

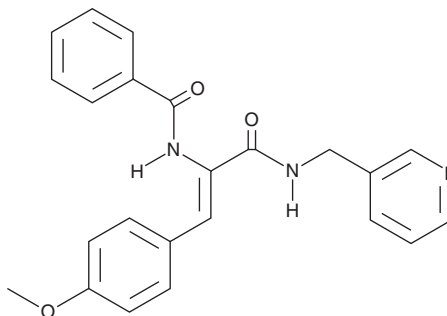
PRODUCT INFORMATION



IS20

Item No. 45324

CAS Registry No.: 1322013-92-2
Formal Name: N-[(1Z)-2-(4-methoxyphenyl)-1-[[[(3-pyridinylmethyl)amino]carbonyl]ethenyl]-benzamide
MF: C₂₃H₂₁N₃O₃
FW: 387.4
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

IS20 is supplied as a crystalline solid. A stock solution may be made by dissolving the IS20 in the solvent of choice, which should be purged with an inert gas. IS20 is sparingly soluble (1-10 mg/ml) in DMSO and ethanol.

Description

IS20 is a nonpeptide agonist of prokineticin receptor 1 (PKR₁) and a positive allosteric modulator (PAM) of PKR₂.¹ It increases intracellular calcium release and Akt phosphorylation in PKR₁-expressing CHO cells (EC₅₀ = 44 nM). It also enhances prokineticin-induced ERK phosphorylation in PKR₂-expressing CHO cells when used at a concentration of 1 nM. IS20 (0.5 mg/kg, i.p.) increases survival, reduces cardiac lesion formation, improves cardiac function, and increases neovasculation in a mouse model of myocardial infarction.

Reference

1. Wang, S., Zhao, L., Shi, X.-J., *et al.* Development of highly potent, selective, and cellular active triazolo [1,5-a]pyrimidine-based inhibitors targeting the DCN1-UBC12 protein-protein interaction. *J. Med Chem.* **62**(5), 2772-2797 (2019).

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, 03/30/2026

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