

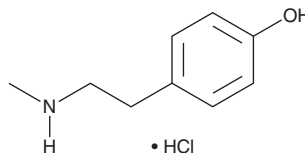
PRODUCT INFORMATION



N-methyl Tyramine (hydrochloride)

Item No. 35233

CAS Registry No.: 13062-76-5
Formal Name: 4-[2-(methylamino)ethyl]-phenol, monohydrochloride
MF: C₉H₁₃NO • HCl
FW: 187.7
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

N-methyl Tyramine (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the N-methyl tyramine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. N-methyl Tyramine (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of N-methyl tyramine (hydrochloride) in these solvents is approximately 1, 3, and 5 mg/ml, respectively.

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 N-methyl tyramine (hydrochloride) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of N-methyl tyramine (hydrochloride) in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

N-methyl Tyramine is a phenethylamine alkaloid that has been found in *C. aurantium* and has antilipolytic activity.¹ It inhibits lipolysis induced by isoprenaline (isoproterenol; Item No. 15592) in isolated rat and human adipocytes when used at a concentration of 5.8 μM.

Reference

1. Mercader, J., Wanecq, E., Chen, J., *et al.* Isopropyl norepinephrine is a stronger lipolytic agent in human adipocytes than synephrine and other amines present in *Citrus aurantium*. *J. Physiol. Biochem.* **67(3)**, 443-452 (2011).

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, 11/17/2022

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