

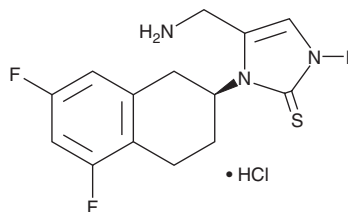
PRODUCT INFORMATION



Nepicastat (hydrochloride)

Item No. 22126

CAS Registry No.: 170151-24-3
Formal Name: 5-(aminomethyl)-1-[(2S)-5,7-difluoro-1,2,3,4-tetrahydro-2-naphthalenyl]-1,3-dihydro-2H-imidazole-2-thione, monohydrochloride
Synonym: RS 25560-197
MF: C₁₄H₁₅F₂N₃S • HCl
FW: 331.8
Purity: ≥98%
UV/Vis.: λ_{max}: 267 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

Nepicastat (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the nepicastat (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Nepicastat (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of nepicastat (hydrochloride) in these solvents is approximately 2 and 10 mg/ml, respectively.

Description

Nepicastat is an inhibitor of dopamine β-hydroxylase (DBH; IC₅₀ = 9 nM for the purified human enzyme).¹ It is selective for DBH over a panel of 12 enzymes and 13 neurotransmitter receptors (IC₅₀s or K_s = >10 μM). Nepicastat dose-dependently reduces norepinephrine content and increases dopamine content in the mesenteric artery, left ventricle, and cerebral cortex in spontaneously hypertensive rats, as well as in the renal artery, left ventricle, and cerebral cortex in beagle dogs. It attenuates increases in diastolic blood pressure and heart rate induced by preganglionic sympathetic nerve stimulation in pithed spontaneously hypertensive rats when administered orally at doses of 10 and 30 mg/kg.² Nepicastat (50 mg/kg) reduces the progressive ratio response for cocaine, but not food or sucrose pellets, in rats.³ It also reduces reinstatement of cocaine-seeking behavior induced by cues, yohimbine (Item No. 19869), or foot-shock in rats.

References

1. Stanley, W.C., Li, B., Bonhaus, D.W., *et al.* Catecholamine modulatory effects of nepicastat (RS-25560-197), a novel, potent and selective inhibitor of dopamine-β-hydroxylase. *Br. J. Pharmacol.* **121**(8), 1803-1809 (1997).
2. Stanley, W.C., Lee, K., Johnson, L.G., *et al.* Cardiovascular effects of nepicastat (RS-25560-197), a novel dopamine β-hydroxylase inhibitor. *J. Cardiovasc. Pharmacol.* **31**(6), 963-970 (1998).
3. Schroeder, J.P., Epps, S.A., Grice, T.W., *et al.* The selective dopamine β-hydroxylase inhibitor nepicastat attenuates multiple aspects of cocaine-seeking behavior. *Neuropsychopharmacology* **38**(6), 1032-1038 (2013).

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, 12/09/2022

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