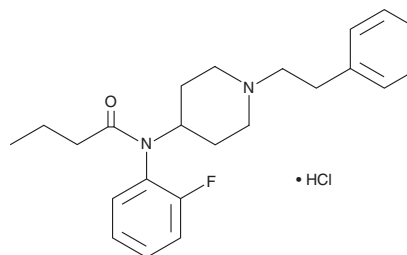


PRODUCT INFORMATION

ortho-Fluorobutyryl fentanyl (hydrochloride) (exempt preparation) Item No. 39596

CAS Registry No.: 2306823-44-7
Formal Name: N-(2-fluorophenyl)-N-[1-(2-phenylethyl)-4-piperidiny]-butanamide, monohydrochloride
Synonyms: 2-FBF, o-FBF, 2-Fluorobutyryl fentanyl, o-Fluorobutyryl fentanyl
MF: C₂₃H₂₉FN₂O • HCl
FW: 405.0
Purity: ≥98%
Supplied as: A 100 µg/ml solution in methanol
Storage: -20°C
Stability: ≥3 years



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

Description

ortho-Fluorobutyryl fentanyl (hydrochloride) (exempt preparation) (Item No. 39596) is an analytical reference standard categorized as an opioid.¹ *ortho*-Fluorobutyryl fentanyl is regulated as a Schedule I compound in the United States. *ortho*-Fluorobutyryl fentanyl (hydrochloride) (exempt preparation) (Item No. 39596) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

Reference

- Hassanien, S.H., Bassman, J.R., Perrien Naccarato, C.M., *et al.* *In vitro* pharmacology of fentanyl analogs at the human mu opioid receptor and their spectroscopic analysis. *Drug Test. Anal.* **12**(8), 1212-1221 (2020).

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, 08/22/2022

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