

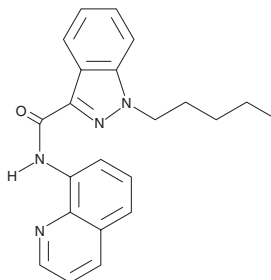
PRODUCT INFORMATION



THJ

Item No. 14889

CAS Registry No.: 2180931-94-4
Formal Name: 1-pentyl-N-(quinolin-8-yl)-1H-indazole-3-carboxamide
MF: C₂₂H₂₂N₄O
FW: 358.4
Purity: ≥98%
UV/Vis.: λ_{max}: 209, 245, 325 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

THJ 018 (Item No. 11962) is an analog of JWH 018 (Item No. 10900) where the core indole structure is substituted with an indazole base. Though no biological activity has been reported for THJ 018, structurally similar analogs bind to the brain central cannabinoid (CB₁) receptors and display agonist properties in functional assays, suggesting that it would have the same *in vivo* effects as Δ⁹-THC (Item No. 12068) and various synthetic CBs.¹ THJ is a derivative of THJ 018 featuring an 8-aminoquinoline replacing the naphthalene group. The physiological and toxicological properties of this compound have not been determined. This product is intended for research and forensic applications.

Reference

1. AKB48 (APINACA) and 5F-AKB48 (5F-APINACA), *Drug Enforcement Administration* (2013).

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