

# PRODUCT INFORMATION



## MJ15

Item No. 11956

**CAS Registry No.:** 944154-76-1  
**Formal Name:** 5-(4-chlorophenyl)-1-(2,4-dichlorophenyl)-4-methyl-N-(3-pyridinylmethyl)-1H-pyrazole-3-carboxamide

**MF:** C<sub>23</sub>H<sub>17</sub>Cl<sub>3</sub>N<sub>4</sub>O

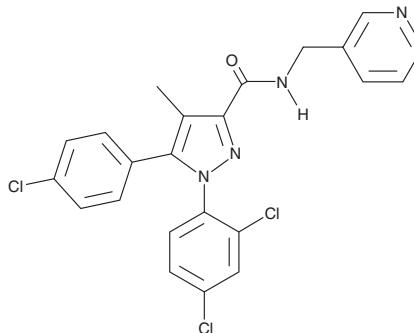
**FW:** 471.8

**Purity:** ≥98%

**Supplied as:** A crystalline solid

**Storage:** -20°C

**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

## Laboratory Procedures

MJ15 is supplied as a crystalline solid. A stock solution may be made by dissolving the MJ15 in the solvent of choice, which should be purged with an inert gas. MJ15 is soluble in the organic solvent DMSO. MJ15 is also slightly soluble in ethanol.

## Description

MJ15 is a potent antagonist of the cannabinoid (CB) receptor CB<sub>1</sub> (IC<sub>50</sub> = 118.9 pM; K<sub>i</sub> = 27.2 pM for rat CB<sub>1</sub>).<sup>1</sup> It is selective for CB<sub>1</sub> over CB<sub>2</sub>, exhibiting only 46% inhibition of CB<sub>2</sub> at a concentration of 10 μM. MJ15 reverses (+)-WIN 55,212-2 (Item No. 10009023) inhibition of forskolin-induced cAMP accumulation in a dose-dependent manner in CHO cells expressing human CB<sub>1</sub>. MJ15 (40 mg/kg) reduces food intake and body weight gain in a rat model of diet-induced obesity. It also reduces body weight gain and white adipose accumulation in a mouse model of early onset diet-induced obesity.

## Reference

1. Chen, W., Tang, H., Liu, H., *et al.* Novel selective antagonist of the cannabinoid CB<sub>1</sub> receptor, MJ15, with prominent anti-obesity effect in rodent models. *Eur. J. Pharmacol.* **637(1-3)**, 178-185 (2010).

### 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, 10/06/2022

## 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