

PRODUCT INFORMATION



AZD 5462

Item No. 40956

CAS Registry No.: 2787501-83-9
Formal Name: cis-4-[2-fluoro-4-methoxy-5-[[[(1S,2R,3S,4R)-3-[[[(1-methylcyclobutyl)methyl]amino]carbonyl]bicyclo[2.2.1]hept-2-yl]amino]carbonyl]phenoxy]-1-methylcyclohexanecarboxylic acid

MF: C₃₀H₄₁FN₂O₆

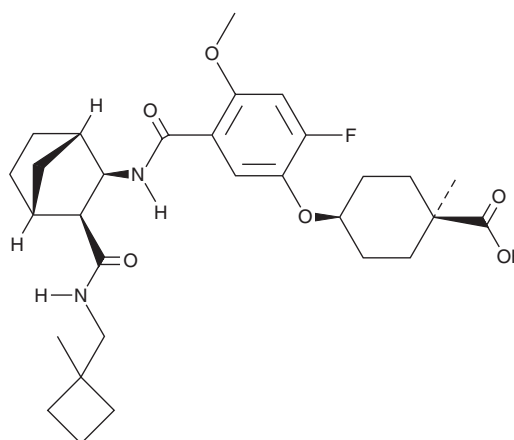
FW: 544.7

Purity: ≥98%

Supplied as: A 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

AZD 5462 is supplied as a solid. A stock solution may be made by dissolving the AZD 5462 in the solvent of choice, which should be purged with an inert gas. AZD 5462 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

AZD 5462 is an agonist of relaxin family peptide receptor 1 (RXFP1).¹ It induces cAMP mobilization in HEK293 cells expressing the human RXFP1 (EC₅₀ = 15.85 nM). AZD 5462 increases heart rate and mean arterial blood pressure (MABP) in rats. It increases left ventricular ejection fraction (LVEF) in a cynomolgus monkey model of heart failure induced by a high-fat diet and age when administered at doses of 1 or 10 mg/kg per day.

Reference

1. Granberg, K.L., Sakamaki, S., Larsson, N., *et al.* Discovery of clinical candidate AZD5462, a selective oral allosteric RXFP1 agonist for treatment of heart failure. *J. Med. Chem.* **67(6)**, 4419-4441 (2024).

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, 06/07/2024

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