

# PRODUCT INFORMATION



## Thiophene fentanyl (hydrochloride)

Item No. 22802

**CAS Registry No.:** 2306823-39-0  
**Formal Name:** N-phenyl-N-[1-(2-phenylethyl)-4-piperidinyl]-2-thiophenecarboxamide, monohydrochloride

**Synonyms:** Thiofuranyl fentanyl, 2-Thiofuranyl fentanyl, 2-Thiophenoyl fentanyl

**MF:** C<sub>24</sub>H<sub>26</sub>N<sub>2</sub>OS • HCl  
**FW:** 427.0

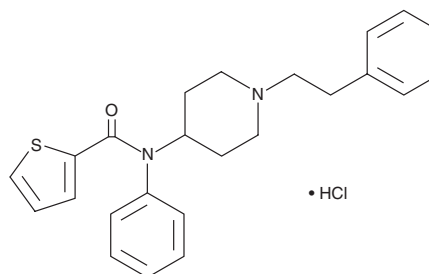
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 246 nm

**Supplied as:** A crystalline solid

**Storage:** -20°C

**Stability:** ≥5 years



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

### Description

Thiophene fentanyl (hydrochloride) (Item No. 22802) is an analytical reference standard categorized as an opioid.<sup>1,2</sup> Thiophene fentanyl induces generalized responses in rats trained to discriminate morphine from saline.<sup>3</sup> Thiophene fentanyl is regulated as a Schedule I compound in the United States. 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. Varshneya, N.B., Walentiny, D.M., Moisa, L.T., *et al.* Opioid-like antinociceptive and locomotor effects of emerging fentanyl-related substances. *Neuropharmacology* **151**, 171-179 (2019).
3. Eneanya, C., Walker, T., Tella, S., *et al.* Morphine drug discrimination: A behavioral model to assess novel synthetic opioid substances. *FASEB J.* **33(S1)**, 663.2-663.2 (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, 06/16/2023

#### 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