

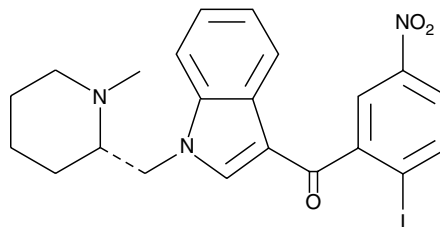
Product Information



(S)-AM1241

Item No. 10490

CAS Registry No.: 444912-53-2
Formal Name: (2-iodo-5-nitrophenyl)[1-[[[(2S)-1-methyl-2-piperidinyl]methyl]-1H-indol-3-yl]-methanone
Synonym: (-)-AM1241
MF: C₂₂H₂₂IN₃O₃
FW: 503.3
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

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

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

(S)-AM1241 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (S)-AM1241 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. (S)-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.

(S)-AM1241 binds the peripheral cannabinoid receptor (CB₂) from human, rat, and mouse with K_i values of 658, 893, and 577 nM, respectively. It exhibits weak binding for central cannabinoid receptor (CB₁) with a K_i value exceeding 10 μM.¹ (S)-1241 is an agonist for CB₂ for all three species, but shows greater activity at human CB₂ (EC₅₀ = 131 nM) than for either rat or mouse CB₂ (EC₅₀ = 785 and 2,000 nM, respectively).¹ Like the racemate AM1241 (Item No. 10010118), (S)-AM1241 produces antinociception in rats to thermal, but not mechanical pain.² The pain-reducing effect of (S)-AM1241 is blocked by the CB₂-specific inhibitor SR 144528, but not by either the CB₁-selective inhibitor rimonabant or the opioid receptor blocker naloxone.²

References

1. Bingham, B., Jones, P.G., Uveges, A.J., *et al.* Species-specific *in vitro* pharmacological effects of the cannabinoid receptor 2 (CB₂) selective ligand AM1241 and its resolved enantiomers. *Br. J. Pharmacol.* **151**, 1061-1070 (2007).
2. Rahn, E.J., Zvonok, A.M., Makriyannis, A., *et al.* Antinociceptive effects of racemic AM1241 and its chirally synthesized enantiomers: Lack of dependence upon opioid receptor activation. *AAPS J.* **12(2)**, 147-157 (2010).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10490

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, 12/02/2011

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