

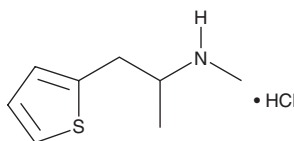
PRODUCT INFORMATION



Methiopropamine (hydrochloride)

Item No. 11161

CAS Registry No.: 7464-94-0
Formal Name: N, α -dimethyl-2-thiopheneethanamine, monohydrochloride
Synonyms: MPA, 2-MPA, NSC 400137
MF: C₈H₁₃NS • HCl
FW: 191.7
Purity: \geq 98%
Stability: \geq 2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{\max} : 233 nm



Laboratory Procedures

For long term storage, we suggest that methiopropamine (MPA) (hydrochloride) be stored as supplied at -20°C. It should be stable for at least two years.

MPA (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the MPA (hydrochloride) in the solvent of choice. MPA (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of MPA (hydrochloride) in ethanol and DMF is approximately 20 mg/ml and approximately 10 mg/ml in DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of MPA (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of MPA (hydrochloride) in PBS, pH 7.2, is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Methiopropamine is a structural analog of methamphetamine in which the phenyl group has been replaced with thiophene. The physiological and toxicological properties of this compound have not been studied. This product is intended for forensic and research applications.

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, 02/26/2016

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