

PRODUCT INFORMATION



Butyryl fentanyl (hydrochloride) (exempt preparation)

Item No. 44604

CAS Registry No.: 1443-52-3

Formal Name: N-phenyl-N-[1-(2-phenylethyl)-4-piperidyl]-butanamide, monohydrochloride

Synonyms: Butyr fentanyl, NIH 10486

MF: $C_{23}H_{30}N_2O \cdot HCl$

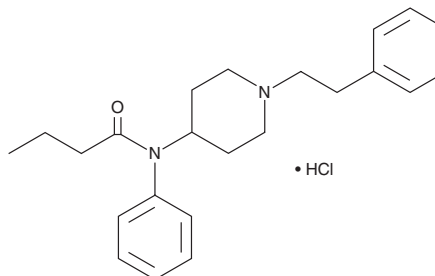
FW: 387.0

Purity: $\geq 98\%$

Supplied as: A 100 $\mu g/ml$ solution in methanol

Storage: $-20^{\circ}C$

Stability: ≥ 3 years



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

Description

Butyryl fentanyl (hydrochloride) (exempt preparation) (Item No. 44604) is an analytical reference standard categorized as an opioid.¹ Butyryl fentanyl is also a potential impurity found in preparations of fentanyl.² Butyryl fentanyl is regulated as a Schedule I compound in the United States. Butyryl fentanyl (hydrochloride) (exempt preparation) (Item No. 44604) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

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. Garg, A., Solas, D.W., Takahashi, L.H., *et al.* Forced degradation of fentanyl: Identification and analysis of impurities and degradants. *J. Pharm. Biomed. Anal.* **53**(3), 325-334 (2010).

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, 09/25/2025

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