

Minoxidil sulfate

K⁺ channel activator

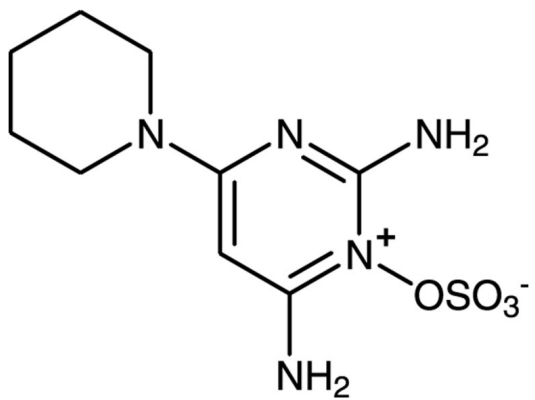
A selective ATP-sensitive K⁺ (K_{ATP}) channel opener. It is the active metabolite of minoxidil (Prod. No. BML-KC125) and is a potent (IC₅₀=0.14 μM) vascular smooth muscle relaxant. Antihypertensive. Stimulates renin secretion.

Citations: 7

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at room temperature. Make solution fresh before use.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	U-58838
Appearance	White solid.
CAS	83701-22-8
Couple Target	Potassium channel
Couple Type	Activator
Formula	$C_9H_{15}N_5O_4S$
MW	289.3
Purity	≥98%
Solubility	Soluble in DMSO (25mg/ml).
Technical Info / Product Notes	Minoxidil sulfate will slowly convert to minoxidil in DMSO solutions.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com