

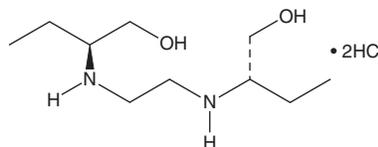
PRODUCT INFORMATION



Ethambutol (hydrochloride)

Item No. 23713

CAS Registry No.: 1070-11-7
Formal Name: 2S,2'S-(1,2-ethanediyldiimino)bis-1-butanol, dihydrochloride
MF: $C_{10}H_{24}N_2O_2 \cdot 2HCl$
FW: 277.2
Purity: $\geq 98\%$
Supplied as: A solid
Storage: 4°C
Stability: ≥ 4 years



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

Laboratory Procedures

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

Description

Ethambutol is an antimycobacterial compound that inhibits the growth of *M. tuberculosis in vitro* at a concentration of 10 $\mu\text{g/ml}$.¹ It is protective against *M. smegmatis* infections in rhesus monkeys (50, 30, and 0% fatalities at doses of 12.5, 25, and 50 mg/kg, respectively) and clears *M. smegmatis* pulmonary infiltrates at a dose of 100 mg/kg.² Formulations containing ethambutol have been used in the treatment of tuberculosis.

References

1. Dickinson, J.M. and Mitchison, D.A. In vitro and in vivo studies to assess the suitability of anti-tuberculous drugs for use in intermittent chemotherapy regimens. *Bull. Int. Union. Tuberc.* **41(S1)**, 309-315 (1968).
2. Schmidt, L.H. Studies on the antituberculous activity of ethambutol in monkey. *Ann. N.Y. Acad. Sci.* **135(2)**, 747-758 (1966).

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, 11/04/2022

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