

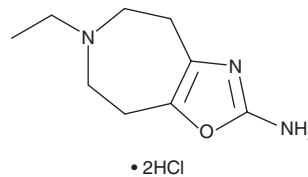
PRODUCT INFORMATION



B-HT 933 (hydrochloride)

Item No. 11941

CAS Registry No.: 36067-72-8
Formal Name: 6-ethyl-5,6,7,8-tetrahydro-4H-oxazolo[4,5-d]azepin-2-amine, dihydrochloride
Synonyms: Azepexole, Oxazoloazepin
MF: C₉H₁₅N₃O • 2HCl
FW: 254.2
Purity: ≥95%
UV/Vis.: λ_{max}: 220 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

B-HT 933 (hydrochloride) is supplied as a crystalline solid. Aqueous solutions of B-HT 933 (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of B-HT 933 (hydrochloride) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

B-HT 933 is an agonist of α₂-adrenergic receptors (α₂-ARs) with an EC₅₀ value of 0.65 μM for contraction of canine saphenous veins.¹ It induces concentration-dependent inhibition of the twitch response in rat vas deferens and guinea pig ileum.² B-HT 933 induces sedation and reduces motor activity in mice, an effect that is reversed by the α₂-AR antagonist yohimbine (Item No. 19869).³ B-HT 933 decreases blood pressure and cardiac output in cats injected with 30 μg/kg intracisternally.⁴ It also exhibits antinociceptive effects in mice with ED₅₀ values of 12.5, 20.5, and 6.1 mg/kg for the tail-immersion, tail-pinch, and acetic acid writhing tests, respectively.⁵

References

1. Ruffolo, R.R., Jr. and Zeid, R.L. Relationship between alpha 2-adrenoceptor occupancy and response for B-HT 933 in canine saphenous vein. *Eur. J. Pharmacol.* **111(2)**, 267-271 (1985).
2. Vargas, M.L. and Brugger, A.J. Mechanism of action of B-HT 933 (azepexole) in rat vas deferens and guinea-pig ileum. *Eur. J. Pharmacol.* **119(3)**, 137-141 (1985).
3. Anden, N.-E. Selective stimulation of dopamine and noradrenaline autoreceptors by B-HT920 and B-HT933, respectively. *Naunyn-Schmiedebergs Arch. Pharmacol.* **321**, 100-104 (1982).
4. Kobinger, W. and Pichler, L. Pharmacological characterization of B-HT 933 (2-amino-6-ethyl-4,5,7,8-tetrahydro-6H-oxazolo-[5,4-d]-azepin-dihydrochloride) as a hypotensive agent of the "clonidine-type". *Naunyn-Schmiedebergs Arch. Pharmacol.* **300**, 39-46 (1977).
5. Vargas, M.L., Bansinath, M., Turndorf, H., et al. Antinociceptive effects of azepexole (BHT 933) in mice. *Pain* **36(1)**, 117-123 (1989).

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, 11/08/2022

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