

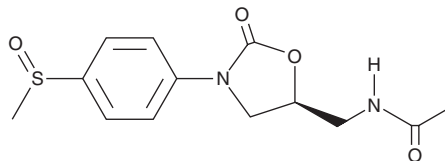
PRODUCT INFORMATION



DuP-105

Item No. 45298

CAS Registry No.: 96800-41-8
Formal Name: N-[[[(5S)-3-[4-(methylsulfinyl)phenyl]-2-oxo-5-oxazolidinyl]methyl]-acetamide
MF: $C_{13}H_{16}N_2O_4S$
FW: 296.3
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

DuP-105 is supplied as a solid. A stock solution may be made by dissolving the DuP-105 in the solvent of choice, which should be purged with an inert gas. DuP-105 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

DuP-105 is an oxazolidinone antibiotic.¹ It is active against *Staphylococcus*, *Streptococcus*, *Bacteroides*, *Fusobacterium*, and *Peptostreptococcus* species ($MIC_{50s} = 0.25-8 \mu g/ml$). *In vivo*, DuP-105 reduces lethality in immunocompetent and immunosuppressed mouse models of *S. aureus* infection ($ED_{50s} = 16.6$ and 32.7 mg/kg, respectively). It also reduces renal bacterial burden in a mouse model of *S. aureus*-induced kidney infection.

Reference

1. Slee, A. M., Wuonola, M. A., McRipley, R. J., *et al.* Oxazolidinones, a new class of synthetic antibacterial agents: in vitro and in vivo activities of DuP 105 and DuP 721. *Antimicrob Agents Chemother.* **31**(11), 1791-7 (1987).

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, 03/06/2026

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