

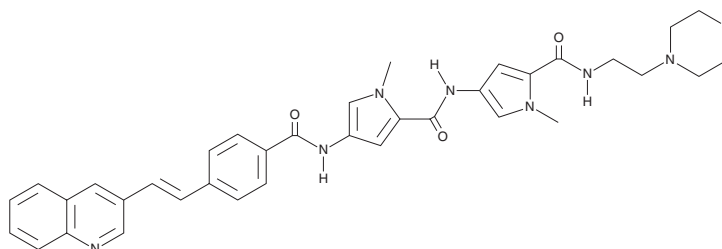
PRODUCT INFORMATION



MGB-BP-3

Item No. 42613

CAS Registry No.: 1000277-08-6
Formal Name: 1-methyl-N-[1-methyl-5-[[[2-(4-morpholinyl)ethyl]amino]carbonyl]-1H-pyrrol-3-yl]-4-[[4-[(1E)-2-(3-quinolinyl)ethenyl]benzoyl]amino]-1H-pyrrole-2-carboxamide
MF: C₃₆H₃₇N₇O₄
FW: 631.7
Purity: ≥98%
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

MGB-BP-3 is supplied as a solid. A stock solution may be made by dissolving the MGB-BP-3 in the solvent of choice, which should be purged with an inert gas. MGB-BP-3 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

MGB-BP-3 is a lexitropsin antibiotic.^{1,2} It binds to bacterial dsDNA and is active against *S. aureus*, *S. faecalis*, and *M. fortuitum*, as well as *C. albicans* (MICs = 3, 25, 50, and 50 µg/ml, respectively). *In vivo*, MGB-BP-3 (40 mg/kg) decreases joint inflammation severity in a mouse model of joint sepsis induced by *S. aureus*.²

References

1. Hind, C., Clifford, M., Woolley, C., *et al.* Insights into the spectrum of activity and mechanism of action of MGB-BP-3. *ACS Infect. Dis.* **8**(12), 2552-2563 (2022).
2. Anthony, N.G., Breen, D., Clarke, J., *et al.* Antimicrobial lexitropsins containing amide, amidine, and alkene linking groups. *J. Med. Chem.* **50**(24), 6116-6125 (2007).

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, 01/15/2025

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