

PRODUCT INFORMATION



Methylfentanyl Isomer Mixture

Item No. 44680

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

Storage: -20°C

Stability: ≥3 years

Fill volume: >1 ml

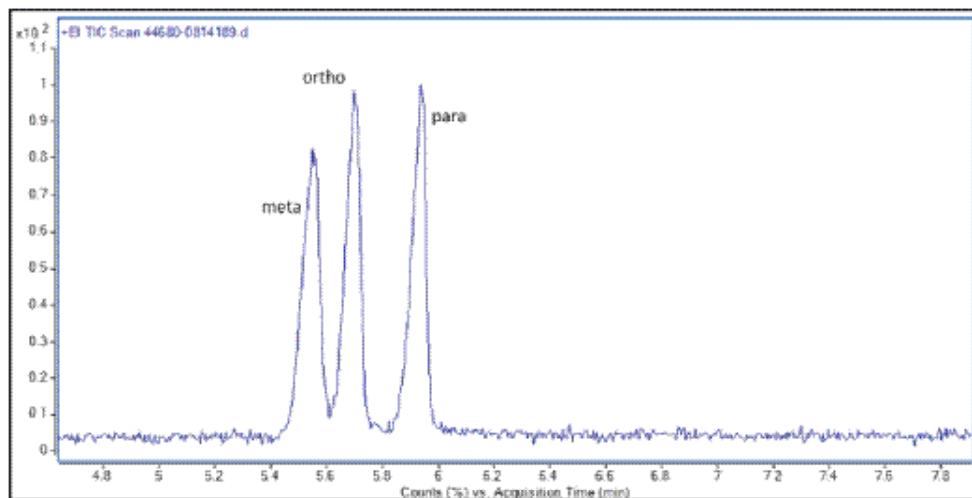
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description and Data

Methylfentanyl Isomer Mixture (Item No. 44680) is a mixture of analytical reference standards containing methylfentanyl isomers. The mixture contains *para*-, *meta*-, and *ortho*-methylfentanyl. The mixture is supplied in an amber ampule in which the headspace has been purged with argon. It is designed for use in method development and as a reference standard for these substances during routine analytical checks using GC-MS, LC-MS, or HPLC-UV instrumentation. Methylfentanyl 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 three methylfentanyl isomers. Included are:

Item No. 44680	Methylfentanyl Isomer Mixture	
Item Name	CAS Number	Concentration (mg/ml)
<i>para</i> -Methylfentanyl (hydrochloride)	1807-12-1	1.0 (as free base)
<i>meta</i> -Methylfentanyl (hydrochloride)	1465-22-1	1.0 (as free base)
<i>ortho</i> -Methylfentanyl (hydrochloride)	1443-53-4	1.0 (as free base)



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, 02/10/2026

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