

PRODUCT INFORMATION



A-834735

Item No. 11163

CAS Registry No.: 895155-57-4

Formal Name: [1-[(tetrahydro-2H-pyran-4-yl)methyl]-1H-indol-3-yl](2,2,3,3-tetramethylcyclopropyl)-methanone

MF: C₂₂H₂₉NO₂

FW: 339.5

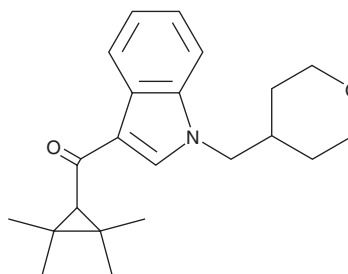
Purity: ≥98%

UV/Vis.: λ_{max}: 217, 246, 304 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

A-834735 is an indole-derived cannabinoid (CB) with a 3-tetramethylcyclopropylmethanone substituent. It acts as a full agonist at both the central CB₁ and peripheral CB₂ receptors in rat with K_i values of 4.6 and 0.31 nM, respectively, and EC₅₀ values of 12 and 0.21 nM, respectively.¹ A-834735 readily crosses the blood-brain barrier, dose-dependently reversing thermal hyperalgesia without adverse side effects in a rat neuropathic pain model.¹

Reference

1. Chin, C.-L., Tovcimak, A.E., Hradil, V.P., *et al.* Differential effects of cannabinoid receptor agonists on regional brain activity using pharmacological MRI. *Br. J. Pharmacol.* **153**(2), 367-379 (2008).

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, 07/07/2023

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