

PRODUCT INFORMATION



GC-MS Drug Standard Mixture 6

Item No. 43673

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

Fill volume: >1 ml

Storage: -20°C

Stability: ≥2 years

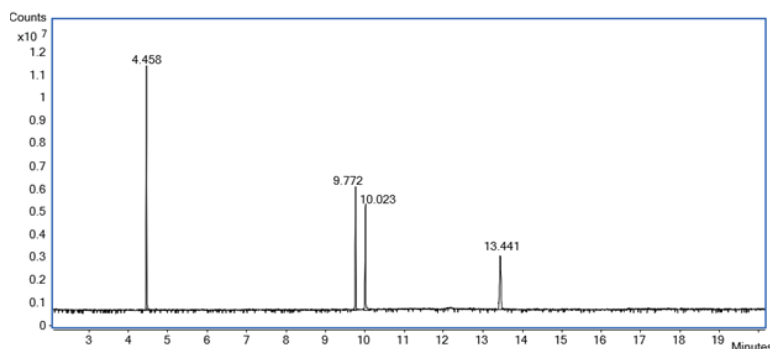
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description and Data

GC-MS drug standard mixture 6 (Item No. 43673) is a mixture of analytical reference standards containing an amphetamine and opioids. The mixture contains D-amphetamine, hydrocodone, oxycodone, and buprenorphine. The mixture is supplied in an amber ampule in which the headspace has been purged with argon. It is designed for use in method development and as a reference standard for these substances during routine analytical checks using GC-MS instrumentation. GC-MS drug standard mixture 6 contains compounds that are regulated as Schedule II and III compounds in the United States. This product is intended for research and forensic applications.

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

Item Number	Item Name	Concentration
40155	D-Amphetamine	1 mg/ml
15461	Hydrocodone	1 mg/ml
15465	Oxycodone	1 mg/ml
14025	Buprenorphine (hydrochloride)	1 mg/ml



Item Number: 43673		Item Name: GC-MS Drug Standard Mixture 6	
Item Number	Item Name	Formula Weight:	RT (min)
40155	D-Amphetamine	149.2	4.46
15461	Hydrocodone	299.4	9.77
15465	Oxycodone	315.4	10.02
14025	Buprenorphine (hydrochloride)	504.1	13.44
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, 08/25/2025

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