

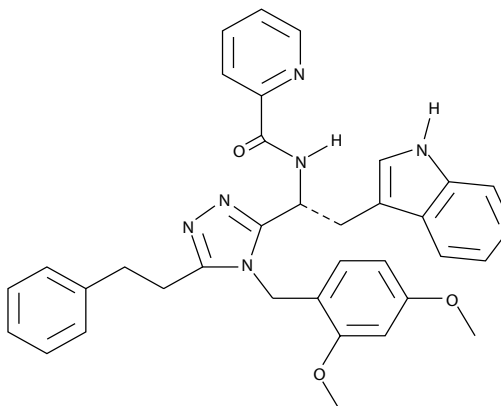
Product Information



JMV3002

Item No. 10012699

CAS Registry No.: 925239-03-8
Formal Name: N-[(1R)-1-[4-[(2,4-dimethoxyphenyl)methyl]-5-(2-phenylethyl)-4H-1,2,4-triazol-3-yl]-2-(1H-indol-3-yl)ethyl]-2-pyridinecarboxamide
MF: C₃₅H₃₄N₆O₃
FW: 586.7
Purity: ≥98%
Stability: ≥1 year at -20°C
Supplied as: A solution in methyl acetate
UV/Vis.: λ_{max}: 270 nm



Laboratory Procedures

For long term storage, we suggest that JMV3002 be stored as supplied at -20°C. It should be stable for at least one year. JMV3002 is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of JMV3002 in these solvents is approximately 20 mg/ml.

JMV3002 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of JMV3002 should be diluted with the aqueous buffer of choice. JMV3002 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Ghrelin is an endogenous ligand for the growth hormone secretagogue receptor that stimulates food intake and transduces signals to hypothalamic regulatory nuclei that control energy homeostasis. JMV3002 is a potent ghrelin receptor antagonist with an IC₅₀ value of 1.1 nM *in vitro*.¹ At 80 µg/kg, JMV3002 inhibits hexarelin-stimulated food intake by as much as 98% in rats.¹ JMV3002 alone does not elicit growth hormone release nor does it inhibit hexarelin-stimulated growth hormone secretion when tested in infant rats at a dose of 160 µg/kg.¹

Reference

1. Moulin, A., Demange, L., Ryan, J., *et al.* New trisubstituted 1,2,4-triazole derivatives as potent ghrelin receptor antagonists. 3. Synthesis and pharmacological *in vitro* and *in vivo* evaluations. *J. Med. Chem.* **51**, 689-693 (2008).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10012699

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY. NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

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 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, 01/07/2014

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