

PRODUCT INFORMATION



GC-MS Opioid Mixture 2

Item No. 23876

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

Fill volume: >1 ml

Storage: -20°C

Stability: ≥7 years

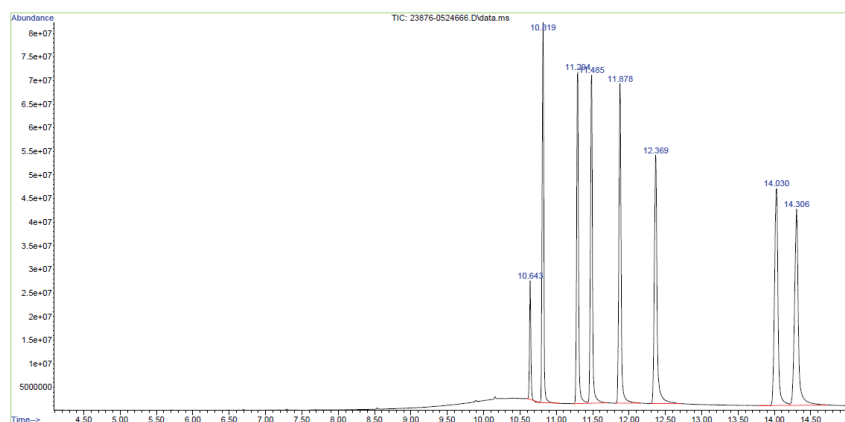
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description and Contents

GC-MS opioid mixture 2 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 research and forensic applications.

This mixture contains compounds that are regulated as Schedule I compounds in the United States as well as non-regulated compounds. Included are:

Item Number	Item Name	Concentration
22278	U-48800 (hydrochloride)	1 mg/ml
21098	U-49900	1 mg/ml
19313	FIBF (hydrochloride)	1 mg/ml
19425	ortho-Fluorofentanyl (hydrochloride)	1 mg/ml
9002747	(±)-cis-3-methyl Fentanyl (hydrochloride)	1 mg/ml
20782	Methoxyacetyl fentanyl (hydrochloride)	1 mg/ml
20439	Cyclopentyl fentanyl (hydrochloride)	1 mg/ml
20859	Tetrahydrofuran fentanyl (hydrochloride)	1 mg/ml



Item Number: 23876		GC-MS Opioid Mixture 2
Item Number	Item Name	RT (min)
22278	U-48800 (hydrochloride)	10.6
21098	U-49900	10.8
19313	FIBF (hydrochloride)	11.3
19425	ortho-Fluorofentanyl (hydrochloride)	11.5
9002747	(±)-cis-3-methyl Fentanyl (hydrochloride)	11.9
20782	Methoxyacetyl fentanyl (hydrochloride)	12.4
20439	Cyclopentyl fentanyl (hydrochloride)	14.0
20859	Tetrahydrofuran fentanyl (hydrochloride)	14.3
LC-MS Conditions:		
Column: 30 m × 0.32 mm, 0.5 µm Rtx-5MS		Flow Rate: 2 ml/min
Inlet Temp: 300°C		Split: 15:1
Carrier gas: He		Transfer Line temp: 300°C
Oven program: hold 50°C for 1 min, 30°C/min to 300°C hold 300°C for 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