

PRODUCT INFORMATION



2-Methoxymethcathinone (hydrochloride)

Item No. 9001186

Formal Name: 1-(2-methoxyphenyl)-2-(methylamino)propan-1-one, monohydrochloride

Synonym: 2-MeOMC

MF: C₁₁H₁₅NO₂ • HCl

FW: 229.7

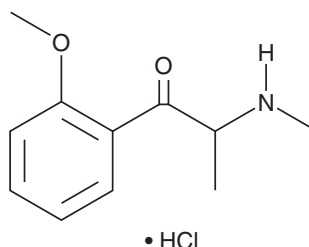
Purity: ≥98%

UV/Vis.: λ_{max}: 212, 254, 317 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥5 years



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

Description

Methedrone (Item No. 10529), also known as *para*-methoxymethcathinone, is a psychoactive compound that has been identified as an adulterant of party pills and bath salts.¹ 2-MeOMC (hydrochloride) is a positional isomer of methedrone, having the methoxy group at the 2 rather than the 4 position. It is also the β-keto derivative of 2-methoxymethamphetamine, a β-adrenergic receptor agonist used as a bronchodilator. The properties of 2-MeOMC have been poorly studied. This product is intended for forensic and research applications.

Reference

1. Kikura-Hanajiri, R., Uchiyama, N., and Goda, Y. Survey of current trends in the abuse of psychotropic substances and plants in Japan. *Leg. Med. (Tokyo)* **13**(3), 109-15 (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, 07/06/2023

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