

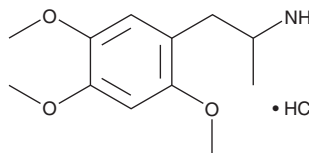
PRODUCT INFORMATION



2,4,5-Trimethoxyamphetamine (hydrochloride)

Item No. 33715

CAS Registry No.: 15995-72-9
Formal Name: 2,4,5-trimethoxy- α -methylbenzeneethanamine, monohydrochloride
Synonyms: TMA-2, 2,4,5-TMA
MF: C₁₂H₁₉NO₃ • HCl
FW: 261.7
Purity: \geq 98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 5 years



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

Description

2,4,5-Trimethoxyamphetamine (hydrochloride) (Item No. 33715) is an analytical reference standard categorized as an amphetamine. 2,4,5-Trimethoxyamphetamine has psychoactive properties.¹ 2,4,5-Trimethoxyamphetamine is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

Reference

1. Ewald, A.H. and Maurer, H.H. 2,5-dimethoxyamphetamine-derived designer drugs: Studies on the identification of cytochrome P450 (CYP) isoenzymes involved in formation of their main metabolites and on their capability to inhibit CYP2D6. *Toxicol. Lett.* **183(1-3)**, 52-57 (2008).

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, 06/23/2023

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