

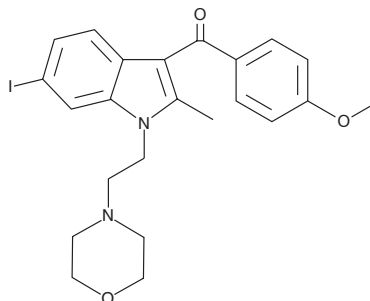
Product Information



AM630

Item No. 10006974

CAS Registry No.: 164178-33-0
Formal Name: [6-iodo-2-methyl-1-[2-(4-morpholinyl)ethyl]-1H-indol-3-yl] (4-methoxyphenyl)-methanone
Synonym: Iodopravadoline
MF: C₂₃H₂₅IN₂O₃
FW: 504.4
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that AM630 be stored as supplied at -20°C. It should be stable for at least two years.

AM630 is supplied as a crystalline solid. A stock solution may be made by dissolving the AM630 in an organic solvent purged with an inert gas. AM630 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of AM630 in these solvents is approximately 3 and 10 mg/ml, respectively.

AM630 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AM630 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. AM630 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Cannabinoids produce biochemical and pharmacological effects by interacting with the central cannabinoid (CB₁) and peripheral cannabinoid (CB₂) G protein-coupled receptors. AM630 is a selective CB₂ receptor antagonist that binds to CB₁ and CB₂ receptors with K_i values of 5.2 μM and 31.2 nM, respectively.¹ AM630 behaves as an inverse agonist at CB₂ receptors, attenuating the antinociceptive effects of a number of cannabinoids, and as a weak partial agonist at CB₁ receptors.¹⁻³

References

1. Ross, R.A., Brockie, H.C., Stevenson, L.A., *et al.* Agonist-inverse agonist characterization at CB₁ and CB₂ cannabinoid receptors of L759633, L759656 and AM630. *Br. J. Pharmacol.* **126**, 665-672 (1999).
2. Malan, T.P., Jr., Ibrahim, M.M., Deng, H., *et al.* CB₂ cannabinoid receptor-mediated peripheral antinociception. *Pain* **93**, 239-245 (2001).
3. Bisogno, T., Ortar, G., Petrosino, S., *et al.* Development of a potent inhibitor of 2-arachidonoylglycerol hydrolysis with antinociceptive activity *in vivo*. *Biochem. Biophys. Acta* **1791**, 53-60 (2009).

Related Products

AM2201-d₅ - Item No. 10706 • AM2201 - Item No. 10707 • AM251 - Item No. 71670 • Pravadoline - Item No. 10006973 • L-759,633 - Item No. 10009280 • AM1241 - Item No. 10010118 • GW 842166X - Item No. 10010372

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent under separate cover to the MSDS supervisor at your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 05/25/2011

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com