

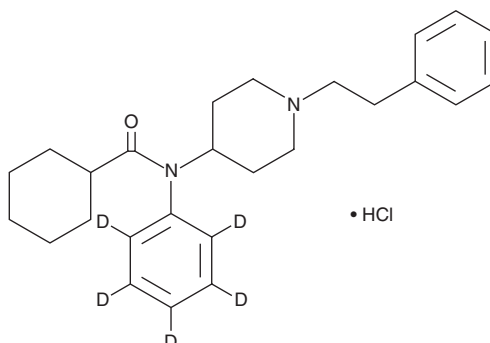
PRODUCT INFORMATION



Cyclohexyl fentanyl-d₅ (hydrochloride) (exempt preparation)

Item No. 35875

CAS Registry No.: 2747915-25-7
Formal Name: N-(phenyl-d₅)-N-[1-(2-phenylethyl)-4-piperidinyl]-cyclohexanecarboxamide, monohydrochloride
MF: C₂₆H₂₉D₅N₂O • HCl
FW: 432.1
Chemical Purity: ≥98% (Cyclohexyl fentanyl-d₅ (hydrochloride))
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀
Supplied as: A 100 µg/ml solution in methanol
Storage: -20°C
Stability: ≥3 years



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

Description

Cyclohexyl fentanyl-d₅ (hydrochloride) (exempt preparation) (Item No. 35875) is intended for use as an internal standard for the quantification of cyclohexyl fentanyl (Item No. 22390) 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).

Cyclohexyl fentanyl is categorized as an opioid.^{1,2} Cyclohexyl fentanyl-d₅ is regulated as a Schedule I compound in the United States. Cyclohexyl fentanyl-d₅ (hydrochloride) (exempt preparation) (Item No. 35875) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

References

1. Åstrand, A., Vikingsson, S., Jakobsen, I., *et al.* Activation of the µ-opioid receptor by alicyclic fentanyls: Changes from high potency full agonists to low potency partial agonists with increasing alicyclic substructure. *Drug Test Anal.* **13**(1), 169-174 (2020).
2. Eshleman, A.J., Nagarajan, S., Wolfrum, K.M., *et al.* Affinity, potency, efficacy, selectivity, and molecular modeling of substituted fentanyls at opioid receptors. *Biochem. Pharmacol.* **182**, 114293 (2020).

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, 12/19/2023

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