

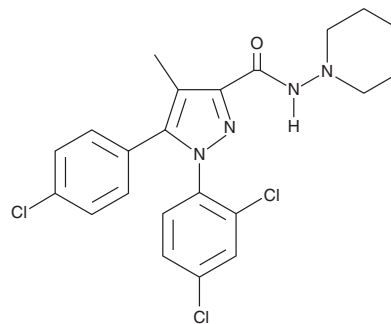
# PRODUCT INFORMATION



## Rimonabant

Item No. 9000484

**CAS Registry No.:** 168273-06-1  
**Formal Name:** 5-(4-chlorophenyl)-1-(2,4-dichlorophenyl)-4-methyl-N-1-piperidinyl-1H-pyrazole-3-carboxamide  
**Synonym:** SR141716  
**MF:** C<sub>22</sub>H<sub>21</sub>Cl<sub>3</sub>N<sub>4</sub>O  
**FW:** 463.8  
**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

Rimonabant is supplied as a crystalline solid. A stock solution may be made by dissolving the rimonabant in the solvent of choice. Rimonabant is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of rimonabant in ethanol is approximately 30 mg/ml and approximately 20 mg/ml in DMSO and DMF.

Rimonabant is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, rimonabant should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Rimonabant has a solubility of approximately 0.3 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Rimonabant is a cannabinoid 1 (CB<sub>1</sub>) receptor antagonist (K<sub>i</sub> = 5.6 nM).<sup>1</sup> It is selective for CB<sub>1</sub> over CB<sub>2</sub> receptors (K<sub>i</sub> = >1,000 nM), as well as a panel of 37 other receptors and channels (IC<sub>50</sub>s = >1,000 nM). Rimonabant (10 μM) inhibits phytohemagglutinin-induced proliferation of isolated human peripheral blood mononuclear cells (PBMCs).<sup>1</sup> Intraperitoneal administration of rimonabant prevents decreases in body temperature and increases in tail-flick latency induced by the CB<sub>1</sub> and CB<sub>2</sub> receptor agonist (+)-WIN 55,212-2 (Item No. 10009023) in mice (ED<sub>50</sub>s = 0.28 and 1.62 mg/kg, respectively) and oral administration reduces body weight in a mouse model of diet-induced obesity when administered at a dose of 10 mg/kg in the drinking water.<sup>2,3</sup> Rimonabant (10 mg/kg) decreases the percentage of time spent in the open arms of the elevated plus maze in mice, indicating anxiety-like activity.<sup>4</sup> Formulations containing rimonabant have previously been used in the treatment of obesity.

### References

1. Malfitano, A.M., Laezza, C., Pisanti, S., et al. Rimonabant (SR141716) exerts anti-proliferative and immunomodulatory effects in human peripheral blood mononuclear cells. *Brit. J Pharmacol.* **153**(5), 1003-1010 (2009).
2. Rinaldi-Carmona, M., Barth, F., Héaulme, M., et al. SR141716A, a potent and selective antagonist of the brain cannabinoid receptor. *FEBS Lett.* **350**(2-3), 240-244 (1994).
3. Lee, S.H., Seo, H.J., Lee, S.H., et al. Biarylpyrazolyl oxadiazole as potent, selective, orally bioavailable cannabinoid-1 receptor antagonists for the treatment of obesity. *J. med. Chem.* **51**, 7216-7233 (2008).
4. Bellocchio, L., Soria-Gómez, E., Quarta, C., et al. Activation of the sympathetic nervous system mediates hypophagic and anxiety-like effects of CB<sub>1</sub> receptor blockade. *Proc. Natl. Acad. Sci. USA* **110**(12), 4786-4791 (2013).

#### 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, 03/21/2023

#### 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