

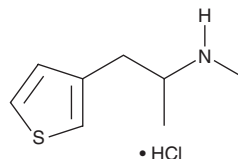
PRODUCT INFORMATION



Methiopropamine 3'-thiophene isomer (hydrochloride)

Item No. 9001980

CAS Registry No.: 857361-90-1
Formal Name: N, α -dimethyl-3-thiopheneethanamine, monohydrochloride
Synonym: 3-MPA
MF: C₈H₁₃NS • HCl
FW: 191.7
Purity: \geq 98%
Supplied as: A neat solid
Storage: -20°C
Stability: \geq 2 years



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

Description

Methiopropamine (Item No. 11161) is a structural analog of methamphetamine (Item Nos. ISO60168 | 13997) in which the phenyl group has been replaced by thiophene with the sulfur atom at the 2 position. Methiopropamine 3'-thiophene isomer is an isomer of methiopropamine that has the sulfur atom of the thiophene group at the 3 position. The physiological and toxicological properties of this compound are not known. This product is intended for forensic and research applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards.

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, 07/11/2019

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