

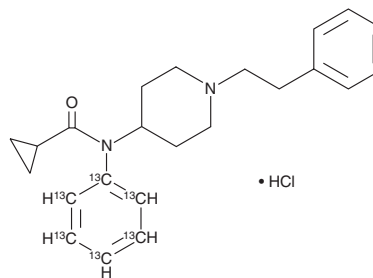
# PRODUCT INFORMATION



## Cyclopropyl fentanyl-<sup>13</sup>C<sub>6</sub> (hydrochloride)

Item No. 25563

**CAS Registry No.:** 2748409-79-0  
**Formal Name:** N-(1-phenethylpiperidin-4-yl)-N-(phenyl-<sup>13</sup>C<sub>6</sub>) cyclopropanecarboxamide, monohydrochloride  
**MF:** C<sub>17</sub>[<sup>13</sup>C]<sub>6</sub>H<sub>28</sub>N<sub>2</sub>O • HCl  
**FW:** 390.9  
**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

Cyclopropyl fentanyl-<sup>13</sup>C<sub>6</sub> (hydrochloride) (Item No. 25563) is intended for use as an internal standard for the quantification of cyclopropyl fentanyl (Item Nos. 23603 | 21739 | 24101) by GC- or LC-MS. Cyclopropyl fentanyl is categorized as an opioid.<sup>1</sup> Cyclopropyl fentanyl-<sup>13</sup>C<sub>6</sub> is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

### Reference

1. Hassanien, S.H., Bassman, J.R., Perrien Naccarato, C.M., *et al.* *In vitro* pharmacology of fentanyl analogs at the human mu opioid receptor and their spectroscopic analysis. *Drug Test. Anal.* **12(8)**, 1212-1221 (2020).

#### 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, 07/14/2023

#### 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