

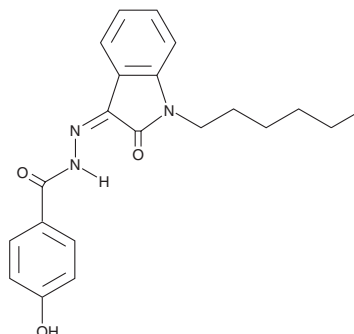
PRODUCT INFORMATION



MDA-19 4-hydroxybenzoyl metabolite

Item No. 36040

Formal Name: (Z)-N'-(1-hexyl-2-oxoindolin-3-ylidene)-4-hydroxybenzohydrazide
Synonym: BZO-HEXOXIZID 4-hydroxybenzoyl metabolite
MF: C₂₁H₂₃N₃O₃
FW: 365.4
Purity: ≥98%
UV/Vis.: λ_{max}: 350 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥5 years



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

Description

MDA-19 4-hydroxybenzoyl metabolite (Item No. 36040) is an analytical reference standard that is structurally similar to known synthetic cannabinoids. MDA-19 4-hydroxybenzoyl metabolite is a potential metabolite of MDA-19 (BZO-HEXOXIZID; Item No. 10563) based on the published metabolism of 5-fluoro CUMYL-PINACA and CUMYL-PeGACLONE (Item No. 22696).^{1,2} At the time MDA-19 4-hydroxybenzoyl metabolite (Item No. 36040) was made available for purchase, specific metabolism data had not been published. Contact us if updated information on this molecule is now available. This product is intended for research and forensic applications.

References

1. Staeheli, S.N., Poetzsch, M., Veloso, V.P., *et al.* In vitro metabolism of the synthetic cannabinoids CUMYL-PINACA, 5F-CUMYL-PINACA, CUMYL-4CN-BINACA, 5F-CUMYL-P7AICA and CUMYL-4CN-B7AICA. *Drug Test Anal.* **10**(1), 148-157 (2018).
2. Mogler, L., Wilde, M., Huppertz, L.M., *et al.* Phase I metabolism of the recently emerged synthetic cannabinoid CUMYL-PEGACLONE and detection in human urine samples. *Drug Test Anal.* **10**(5), 886-891 (2018).

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, 05/22/2024

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