

PRODUCT INFORMATION



Polymyxin E₁ (sulfate)

Item No. 39718

CAS Registry No.: 207990-63-4

Formal Name: N²-(6-methyl-1-oxooctyl)-L-2,4-diaminobutanoyl-L-threonyl-L-2,4-diaminobutanoyl-L-2,4-diaminobutanoyl-D-leucyl-L-leucyl-L-2,4-diaminobutanoyl-L-threonine, cyclic (10→4)-peptide, sulfate

Synonym: Colistin A sulfate

MF: C₅₃H₁₀₀N₁₆O₁₃ • XH₂SO₄

FW: 1,169.5

Purity: ≥90%

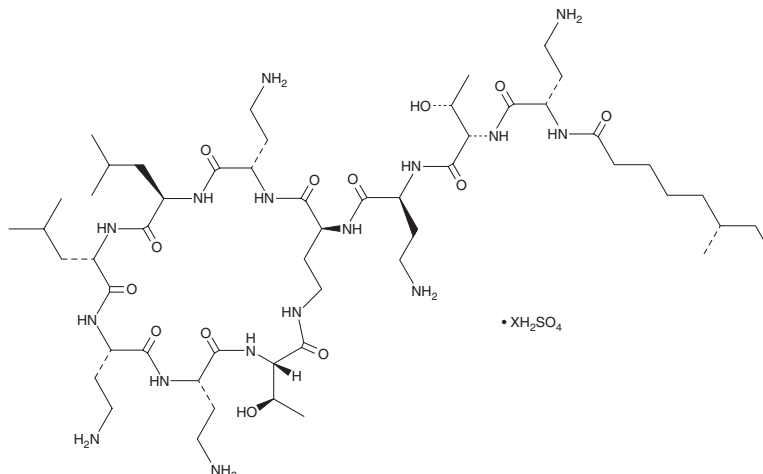
Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Bacterium/*Bacillus polymyxa* var. *colistinus*

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



Laboratory Procedures

Polymyxin E₁ (sulfate) is supplied as a solid. Aqueous solutions of polymyxin E₁ (sulfate) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of polymyxin E₁ (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

Polymyxin E₁ is a peptide antibiotic and a major constituent of the cationic lipopeptide antibiotic colistin (Item No. 17584) originally isolated from *B. polymyxa*.¹ It is active against *P. aeruginosa* and *E. coli* (MICs = 2 and 0.8 µg/ml, respectively).

Reference

1. Kline, T., Holub, D., Therrien, J., et al. Synthesis and characterization of the colistin peptide polymyxin E1 and related antimicrobial peptides. *J. Pept. Res.* **57**(3), 175-187 (2001).

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, 10/05/2023

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