

# PRODUCT INFORMATION



## $\alpha$ -methyl Thiofentanyl (hydrochloride)

Item No. 20821

CAS Registry No.: 117332-94-2

Formal Name: N-[1-[1-methyl-2-(2-thienyl)ethyl]-4-piperidyl]-N-phenylpropanamide, monohydrochloride

Synonym: NIH 10538

MF:  $C_{21}H_{28}N_2OS \cdot HCl$

FW: 393.0

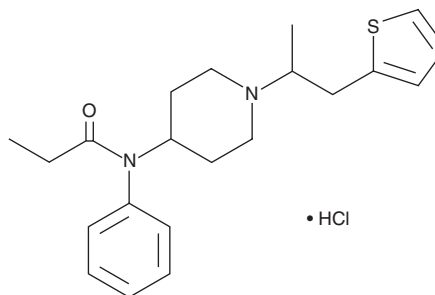
Purity:  $\geq 98\%$

UV/Vis.:  $\lambda_{max}$ : 229 nm

Supplied as: A crystalline solid

Storage:  $-20^{\circ}C$

Stability:  $\geq 5$  years



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

### Description

$\alpha$ -methyl Thiofentanyl (hydrochloride) (Item No. 20821) is an analytical reference standard that is structurally categorized as an opioid. It is reported to be a selective  $\mu$ -opioid receptor agonist, although with 100-fold increased potency compared to morphine (Item No. ISO60147).<sup>1</sup> This product is intended for research and forensic applications.

### Reference

1. J. Woods, F. Medzihradsky, C. Smith, *et al.* Evaluation of new compounds for opioid activity: 1987 Annual report. **81** 543-590.

#### 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, 06/22/2023

#### 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