

Bio-Scale™ Mini Nuvia™ IMAC Ni-Charged, Cartridges 1 and 5 ml

Quick Start Guide

Overview

Bio-Scale Mini Nuvia IMAC Ni-Charged Cartridges are convenient, disposable prepacked low-pressure chromatography columns. Bio-Scale Mini Cartridges offer both increased run-to-run uniformity and high purity of protein through a patent pending column design and novel resin technology. This quick start guide will guide you in the basic use of the cartridges. For additional information, please consult the instruction manual, available online at bio-rad.com.

Specifications

Dimensions	1 ml: 40 mm length x 5.6 mm inner diameter 5 ml: 40 mm length x 12.6 mm inner diameter
Maximum pressure tolerance	45 psi
Recommended flow rates	1 ml: 1.5 ml/min; 5 ml: 7 ml/min
Maximum recommended flow rate	1 ml: 6 ml/min; 5 ml: 20 ml/min
Fittings	Female luer inlet and male luer outlet
Material of construction	Polypropylene column body, polyethylene (HDPE) frit
Shipping and storage conditions	20% ethanol
Autoclavability	Not autoclavable

Recommended Protocol

Step	Volume, CV*	Buffer/Lysate Solution
Equilibration/binding	5	50 mM sodium phosphate, 0.3 M NaCl, pH 8.0
Sample loading	Varies	25–40 mg/ml resin of histidine-tagged protein in 10x volume binding buffer
Wash 1	6	50 mM sodium phosphate, 0.3 M NaCl, pH 8.0
Wash 2	6	50 mM sodium phosphate, 0.3 M NaCl, pH 8.0, 15 or 30 mM imidazole
Elution	5	50 mM sodium phosphate, 0.3 M NaCl, 500 mM imidazole
Re-equilibration	5	50 mM sodium phosphate, 0.3 M NaCl, pH 8.0
Storage	5 and 5	deionized H ₂ O followed by 20% ethanol in deionized H ₂ O

* CV = column volume

Ordering Information

Bio-Scale Mini Nuvia IMAC Ni-Charged Cartridges and Resin

Catalog # Description

780-0811	Bio-Scale Mini Nuvia IMAC Ni-Charged Cartridge , 1 x 5 ml columns
780-0812	Bio-Scale Mini Nuvia IMAC Ni-Charged Cartridge , 5 x 5 ml columns
780-0800	Nuvia IMAC Ni-Charged Resin , 25 ml bottle
780-0801	Nuvia IMAC Ni-Charged Resin , 100 ml bottle
780-0802	Nuvia IMAC Ni-Charged Resin , 500 ml bottle

Fittings, Tubing, and Fittings Kits

Catalog # Description

731-8225	1.6 mm Barb to Male Luer
731-8222	1.6 mm Barb to Female Luer
732-0112	Luer to 10-32 Adaptor Fittings Kit , includes 10-32 F to M luer and 10-32 F to F luer
732-0113	Luer to BioLogic System Fittings Kit , includes 1/4-28 F to M luer and 1/4-28 F to F luer

Please visit us at bio-rad.com for the most up-to-date list of bottled resin and cartridge offerings. Larger package sizes of media are available for process scale chromatography.