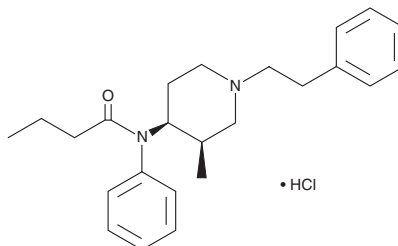


PRODUCT INFORMATION

(±)-cis-3-methyl Butyryl fentanyl (hydrochloride) (exempt preparation)
Item No. 39594

CAS Registry No.: 88641-20-7
Formal Name: cis-N-[3-methyl-1-(2-phenylethyl)-4-piperidiny]-N-phenyl-butanamide, monohydrochloride
MF: C₂₄H₃₂N₂O • HCl
FW: 401.0
Purity: ≥98%
Supplied as: A 100 µg/ml solution in methanol
Storage: -20°C
Stability: ≥3 years



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

Description

(±)-cis-3-methyl Butyryl fentanyl (hydrochloride) (exempt preparation) (Item No. 39594) is an analytical reference standard that is structurally similar to known opioids. (±)-cis-3-methyl Butyryl fentanyl is regulated as a Schedule I compound in the United States. (±)-cis-3-methyl Butyryl fentanyl (hydrochloride) (exempt preparation) (Item No. 39594) is provided as a DEA exempt preparation. This product is intended for research and forensic applications.

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, 08/21/2023

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