

# Minoxidil

K<sup>+</sup> channel activator

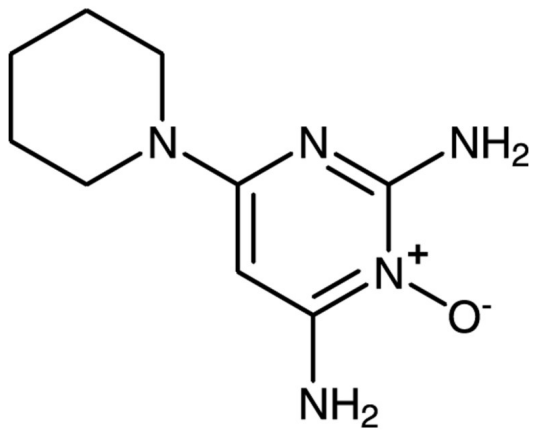
A selective ATP-sensitive K<sup>+</sup> (K<sub>ATP</sub>) channel opener. It is metabolized to the more active sulfate *in vivo*. A clinically useful anti-alopecia. Antihypertensive. Stimulates renin secretion. Potent vascular smooth muscle relaxant.

Citations: 7

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	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. Stock solutions are stable for up to 3 months at -20°C.
<b>Long Term Storage</b>	Ambient
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	U-10858
<b>Appearance</b>	White solid.
<b>CAS</b>	38304-91-5
<b>Couple Target</b>	Potassium channel
<b>Couple Type</b>	Activator
<b>Formula</b>	$C_9H_{15}N_5O$
<b>MI</b>	<b>14:</b> 6203
<b>MW</b>	209.3
<b>Purity</b>	≥98% (HPLC)
<b>Solubility</b>	Soluble in 100% ethanol (29mg/ml), DMSO (6.5mg/ml) or water (2.2mg/ml).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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