

PRODUCT INFORMATION

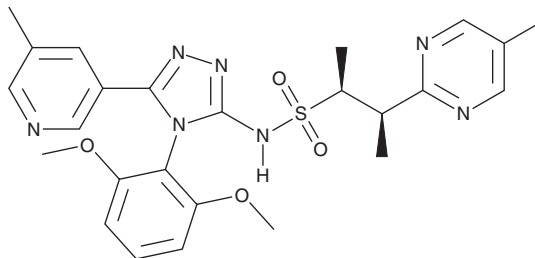


Azelaprag

Item No. 44644

CAS Registry No.: 2049980-18-7
Formal Name: (α S, β R)-N-[4-(2,6-dimethoxyphenyl)-5-(5-methyl-3-pyridinyl)-4H-1,2,4-triazol-3-yl]- α , β ,5-trimethyl-2-pyrimidineethanesulfonamide

Synonyms: AMG 986, BGE-105
MF: C₂₅H₂₉N₇O₄S
FW: 523.6
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

Azelaprag is supplied as a solid. A stock solution may be made by dissolving the azelaprag in the solvent of choice, which should be purged with an inert gas. Azelaprag is slightly soluble (0.1-1 mg/ml) in DMSO and a 1:1 solution of acetonitrile:water.

Description

Azelaprag is an agonist of the apelin receptor (APJ).¹ It inhibits forskolin-induced cAMP accumulation in CHO cells expressing human APJ (EC₅₀ = 0.23 nM). Azelaprag (1 mg/kg, i.v.) increases stroke volume and heart rate and decreases systemic vascular resistance, indicating increased cardiac reserve, in ZSF1 obese rats in a model of diastolic dysfunction driven by impaired metabolism, endothelial dysfunction, and hypertension.

Reference

1. Ason, B., Chen, Y., Guo, Q., *et al.* Cardiovascular response to small-molecule APJ activation. *JCI Insight* 5(8), e132898 (2020).

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, 02/05/2026

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