

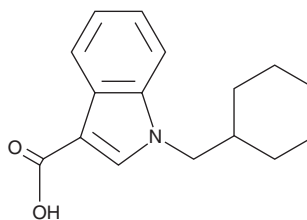
# PRODUCT INFORMATION



## BB-22 3-carboxyindole metabolite

Item No. 14378

**CAS Registry No.:** 858515-71-6  
**Formal Name:** 1-(cyclohexylmethyl)-1H-indole-3-carboxylic acid  
**MF:** C<sub>16</sub>H<sub>19</sub>NO<sub>2</sub>  
**FW:** 257.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 217, 289 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 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).<sup>1,2</sup> Both BB-22 and PB-22 are distinctive in that an 8-hydroxyquinoline replaces the naphthalene group of JWH 018. BB-22 3-carboxyindole metabolite is expected to be a metabolite of BB-22 that would be detectable both in serum and urine. 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**, 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, 07/13/2017

#### 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