

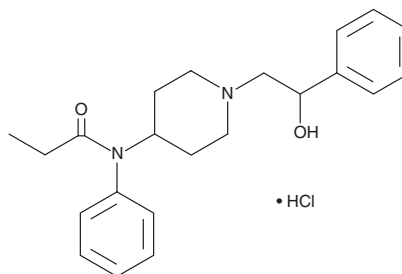
PRODUCT INFORMATION



β -hydroxy Fentanyl (hydrochloride)

Item No. 20789

CAS Registry No.: 1473-95-6
Formal Name: N-[1-(2-hydroxy-2-phenylethyl)-4-piperidiny]-N-phenylpropanamide, monohydrochloride
Synonyms: NIH 10506, R-4333
MF: C₂₂H₂₈N₂O₂ • HCl
FW: 388.9
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years



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

Description

β -hydroxy Fentanyl (hydrochloride) (Item No. 20789) is an analytical reference standard that is structurally categorized as an opioid. It is more potent than fentanyl (Item Nos. ISO60197 | 14719) at the wild-type μ -opioid receptor *in vitro* (EC_{50} s = 38.1 and 64.9 nM, respectively).¹ This compound differs from the hydroxypropionyl compound that has been described as a metabolite of fentanyl.² The physiological and toxicological properties of β -hydroxy fentanyl are not known. This product is intended for research and forensic applications.

References

1. Pil, J. and Tytgat, J. Serine 329 of the μ -opioid receptor interacts differently with agonists. *J. Pharm. Exp. Ther.* **304**(3), 924-930 (2003).
2. 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).

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, 06/22/2023

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