

Nicotinic acid

Nicotinic acid (vitamin B3, niacin) is a naturally occurring organic compound and essential nutrient. It is a precursor of the coenzymes NAD and NADP, which play a role in hydrogen transfer processes, such as fat, carbohydrate, protein, and alcohol catabolism, DNA repair, cell signaling, etc. Insufficiency of nicotinic acid leads to nausea, skin and mouth lesions, anemia, headaches, tiredness, and in severe cases the disease pellagra. Medically used in the treatment of pellagra.

Materials Provided

Catalog Number:	92-1168
Quantity Supplied:	1 vial(s), 10 mg (re-packaged) per vial

Description

DiscoverX control ligands have been functionally tested and validated for optimal use with all cell line targets and Assay Ready kits.

Product Information

Molecular Weight:	123.11 g/mol
Source:	Synthetic
Purity:	≥ 98%
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	23-25°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

Stock Concentration:	100 mM
Reconstitution Volume:	812 µL with 10 mg vial
Reconstitution Solvent:	1N NaOH

For Research Use Only

Ordering: +1.510.979.1415 option 4 or e-mail CustomerServiceDRX@eurofins.com
Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com
General product information: www.discoverx.com