

PRODUCT INFORMATION

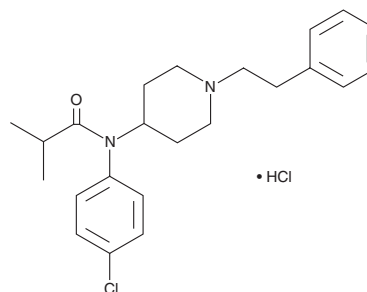


para-Chloroisobutyryl fentanyl (hydrochloride)

Item No. 20018

CAS Registry No.: 2306827-85-8
Formal Name: N-(4-chlorophenyl)-2-methyl-N-[1-(2-phenylethyl)-4-piperidinyl]-propanamide, monohydrochloride
Synonyms: 4-Chloride Isobutyryl fentanyl, 4-chloro iBF, 4-Chloroisobutyryl fentanyl, *p*-Chloride Isobutyryl fentanyl, *p*-chloro iBF, *p*-Chloroisobutyryl fentanyl, *para*-Chloride Isobutyryl fentanyl, *para*-chloro iBF

MF: C₂₃H₂₉ClN₂O • HCl
FW: 421.4
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

para-Chloroisobutyryl fentanyl (hydrochloride) (Item No. 20018) is an analytical reference material categorized as an opioid.¹ *para*-Chloroisobutyryl fentanyl 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.

Reference

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).

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, 08/19/2022

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