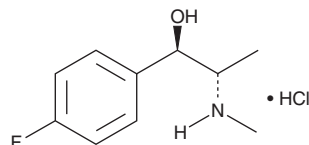


PRODUCT INFORMATION

4-Fluoromethcathinone metabolite (hydrochloride) (±)-Ephedrine stereochemistry

Item No. 14757

CAS Registry No.: 7799-72-6
Formal Name: 4-fluoro-α-[(1R)-1-(methylamino)ethyl]-benzenemethanol, monohydrochloride
Synonyms: Flephedrone metabolite, 4-FMC metabolite
MF: C₁₀H₁₄FNO • HCl
FW: 219.7
Purity: ≥98%
UV/Vis.: λ_{max}: 264, 270 nm
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

4-Fluoromethcathinone (4-FMC) (Item No. 10859) is a fluorinated cathinone derivative identified in several designer drugs that are marketed as plant feeders.^{1,2} This metabolite of 4-FMC features conversion of the β-keto group to β-hydroxy and is an enantiomeric mixture of the R,S and S,R orientations at carbons one and two, as in ephedrine. The physiological and toxicological properties of this compound have not been evaluated. This product is intended for forensic and research purposes.

References

1. Brandt, S.D., Sumnall, H.R., Measham, F., et al. Analyses of second-generation 'legal-highs' in the UK: Initial findings. *Drug Test. Anal.* **2**(8), 377-382 (2010).
2. Gibbons, S. and Zloh, M. An analysis of the 'legal high' mephedrone. *Bioorg. Med. Chem. Lett.* **20**(14), 4135-4139 (2010).

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, 06/22/2023

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