

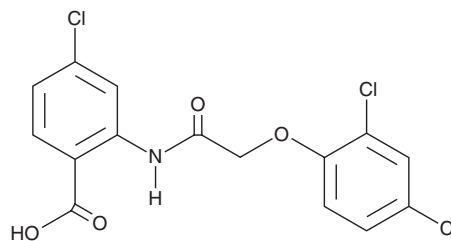
PRODUCT INFORMATION



MRSA Antibiotic 1

Item No. 40432

CAS Registry No.: 667410-81-3
Formal Name: 4-chloro-2-[[2-(2,4-dichlorophenoxy)acetyl]amino]-benzoic acid
MF: C₁₅H₁₀Cl₃NO₄
FW: 374.6
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 1 is supplied as a solid. A stock solution may be made by dissolving the MRSA antibiotic 1 in the solvent of choice, which should be purged with an inert gas. MRSA antibiotic 1 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

MRSA antibiotic 1 is an antibiotic.¹ It is active against *S. aureus*, methicillin-resistant *S. aureus* (MRSA), and *B. subtilis* (MICs = 2, 4, and 2 µg/ml, respectively) without inducing cytotoxicity in HepG2, HSkMC, or IMR-90 cells at 10 µM. MRSA antibiotic 1 is also active against 40 antibiotic-resistant clinical isolates (median MICs = 2-16 µg/ml). It decreases MRSA viral titers in the skin in a neutropenic mouse wound infection model when applied topically at 1% (w/w), as well as reduces MRSA viral titers in the thigh in a neutropenic mouse thigh infection model when administered intraperitoneally at 80 mg/kg.

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, 07/03/2024

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