

Nicotinamide

Nicotinamide, the water-soluble amide of nicotinic acid, is a component of the two most important coenzymes - nicotinamide adenine dinucleotide and nicotinamide adenine dinucleotide phosphate. Thus nicotinamide is involved in numerous oxidation-reduction reactions in mammalian biological systems acting as an antioxidant.

Citations: 4

[View Online »](#)

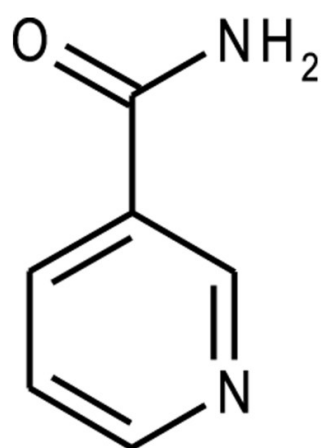
Ordering Information

[Order Online »](#)

BML-KI283-0500	500µl
----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -70°C for up to 1 year.
Long Term Storage	-80°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Clear solution.
CAS	98-92-0
Formula	$C_6H_6N_2O$
Formulation	50 mM Nicotinamide in 50 mM Tris/Cl, pH 8.0, 137 mM NaCl, 2.7 mM KCl, 1 mM $MgCl_2$.
MI	14: 6523
MW	122.1
Solubility	Soluble in water, 100% ethanol or glycerol. Slightly soluble in chloroform, almost insoluble in ether or acetone.

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)