

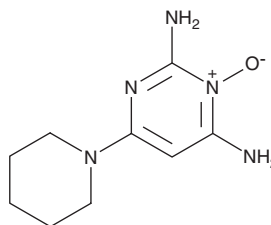
PRODUCT INFORMATION



Minoxidil

Item No. 15302

CAS Registry No.: 38304-91-5
Formal Name: 6-(1-piperidiny)-2,4-pyrimidinediamine 3-oxide
Synonyms: Loniten, U-10858
MF: C₉H₁₅N₅O
FW: 209.3
Purity: ≥98%
UV/Vis.: λ_{max}: 231, 262, 284 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

Minoxidil is supplied as a crystalline solid. A stock solution may be made by dissolving the minoxidil in the solvent of choice. Minoxidil is soluble in ethanol at a concentration of approximately 1 mg/ml.

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 minoxidil can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of minoxidil in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Minoxidil is a pyrimidine derivative that was originally developed as an anti-hypertensive agent and unexpectedly found to stimulate hair growth. Minoxidil directly affects arteriolar smooth muscle to decrease vascular resistance in part by opening ATP-sensitive potassium channels.^{1,2} It can activate cyclooxygenase 1 (AC₅₀ = 80 μM), which is expressed in the dermal papilla of hair follicles, increasing prostaglandin E₂ production in BALB/c 3T3 fibroblasts and human dermal papilla fibroblasts.³

References

1. Limas, C.J. and Cohn, J.N. Stimulation of vascular smooth muscle sodium, potassium--adenosinetriphosphatase by vasodilators. *Circ. Res.* **35(4)**, 601-607 (1974).
2. Imenshahidi, M., Hadizadeh, F., Firoozeh-Moghadam, A., *et al.* Synthesis and vasorelaxant effect of 9-aryl-1,8-acridinediones as potassium channel openers in isolated rat aorta. *Iran. J. Pharm. Res.* **11(1)**, 229-233 (2012).
3. Michelet, J.-F., Commo, S., Billoni, N., *et al.* Activation of cytoprotective prostaglandin synthase-1 by minoxidil as a possible explanation for its hair growth-stimulating effect. *J. Invest. Dermatol.* **108**, 205-209 (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, 11/02/2022

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