

PRODUCT INFORMATION



GC-MS Drug Standard Mixture 2

Item No. 21724

Supplied as: A solution in methanol (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 Contents

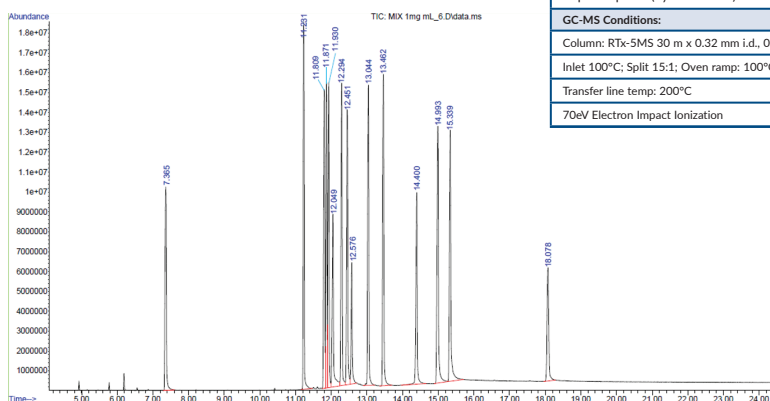
GC-MS drug standard mixture 2 (Item No. 21724) is a mixture of analytical reference standards and analytical reference materials. It contains common opioids, benzodiazepines, and cannabinoids as well as an androgenic anabolic steroid and a cathinone. The mixture is supplied in an amber ampule in which 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-IV compounds in the United States. Included are:

Item Name	Concentration
bk-MDEA (hydrochloride)	1 mg/ml
(±)-Cannabichromene	1 mg/ml
Diazepam	1 mg/ml
Δ ⁹ -THC	1 mg/ml
Hydrocodone	1 mg/ml
Hydromorphone	1 mg/ml
Cannabinol	1 mg/ml
Oxycodone	1 mg/ml
Oxymorphone	1 mg/ml
Testosterone	1 mg/ml
Fentanyl (hydrochloride)	1 mg/ml
Clonazepam	1 mg/ml
Alprazolam	1 mg/ml
AB-FUBINACA	1 mg/ml
Buprenorphine (hydrochloride)	1 mg/ml

Item Number: 21724	GC-MS Drug Standard Mixture 2	
Item Name	Retention Time (min)	GC-MS Library (Pass/Fail)
bk-MDEA (hydrochloride)	7.4	Pass
(±)-Cannabichromene	11.2	Pass
Diazepam	11.8	Pass
Δ ⁹ -THC	11.9	Pass
Hydrocodone	11.9	Pass
Hydromorphone	12.0	Pass
Cannabinol	12.3	Pass
Oxycodone	12.5	Pass
Oxymorphone	12.6	Pass
Testosterone	13.0	Pass
Fentanyl (hydrochloride)	13.5	Pass
Clonazepam	14.4	Pass
Alprazolam	15.0	Pass
AB-FUBINACA	15.3	Pass
Buprenorphine (hydrochloride)	18.1	Pass

GC-MS Conditions:	
Column: RTX-5MS 30 m x 0.32 mm i.d., 0.5 μm	
Inlet 100°C; Split 15:1; Oven ramp: 100°C hold for 2 min, 30°C/min to 325°C, hold 10 min	
Transfer line temp: 200°C	Flow Rate: 2 ml/min
70eV Electron Impact Ionization	



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, 07/17/2023

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