

PRODUCT INFORMATION



MRSA Antibiotic 2

Item No. 40433

CAS Registry No.: 667410-02-8
Formal Name: 5-bromo-2-[[2-(2,4-dichlorophenoxy)acetyl]amino]-benzoic acid

Synonyms: BRD-K80450985, Methicillin-resistant *S. aureus* Antibiotic 2, Methicillin-resistant *Staphylococcus aureus* Antibiotic 2

MF: C₁₅H₁₀BrCl₂NO₄

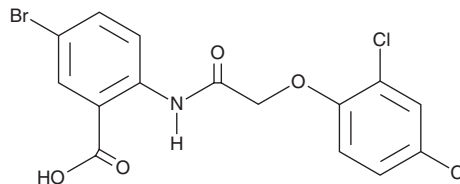
FW: 419.1

Purity: ≥95%

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

MRSA antibiotic 2 is supplied as a solid. A stock solution may be made by dissolving the MRSA antibiotic 2 in the solvent of choice, which should be purged with an inert gas. MRSA antibiotic 2 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

MRSA antibiotic 2 is an antibiotic.¹ It is active against the Gram-positive bacteria *B. subtilis*, *S. aureus*, and methicillin-resistant *S. aureus* (MRSA; MICs = 2, 2, and 4 µg/ml, respectively) but not the Gram-negative bacteria *E. coli*, *A. baumannii*, and *P. aeruginosa* (MICs = >128 µg/ml for all). MRSA antibiotic 2 (32 µg/ml) disrupts the pH gradient of *S. aureus* and *B. subtilis* in a proton motive force assay using DiSC₃(5) (3,3'-dipropylthiadicarbocyanine; Item No. 36308).

Reference

1. Wong, F., Zheng, E.J., Valeri, J.A., *et al.* Discovery of a structural class of antibiotics with explainable deep learning. *Nature* **626(7997)**, 177-185 (2024).

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, 12/10/2024

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