

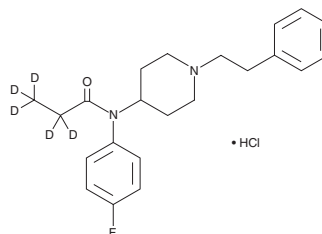
PRODUCT INFORMATION



para-Fluorofentanyl-d₅ (hydrochloride)

Item No. 19597

CAS Registry No.: 2747915-22-4
Formal Name: N-(4-fluorophenyl)-N-(1-phenethylpiperidin-4-yl)propanamide-d₅, monohydrochloride
Synonyms: *p*-FF-d₅, *p*-Fluorofentanyl-d₅
MF: C₂₂H₂₂D₅FN₂O • HCl
FW: 396.0
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥3 years



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

Description

para-Fluorofentanyl-d₅ (hydrochloride) (Item No. 19597) is intended for use as an internal standard for the quantification of *para*-fluorofentanyl (Item Nos. 19735 | 15496 | 39713) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

para-Fluorofentanyl is categorized as an opioid.¹ *para*-Fluorofentanyl has antinociceptive activity in rats.² It has been found in samples seized by law enforcement.³ *para*-Fluorofentanyl-d₅ 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. Ulens, C., Van Boven, M., Daenens, P., *et al.* Interaction of *p*-fluorofentanyl on cloned human opioid receptors and exploration of the role of Trp-318 and His-319 in μ -opioid receptor selectivity. *J. Pharmacol. Exp. Ther.* **294**(3), 1024-1033 (2000).
2. Casy, A.F. and Huckstep, M.R. Structure-activity studies of fentanyl. *J. Pharm. Pharmacol.* **40**(9), 605-608 (1988).
3. de Boer, D., Goemans, W.P., Ghezavat, V.R., *et al.* Seizure of illicitly produced *para*-fluorofentanyl: Quantitative analysis of the content of capsules and tablets. *J. Pharm. Biomed. Anal.* **31**(3), 557-562 (2003).

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, 04/17/2024

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