

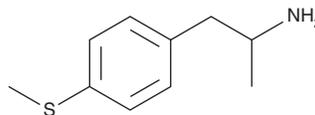
PRODUCT INFORMATION



4-MTA

Item No. 11555

CAS Registry No.: 14116-06-4
Formal Name: α -methyl-4-(methylthio)-benzeneethanamine
Synonyms: 4-Methylthioamphetamine, P 1882
MF: C₁₀H₁₅NS
FW: 181.3
Purity: \geq 98%
UV/Vis.: λ_{max} : 205, 257 nm
Supplied as: A solution in ethanol
Storage: -20°C
Stability: \geq 4 years



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

Description

4-MTA is a methylthio analog of para-methoxyamphetamine (PMA), a hallucinogenic drug that has been scheduled in many countries.¹ In animal models, the effects of 4-MTA are qualitatively and quantitatively indistinguishable from those of PMA.¹ The metabolism of this compound has recently been described.¹ This product is intended for forensic applications.

Reference

1. Staack, R.F. and Maurer, H.H. Metabolism of designer drugs of abuse. *Curr. Drug Metab.* **6(3)**, 259-274 (2005).

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, 06/13/2023

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