

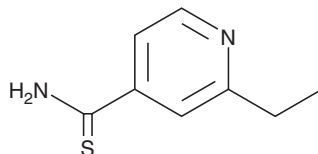
PRODUCT INFORMATION



Ethionamide

Item No. 23714

CAS Registry No.: 536-33-4
Formal Name: 2-ethyl-4-pyridinecarbothioamide
Synonym: NSC 255115
MF: C₈H₁₀N₂S
FW: 166.2
Purity: ≥98%
UV/Vis.: λ_{max}: 290 nm
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

Ethionamide is supplied as a solid. A stock solution may be made by dissolving the ethionamide in the solvent of choice. Ethionamide is soluble in the organic solvent DMSO, which should be purged with an inert gas. Ethionamide is also slightly soluble in methanol.

Description

Ethionamide is an antimycobacterial compound that is active against *M. tuberculosis* (MICs = 0.3-1.25 µg/ml).^{1,2} It is activated *via* oxidation by flavin monooxygenase and inhibits the InhA enzyme involved in mycolic acid biosynthesis.¹ Formulations containing ethionamide have been used in the second-line treatment of multi-drug resistant tuberculosis.

References

1. Laborde, J., Deraeve, C., and Bernardes-Génisson, V. Update of antitubercular prodrugs from a molecular perspective: Mechanisms of action, bioactivation pathways, and associated resistance. *ChemMedChem* **12(20)**, 1657-1676 (2017).
2. Heifets, L.B., Lindholm-Levy, P.J., and Flory, M. Comparison of bacteriostatic and bactericidal activity of isoniazid and ethionamide against *Mycobacterium avium* and *Mycobacterium tuberculosis*. *Am. Rev. Respir. Dis.* **143(2)**, 268-270 (1991).

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, 10/21/2022

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