

PRODUCT INFORMATION



Synthetic Cannabinoid Custom Mix 1

Item No. 9002566

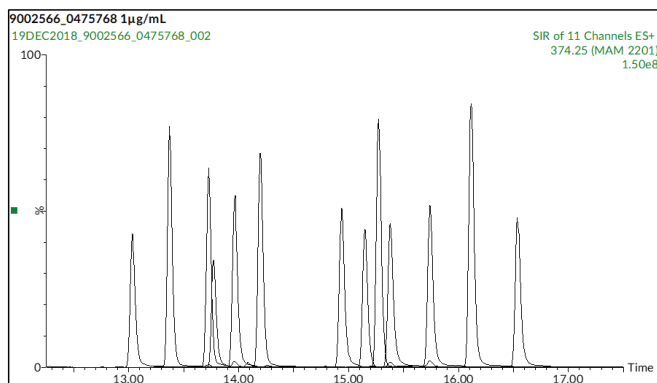
Supplied as: A solution in methanol (100 µg/ml of each compound)
Fill Volume: >1 ml
Storage: -20°C
Stability: ≥7 years

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

Description and Data

Synthetic cannabinoid mixture 1 (Item No. 9002566) is a mixture of analytical reference standards and analytical reference materials. It contains common synthetic cannabinoids. 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 LC-MS instrumentation. This product is intended for forensic and research applications.

This mixture contains compounds that are regulated as Schedule I compounds in the United States. Included are UR-144 N-(5-chloropentyl) analog (Item No. 11951), AM2201 (Item No. 10707), JWH 018 (Item No. 10900), JWH 018-quinolinyl carboxamide (Item No. 14891), JWH 073 (Item No. 10904), JWH 122 (Item No. 10591), JWH 210 (Item No. 10644), JWH 250 (Item No. 13634), MAM2201 (Item No. 9001219), UR-144 (Item No. ISO00055), UR-144 degradant (Item No. 11928), XLR11 (Item No. 11565), and XLR11 degradant (Item No. 14055).



Item Number: 9002566		Synthetic Cannabinoid Custom Mix 1	
Item Number	Item Name	Formula Weight:	RT (min)
11951	UR-144-N-(5-chloropentyl) analog	345.9	15.15
10707	AM2201	359.4	13.03
10900	JWH 018	341.5	14.94
14891	JWH 018 8-quinolinyl carboxamide	357.5	15.38
10904	JWH 073	327.4	13.97
10591	JWH 122	355.5	15.74
10644	JWH 210	369.5	16.53
13634	JWH 250	335.4	13.73
9001219	MAM2201	373.5	13.77
ISO00055	UR-144	311.5	15.27/16.11*
11928	UR-144 degradant	311.5	15.27/16.11*
11565	XLR11	329.5	13.37/14.20*
14055	XLR11 degradant	329.5	13.37/14.20*

*individual retention times not confirmed

LC-MS Conditions: Waters Acquity UPLC – Xevo TQD

Mobile Phase A: Water with 0.1% formic acid

Mobile Phase B: Acetonitrile with 0.1% formic acid

Column: Waters Cortecs C18, 2.1x100mm, 1.6 µm Flow Rate: 400 µl/min

LC Gradient: 0min 10%B, 20min 95%B, 21.5min 95%B, 21.6-25min 10%B

Positive Electrospray Ionization Full Scan: SIR

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, 09/22/2023

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