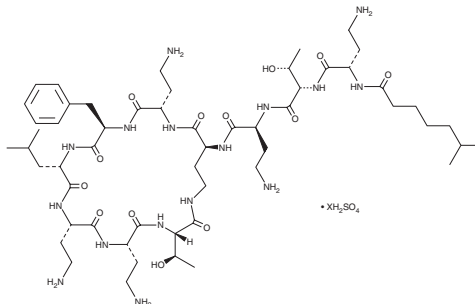


# PRODUCT INFORMATION

## Polymyxin B<sub>2</sub> (sulfate)

Item No. 39720

CAS Registry No.: 108965-69-1  
 MF: C<sub>55</sub>H<sub>96</sub>N<sub>16</sub>O<sub>13</sub> • XH<sub>2</sub>SO<sub>4</sub>  
 FW: 1,189.5  
 Purity: ≥90%  
 Supplied as: A solid  
 Storage: -20°C  
 Stability: ≥4 years  
 Item Origin: Bacterium/*Pseudomonas* sp.



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

### Laboratory Procedures

Polymyxin B<sub>2</sub> (sulfate) is supplied as a solid. A stock solution may be made by dissolving the polymyxin B<sub>2</sub> (sulfate) in water. The solubility of polymyxin B<sub>2</sub> (sulfate) in water is approximately 50 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Polymyxin B<sub>2</sub> is an antibiotic and a major constituent of the cationic lipopeptide antibiotic polymyxin B (Item No. 14157) originally isolated from *B. polymyxa*.<sup>1</sup> It is active against *P. aeruginosa*, *A. baumannii*, *K. pneumoniae*, and *E. cloacae* strains (MICs = 1-2, 0.5-1, 0.25-0.5, and 0.25-1 µg/ml, respectively). *In vivo*, polymyxin B<sub>2</sub> (4 mg/kg) reduces the number of blood colony-forming units (CFUs) in a mouse model of *P. aeruginosa* blood infection.

### Reference

1. Roberts, K.D., Azad, M.A.K., Wang, J., *et al.* Antimicrobial activity and toxicity of the major lipopeptide components of polymyxin B and colistin: Last-line antibiotics against multidrug-resistant Gram-negative bacteria. *ACS Infect. Dis.* **1**(11), 568-575 (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, 10/03/2023

#### 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