

PRODUCT INFORMATION



AZD 8329

Item No. 45334

CAS Registry No.: 1048668-70-7

Formal Name: 4-[5-(1,1-dimethylethyl)-4-
[[tricyclo[3.3.1.1^{3,7}]dec-2-ylamino]
carbonyl]-1H-pyrazol-1-yl]-benzoic acid

MF: C₂₅H₃₁N₃O₃

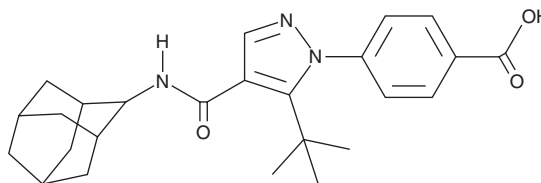
FW: 421.5

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

AZD 8329 is supplied as a solid. A stock solution may be made by dissolving the AZD 8329 in the solvent of choice, which should be purged with an inert gas. AZD 8329 is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

AZD 8329 is an inhibitor of 11β-hydroxysteroid dehydrogenase 1 (11β-HSD1; IC₅₀ = 9 nM for the human enzyme).¹ It is selective for 11β-HSD1 over 11β-HSD2 and 17β-HSD1 (IC₅₀s = >30 μM for both), and over 17β-HSD3 with less than 5% inhibition at 10 μM. AZD 8329 (50 mg/kg per day) prevents impairments in cardiac function in a minipig model of myocardial infarction.² Repeated oral administration of AZD 8329 over seven days induces tachyphylaxis in rats.³

References

1. Scott, J.S., de Schoolmeester, J., Kilgour, E., *et al.* Novel acidic 11β-hydroxysteroid dehydrogenase type 1 (11β-HSD1) inhibitor with reduced acylglucuronide liability: The discovery of 4-[4-(2-adamantylcarbamoyl)-5-tert-butyl-pyrazol-1-yl]benzoic acid (AZD8329). *J. Med. Chem.* **55**(22), 10136-10147 (2012).
2. Al Disi, S., Ascione, R., Khan, S., *et al.* Pharmacological inhibition of hydroxysteroid 11-beta dehydrogenase type 1 (11-βHSD1) after myocardial infarction preserves cardiac function in a translational mini-pig model. *Br. J. Pharmacol.* (2025).
3. Morentin Gutierrez, P., Gyte, A., de Schoolmeester, J., *et al.* Continuous inhibition of 11β-hydroxysteroid dehydrogenase type I in adipose tissue leads to tachyphylaxis in humans and rats but not in mice. *Br. J. Pharmacol.* **172**(20), 4806-4816 (2015).

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, 04/03/2026

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