

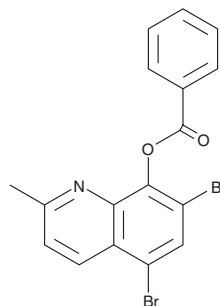
PRODUCT INFORMATION



Broxaldine

Item No. 31113

CAS Registry No.: 3684-46-6
Formal Name: 5,7-dibromo-2-methyl-8-quinolinol,
8-benzoate
Synonym: Brobenzoxaldine
MF: C₁₇H₁₁Br₂NO₂
FW: 421.1
Purity: ≥98%
UV/Vis.: λ_{max}: 239 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

Broxaldine is supplied as a solid. A stock solution may be made by dissolving the broxaldine in the solvent of choice, which should be purged with an inert gas. Broxaldine is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of broxaldine in these solvents is approximately 1 and 5 mg/ml, respectively.

Broxaldine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, broxaldine should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Broxaldine has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Broxaldine is an antimicrobial agent.^{1,2} It is active against the bacterium *C. difficile* (MIC = 4 μM) and decreases the viability of *S. mansoni* larvae at 10 μM.

References

1. AbdelKhalek, A., Mohammad, H., Mayhoub, A.S., *et al.* Screening for potent and selective anticlostridial leads among FDA-approved drugs. *J. Antibiot. (Tokyo)* **73(6)**, 392-409 (2020).
2. Panic, G., Vargas, M., Scandale, I., *et al.* Activity profile of an FDA-approved compound library against *Schistosoma mansoni*. *PLoS Negl. Trop. Dis.* **9(7)**, e0003962 (2015).

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, 12/09/2022

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