

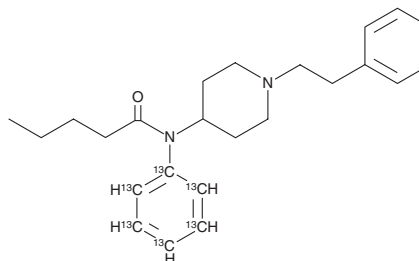
PRODUCT INFORMATION



Valeryl fentanyl-¹³C₆

Item No. 26220

CAS Registry No.: 2748625-09-2
Formal Name: N-phenyl-¹³C₆-N-[1-(2-phenylethyl)-4-piperidinyl]-pentanamide
Synonyms: Pentanoyl fentanyl-¹³C₆, Pentanyl fentanyl-¹³C₆
MF: C₁₈[¹³C]₆H₃₂N₂O
FW: 370.5
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

Valeryl fentanyl-¹³C₆ (Item No. 26220) is an analytical reference material intended for use as an internal standard for the quantification of valeryl fentanyl (Item Nos. 19798 | 18934) by GC- or LC-MS. Valeryl fentanyl is categorized as an opioid.¹ Valeryl fentanyl-¹³C₆ 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.

Reference

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).

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, 01/26/2022

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