

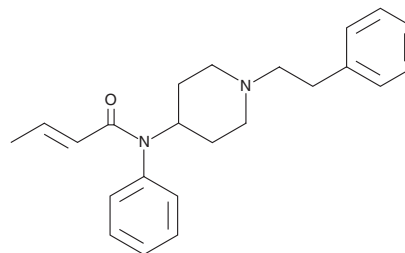
# PRODUCT INFORMATION



## Crotonyl fentanyl

Item No. 22801

**CAS Registry No.:** 760930-59-4  
**Formal Name:** (2E)-N-phenyl-N-[1-(2-phenylethyl)-4-piperidinyl]-2-butenamide  
**MF:** C<sub>23</sub>H<sub>28</sub>N<sub>2</sub>O  
**FW:** 348.5  
**Purity:** ≥98%  
**Supplied as:** A neat solid  
**Storage:** -20°C  
**Stability:** ≥6 years



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

### Description

Crotonyl fentanyl (Item No. 22801) is an analytical reference material categorized as an opioid.<sup>1-3</sup> Crotonyl fentanyl induces generalized responses in mice trained to discriminate oxycodone from saline.<sup>4</sup> Crotonyl fentanyl 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 for reference materials.

### 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, 10/08/2021

#### 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