

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

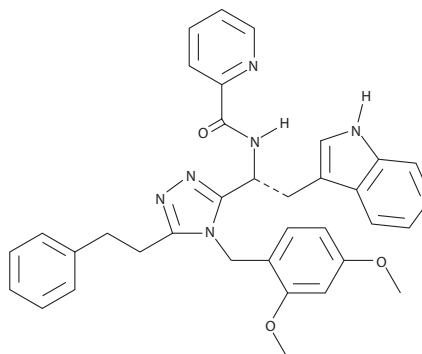
Purity: ≥98%

UV/Vis.: λ_{max}: 270 nm

Supplied as: A solution in methyl acetate

Storage: -20°C

Stability: ≥2 years



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

Laboratory Procedures

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 (DMF) purged with an inert gas can be used. The solubility of JMV3002 in ethanol and DMF is approximately 30 mg/ml and approximately 20 mg/ml in DMSO.

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.

Description

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).

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, 03/01/2024

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