

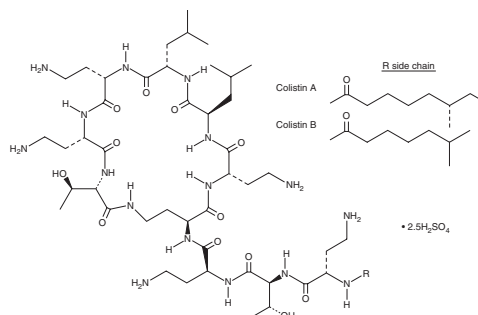
PRODUCT INFORMATION



Colistin (sulfate)

Item No. 17584

Formal Name: Colistin, sulfate salt (2:5)
Synonym: Polymyxin E Complex
MF: $[C_{52}H_{98}N_{16}O_{13}]_2 \cdot 5H_2SO_4$
FW: 2,801.3
Purity: $\geq 95\%$
Supplied as: A crystalline solid
Storage: Room temperature
Stability: ≥ 4 years



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

Laboratory Procedures

Colistin (sulfate) is supplied as a crystalline solid. Aqueous solutions of colistin (sulfate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of colistin (sulfate) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Colistin is a complex antibiotic containing greater than 30 components, with the cyclic polypeptide antibiotics polymyxins E₁ (colistin A) and E₂ (colistin B) as the major components.^{1,2} It was originally isolated from *B. polymyxa*.¹ Colistin inhibits the growth of the Gram-negative bacteria *E. coli*, *P. aeruginosa*, *P. fluorescens*, and *S. enterica* (MICs = 0.04-2.08 µg/ml) and Gram-positive *L. ivanovii* and *L. monocytogenes* (MICs = 2.5-10 µg/ml) but is not active against Gram-positive *L. lactis*, *P. polymyxa*, *P. acidilactici*, or *S. aureus* at concentrations up to 5 µg/ml.³ It also inhibits the growth of clinical isolates of both susceptible and multidrug-resistant *P. aeruginosa* (MICs = 1-2 mg/l).⁴ Colistin binds selectively to LPS from susceptible strains of *K. pneumoniae* compared to resistant strains (K_D s = 0.56 and 2.83 µM, respectively), which may contribute to its mechanism of action against Gram-negative bacteria.⁵ *In vivo*, colistin slows the growth of *P. aeruginosa* and *A. baumannii* in a neutropenic mouse model of thigh infection.⁶

References

1. Suzuki, T., Inouye, H., Fujikawa, K., *et al.* Studies on the chemical structure of colistin. I. Fractionation, molecular weight determination, amino acid and fatty acid composition. *J. Biochem.* **54**(1), 25-33 (1963).
2. Dotsikas, Y., Markopoulou, C.K., Koundourellis, J.E., *et al.* Validation of a novel LC-MS/MS method for the quantitation of colistin A and B in human plasma. *J. Sep. Sci.* **34**(1), 37-45 (2011).
3. Naghmouchi, K., Hammami, R., Fliss, I., *et al.* Colistin A and colistin B among inhibitory substances of *Paenibacillus polymyxa* JB05-01-1. *Arch. Microbiol.* **194**(5), 363-370 (2012).
4. Mitsugui, C.S., Tognim, M.C.B., Cardoso, C.L., *et al.* *In vitro* activity of polymyxins in combination with β -lactams against clinical strains of *Pseudomonas aeruginosa*. *Int. J. Antimicrob. Agents* **38**(5), 447-450 (2011).
5. Velkov, T., Deris, Z.Z., Huang, J.X., *et al.* Surface changes and polymyxin interactions with a resistant strain of *Klebsiella pneumoniae*. *Innate Immun.* **20**(4), 350-363 (2014).
6. Cheah, S.-E., Wang, J., Nguyen, V.T.T., *et al.* New pharmacokinetic/pharmacodynamic studies of systemically administered colistin against *Pseudomonas aeruginosa* and *Acinetobacter baumannii* in mouse thigh and lung infection models: Smaller response in lung infection. *J. Antimicrob. Chemother.* **70**(12), 3291-3297 (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, 11/30/2022

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