

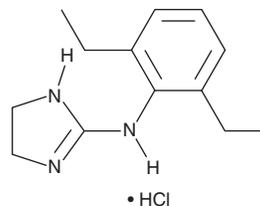
PRODUCT INFORMATION



ST 91

Item No. 43562

CAS Registry No.: 4749-61-5
Formal Name: N-(2,6-diethylphenyl)-4,5-dihydro-1H-imidazol-2-amine, monohydrochloride
MF: C₁₃H₁₉N₃ • HCl
FW: 253.8
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

ST 91 is supplied as a solid. A stock solution may be made by dissolving the ST 91 in the solvent of choice, which should be purged with an inert gas. ST 91 is soluble (≥10 mg/ml) in ethanol and DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of ST 91 can be prepared by directly dissolving the solid in aqueous buffers. ST 91 is soluble (≥10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

ST 91 is an agonist of α_2 -adrenergic receptors (α_2 -ARs).¹ It induces cAMP accumulation in HEK293 cells expressing α_{2A} , α_{2B} , or α_{2C} -ARs (EC_{50s} = 155, 371, and 851 nM, respectively). ST 91 (100 nM) inhibits isoproterenol-induced relaxation in phenylephrine-precontracted isolated rat mesenteric arterial rings.² Intrathecal administration of ST 91 (10 and 30 μ g/animal) reduces paw flinching in the first and second phases of the formalin test in rats.³ It increases plasma atrial natriuretic peptide (ANP) levels, urinary output, and sodium, potassium, and cGMP excretion in conscious rats when administered at a dose of 250 μ g/animal.⁴

References

1. Jasper, J.R., Lesnick, J.D., Chang, L.K., *et al.* Ligand efficacy and potency at recombinant α_2 adrenergic receptors: Agonist-mediated [³⁵S]GTP γ S binding. *Biochem. Pharmacol.* **55(7)**, 1035-1043 (1998).
2. Kató, E., Lipták, L., Shujaa, N., *et al.* α_{2B} -adrenoceptor agonist ST-91 antagonizes β_2 -adrenoceptor-mediated relaxation in rat mesenteric artery rings. *Eur. J. Pharmacol.* **580(3)**, 361-365 (2008).
3. Nazarian, A., Christianson, C.A., Hua, X.-Y., *et al.* Dexmedetomidine and ST-91 analgesia in the formalin model is mediated by α_{2A} -adrenoceptors: A mechanism of action distinct from morphine. *Br. J. Pharmacol.* **155(7)**, 1117-1126 (2008).
4. Gutkowska, J., Mukaddam-Daher, S., and Tremblay, J. The peripheral action of clonidine analog ST-91: Involvement of atrial natriuretic factor. *J. Pharmacol. Exp. Ther.* **281(2)**, 670-676 (1997).

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, 08/07/2025

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