

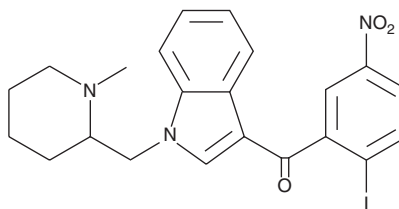
# Product Information



## AM1241

Item No. 10010118

**CAS Registry No.:** 44912-48-5  
**Formal Name:** (2-iodo-5-nitrophenyl)-(1-(1-methylpiperidin-2-ylmethyl)-1H-indol-3-yl)methanone  
**MF:** C<sub>22</sub>H<sub>22</sub>IN<sub>3</sub>O<sub>3</sub>  
**FW:** 503.3  
**Purity:** ≥97%  
**Stability:** ≥2 years at -20°C  
**Supplied as:** A crystalline solid



### Laboratory Procedures

For long term storage, we suggest that AM1241 be stored as supplied at -20°C. It will be stable for at least two years.

AM1241 is supplied as a crystalline solid. A stock solution may be made by dissolving the AM1241 in an organic solvent purged with an inert gas. AM1241 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of AM1241 in these solvents is approximately 5, 10, and 25 mg/ml, respectively.

AM1241 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AM1241 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. AM1241 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

AM1241 is a peripheral cannabinoid (CB<sub>2</sub>) receptor agonist with a K<sub>i</sub> value of 2 nM and greater than 100-fold selectivity over the central cannabinoid (CB<sub>1</sub>) receptor *in vitro*.<sup>1</sup> AM1241 produces antinociception to thermal stimuli in rat hindpaw. The antinociceptive actions of AM1241 were blocked by the CB<sub>2</sub> receptor-selective antagonist AM630 but not by the CB<sub>1</sub> receptor-selective antagonist AM251 (Catalog No. 71670).<sup>1</sup>

### Reference

1. Malan, T.P., Jr., Ibrahim, M.M., Deng, H., *et al.* CB<sub>2</sub> cannabinoid receptor-mediated peripheral antinociception. *Pain* **93**, 239-245 (2001).

### Related Products

AM2201-d<sub>5</sub> - Item No. 10706 • AM2201 - Item No. 10707 • AM251 - Item No. 71670 • CP 55,940 - Item No. 90084 • JWH 015 - Item No. 10009018 • WIN 55212-2 (mesylate) - Item No. 10009023 • (±)-SLV 319 - Item No. 10009226

**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

## 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