



Certificate of Analysis

GC-MS Drug Standard Mixture 2

Item #: 21724

Batch #: 0660536

Expiry: 03OCT2027 (valid from date of certification)

Contents: 1 mg/mL of each compound in methanol

Storage and Handling: Temperature sensitive. Product should be stored as supplied at -20°C.

Results:

Method	Specification	Specification Met
GC-MS	All analytes identified	Yes

Catalog Number: 21724 Batch Number: 0660536		GC-MS Drug Test Mixture 2	
Item #	Name	Retention Time (min)	GC-MS Library (Pass/Fail)
9001123	Ethylone (aka bk-MDEA) HCl	6.8	Pass
21721	Cannabichromene	10.4	Pass
15887	Diazepam	10.9	Pass
ISO00144	Hydrocodone	11.0	Pass
ISO00157	Δ^9 -THC	11.03	Pass
ISO00146	Hydromorphone	11.1	Pass
12066	Cannabinol	11.4	Pass
15465	Oxycodone	11.5	Pass
15911	Oxymorphone	11.6	Pass
ISO00154	Testosterone	12.1	Pass
14719	Fentanyl HCl	12.6	Pass
14263	Clonazepam	13.4	Pass
14255	Alprazolam	14.0	Pass
14039	AB-FUBINACA	14.5	Pass
14025	Buprenorphine HCl	16.6	Pass
GC-MS Conditions:			
Column: HP-5MS 30m x 0.25mm i.d., 0.25uM			
Inlet: 200°C; Split 15:1			
Oven Ramp: 100°C hold for 2 min., 30°C/min. to 210°C, 12°C/min to 325°C. Hold 325°C for 10 min.			
MSD Transfer Line: 300°C		Flow Rate: 2mL/min.	
70eV Electron Impact Ionization			

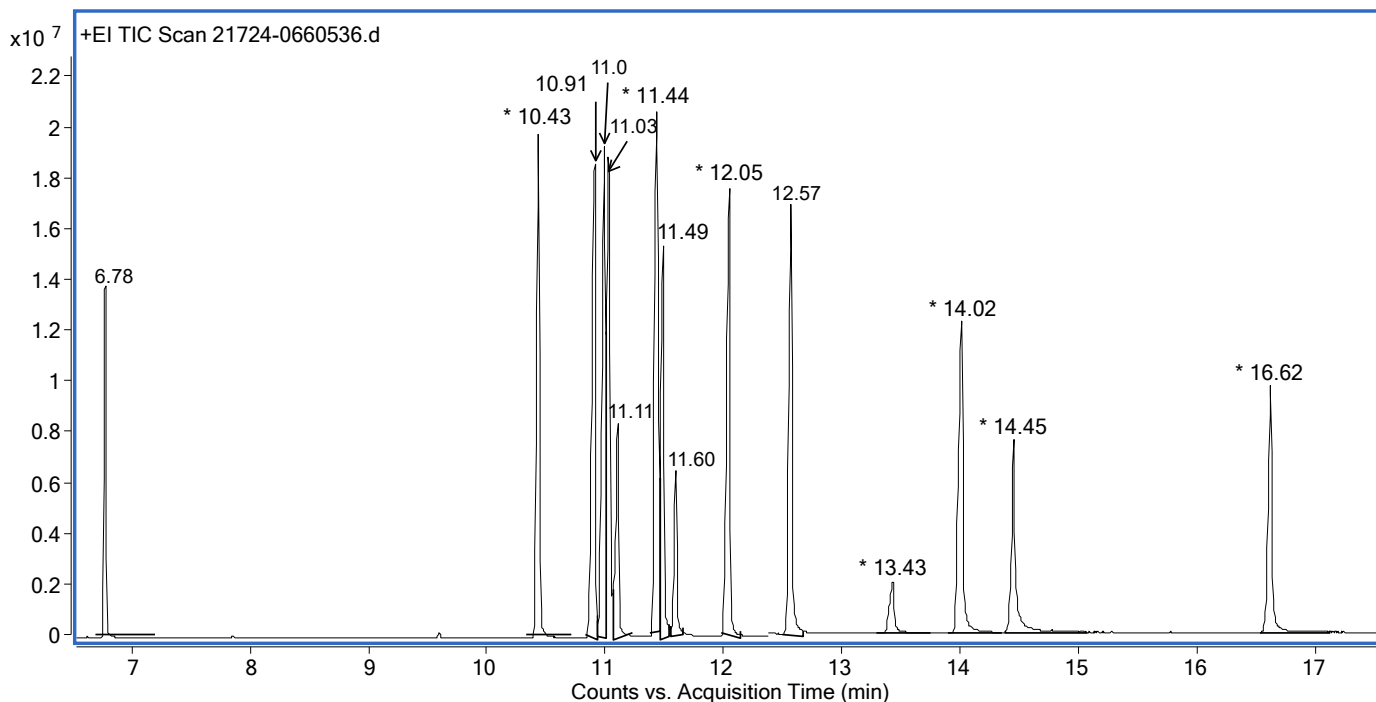
Cayman Chemical Co.

1180 E. Ellsworth Rd
Ann Arbor, MI 48108
www.caymanchem.com

Contact Us:

1-800-364-9897 or (734)971-3335
Fax: (734)971-3640
Email: techserv@caymanchem.com

For R&D purposes only. Not intended for human or animal use.



Mixture was prepared using gravimetric and volumetric solution preparation procedures from a combination of neat and formulated pure standards. The concentration of this mixture was prepared with an estimated error of $\pm 10\%$.

Cayman Chemical certifies that this product meets the specifications stated in this certificate and warrants this product to meet the stated acceptance criteria through the expiration date when stored unopened as recommended.

Approval:  Title: Director of Analytical Chemistry

Certification Date: 03OCT2022