

PRODUCT INFORMATION



Fluorofentanyl Isomer Mixture

Item No. 37028

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

Fill Volume: >1 ml

Storage: -20°C

Stability: ≥5 years

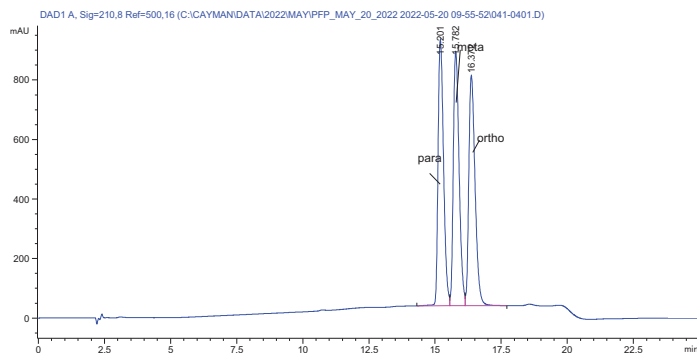
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description and Data

Fluorofentanyl isomer mixture (Item No. 37028) is a mixture of analytical reference standards and analytical reference materials containing *para*-, *meta*-, and *ortho*-fluorofentanyl. 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 HPLC-UV or LC- or GC-MS instrumentation. Fluorofentanyl isomer 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 fluorofentanyl isomers. Included are:

Item Number	Item Name	Concentration
15496	<i>para</i> -Fluorofentanyl (hydrochloride)	1 mg/ml
19424	<i>meta</i> -Fluorofentanyl (hydrochloride)	1 mg/ml
19425	<i>ortho</i> -Fluorofentanyl (hydrochloride)	1 mg/ml



Item Number: 37028		Fluorofentanyl Isomer Mixture
Item Number	Item Name	RT (min)
15496	<i>para</i> -Fluorofentanyl (hydrochloride)	15.20
19424	<i>meta</i> -Fluorofentanyl (hydrochloride)	15.78
19425	<i>ortho</i> -Fluorofentanyl (hydrochloride)	16.37
LC-MS Conditions:		
Column: Phenomenex Luna biphenyl 250x4.6mm		Flow Rate: 1 ml/min
Wavelength: 210 nm		
Mobile Phase A: 10:90:0.1 MeOH:H ₂ O 10mM Amm. Acetate		
Mobile Phase B: 90:10:0.1 MeOH:H ₂ O 10mM Amm. Acetate		
Gradient: 70%B to 85%B over 15.1 minutes		

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, 07/18/2023

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