

CERTIFICATE of ANALYSIS



meta-methoxy Furanyl fentanyl

N-(3-methoxyphenyl)-N-(1-phenethylpiperidin-4-yl)furan-2-carboxamide

Item No. 26050 • Batch No. 0539202

Purity Specification: $\geq 98\%$

Molecular Formula : C₂₅H₂₈N₂O₃

CAS Number: (unspecified)

Formula Weight : 404.5015

Expiry date: 03DEC2022

Tests	Results
HPLC	Purity: 100.0 %
IR	Conforms
Mass spec	MH ⁺ : 405.2
Melting Point	133 - 136 °C
TLC	Purity: 100 %
UV	λ max: 260 nm
NMR	Conforms

Reviewed and approved by: Libby Perry

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

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