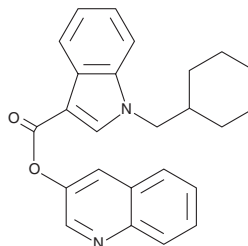


PRODUCT INFORMATION



BB-22 3-hydroxyquinoline isomer Item No. 14505

CAS Registry No.: 2365471-57-2
Formal Name: 1-(cyclohexylmethyl)-1H-indole-3-carboxylic acid, 3-quinolinyl ester
MF: C₂₅H₂₄N₂O₂
FW: 384.5
Purity: ≥98%
UV/Vis.: λ_{max}: 233, 294 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

BB-22 (Item No. 14099) is an analog of the potent synthetic cannabinoid, JWH 018 (Item No. 10900) that is structurally similar to PB-22 (Item No. 14096) and its derivative, 5-fluoro PB-22 (Item No. 14095).^{1,2} Both BB-22 and PB-22 are distinctive in that an 8-hydroxyquinoline replaces the naphthalene group of JWH 018. BB-22 3-hydroxyquinoline isomer differs from BB-22 structurally by having the hydroxyquinoline group attached at the three carbon instead of through the eight position of the ring. The physiological and toxicological properties of this compound are not known. This product is intended for forensic and research applications.

References

1. Sobolevsky, T., Prasolov, I., and Rodchenkov, G. Detection of JWH-018 metabolites in smoking mixture post-administration urine. *Forensic Sci. Int.* **200(1-3)**, 141-147 (2010).
2. Moran, C.L., Le, V.H., Chimalakonda, K.C., *et al.* Quantitative measurement of JWH-018 and JWH-073 metabolites excreted in human urine. *Anal. Chem.* **83(11)**, 4228-4236 (2011).

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, 06/19/2023

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