

# PRODUCT INFORMATION



## GC-MS Opioid Mixture 1

Item No. 23629

**Supplied as:** A solution in methanol  
**Fill volume:** >1 ml  
**Storage:** -20°C  
**Stability:** ≥5 years

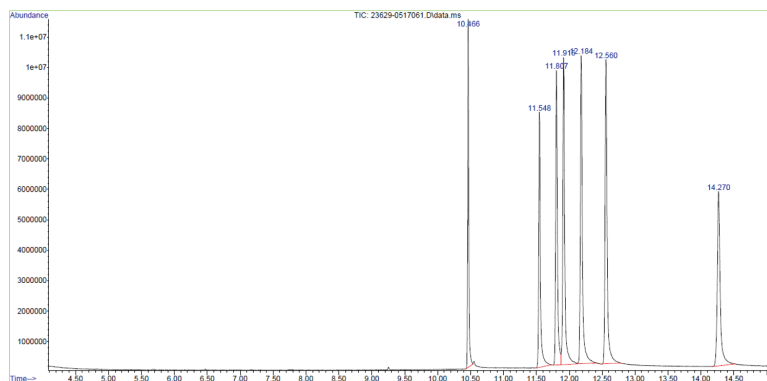
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Description and Data

GC-MS opioid mixture 1 (Item No. 23629) is a mixture of analytical reference standards and analytical reference materials. It contains common opioids. 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 and Schedule II compounds in the United States. Included are:

Item Number	Item Name	Concentration
18596	U-47700	1 mg/ml
14719	Fentanyl (hydrochloride)	1 mg/ml
ISO00128	Acetyl fentanyl (hydrochloride)	1 mg/ml
19312	Acrylfentanyl (hydrochloride)	1 mg/ml
14728	Butyryl fentanyl (hydrochloride)	1 mg/ml
21739	Cyclopropyl fentanyl (hydrochloride)	1 mg/ml
18705	Furanyl fentanyl (hydrochloride)	1 mg/ml



Item Number: 23629		GC-MS Opioid Mixture 1
Item Number	Item Name	Retention Time (min)
18596	U-47700	10.5
14719	Fentanyl (hydrochloride)	11.5
ISO00128	Acetyl fentanyl (hydrochloride)	11.8
19312	Acrylfentanyl (hydrochloride)	11.9
14728	Butyryl fentanyl (hydrochloride)	12.2
21739	Cyclopropyl fentanyl (hydrochloride)	12.6
18705	Furanyl fentanyl (hydrochloride)	14.3
<b>GC-MS Conditions:</b>		
Initial Temp: 50°C		
Initial Time: 1 min		
Ramp: 30°C/min to 300°C		
Final Time: 5.67 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.

Copyright Cayman Chemical Company, 07/14/2023

### 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