

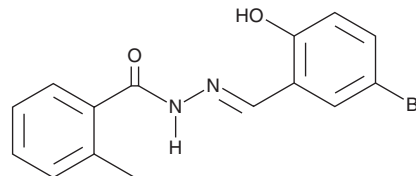
# PRODUCT INFORMATION



## BHBM

Item No. 29817

**CAS Registry No.:** 302807-84-7  
**Formal Name:** 2-methyl-benzoic acid 2-[(5-bromo-2-hydroxyphenyl)methylene]hydrazide  
**MF:** C<sub>15</sub>H<sub>13</sub>BrN<sub>2</sub>O<sub>2</sub>  
**FW:** 333.2  
**Purity:** ≥95%  
**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

BHBM is supplied as a crystalline solid. A stock solution may be made by dissolving the BHBM in the solvent of choice, which should be purged with an inert gas. BHBM is soluble in DMSO.

### Description

BHBM is an acylhydrazone with antifungal activity.<sup>1,2</sup> It is active against *C. neoformans* *in vitro* (MIC<sub>80</sub> = 1 µg/ml).<sup>2</sup> BHBM (0.25, 1, and 4 µg/ml) inhibits the synthesis of glucosylceramide, which is essential to fungal cell division, in *C. neoformans* but not J774 murine macrophages.<sup>1</sup> *In vivo*, BHBM (1.2 mg/kg per day) increases survival in a mouse model of *C. neoformans* infection.

### References

1. Haranahalli, K., Lazzarini, C., Sun, Y., et al. SAR studies on aromatic acylhydrazone-based inhibitors of fungal sphingolipid synthesis as next-generation antifungal agents. *J. Med. Chem.* **62(17)**, 8249-8273 (2019).
2. 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