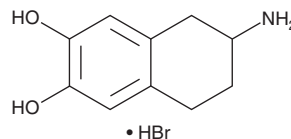


PRODUCT INFORMATION



(±)-2-Amino-6,7-dihydroxy-1,2,3,4-tetrahydronaphthalene (hydrobromide)
Item No. 23866

CAS Registry No.: 13575-86-5
Formal Name: 6-amino-5,6,7,8-tetrahydro-2,3-naphthalenediol, monohydrobromide
Synonyms: (±)-ADTN, 6,7-ADTN, 6,7-diOHATN, NSC 287353
MF: C₁₀H₁₃NO₂ • HBr
FW: 260.1
Purity: ≥95%
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

(±)-2-Amino-6,7-dihydroxy-1,2,3,4-tetrahydronaphthalene (6,7-ADTN) is supplied as a solid. A stock solution may be made by dissolving the 6,7-ADTN in water. The solubility of 6,7-ADTN in water is approximately 100 mM. We do not recommend storing the aqueous solution for more than one day.

Description

6,7-ADTN is a dopamine receptor agonist (EC₅₀s = 3.5 and 0.65 μM in rat striatal and nucleus accumbens homogenates, respectively).¹ 6,7-ADTN stimulates production of cAMP in rat striatal homogenates, a process that is reduced by 70% in the presence of the dopamine antagonist fluphenazine (Item No. 23555) at a concentration of 1 μM. *In vivo*, 6,7-ADTN induces hyperlocomotion in rats when injected bilaterally at a dose of 100 nmol per side into the nucleus accumbens. It increases cAMP production and glycoprotein secretion in the albumen gland of *H. duryi* (snail).² It also reduces sucrose feeding in *R. maderae* (cockroach) nymphs.³

References

1. Woodruff, G.N., Watling, K.J., Andrews, C.D., *et al.* Dopamine receptors in rat striatum and nucleus accumbens; conformational studies using rigid analogues of dopamine. *J. Pharm. Pharmacol.* **29(7)**, 422-427 (1977).
2. Mukai, S.T., Kiehn, L., and Saleuddin, A.S.M. Dopamine stimulates snail albumen gland glycoprotein secretion through the activation of a D1-like receptor. *J. Exp. Biol.* **207(Pt 14)**, 2507-2518 (2004).
3. Allen, J.M., Van Kummer, B.H., and Cohen, R.W. Dopamine as an anorectic neuromodulator in the cockroach *Rhyarobia maderae*. *J. Exp. Biol.* **214(Pt 22)**, 3843-3849 (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, 02/01/2023

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