

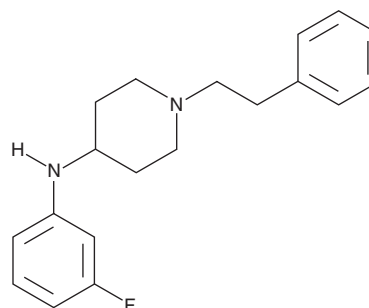
PRODUCT INFORMATION



Despropionyl *meta*-Fluorofentanyl

Item No. 20125

CAS Registry No.: 416881-38-4
Formal Name: N-(3-fluorophenyl)-1-(2-phenylethyl)-4-piperidinamine
Synonyms: *m*-fluoro 4-ANPP, *meta*-fluoro 4-ANPP, Despropionyl *m*-FF, Despropionyl *meta*-FF, Despropionyl *m*-Fluorofentanyl
MF: C₁₉H₂₃FN₂
FW: 298.4
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥6 years



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

Description

Despropionyl *meta*-fluorofentanyl (Item No. 20125) is an analytical reference material that is structurally similar to known opioids. Despropionyl *meta*-fluorofentanyl is a precursor to *meta*-fluorofentanyl and *meta*-fluorobutyryl fentanyl and a potential impurity in illicit preparations of this class of compounds. It is also a potential metabolite of *meta*-fluorofentanyl and *meta*-fluorobutyryl fentanyl based on the published metabolism of fentanyl.¹⁻⁵ At the time despropionyl *meta*-fluorofentanyl (Item No. 20125) was made available for purchase, specific metabolism data had not been published. Contact us if updated information on this molecule is now available. This product is intended for research and forensic applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards for reference materials.

References

1. Pease, J.P., LePine, A.J., and Smith, C.M. Methods for preparing fentanyl and fentanyl intermediates. *Patent Application Publication US20130281702A1*, (2013).
2. Watanabe, S., Vikingsson, S., Roman, M., *et al.* *In vitro* and *in vivo* metabolite identification studies for the new synthetic opioids acetylfentanyl, acrylfentanyl, furanylfentanyl, and 4-fluoro-isobutyrylfentanyl. *AAPS J.* **19(4)**, 1102-1122 (2017).
3. Labroo, R.B., Paine, M.F., Thummel, K.E., *et al.* Fentanyl metabolism by human hepatic and intestinal cytochrome P450 3A4: Implications for interindividual variability in disposition, efficacy, and drug interactions. *Drug Metab. Dispos.* **25(9)**, 1072-1079 (1997).
4. Melent'ev, A.B., Kataev, S.S., and Dvorskaya, O.N. Identification and analytical properties of acetyl fentanyl metabolites. *J. Anal. Chem.* **70(2)**, 216-224 (2015).
5. Steuer, A.E., Williner, E., Staeheli, S.N., *et al.* Studies on the metabolism of the fentanyl-derived designer drug butyrylfentanyl in human *in vitro* liver preparations and authentic human samples using liquid chromatography-high resolution mass spectrometry (LC-HRMS). *Drug Test Anal.* **9(7)**, 1085-1092 (2017).

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, 12/02/2020

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