

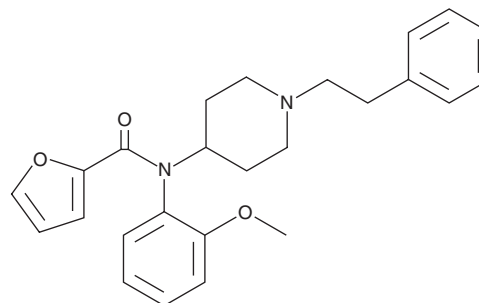
PRODUCT INFORMATION



ortho-methoxy Furanyl fentanyl

Item No. 22805

CAS Registry No.: 101343-50-4
Formal Name: N-(2-methoxyphenyl)-N-[1-(2-phenylethyl)-4-piperidinyl]-2-furancarboxamide
Synonyms: *o*-methoxy Fu-F, *ortho*-methoxy Fu-F, *o*-methoxy Furanyl fentanyl
MF: C₂₅H₂₈N₂O₃
FW: 404.5
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥5 years



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

Description

ortho-methoxy Furanyl fentanyl (Item No. 22805) is an analytical reference material that is structurally similar to known opioids. *ortho*-methoxy Furanyl fentanyl is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

This product is a qualified Reference Material (RM) that has been manufactured and tested to meet ISO/IEC 17025 and ISO Guide 34 guidelines. These materials are tested using validated analytical methods on qualified instrumentation to ensure traceability of measurements. All traceable RMs may be distinguished by their CofAs and can be downloaded using the batch number located on the product label. For a representative CofA please contact our technical support.

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, 12/03/2020

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