

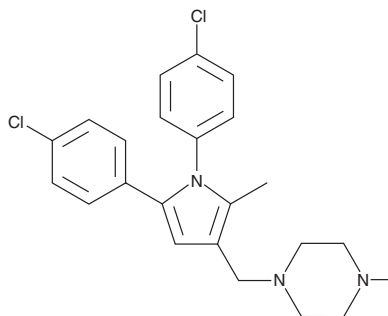
PRODUCT INFORMATION



BM212

Item No. 22986

CAS Registry No.: 146204-42-4
Formal Name: 1-[[1,5-bis(4-chlorophenyl)-2-methyl-1H-pyrrol-3-yl]methyl]-4-methyl-piperazine
MF: C₂₃H₂₅Cl₂N₃
FW: 414.4
Purity: ≥98%
UV/Vis.: λ_{max}: 247, 297 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

BM212 is supplied as a crystalline solid. A stock solution may be made by dissolving the BM212 in the solvent of choice, which should be purged with an inert gas. BM212 is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of BM212 in these solvents is approximately 0.5 mg/ml.

Description

BM212 is an antimycobacterial compound.¹ It inhibits the growth of laboratory and clinical isolate *M. tuberculosis* strains (MICs = 0.7-1.5 µg/ml), including strains resistant to ethambutol (Item No. 23713), isoniazid (Item No. 20378), rifampicin (Item No. 14423), and rifabutin (Item No. 16468). BM212 also inhibits the growth of *M. fortuitum*, *M. smegmatis*, *M. kansasii*, and *M. avium* strains (MICs = 3.1-12.5, 3.1-25, 3.1-6.2, and 0.4-3.1 µg/ml, respectively). It inhibits replication of intracellular *M. tuberculosis* in infected U937 macrophages in a dose-dependent manner at concentrations ranging from 0.5-10 µg/ml. Some *M. smegmatis*, *M. bovis* BCG, and *M. tuberculosis* H37Rv mutant strains carry mutations in the mycobacterium membrane protein large 3 (mmpL3) gene and are resistant to BM212, indicating mmpL3 is likely a cellular target of BM212.²

References

1. Deidda, D., Lampis, G., Fioravanti, R., *et al.* Bactericidal activities of the pyrrole derivative BM212 against multidrug-resistant and intramacrophagic *Mycobacterium tuberculosis* strains. *Antimicrob. Agents Chemother.* **42(11)**, 3035-3037 (1998).
2. LaRosa, V., Poce, G., Canseco, J.O., *et al.* MmpL3 is the cellular target of the antitubercular pyrrole derivative BM212. *Antimicrob. Agents Chemother.* **56(1)**, 324-331 (2012).

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, 12/21/2022

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