

PRODUCT INFORMATION



meta-methoxy Furanyl fentanyl

Item No. 26050

CAS Registry No.: 2749434-68-0
Formal Name: N-(3-methoxyphenyl)-N-[1-(2-phenylethyl)-4-piperidinyl]-2-furancarboxamide

Synonyms: *m*-methoxy Fu-F, *meta*-methoxy Fu-F, *m*-methoxy Furanyl fentanyl

MF: C₂₅H₂₈N₂O₃

FW: 404.5

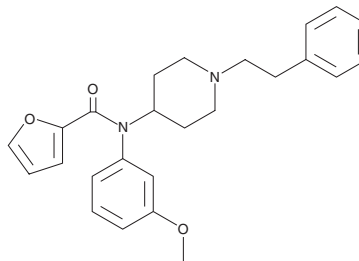
Purity: ≥98%

UV/Vis.: λ_{max}: 260 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥5 years



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

Description

meta-methoxy Furanyl fentanyl (Item No. 26050) is an analytical reference standard that is structurally similar to known opioids. *meta*-methoxy Furanyl fentanyl is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

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, 07/19/2023

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