

PRODUCT INFORMATION



Synthetic Cannabinoid Mixture 3

Item No. 9002568

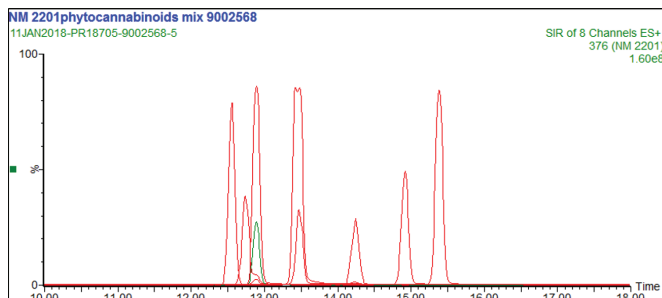
Supplied as: A solution in acetonitrile (100 µg/ml of each compound)
Fill Volume: >1 ml
Storage: -20°C
Stability: ≥7 years

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

Description and Data

Synthetic cannabinoid mixture 3 (Item No. 9002568) 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 5-fluoro NNEI (Item No. 14147), 5-fluoro PB-22 (Item No. 14095), 5-fluoro PB-22 3-hydroxyquinoline isomer (Item No. 14511), FUB-144 (Item No. 15537), MMB-FUBINACA (Item No. 9001960), FUB-PB-22 (Item No. 14949), NM2201 (Item No. 15334), PB-22 (Item No. ISO00122), and THJ2201 (Item No. 14789).



Item Number: 9002568		Synthetic Cannabinoid Mixture 3	
Item Number	Item Name	Formula Weight	RT (min)
14147	5-fluoro NNEI	374.5	12.9
14095	5-fluoro PB-22	376.4	12.7/13.4
14511	5-fluoro PB-22 3-hydroxyquinoline isomer	376.4	12.7/13.4
15537	FUB-144	349.5	15.4
9001960	MMB-FUBINACA	383.4	12.5
14949	FUB-PB-22	396.4	13.5
15334	NM2201	375.4	12.9
ISO00122	PB-22	358.4	14.2
14789	THJ2201	360.4	14.9

LC-MS Conditions:

Mobile Phase A: Water + 0.1% Formic Acid	
Mobile Phase B: Acetonitrile:Water (90:10)	
Column: Waters Cortecs C18, 2.1 x 100 mm, 1.6 µm	Flow Rate: 400 µl/min
LC Gradient: 0 minutes 10% B, 20 minutes 100% B, 21.5 minutes 100% B, 21.6 - 25 minutes 10% B	
Positive Electrospray Ionization	SIR Scan

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