

PRODUCT INFORMATION



GC-MS Drug Standard Mixture 5

Item No. 24639

Supplied as: A solution in methanol (1 mg/ml of each compound)

Fill volume: >1 ml

Storage: -20°C

Stability: ≥6 years

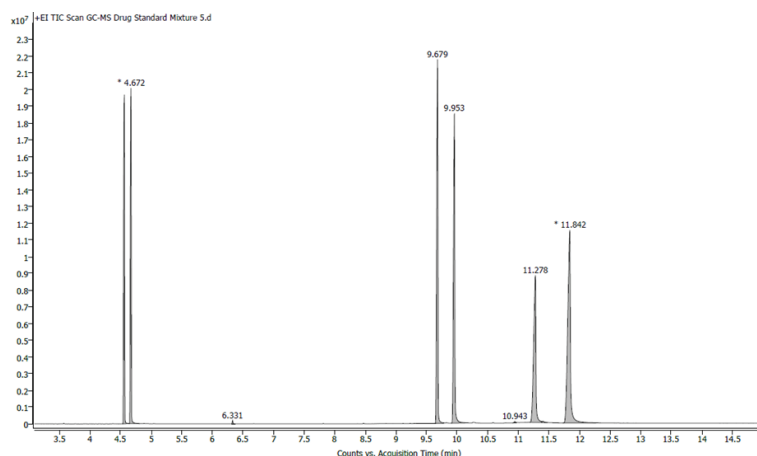
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description and Data

GC-MS drug standard mixture 5 (Item No. 24639) is a mixture of analytical reference standards and analytical reference materials. It contains common amphetamines, opioids, and benzodiazepines. 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. This product is intended for research and forensic applications.

This mixture contains compounds that are regulated as Schedule I compounds in the United States. Included are:

Item Number	Item Name	Concentration
38010	Phentermine	1 mg/ml
36566	(±)-Methamphetamine	1 mg/ml
15465	Oxycodone	1 mg/ml
15461	Hydrocodone	1 mg/ml
14263	Clonazepam	1 mg/ml
14255	Alprazolam	1 mg/ml



Item Number: 24639		Item Name: GC-MS Drug Standard Mixture 5	
Item Number	Item Name	Formula Weight:	RT (min)
38010	Phentermine	149.2	4.56
36566	(±)-Methamphetamine	149.2	4.67
15465	Oxycodone	315.4	9.95
15461	Hydrocodone	146.0	9.67
14263	Clonazepam	315.7	11.27
14255	Alprazolam	308.8	11.84
GC-MS Conditions:			
Column: 30 m x 0.25 mm, 0.25 µm HP-5MS			
Flow Rate: 2 ml/min			
Inlet Temp: 300°C			
Split: 15:1			
Carrier gas: He			
Oven Program: Hold 50°C for 1 min, 30°C/min to 300°C hold 300°C for 5.67 min.			
Negative Electrospray Ionization		Full MS/Survey Scan	

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, 02/10/2026

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