

PRODUCT INFORMATION



WIN 54,461

Item No. 9001204

CAS Registry No.: 166599-63-9

Formal Name: [6-bromo-2-methyl-1-[2-(4-morpholinyl)ethyl]-1H-indol-3-yl](4-methoxyphenyl)-methanone

Synonym: 6-Bromopravadoline

MF: C₂₃H₂₅BrN₂O₃

FW: 457.43

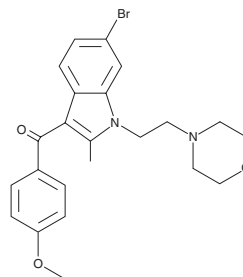
Purity: ≥98%

UV/Vis.: λ_{max}: 226, 283, 319 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

WIN 54,461 is an aminoalkylindole which acts as a cannabinoid (CB) mimetic, displacing WIN 55,212-2 from CB receptors on rat cerebellum membranes (IC₅₀ = 515 nM).¹ As it antagonizes the inhibition of electrically-induced contractions in isolated mouse vas deferens preparations by tetrahydrocannabinol or pravadoline, it may act as a receptor antagonist.¹ The physiological effects of WIN 54,461 *in vivo* have not been examined. This product is intended for research and forensic applications.

Reference

1. Eissenstat, M.A., Bell, M.R., D'Ambra, T.E., *et al.* Aminoalkylindoles: Structure-activity relationships of novel cannabinoid mimetics. *J. Med. Chem.* **38**(16), 3094-3105 (1995).

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