

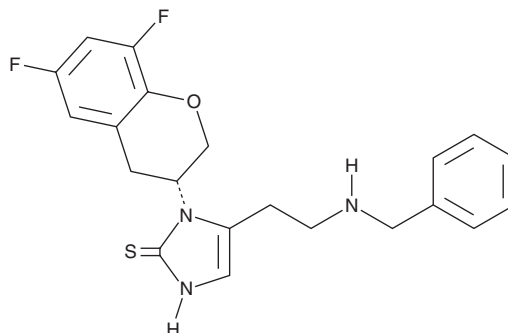
PRODUCT INFORMATION



Zamicastat

Item No. 42101

CAS Registry No.: 1080028-80-3
Formal Name: 1-[(3R)-6,8-difluoro-3,4-dihydro-2H-1-benzopyran-3-yl]-1,3-dihydro-5-[2-[(phenylmethyl)amino]ethyl]-2H-imidazole-2-thione
Synonym: BIA 5-1058
MF: C₂₁H₂₁F₂N₃OS
FW: 401.5
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

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

Description

Zamicastat is an inhibitor of dopamine β-hydroxylase (DβH).¹ It increases levels of norepinephrine and decreases levels of dopamine in the hearts of aged spontaneously hypertensive rats (SHRs) when administered at a dose of 30 mg/kg per day. Zamicastat decreases diastolic and systolic blood pressure in the same model. It increases serum levels of erythropoietin (EPO), decreases the heart-to-body weight ratio, and increases survival in a mouse model of hypertension using aged rats fed a high-salt diet.²

References

1. Igreja, B., Pires, N., Loureiro, A., *et al.* Cardiometabolic and inflammatory benefits of sympathetic down-regulation with zamicastat in aged spontaneously hypertensive rats. *ACS Pharmacol. Transl. Sci.* **2**(5), 353-360 (2019).
2. Igreja, B., Pires, N.M., Wright, L.C., *et al.* Effects of zamicastat treatment in a genetic model of salt-sensitive hypertension and heart failure. *Eur. J. Pharmacol.* **842**, 125-132 (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, 12/03/2025

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