broad spectrum antifungal azole, ravuconazole: synthesis and biological properties. *Bioorg. Med. Chem. Lett.* **13(21)**, 3669-3672 (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
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, 10/29/2024

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