

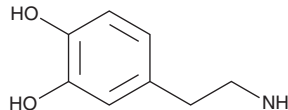
PRODUCT INFORMATION



Dopamine

Item No. 36532

CAS Registry No.: 51-61-6
Formal Name: 4-(2-aminoethyl)-1,2-benzenediol
Synonyms: DA, 3-hydroxy Tyramine, KW 3160, NSC 173182
MF: C₈H₁₁NO₂
FW: 153.2
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

Dopamine is supplied as a solid. A stock solution may be made by dissolving the dopamine in the solvent of choice, which should be purged with an inert gas. Dopamine is soluble in the organic solvent DMSO (heated) at a concentration of approximately 1 mg/ml.

Description

Dopamine is an endogenous catecholamine neurotransmitter synthesized from the amino acid L-tyrosine that acts as an agonist at dopamine receptors (D₁₋₅).¹ Dopamine is mainly synthesized in the substantia nigra and ventral tegmental area and is a precursor in norepinephrine and epinephrine biosynthesis. Dopamine-containing neurons in the brain are involved in reward-motivated behavior, motor control, and hormone release. Dopamine is also synthesized in the adrenal glands where it exerts peripheral paracrine functions including control of vasodilation, sodium excretion, insulin production, gastrointestinal motility, and the activity of lymphocytes.^{2,3} Loss or damage of dopaminergic neurons in the substantia nigra is associated with Parkinson's disease.⁴

References

1. Missale, C., Nash, S.R., Robinson, S.W., *et al.* Dopamine receptors: From structure to function. *Physiol. Rev.* **78(1)**, 190-225 (1998).
2. Hayaishi, O. Molecular genetic studies on sleep-wake regulation, with special emphases on the prostaglandin D₂ system. *J. Appl. Physiol.* **92(2)**, 863-868 (2015).
3. Garza, J.H.H. and Carr, D.J.J. Neuroendocrine peptide receptors on cells of the immune system. *Chem. Immunol.* **69**, 132-154 (1997).
4. Angeles, D.C., Ho, P., Dymock, B.W., *et al.* Antioxidants inhibit neuronal toxicity in Parkinson's disease-linked LRRK2. *Ann. Clin. Transl. Neurol.* **3(4)**, 288-294 (2016).

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, 10/21/2022

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