

PRODUCT INFORMATION



Azelnidipine

Item No. 21249

CAS Registry No.: 123524-52-7
Formal Name: 2-amino-1,4-dihydro-6-methyl-4-(3-nitrophenyl)-3,5-pyridinedicarboxylic acid, 3-[1-(diphenylmethyl)-3-azetidiny] 5-(1-methylethyl) ester

Synonyms: CS 905, RS 9054

MF: C₃₃H₃₄N₄O₆

FW: 582.7

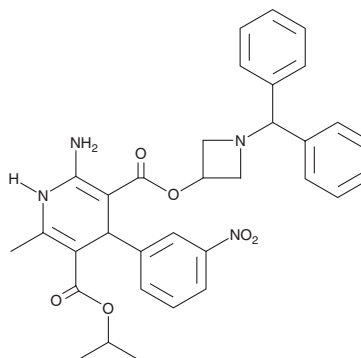
Purity: ≥98%

UV/Vis.: λ_{max}: 255, 343 nm

Supplied as: A crystalline 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

Azelnidipine is supplied as a crystalline solid. A stock solution may be made by dissolving the azelnidipine in the solvent of choice, which should be purged with an inert gas. Azelnidipine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of azelnidipine in ethanol is approximately 15 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Azelnidipine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, azelnidipine should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Azelnidipine has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Azelnidipine is a third generation, dihydropyridine L-type calcium channel blocker with a slow onset profile that displays long-lasting antihypertensive activity in hypertensive rat models compared to other L-type calcium channel blockers.^{1,2}

References

1. Oizumi, K., Nishino, H., Koike, H., *et al.* Antihypertensive effects of CS-905, a novel dihydropyridine Ca⁺⁺ channel blocker. *Jpn. J. Pharmacol.* **51(1)**, 57-64 (1989).
2. Chen, B.-L., Zhang, Y.-Z., Luo, J.-Q., *et al.* Clinical use of azelnidipine in the treatment of hypertension in Chinese patients. *Ther. Clin. Risk Manag.* **11**, 309-318 (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, 11/11/2022

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