

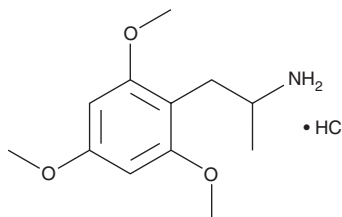
# PRODUCT INFORMATION



## 2,4,6-Trimethoxyamphetamine (hydrochloride)

Item No. 12099

**CAS Registry No.:** 23815-74-9  
**Formal Name:** 2,4,6-trimethoxy- $\alpha$ -methylbenzeneethanamine, monohydrochloride  
**Synonym:** TMA-6  
**MF:**  $C_{12}H_{19}NO_3 \cdot HCl$   
**FW:** 261.7  
**Purity:**  $\geq 98\%$   
**Supplied as:** A crystalline solid  
**Storage:**  $-20^{\circ}C$   
**Stability:**  $\geq 5$  years



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

### Description

The trimethoxyamphetamines (TMA) are a family of designer drugs developed on an amphetamine base. TMA-6, known more formally as 2,4,6-trimethoxyamphetamine, is a psychedelic hallucinogen which has appeared on the illicit drug market.<sup>1,2</sup> Unlike amphetamines and cathinones, TMA-6 has no effect on the release or re-uptake of monoamines (dopamine, serotonin, and norepinephrine).<sup>3</sup> This product is intended for research and forensic applications.

### References

1. Nieddu, M., Boatto, G., Pirisi, M.A., *et al.* LC-MS analysis of trimethoxyamphetamine designer drugs (TMA series) from urine samples. *J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.* **867(1)**, 126-130 (2008).
2. Zhang, Z. and Simpkins, J.W. Okadaic acid induces tau phosphorylation in SH-SY5Y cells in an estrogen-preventable manner. *Brain Res.* **1345**, 176-181 (2010).
3. Nagai, F., Nonaka, R., and Satoh Hisashi Kamimura, K. The effects of non-medically used psychoactive drugs on monoamine neurotransmission in rat brain. *Eur. J. Pharmacol.* **559(2-3)**, 132-137 (2007).

#### 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, 07/25/2023

#### 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