

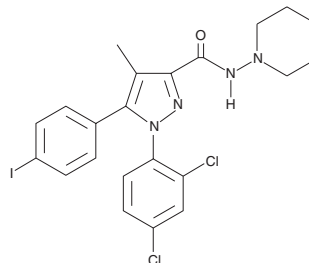
Product Information



AM251

Item No. 71670

CAS Registry No.: 183232-66-8
Formal Name: 1-(2,4-dichlorophenyl)-5-(4-iodophenyl)-4-methyl-N-1-piperidinyl-1H-pyrazole-3-carboxamide
MF: C₂₂H₂₁Cl₂IN₄O
FW: 555.2
Purity: ≥98%
Stability: ≥1 year at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that AM251 be stored as supplied at -20°C. It will be stable for at least one year.

AM251 is supplied as a crystalline solid. A stock solution may be made by dissolving the AM251 in an organic solvent purged with an inert gas. AM251 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of AM251 in these solvents is approximately 10 mg/ml. AM251 will be stable for at least six months in these solvents if stored at -20°C.

AM251 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AM251 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. AM251 has a solubility of approximately 14 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

SR 141716A is a biarylpyrazole compound which has been considered for many years as the prototypical cannabinoid receptor antagonist.¹ Some very close structural relatives of SR 141716A were recently prepared and tested for their ability to act as antagonists of the CB₁ receptor.² AM251 is one of these analogs, wherein the *p*-chloro group attached to the phenyl substituent at C-5 of the pyrazole ring is replaced with a *p*-iodo group. The resulting compound exhibits slightly better binding affinity for the CB₁ receptor with a K_i value of 7.5 nM compared to SR 141716A, which has a K_i value of 11.5 nM. However AM251 is about two-fold more selective for the CB₁ receptor as compared with SR 141716A.²

References

1. Rinaldi-Carmona, M., Barth, F., Héaulme, M., *et al.* SR141716A, a potent and selective antagonist of the brain cannabinoid receptor. *FEBS Lett.* **350**, 240-244 (1994).
2. Lan, R., Liu, Q., Fan, P., *et al.* Structure-activity relationships of pyrazole derivatives as cannabinoid receptor antagonists. *J. Med. Chem.* **42**, 769-776 (1999).

Related Products

AM2201-d₅ - Item No. 10706 • AM2201 - Item No. 10707 • Arachidonic Acid-(2-aminoethyl)-ester - Item No. 91050 • Pravadoline - Item No. 10006973

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

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 *via* email to 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