

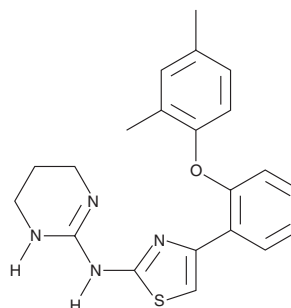
PRODUCT INFORMATION



Abafungin

Item No. 35465

CAS Registry No.: 129639-79-8
Formal Name: N-[4-[2-(2,4-dimethylphenoxy)phenyl]-2-thiazolyl]-1,4,5,6-tetrahydro-2-pyrimidinamine
Synonym: BAY-W 6341
MF: C₂₁H₂₂N₄OS
FW: 378.5
Purity: ≥98%
UV/Vis.: λ_{max}: 244, 301 nm
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

Abafungin is supplied as a solid. A stock solution may be made by dissolving the abafungin in the solvent of choice, which should be purged with an inert gas. Abafungin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of abafungin in these solvents is approximately 0.3, 2.5, and 10 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of abafungin can be prepared by directly dissolving the solid in aqueous buffers. The solubility of abafungin in PBS (pH 7.2) is approximately 0.3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Abafungin is an antifungal agent.¹ It is active against a variety of dermatophytes, yeast, molds, and other fungi (MICs = ≤0.06-2, 0.5-16, ≤0.06-1, and 0.25-8 µg/ml, respectively). Abafungin inhibits ergosterol biosynthesis, as well as acts directly on the fungal cell membrane to induce cell death.

Reference

1. Borelli, C., Schaller, M., Niewerth, M., *et al.* Modes of action of the new arylguanidine abafungin beyond interference with ergosterol biosynthesis and in vitro activity against medically important fungi. *Chemotherapy* **54**(4), 245-259 (2022).

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, 12/12/2022

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