

# PRODUCT INFORMATION



## FIBF (hydrochloride)

Item No. 19313

**CAS Registry No.:** 2309383-06-8  
**Formal Name:** N-(4-fluorophenyl)-2-methyl-N-[1-(2-phenylethyl)-4-piperidinyl]-propanamide, monohydrochloride

**Synonyms:** *p*-FIBF, *para*-FIBF, *p*-Fluoroisobutyryl fentanyl, *para*-Fluoroisobutyryl fentanyl, *p*-fluoro iBF, *para*-fluoro iBF, *p*-fluoro Isobutyryl fentanyl, *para*-fluoro Isobutyryl fentanyl

**MF:** C<sub>23</sub>H<sub>29</sub>FN<sub>2</sub>O • HCl

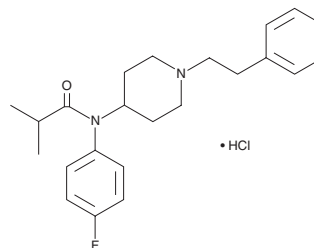
**FW:** 405.0

**Purity:** ≥98%

**Supplied as:** A neat solid

**Storage:** -20°C

**Stability:** ≥4 years



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

### Description

FIBF (hydrochloride) (Item No. 19313) is an analytical reference material categorized as an opioid.<sup>1</sup> FIBF has been associated with fatal multidrug overdoses, as well as overdoses where it was the sole intoxicant responsible for death.<sup>2</sup> FIBF is regulated as a Schedule I compound in the United States. 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. 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).
2. Pardi, J., Toriello, A., and Cooper, G. Evaluation of 4-fluoroisobutyrylfentanyl in blood samples from 247 authentic cases submitted to the New York City Office of Chief Medical Examiner in 2017–2018. *Forensic Toxicol.* **38**, 340-351 (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, 05/04/2021

#### 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