

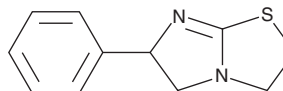
PRODUCT INFORMATION



Tetramisole

Item No. 39242

CAS Registry No.: 5036-02-2
Formal Name: 2,3,5,6-tetrahydro-6-phenyl-imidazo[2,1-b]thiazole
Synonyms: Phenyltetrahydroimidazothiazole, PTHIT
MF: C₁₁H₁₂N₂S
FW: 204.3
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.

Description

Tetramisole (Item No. 39242) is an analytical reference standard categorized as an imidazothiazole. Tetramisole is a racemic form of (-)-levamisole (Item Nos. 21321 | 14874 | 33972) and has been found as an adulterant in cocaine seized by law enforcement.¹ This product is intended for research and forensic applications.

Reference

1. Hundertmark, M., Hess, C., Röhrich, J., *et al.* Analysis of tetramisole metabolites - Is "Aminorex" found in forensic samples of cocaine users actually 4-phenyl-2-imidazolidinone? *Drug Test Anal.* **15**(8), 865-878 (2023).

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, 03/10/2025

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