

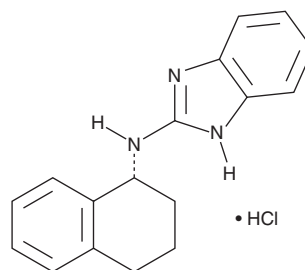
PRODUCT INFORMATION



NS 8593 (hydrochloride)

Item No. 29774

CAS Registry No.: 875755-24-1
Formal Name: N-[(1R)-1,2,3,4-tetrahydro-1-naphthalenyl]-1H-benzimidazol-2-amine, monohydrochloride
MF: C₁₇H₁₇N₃ • HCl
FW: 299.8
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

NS 8593 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the NS 8593 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. NS 8593 (hydrochloride) is soluble in organic solvents such as ethanol and DMSO. The solubility of NS 8593 (hydrochloride) in these solvents is approximately 20 and 100 mM, respectively.

Description

NS 8593 is an inhibitory gating modifier of small-conductance calcium-activated potassium (K_{Ca}2/SK) channels (K_ds = 0.42, 0.6, and 0.73 μM for K_{Ca}2.1/SK1, K_{Ca}2.2/SK2, and K_{Ca}2.3/SK3, respectively, in the presence of calcium) that decreases the calcium sensitivity of SK channels.¹ It is selective for K_{Ca}2/SK channels over intermediate- (K_{Ca}3.1/IK_{Ca}1) and large-conductance (BK) potassium channels at 10 μM. NS 8593 induces relaxation of potassium- or acetylcholine chloride-precontracted isolated tracheal rings from wild-type mice (IC₅₀s = 8.9 and 39.8 μM, respectively) or from mice in an ovalbumin-induced model of allergic asthma (IC₅₀s = 16.4 and 32.2 μM, respectively).² It inhibits aerosolized acetylcholine chloride-induced increases in respiratory system resistance in mice when administered as an aerosol at a dose of 500 μM. NS 8593 (5 mg/kg) decreases the duration of burst-pacing-induced atrial fibrillation in normotensive and spontaneously hypertensive rats.³

References

1. Strøbaek, D., Hougaard, C., Johansen, T.H., *et al.* Inhibitory gating modulation of small conductance Ca²⁺-activated K⁺ channels by the synthetic compound (R)-N-(benzimidazol-2-yl)-1,2,3,4-tetrahydro-1-naphthylamine (NS8593) reduces afterhyperpolarizing current in hippocampal CA1 neurons. *Mol. Pharmacol.* **70**(5), 1771-1782 (2006).
2. Liu, B.-B., Peng, Y.-B., Zhang, W.-J., *et al.* NS8593 inhibits Ca²⁺ permeant channels reversing mouse airway smooth muscle contraction. *Life Sci.* **238**, 116953 (2019).
3. Diness, J.G., Skibsbay, L., Jespersen, T., *et al.* Effects on atrial fibrillation in aged hypertensive rats by Ca²⁺-activated K⁺ channel inhibition. *Hypertension* **57**(6), 1129-1135 (2011).

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