

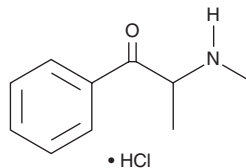
# PRODUCT INFORMATION



## Methcathinone (hydrochloride) (exempt preparation)

Item No. 15656

**CAS Registry No.:** 49656-78-2  
**Formal Name:** 2-(methylamino)-1-phenyl-1-propanone, monohydrochloride  
**Synonym:** DL-Ephedrone  
**MF:** C<sub>10</sub>H<sub>13</sub>NO • HCl  
**FW:** 199.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 249 nm  
**Supplied as:** A solution in methanol  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

Methcathinone (hydrochloride) (exempt preparation) is supplied as a solution in methanol. To change the solvent, simply evaporate the methanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of methcathinone (hydrochloride) (exempt preparation) in these solvents is approximately 5, 2, and 1 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. If an organic solvent-free solution of methcathinone (hydrochloride) (exempt preparation) is needed, it can be prepared by evaporating the methanol and directly dissolving the neat oil in aqueous buffers. The solubility of methcathinone (hydrochloride) (exempt preparation) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Methcathinone (hydrochloride) (exempt preparation) (Item No. 15656) is an analytical reference standard categorized as a psychoactive cathinone. It is a β-ketone analog of methamphetamine and has comparable effects on monoamine uptake.<sup>1</sup> Methcathinone is regulated as a Schedule I controlled substance in the United States. This product is intended for forensic and research applications only.

### Reference

1. Cozzi, N.V., Sievert, M.K., Shulgin, A.T., *et al.* Inhibition of plasma membrane monoamine transporters by β-ketoamphetamines. *Eur. J. Pharmacol.* **381(1)**, 63-69 (1999).

#### 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, 04/17/2018

#### 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