

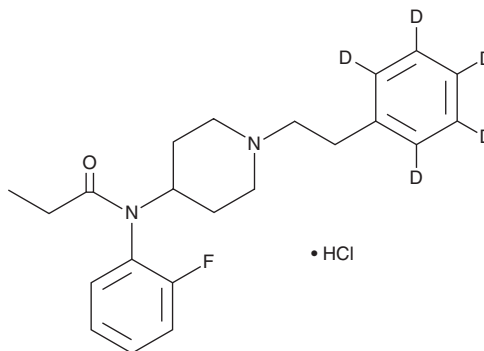
PRODUCT INFORMATION



ortho-Fluorofentanyl-d₅ (hydrochloride)

Item No. 42428

Formal Name: N-(2-fluorophenyl)-N-[1-(2-phenylethyl-d₅)-4-piperidinyl]-propanamide, monohydrochloride
Synonyms: *o*-FF-d₅, *ortho*-FF-d₅, *o*-Fluorofentanyl-d₅
MF: C₂₂H₂₂D₅FN₂O • HCl
FW: 396.0
Chemical Purity: ≥98% (*ortho*-Fluorofentanyl)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Description

ortho-Fluorofentanyl-d₅ (hydrochloride) (Item No. 42428) is an analytical reference material intended for use as an internal standard for the quantification of *ortho*-fluorofentanyl (Item Nos. 19425 | 38651) 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).

ortho-Fluorofentanyl is categorized as an opioid.¹ It has been associated with fatal overdoses.² *ortho*-Fluorofentanyl-d₅ is regulated as a Schedule I compound in the United States. 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. Helland, A., Brede, W.R., Michelsen, L.S., *et al.* Two hospitalizations and one death after exposure to *ortho*-fluorofentanyl. *J. Anal. Toxicol.* **41**(8), 708-709 (2017).

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, 02/12/2024

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