

PRODUCT INFORMATION



Ocfentanil-d₅ (hydrochloride)

Item No. 19599

CAS Registry No.: 2747917-28-6
Formal Name: N-(2-fluorophenyl)-2-methoxy-N-[1-(2-phenylethyl-d₅)-4-piperidiny]-acetamide, monohydrochloride

Synonyms: o-fluoro MAF-d₅, ortho-fluoro MAF-d₅, o-fluoro Methoxyacetyl fentanyl-d₅, ortho-fluoro Methoxyacetyl fentanyl-d₅

MF: C₂₂H₂₂D₅FN₂O₂ • HCl

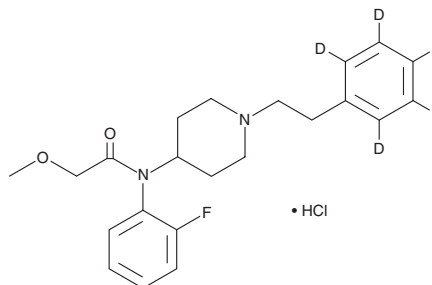
FW: 412.0

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

Ocfentanil-d₅ (hydrochloride) (Item No. 19599) is an analytical reference material intended for use as an internal standard for the quantification of ocfentanil (Item No. 18583) 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).

Ocfentanil is categorized as an opioid.¹ It induces generalized responses in mice trained to discriminate oxycodone from saline.² Ocfentanil-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.

References

1. Vasudevan, L., Vandeputte, M.M., Deventer, M., *et al.* Assessment of structure-activity relationships and biased agonism at the mu opioid receptor of novel synthetic opioids using a novel, stable bio-assay platform. *Biochem. Pharmacol.* **177**, 113910 (2020).
2. Walentiny, D.M., Moisa, L.T., and Beardsley, P.M. Oxycodone-like discriminative stimulus effects of fentanyl-related emerging drugs of abuse in mice. *Neuropharmacology* **150**, 210-216 (2019).

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, 08/18/2022

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