

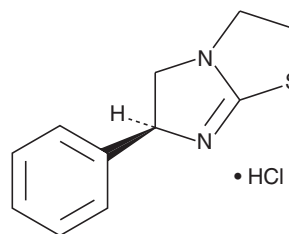
PRODUCT INFORMATION



(-)-Levamisole (hydrochloride)

Item No. 14874

CAS Registry No.: 16595-80-5
Formal Name: (6S)-2,3,5,6-tetrahydro-6-phenyl-imidazo[2,1-b]thiazole, monohydrochloride
Synonyms: NSC 177023, R 12654, L-Tetramisole
MF: C₁₁H₁₂N₂S • HCl
FW: 240.8
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥1 year



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

Description

(-)-Levamisole (hydrochloride) (Item No. 14874) is an analytical reference material categorized as an imidazothiazole. (-)-Levamisole is used as a cutting agent with cocaine, and levamisole-adulterated cocaine has led to cutaneous vasculopathy in some individuals.¹⁻³ This product is intended for research and forensic applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards for reference materials.

References

1. Broséus, J., Gentile, N., and Esseiva, P. The cutting of cocaine and heroin: A critical review. *Forensic Sci. Int.* **262**, 73-83 (2016).
2. Pillow, M.T. and Hughes, A. Levamisole-adulterated cocaine induced vasculitis with skin ulcerations. *West J. Emerg. Med.* **14(2)**, 149-150 (2013).
3. Tran, H., Tan, D., and Marnejon, T.P. Cutaneous vasculopathy associated with levamisole-adulterated cocaine. *Clin. Med. Res.* **11(1)**, 26-30 (2013).

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, 09/26/2023

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