

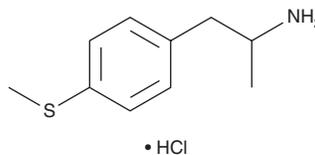
PRODUCT INFORMATION



4-MTA (hydrochloride)

Item No. 11862

CAS Registry No.: 94784-92-6
Formal Name: α -methyl-4-(methylthio)-benzeneethanamine, monohydrochloride
Synonyms: 4-Methylthioamphetamine, P 1882
MF: $C_{10}H_{15}NS \cdot HCl$
FW: 217.8
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 258 nm
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 5 years



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

Description

4-MTA (Item No. 11862), is a selective serotonergic agent which increases serotonin release, blocks its re-uptake ($IC_{50} = 102$ nM), and inhibits monoamine oxidase A.¹⁻³ The detection of this designer drug in serum and urine has been described, as has its metabolism.³⁻⁶ This product is intended for research and forensic applications.

References

- Huang, X., Marona-Lewicka, D., and Nichols, D.E. p-Methylthioamphetamine is a potent new non-neurotoxic serotonin-releasing agent. *Eur. J. Pharmacol.* **229(1)**, 31-38 (1992).
- Gobbi, M., Moia, M., Pirona, L., et al. p-Methylthioamphetamine and 1-(m-chlorophenyl)piperazine, two non-neurotoxic 5-HT releasers *in vivo*, differ from neurotoxic amphetamine derivatives in their mode of action at 5-HT nerve endings *in vitro*. *J. Neurochem.* **82(6)**, 1435-1443 (2002).
- Staack, R.F. and Maurer, H.H. Metabolism of designer drugs of abuse. *Curr. Drug Metab.* **6(3)**, 259-274 (2005).
- Wohlfarth, A., Weinmann, W., and Dresen, S. LC-MS/MS screening method for designer amphetamines, tryptamines, and piperazines in serum. *Anal. Bioanal. Chem.* **396(7)**, 2403-2414 (2010).
- Kerrigan, S., Banuelos, S., Perrella, L., et al. Simultaneous detection of ten psychedelic phenethylamines in urine by gas chromatography-mass spectrometry. *J. Anal. Toxicol.* **35(7)**, 459-469 (2011).
- Ewald, A.H., Peters, F.T., Weise, M., et al. Studies on the metabolism and toxicological detection of the designer drug 4-methylthioamphetamine (4-MTA) in human urine using gas chromatography-mass spectrometry. *J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.* **824(1-2)**, 123-131 (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, 06/15/2023

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