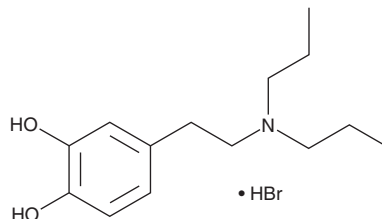


PRODUCT INFORMATION

N,N-Dipropyldopamine (hydrobromide)

Item No. 28937

CAS Registry No.: 65273-66-7
Formal Name: 4-[2-(dipropylamino)ethyl]-1,2-benzenediol, monohydrobromide
MF: $C_{14}H_{23}NO_2 \cdot HBr$
FW: 318.2
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 283 nm
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

N,N-Dipropyldopamine (hydrobromide) is supplied as a solid. A stock solution may be made by dissolving the N,N-dipropyldopamine (hydrobromide) in the solvent of choice, which should be purged with an inert gas. N,N-Dipropyldopamine (hydrobromide) is soluble in the organic solvent DMSO.

Description

N,N-Dipropyldopamine is a dopamine receptor agonist.¹⁻³ It decreases dihydrophenylalanine (DOPA) levels in the limbic forebrain and striatum of reserpinized rats (ED_{50} s = 25 and 20 μ mol/kg, respectively), as well as reduces homovanillic acid (HVA; Item No. 27307) levels in rat striatum when administered at a dose of 80 μ mol/kg.¹ N,N-Dipropyldopamine (0.5-16 mg/kg) reduces spontaneous locomotor activity in mice, an effect that can be reversed by the antipsychotic spiroperidol.^{2,3}

References

1. Wikström, H., Lindberg, P., Martinson, P., *et al.* Pivaloyl esters of N,N-dialkylated dopamine congeners. Central dopamine-receptor stimulating activity. *J. Med. Chem.* **21**(9), 864-867 (1978).
2. Feenstra, M.G., Rollema, H., Horn, A.S., *et al.* Effect of dihydroxy-2-aminotetralin derivatives on dopamine metabolism in the rat striatum. *Naunyn Schmiedebergs Arch. Pharmacol.* **310**(3), 219-225 (1980).
3. Costall, B., Lim, S.K., and Naylor, R.J. Characterisation of the mechanisms by which purported dopamine agonists reduce spontaneous locomotor activity of mice. *Eur. J. Pharmacol.* **73**(2-3), 175-188 (1981).

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, 11/11/2022

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