

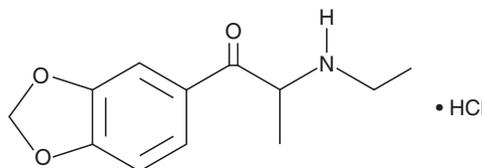
PRODUCT INFORMATION



Ethylone (hydrochloride) polymorph B

Item No. 9002749

CAS Registry No.: 1454266-19-3
Formal Name: 1-(1,3-benzodioxol-5-yl)-2-(ethylamino)-1-propanone, monohydrochloride
MF: C₁₂H₁₅NO₃ • HCl
FW: 257.7
Purity: ≥98%
UV/Vis.: λ_{max}: 234, 280, 317 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

Ethylone (hydrochloride) polymorph B (Item No. 9002749) is an analytical reference standard that is structurally categorized as a cathinone. Ethylone can exist in two conformations (polymorphs) at the C-C bond linking the side chain to the aromatic ring.¹ These polymorphs are identical by mass spectroscopy and nuclear magnetic resonance (NMR), but differ when analyzed by Fourier transform infrared spectroscopy (FTIR), FT-Raman, X-ray diffraction, and solid-state NMR spectroscopy.² The physiological and toxicological properties of this compound are not known. This product is intended for research and forensic applications.

References

1. Cameron, T.S., Grossert, J.S., Maheux, C.R., *et al.* Inversion twinning in a second polymorph of the hydrochloride salt of the recreational drug ethylone. *Acta. Crystallogr. C. Struct. Chem.* **71(Pt.4)**, 266-270 (2015).
2. Maheux, C.R., Alarcon, I.Q., Copeland, C.R., *et al.* Identification of polymorphism in ethylone hydrochloride: Synthesis and characterization. *Drug Test. Anal.* **092015**, (2015).

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, 07/17/2023

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