

PRODUCT INFORMATION



Methylmethcathinone Isomer Mixture

Item No. 44681

Synonym: MMC Isomer Mixture
Supplied as: A solution in methanol (1 mg/ml of each compound)
Fill Volume: >1 ml
Storage: -20°C
Stability: ≥3 years

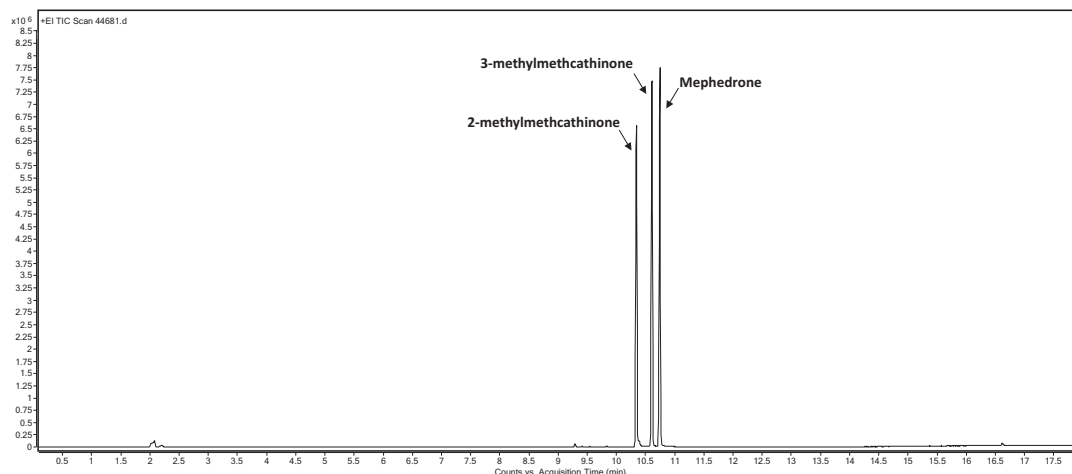
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description and Data

Methylmethcathinone Isomer Mixture (Item No. 44681) is an analytical reference standard mixture containing methylmethcathinone isomers. The mixture contains 2-methylmethcathinone, 3-methylmethcathinone, and mephedrone. The mixture is supplied in an amber ampule in which the headspace has been purged with argon. It is designed for use as a reference standard for these major controlled substances during routine analytical checks using GC-MS instrumentation. Methylmethcathinone Isomer Mixture contains compounds that are regulated as Schedule I compounds in the United States. This product is intended for research and forensic applications.

This mixture contains methylmethcathinone. Included are:

Item No. 44681	Methylmethcathinone Isomer Mixture	
Item Name	CAS Number	Concentration (mg/ml)
2-Methylmethcathinone (hydrochloride)	1246815-51-9	1.0 (as free base)
3-Methylmethcathinone (hydrochloride)	1246816-62-5	1.0 (as free base)
Mephedrone (hydrochloride)	1189726-22-4	1.0 (as free base)



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, 03/05/2026

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