

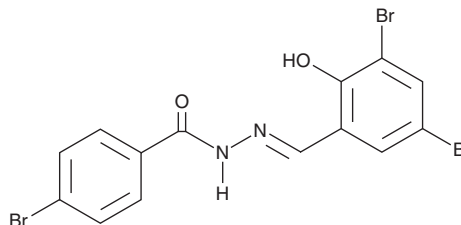
PRODUCT INFORMATION



D13

Item No. 29818

CAS Registry No.: 321945-27-1
Formal Name: 4-bromo-benzoic acid,
2-[(3,5-dibromo-2-hydroxyphenyl)
methylene]hydrazide
MF: C₁₄H₉Br₃N₂O₂
FW: 476.9
Purity: ≥98%
UV/Vis.: λ_{max}: 244, 295, 307, 341 nm
Supplied as: A crystalline 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

D13 is supplied as a crystalline solid. A stock solution may be made by dissolving the D13 in the solvent of choice, which should be purged with an inert gas. D13 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of D13 in these solvents is approximately 10 and 30 mg/ml, respectively.

Description

D13 is an acylhydrazone antifungal.¹ It is active against *C. neoformans* *in vitro* (MIC₈₀ = 0.06 µg/ml). D13 (20 mg/kg per day, p.o.) increases survival in mouse models of *C. neoformans*, *C. albicans*, or *A. fumigatus* infection.

Reference

1. Lazzarini, C., Haranahalli, K., Rieger, R., *et al.* Acylhydrazones as antifungal agents targeting the synthesis of fungal sphingolipids. *Antimicrob. Agents Chemother.* **62(5)**, e00156-18 (2018).

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, 03/20/2024

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