

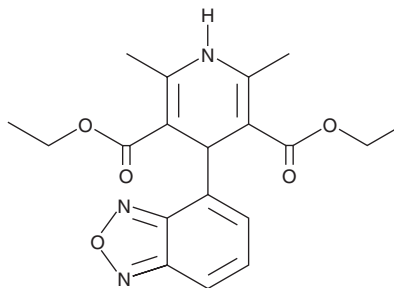
PRODUCT INFORMATION



Darodipine

Item No. 41667

CAS Registry No.: 72803-02-2
Formal Name: 4-(2,1,3-benzoxadiazol-4-yl)-1,4-dihydro-2,6-dimethyl-3,5-pyridinedicarboxylic acid, 3,5-diethyl ester
Synonym: PY 108-068
MF: C₁₉H₂₁N₃O₅
FW: 371.4
Purity: ≥95%
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

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

Description

Darodipine is a calcium channel inhibitor.¹ It selectively inhibits calcium-induced contractions of isolated rabbit aorta (pD₂ = 8.8) over norepinephrine-, serotonin-, or angiotensin II-induced contractions at 10 μM. *In vivo*, continuous infusion of darodipine (2 μg/kg per minute) increases cerebral tissue oxygenation in a cat model of focal ischemia induced by cerebral artery occlusion (CAO).² Darodipine (5 mg/kg) reduces mean arterial blood pressure and increases arterial luminal area in spontaneously hypertensive rats (SHRs).³

References

1. Hof, R.P., Vuorela, H.J., and Neumann, P. PY 108-068, a new, potent, and selective inhibitor of calcium-induced contraction of rabbit aortic rings. *J. Cardiovasc. Pharmacol.* **4**(3), 344-351 (1982).
2. Wiernsperger, N., Gyax, P., and Hofmann, A. Calcium antagonist PY 108-068: Demonstration of its efficacy in various types of experimental brain ischemia. *Stroke* **15**(4), 679-685 (1984).
3. Ferrante, F., Ricci, A., Rossodivita, I., et al. Influence of treatment with the calcium channel blocker darodipine (PY 108-068) on the morphology of pial and coronary arteries in spontaneously hypertensive rats. *Clin. Exp. Hypertens.* **16**(3), 341-357 (1994).

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, 09/19/2024

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