

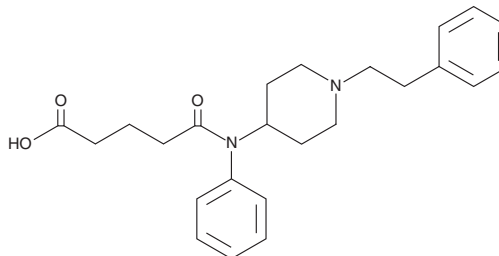
PRODUCT INFORMATION



Valeryl fentanyl carboxy metabolite

Item No. 20497

CAS Registry No.: 526191-19-5
Formal Name: 5-oxo-5-[phenyl[1-(2-phenylethyl)-4-piperidiny]amino]-pentanoic acid
Synonyms: Pentanoyl fentanyl carboxy metabolite, Pentanyl fentanyl carboxy metabolite
MF: C₂₄H₃₀N₂O₃
FW: 394.5
Purity: ≥98%
UV/Vis.: λ_{max}: 206 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Description

Valeryl fentanyl carboxy metabolite (Item No. 20497) is an analytical reference standard that is structurally similar to known opioids. It is a metabolite of valeryl fentanyl (Item Nos. 18934 | 19798).^{1,2} Valeryl fentanyl carboxy metabolite is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

References

1. Labroo, R.B., Paine, M.F., Thummel, K.E., *et al.* Fentanyl metabolism by human hepatic and intestinal cytochrome P450 3A4: Implications for interindividual variability in disposition, efficacy, and drug interactions. *Drug Metab. Dispos.* **25(9)**, 1072-1079 (1997).
2. Staeheli, S.N., Baumgartner, M.R., Gauthier, S., *et al.* Time-dependent postmortem redistribution of butyrfentanyl and its metabolites in blood and alternative matrices in a case of butyrfentanyl intoxication. *Forensic Sci. Int.* **266**, 170-177 (2016).

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, 05/09/2018

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