

CERTIFICATE of ANALYSIS



Methylmethcathinone Isomer Mixture

Item No.: 44681
Batch No.: 0814188

Expiry Date: 23FEB2029 (valid from date of certification)
Supplied as: Each analyte at 1 mg/ml solution (as free base) in methanol; in a deactivated glass ampule

Volume per Ampule: Not less than 1 ml. Ampules are overfilled. It is advised that labs use measured volumes.

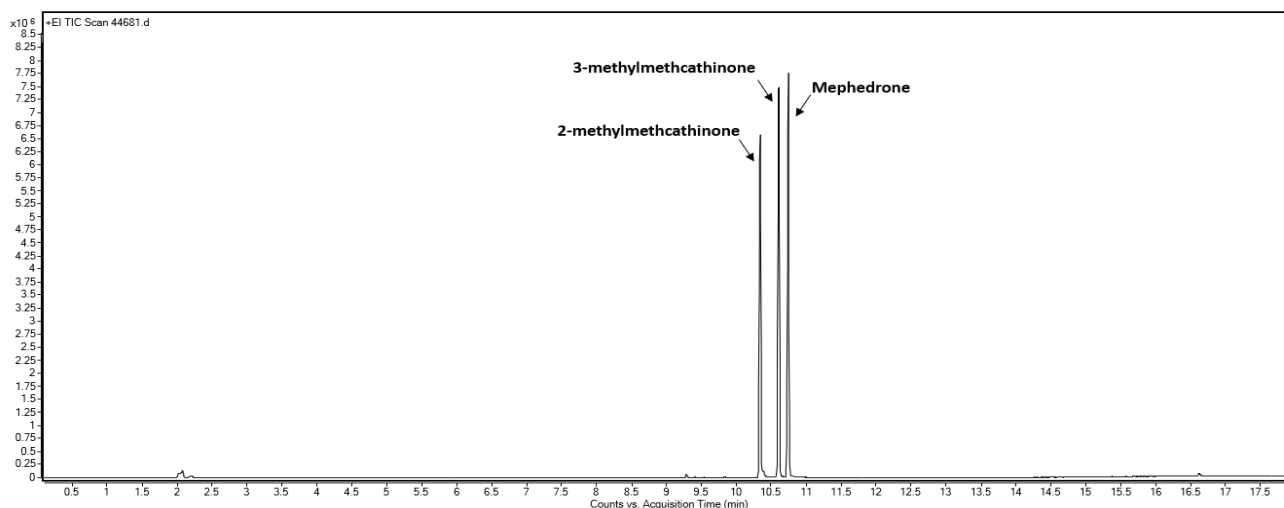
Storage & Handling: Flammable. Store upright and unopened at -20°C. Warm to room temperature prior to opening.

Description

A mixture of methylmethcathinone isomers in methanol solution intended for GC-MS, LC-MS, and/or HPLC-UV applications.

Quality Information

Test	Limit	Result	Meets Specification
GC-MS	All analytes identified	Pass	✓



Approval:

Title: Manager, Analytical Chemistry

Certification Date: 23FEB2026

Cayman Chemical certifies that this product meets the specifications stated in this certificate and warrants this product to meet the stated acceptance criteria through the expiry date when stored unopened as recommended.

CERTIFICATE of ANALYSIS



Quality Control Table

Item Number: 44681		Item Name: Methylmethcathinone Isomer Mixture	Batch Number: 0814188
Item No.	Item Name	Formula	RT (min)
11223	2-Methylmethcathinone (hydrochloride)	$C_{11}H_{15}NO \cdot HCl$	10.34
11224	3-Methylmethcathinone (hydrochloride)	$C_{11}H_{15}NO \cdot HCl$	10.61
10801	Mephedrone (hydrochloride)	$C_{11}H_{15}NO \cdot HCl$	10.74
Conditions:			
Column: HP-5ms Ultra Inert (30 m x 250 μ m x 0.25 μ m)			
Flow Rate: 1.0 ml/min			
Inlet Temp: 250°C			
Split: 20:1			
Carrier gas: He			
Oven Program: 60°C hold for 3 min; 20°C/min to 300°C; 300°C hold for 18 min			

CERTIFICATE of ANALYSIS



Revision History

Revision No.	Date	Reason for Revision
01	23FEB2026	Initial version

Disclaimers

Material Safety Data

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some but not all of the information required for the safe and proper use of this material. Before use, review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

Warranty and Limitation of Remedy

Cayman Chemical Company makes no warranty or guarantee of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specification.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver of Buyer of all claims hereunder with respect to said material. For further details, please refer to our Warranty and Limitations of Remedy located on our website and in our catalog.

This Certificate shall not be reproduced except in full, without written approval from the Cayman Chemical Company.

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 • USA

PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640

techserv@caymanchem.com
www.caymanchem.com