

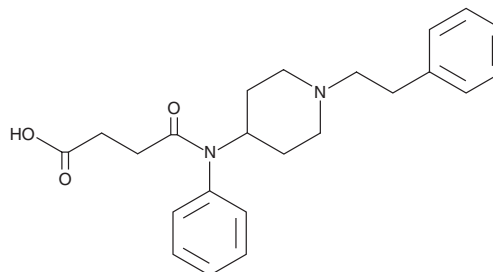
PRODUCT INFORMATION



Butyryl fentanyl carboxy metabolite

Item No. 20494

CAS Registry No.: 52994-23-7
Formal Name: 4-oxo-4-[phenyl[1-(2-phenylethyl)-4-piperidinyl]amino]-butanoic acid
Synonyms: Carboxybutyrfentanyl, Carboxyfentanyl
MF: C₂₃H₂₈N₂O₃
FW: 380.5
Purity: ≥98%
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

Butyryl fentanyl carboxy metabolite (Item No. 20494) is an analytical reference standard that is structurally similar to known opioids. It is a metabolite of butyryl fentanyl (Item Nos. 14728 | 19734).¹ Butyryl fentanyl carboxy metabolite is regulated as a Schedule I compound in the United States. This product is intended for research and forensic applications.

Reference

1. 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, 06/21/2023

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