

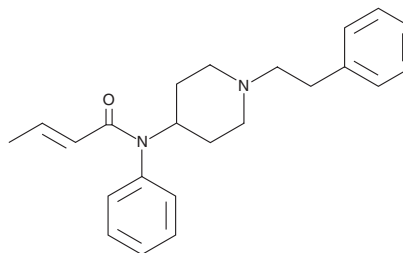
PRODUCT INFORMATION



Crotonyl fentanyl (exempt preparation)

Item No. 39889

CAS Registry No.: 760930-59-4
Formal Name: (2E)-N-phenyl-N-[1-(2-phenylethyl)-4-piperidinyl]-2-butenamide
MF: $C_{23}H_{28}N_2O$
FW: 348.5
Purity: $\geq 98\%$
Supplied as: A 100 $\mu\text{g/ml}$ 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

Crotonyl fentanyl (exempt preparation) (Item No. 39889) is an analytical reference standard categorized as an opioid.¹⁻³ Crotonyl fentanyl induces generalized responses in mice trained to discriminate oxycodone from saline.⁴ Crotonyl fentanyl is regulated as a Schedule I compound in the United States. Crotonyl fentanyl (exempt preparation) (Item No. 39889) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

References

1. 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).
2. 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).
3. Varshneya, N.B., Valentiny, D.M., Moisa, L.T., et al. Opioid-like antinociceptive and locomotor effects of emerging fentanyl-related substances. *Neuropharmacology* **151**, 171-179 (2019).
4. Valentiny, 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, 10/06/2023

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