

PRODUCT INFORMATION



GC-MS Heroin Adulterant Mixture

Item No. 37356

Supplied as: A solution in acetonitrile (1 mg/ml of each compound)

Fill Volume: >1 ml

Storage: -20°C

Stability: ≥5 years

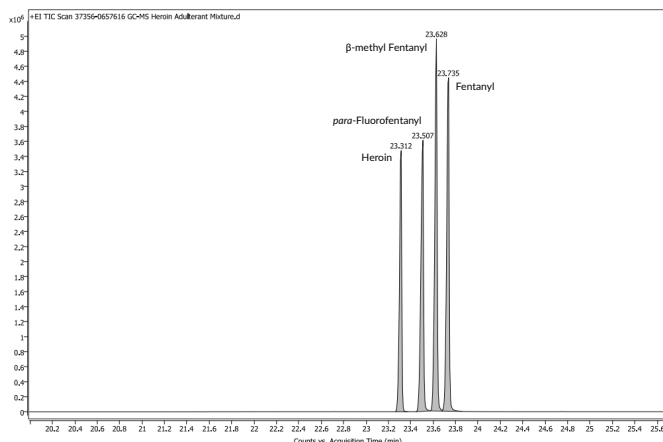
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description and Data

GC-MS heroin adulterant mixture (Item No. 37356) is an analytical reference standard mixture that contains heroin, *para*-fluorofentanyl, β -methyl fentanyl, and fentanyl. 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. GC-MS heroin adulterant mixture contains compounds that are regulated as Schedule I compounds in the United States. This product is intended for research and forensic applications.

This mixture contains opioids and opioid adulterants. Included are:

Item Number	Item Name	Concentration
9001543	Heroin	1 mg/ml
15496	<i>para</i> -Fluorofentanyl (hydrochloride)	1 mg/ml (as free base)
9002860	β -methyl Fentanyl (hydrochloride)	1 mg/ml (as free base)
23580	Fentanyl	1 mg/ml



Item No.	Item Name	RT (min)	NIST Match
9001543	Heroin	23.3	Pass
15496	<i>para</i> -Fluorofentanyl (hydrochloride)	23.5	Pass
9002860	β -methyl Fentanyl (hydrochloride)	23.6	Pass
23580	Fentanyl	23.7	Pass
GC-MS Conditions			
Column: 30 m x 250 μ m x 0.25 μ m		Flow Rate: 1 ml/min	
Inlet Temperature: 250°C			
Injection Volume: 1 μ l			
Split: 20:1			
Carrier Gas: He			
Over Program: Hold 60°C for 10 min. 20°C/min. to 300°C; hold 300°C for 10 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.

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