

PRODUCT INFORMATION

Furanyl fentanyl 3-furancarboxamide isomer (hydrochloride) (exempt preparation)

Item No. 37947

CAS Registry No.: 2306823-47-0

Formal Name: N-phenyl-N-[1-(2-phenylethyl)-4-piperidinyl]-3-furancarboxamide, monohydrochloride

Synonym: Fu-F 3-furancarboxamide isomer

MF: C₂₄H₂₆N₂O₂ • HCl

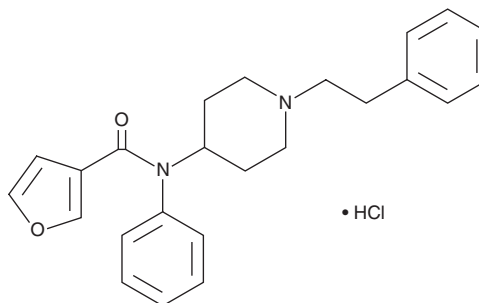
FW: 410.9

Purity: ≥98%

Supplied as: A 100 µg/ml solution in methanol

Storage: -20°C

Stability: ≥5 years



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

Description

Furanyl fentanyl 3-furancarboxamide isomer (hydrochloride) (exempt preparation) (Item No. 37947) is an analytical reference standard that is structurally similar to known opioids. Furanyl fentanyl 3-furancarboxamide isomer induces generalized responses in mice trained to discriminate oxycodone from saline.¹ Furanyl fentanyl 3-furancarboxamide isomer is regulated as a Schedule I compound in the United States. Furanyl fentanyl 3-furancarboxamide isomer (hydrochloride) (exempt preparation) (Item No. 37947) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

Reference

1. 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, 11/15/2022

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