

## **PureLink<sup>™</sup> Nucleic Acid Purification Rack**

Catalog no. K2100-13

**Quantity: 3** 

Store at room temperature

## Instructions for Use

**Description** The PureLink<sup>™</sup> Nucleic Acid Purification Rack is designed for use with PureLink<sup>™</sup> HiPure Plasmid DNA Miniprep, Midiprep, and Maxiprep Kits (see next page for ordering information). The dimensions of the rack are 10.25 inches (l) x 4 inches (w) x 7.5 inches (h). The PureLink<sup>™</sup> Nucleic Acid Purification Rack consists of a:

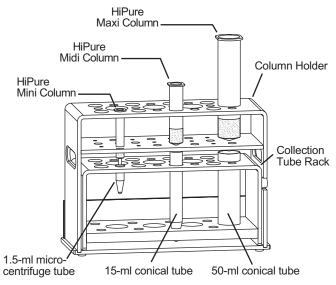
- Column holder rack (for processing 12 miniprep, 8 midiprep, and 4 maxiprep columns)
- Collection Tube Rack (capable of accommodating various types and sizes of recovery tubes)
- Large capacity Waste Tray for collecting waste

The PureLink<sup>™</sup> Nucleic Acid Purification Rack is constructed of clear acrylic for easy monitoring of samples. The rack is compatible with alcohol, but not compatible with chloroform, toluene, benzene, phenol, or acetone.

## Using the Rack

To use the rack with PureLink<sup>™</sup> HiPure Plasmid DNA Kits:

- 1. Separate the nested rack system components and place the Waste Tray into the large column holder. Set aside the Collection Tube Rack for use in the nucleic acid elution step.
- 2. Place the desired columns onto the column holder. Follow the directions outlined in the mini, midi, or maxi prep manual for the purification of nucleic acid. After the last wash step, remove the Waste Tray and insert appropriate Collection Tubes (1.5 ml, 15 ml, or 50 ml tubes) into the Collection Tube Rack corresponding to the placement of the columns (see figure below).



3. Place the loaded Collection Tube Rack into the column holder, lifting up the columns to slightly allow the columns to sit inside the collection tubes. Elute the nucleic acid as described in the mini, midi, or maxi prep manual.

Part No: 25-0792 version A

Rev. Date: 10 Dec 2004

This product is distributed for laboratory research use only. CAUTION: Not for diagnostic use. The safety and efficacy of this product in diagnostic or other clinical uses has not been established.

For technical questions about this product, call the Invitrogen TECH-LINE<sup>SM</sup> 800 955 6288

## Instructions for Use, Continued

Cleaning	Clean the PureLink <sup>™</sup> Nucleic Acid Purification Rack with a mild laboratory detergent and rinse with deionized water after use.		
Accessory Products	Ordering information for PureLink <sup>™</sup> HiPure Plasmid DNA Purification Kits is listed below. For more information, visit our web site at www.invitrogen.com or contact Technical Service.		
	Product	Quantity	Catalog no.
	PureLink <sup>™</sup> HiPure Plasmid DNA Miniprep	25 reactions	K2100-02
		100 reactions	K2100-03
	PureLink <sup>™</sup> HiPure Plasmid DNA Midiprep	25 reactions	K2100-04
		50 reactions	K2100-05
	PureLink <sup>™</sup> HiPure Plasmid DNA Maxiprep	10 reactions	K2100-06
		25 reactions	K2100-07
	should have any questions or concerns about an Invitrogen product or service, contact our Technical Service Representatives. Invitrogen warrants that all of its products will perform according to specifications stated on the certificate of analysis. The company will replace, free of charge, any product that does not meet those specifications. <u>This warranty limits Invitrogen</u> <u>Corporation's liability only to the cost of the product</u> . No warranty is granted for products beyond their listed expiration date. No warranty is applicable unless all product components are stored in accordance with instructions. Invitrogen reserves the right to select the method(s) used to analyze a product unless Invitrogen agrees to a specified method in writing prior to acceptance of the order. Invitrogen makes every effort to ensure the accuracy of its publications, but realizes that the occasional typographical or other error is inevitable. Therefore Invitrogen makes no warranty of any kind regarding the contents of any publication or documentation. If you discover an error in any of our publications, please report it to our Technical Service Representatives. <b>Invitrogen assumes no responsibility or liability for any special, incidental, indirect or consequential loss or damage whatsoever. The above limited warranty is sole and exclusive.</b> <b>No other warranty is made, whether expressed or implied, including any warranty of merchantability or fitness for a particular purpose.</b>		