

# PRODUCT INFORMATION



## GC-MS Drug Standard Mixture 3

Item No. 23539

**Supplied as:** A solution in acetonitrile (1 mg/ml of each compound)  
**Fill volume:** >1 ml  
**Storage:** -20°C  
**Stability:** ≥5 years

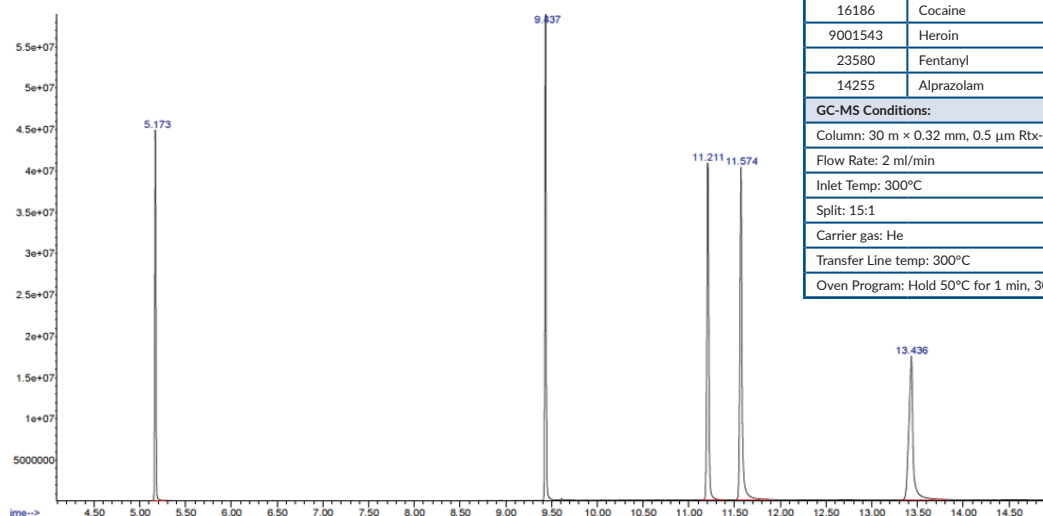
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Description and Data

GC-MS drug standard mixture 3 (Item No. 23539) is a mixture of analytical reference materials and an analytical reference standard. It contains common opioids and stimulants 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, and IV compounds in the United States. Included are:

Item Number	Item Name	Concentration
	(±)-Methamphetamine	1 mg/ml
16186	Cocaine	1 mg/ml
9001543	Heroin	1 mg/ml
23580	Fentanyl	1 mg/ml
14255	Alprazolam	1 mg/ml



Item Number: 23539		GC-MS Drug Standard Mixture 3
Item Number	Item Name	RT (min)
	(±)-Methamphetamine	5.2
16186	Cocaine	9.4
9001543	Heroin	11.2
23580	Fentanyl	11.6
14255	Alprazolam	13.4

**GC-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 5.67 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, 01/04/2024

#### 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