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

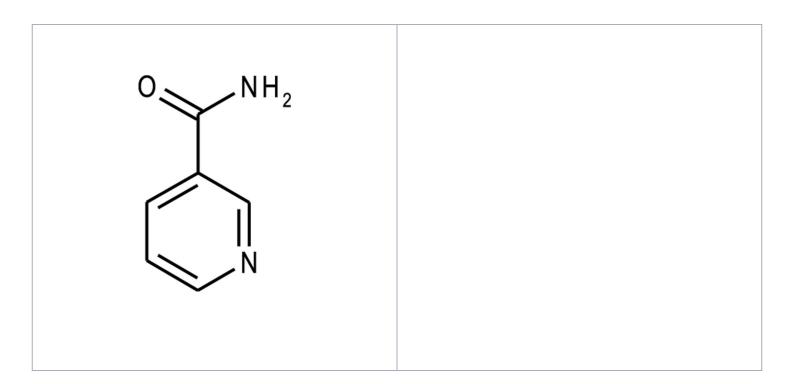
Order Online »

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 MgCl₂.

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

