

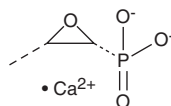
PRODUCT INFORMATION



Fosfomycin (calcium salt)

Item No. 35535

CAS Registry No.: 26016-98-8
Formal Name: P-[(2R,3S)-3-methyl-2-oxiranyl]-
phosphonic acid, monocalcium salt
Synonym: Phosphomycin Calcium
MF: C₃H₅O₄P • Ca
FW: 176.1
Purity: ≥95%
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

Fosfomycin (calcium salt) is supplied as a solid. A stock solution may be made by dissolving the fosfomycin (calcium salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Fosfomycin is a broad-spectrum bactericidal antibiotic.^{1,2} It is active against *Salmonella* and *P. aeruginosa* at 1.56 and 6.25 µg/ml, respectively, as well as methicillin-sensitive *S. aureus* (MSSA), cephalosporin- and penicillin-resistant *S. pneumoniae*, methicillin-resistant *S. aureus* (MRSA), and vancomycin-resistant *Enterococcus*. Formulations containing fosfomycin have been used in the treatment of urinary tract infections.

References

1. Goto, S. Fosfomycin, antimicrobial activity *in vitro* and *in vivo*. *Chemotherapy* **23(Suppl. 1)**, 63-74 (1977).
2. Michalopoulos, A.S., Livaditis, I.G., and Gougoutas, V. The revival of fosfomycin. *Int. J. Infect. Dis.* **15(11)**, e732-e739 (2011).

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, 10/25/2022

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