

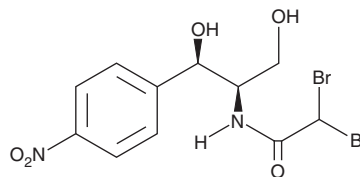
PRODUCT INFORMATION



Bromamphenicol

Item No. 28016

CAS Registry No.: 17371-30-1
Formal Name: [R-(R*,R*)]-2,2-dibromo-N-[2-hydroxy-1-(hydroxymethyl)-2-(4-nitrophenyl)ethyl]-acetamide
MF: C₁₁H₁₂Br₂N₂O₅
FW: 412.0
Purity: ≥98%
UV/Vis.: λ_{max}: 273 nm
Supplied as: A crystalline 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

Bromamphenicol is supplied as a crystalline solid. A stock solution may be made by dissolving the bromamphenicol in the solvent of choice, which should be purged with an inert gas. Bromamphenicol is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

Bromamphenicol is a dibrominated derivative of the antibiotic chloramphenicol.¹ It inhibits rat liver mitochondrial and *E. coli* protein synthesis by 90.6 and 98.8%, respectively, when used at a concentration of 93 μM, and inhibits DNA synthesis in human lymphoblastoid cells by 83% at 1 mM. Bromamphenicol can also bind to the major adhesin subunit DraE from *E. coli*.²

References

1. Manyan, D.R., Arimura, G.K., and Yunis, A.A. Comparative metabolic effects of chloramphenicol analogues. *Mol. Pharmacol.* **11(5)**, 520-527 (1975).
2. Pettigrew, D.M., Roversi, P., Davies, S.G., *et al.* A structural study of the interaction between the Dr haemagglutinin DraE and derivatives of chloramphenicol. *Acta Crystallogr. D Biol. Crystallogr.* **65(Pt 6)**, 513-522 (2009).

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, 12/02/2022

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