

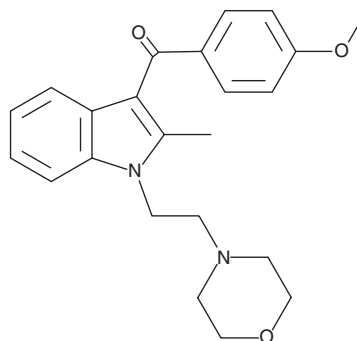
PRODUCT INFORMATION



Pravadoline

Item No. 10006973

CAS Registry No.: 92623-83-1
Formal Name: (4-methoxyphenyl)[2-methyl]-1-[2-(4-morpholinyl)ethyl]-1H-indol-3-yl]-methanone
Synonym: WIN 48,098
MF: C₂₃H₂₆N₂O₃
FW: 378.5
Purity: ≥98%
UV/Vis.: λ_{max}: 219, 272, 322 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

Pravadoline is a nonacidic, aminoalkylindole analgesic agent that inhibits cyclooxygenase.¹ Like NSAIDs, pravadoline inhibits prostaglandin synthesis (IC₅₀ = 5 μM in mouse brain microsomes) and displays antinociceptive activity (ED₅₀ = 26 mg/kg in an acetylcholine-induced writhing assay in mice).¹ In contrast to NSAIDs, pravadoline inhibits neuronally stimulated contractions in mouse vas deferens preparations (IC₅₀ = 0.45 μM), an effect similar to that which is produced by opioid analgesics, yet is not attenuated by the opioid receptor antagonist, naloxone.¹ Pravadoline does not produce gastrointestinal irritation following acute or chronic administration to rodents.¹

Reference

1. D'Ambra, T.E., Estep, K.G., Bell, M.R., *et al.* Conformationally restrained analogues of pravadoline: Nanomolar potent, enantioselective, (aminoalkyl)indole agonists of the cannabinoid receptor. *J. Med. Chem.* **35**(1), 124-135 (1992).

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, 07/17/2023

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