

PRODUCT INFORMATION



GC-MS Drug Standard Mixture 4

Item No. 24632

Supplied as: A solution in methanol (250 µg/ml of each compound)
Fill volume: >1 ml
Storage: -20°C
Stability: ≥7 years

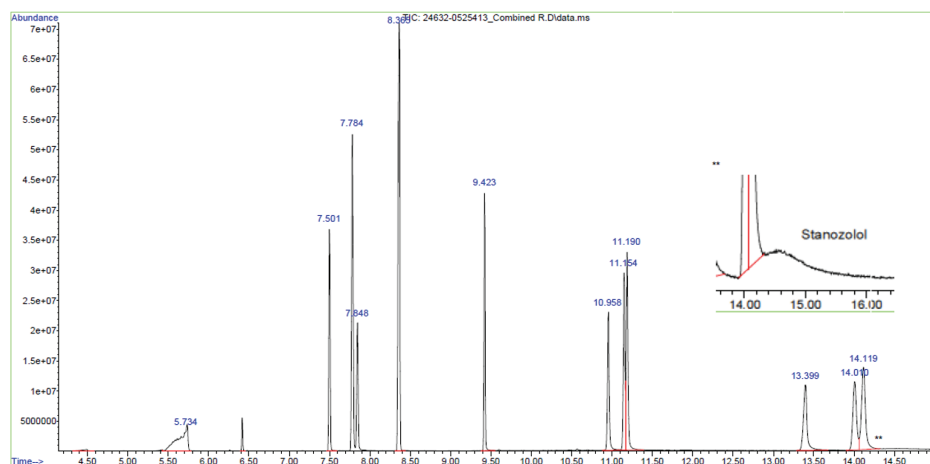
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description and Data

GC-MS drug standard mixture 4 (Item No. 24632) is a mixture of analytical reference standards and analytical reference materials. It contains common opioids, stimulants, and steroids as well as a synthetic cannabinoid and a benzodiazepine. 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 forensic and research applications.

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

Item Number	Item Name	Concentration
9001195	α-Pyrrolidinobutophenone (hydrochloride)	250 µg/ml
9001083	α-Pyrrolidinopentiophenone (hydrochloride)	250 µg/ml
9001123	bk-MDEA (hydrochloride)	250 µg/ml
10393	Butylone (hydrochloride)	250 µg/ml
17014	Tenocyclidine (hydrochloride)	250 µg/ml
14276	PCP (hydrochloride)	250 µg/ml
16186	Cocaine	250 µg/ml
21173	Nandrolone	250 µg/ml
23631	(S)-5-fluoro ADB	250 µg/ml
9001543	Heroin	250 µg/ml
14255	Alprazolam	250 µg/ml
18705	Furanyl fentanyl (hydrochloride)	250 µg/ml
21213	Furanyl fentanyl 3-furancarboxamide isomer (hydrochloride)	250 µg/ml
21167	Stanozolol	250 µg/ml



Item Number: 24632		GC-MS Drug Standard Mixture 4
Item Number	Item Name	RT (min)
9001195	α-Pyrrolidinobutophenone (hydrochloride)	7.50
9001083	α-Pyrrolidinopentiophenone (hydrochloride)	7.78
9001123	bk-MDEA (hydrochloride)	7.78
10393	Butylone (hydrochloride)	7.84
17014	Tenocyclidine (hydrochloride)	8.36
14276	PCP (hydrochloride)	8.36
16186	Cocaine	9.42
21173	Nandrolone	10.96
23631	(S)-5-fluoro ADB	10.15
9001543	Heroin	11.19
14255	Alprazolam	13.40
18705	Furanyl fentanyl (hydrochloride)	11.01
21213	Furanyl fentanyl 3-furancarboxamide isomer (hydrochloride)	14.12
21167	Stanozolol	14.50

LC-MS Conditions:

Column: 30 m × 0.32 mm, 0.5 µm Rtx-5MS
 Flow Rate: 2 ml/min
 Inlet Temp: 300° C
 Split: 15:1
 Carrier Gas: He
 Transfer Line Temp: 300° C
 Oven Program: hold 50° C for 1 min, 30° C/min to 300° C hold 300° C for 20 min

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, 07/14/2023

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