

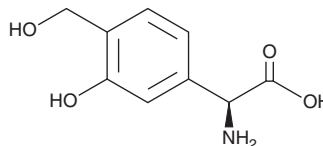
PRODUCT INFORMATION



Forphenicol

Item No. 36382

CAS Registry No.: 71522-58-2
Formal Name: αS-amino-3-hydroxy-4-(hydroxymethyl)-benzeneacetic acid
Synonym: Forfenimex
MF: C₉H₁₁NO₄
FW: 197.2
Purity: ≥90%
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

Forphenicol is supplied as a solid. A stock solution may be made by dissolving the forphenicol in the solvent of choice. Forphenicol is soluble in organic solvents such as methanol and DMSO, which should be purged with an inert gas. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Forphenicol is an immunomodulator and a derivative of the bacterial metabolite forphenicine (Item No. 33504).¹ It increases the phagocytosis of yeast by peritoneal macrophages isolated from thioglycolate-stimulated mice. Forphenicol (100 µg/animal) prevents cyclophosphamide-induced suppression of delayed-type hypersensitivity (DTH), as well as enhances DTH in response to the hapten oxazolone or sheep red blood cells in mice. It enhances the bactericidal activity of macrophages against *P. aeruginosa* in mice when administered at a dose of 0.5 mg/kg.³ Forphenicol (15.6-1,000 µg/animal) increases survival in a mouse model of *P. aeruginosa* infection.² It also inhibits tumor growth in S180 sarcoma and IMC carcinoma mouse xenograft models when administered at doses ranging from 0.05 to 5 mg/kg per day.

References

1. Ishizuka, M., Ishizeki, S., Masuda, T., *et al.* Studies on effects of forphenicol on immune responses. *J. Antibiot. (Tokyo)* **35(8)**, 1042-1048 (1982).
2. Ishizuka, M., Masuda, T., Kanbayashi, N., *et al.* Antitumor effect of forphenicol, a low molecular weight immunomodifier, on murine transplantable tumors and microbial infections. *J. Antibiot. (Tokyo)* **35(8)**, 1049-1054 (1982).
3. Naito, K., Arakawa, H., Watanabe, Y., *et al.* Protective effect of forphenicol, a low molecular weight immunomodifier, against infection with *Pseudomonas aeruginosa* in mice and its mechanisms. *J. Antibiot. (Tokyo)* **40(4)**, 547-545 (1987).

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, 12/19/2022

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