

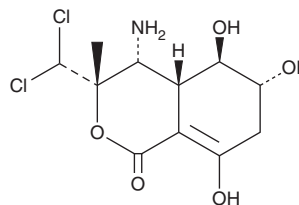
PRODUCT INFORMATION



Bactobolamine

Item No. 38896

CAS Registry No.: 78013-07-7
Formal Name: (3S,4R,4aR,5R,6R)-4-amino-3-(dichloromethyl)-3,4,4a,5,6,7-hexahydro-5,6,8-trihydroxy-3-methyl-1H-2-benzopyran-1-one
MF: C₁₁H₁₅Cl₂NO₅
FW: 312.1
Purity: ≥98%
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

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

Description

Bactobolamine is an immunosuppressant.¹ It inhibits IL-2 production by CTLL cells when used at a concentration of 200 µg/ml. *In vivo*, bactobolamine (200 mg/kg) prolongs heart allograft survival in MHC-compatible and MHC-incompatible rat pairings. It also prolongs survival in a rat liver allograft model and a canine kidney allograft model.

Reference

1. Tomikawa, S., Beck, Y., Watanabe, K., *et al.* Effectiveness of bactobolamine as a new immunosuppressive agent. *Transplant Proc.* **29(1-2)**, 1276-1278 (1997).

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, 10/10/2023

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