

Metformin . hydrochloride

AMPK activator.

Ordering Information

[Order Online »](#)

LKT-M2076-G100	100g
----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Dimethylbiguanide hydrochloride, Imidodicaronimidic diamide, N,N-dimethyl-, monohydrochloride, N,N-Dimethylbiguanide hydrochloride

CAS 1115-70-4

Formula $C_4H_{11}N_5 \cdot HCl$

MW 165.62

Purity $\geq 98\%$

Solubility 100 mM water, 50 mM DMSO



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