

PRODUCT INFORMATION



para-Fluorobutyryl fentanyl-¹³C₆ (hydrochloride)

Item No. 26216

Formal Name: N-(4-fluorophenyl)-N-[1-(2-(phenyl-¹³C₆)ethyl)-4-piperidinyl]-butanamide, monohydrochloride

Synonyms: *p*-FBF-¹³C₆, *para*-FBF-¹³C₆, *p*-Fluorobutyryl fentanyl-¹³C₆, 4-FBF-¹³C₆, 4-Fluorobutyryl fentanyl-¹³C₆

MF: C₁₇[¹³C]₆H₂₉FN₂O • HCl

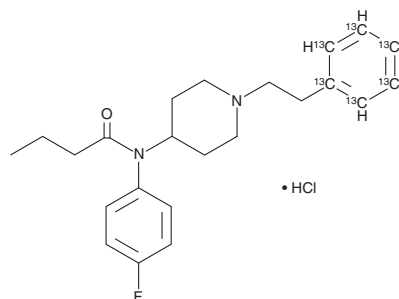
FW: 410.9

Purity: ≥95%

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

para-Fluorobutyryl fentanyl-¹³C₆ (hydrochloride) (Item No. 26216) is intended for use as an internal standard for the quantification of *para*-fluorobutyryl fentanyl (Item Nos. 19826 | 17049) by GC- or LC-MS. *para*-Fluorobutyryl fentanyl is structurally similar to known opioids. It has been associated with fatal overdoses.¹ *para*-Fluorobutyryl fentanyl-¹³C₆ (hydrochloride) 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. Danaceau, J.P., Wood, M., Ehlers, M.D., *et al.* Analysis of 17 fentanyl derivatives in plasma and blood by UPLC-MS/MS with interpretation of findings in surgical and postmortem casework. *Clin. Mass Spec.* **18**, 38-47 (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, 10/11/2021

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