

Metformin

Antidiabetic agent

Metformin is an antidiabetic agent. Reduces the cardiovascular complications of diabetes. Does not induce hypoglycemia. Inhibits hepatic gluconeogenesis via activation of the LKB1/AMPK pathway. Has antiproliferative effects in cancer cell lines.

Citations: 26

[View Online »](#)

Ordering Information

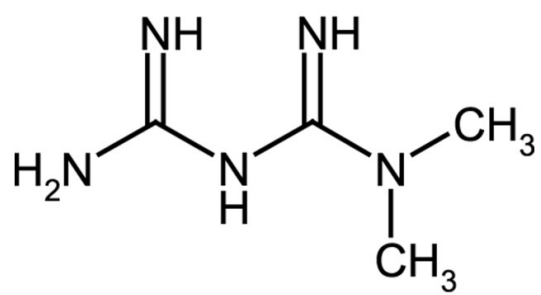
[Order Online »](#)

ALX-270-432-G005	5g
------------------	----

Manuals, SDS & CofA

[View Online »](#)

HCl



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions are stable for up to 6 months when stored at -20°C.
Handling	Hygroscopic.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	1,1-Dimethylbiguanide . hydrochloride
Appearance	White crystalline powder.
CAS	1115-70-4
Formula	$C_4H_{11}N_5 \cdot HCl$
MI	14: 5938
MW	165.6
Purity	≥97% (Assay)
RTECS	DU1800000
Solubility	Soluble in water (50mg/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-
fr@enzolifesciences.com](mailto:info-fr@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)