



TECHNICAL BULLETIN

Vac-Man[®] Jr. Laboratory Vacuum Manifold

Instructions for Use of Products
A7660 and A3110

Vac-Man[®] Jr. Laboratory Vacuum Manifold

All technical literature is available at: www.promega.com/protocols/
 Visit the web site to verify that you are using the most current version of this Technical Bulletin.
 E-mail Promega Technical Services if you have questions on use of this system: techserv@promega.com

1. Description.....		1
2. Product Components		1
3. Setup and Operation.....		2
4. Cleanup		2
5. Related Products		2
6. Summary of Changes		4

1. Description

The Vac-Man[®] Jr. Laboratory Vacuum Manifold (Cat.# A7660) is composed of a #8 neoprene stopper and two, one-way Luer-Lok[®] stopcocks as a single unit that inserts easily into a 1- or 2-liter side-arm flask with a 34–43mm opening. The Vac-Man[®] Jr. Laboratory Vacuum Manifold (Cat.# A3110) is composed of a #10 neoprene stopper and two, one-way Luer-Lok[®] stopcocks as a single unit that inserts easily into a 1- or 2-liter side-arm flask with a 41–48.5mm opening. These small manifolds are ideal for rapid and efficient nucleic acid purification when used in conjunction with a water aspirator-type vacuum source and the columns in any of the PureYield[™], Wizard[®] or Wizard[®] Plus Purification Systems. The Vac-Man[®] Jr. Manifolds are an inexpensive alternative to the popular 20-sample capacity Vac-Man[®] Laboratory Vacuum Manifold (Cat.# A7231).

The Vac-Man[®] Jr. Manifolds are designed for use with a water aspirator vacuum source and **not** a high-pressure vacuum source such as a building-integrated, benchtop-accessible vacuum source. For safe operation, carefully read the operating instructions and cautionary notes before using the Vac-Man[®] Jr. Laboratory Vacuum Manifold.

2. Product Components

PRODUCT	CAT.#
Vac-Man[®] Jr. Laboratory Vacuum Manifold, #8	A7660
Vac-Man[®] Jr. Laboratory Vacuum Manifold, #10	A3110

3 Setup and Operation

Use only PureYield™ or Wizard® Purification Systems with the Vac-Man® Jr. Manifold. Use of other chemicals or purification systems with the manifold may create a safety hazard. Wear a lab coat and safety glasses when using PureYield™ or Wizard® Purification Systems with the Vac-Man® Jr. Manifold. Before beginning a purification procedure, inspect the components of the vacuum manifold. Do not use if cracks or other damage are visible. Ensure that the one-way Luer-Lok® stopcocks are attached firmly to the neoprene stopper.

1. Insert the neoprene stopper into the mouth of a 1- or 2-liter side-arm flask. Do not restrict the stopper in any way (e.g., wire or tape). The stopper should be allowed to come loose, thereby serving as a safety valve in the event of excessive pressure.
2. Insert columns into the stopcocks. Apply the lysate or resin/DNA mix according to the protocol provided in the PureYield™ or Wizard® Purification System.
3. Connect the side-arm flask to the water aspirator vacuum source with a sturdy vacuum-grade hose.
4. Open the stopcocks, and apply vacuum.

Note: Additional DNA purifications can be performed by using multiple manifolds in series or by sequentially using a single manifold and replacing with fresh columns each time.

4. Cleanup

1. With the vacuum off, disconnect the vacuum hose from the side-arm flask.
2. Remove the neoprene stopper, and drain the contents of the flask over a sink or other appropriate liquid disposal area.
3. Replace the neoprene stopper for the next use.

5. Related Products

Product	Size	Cat.#
Vac-Man® Laboratory Vacuum Manifold, 20-sample capacity	1 each	A7231
Vacuum Adapters	20 each	A1331

Nucleic Acid Purification Systems

Product	Size	Cat. #
PureYield™ Plasmid Miniprep System	10 preps	A1220
	100 preps	A1223
	250 preps	A1222
PureYield™ Plasmid Midiprep System	25 preps	A2492
	100 preps	A2495
PureYield™ Plasmid Maxiprep System	10 preps	A2392
	25 preps	A2393
Wizard® SV Gel and PCR Clean-Up System*	50 preps	A9281
	250 preps	A9282
Wizard® SV Genomic DNA Purification System	50 preps	A2360
	250 preps	A2361
Wizard® SV 96 Genomic DNA Purification System	1 × 96 preps	A2370
	4 × 96 preps	A2371
SV Total RNA Isolation System*	50 preps	Z3100
	250 preps	Z3105
SV Total RNA Isolation System, Trial Size*	10 preps	Z3101
Wizard® <i>Plus</i> SV Minipreps DNA Purification System*	50 preps	A1330
	250 preps	A1460
Wizard® <i>Plus</i> SV Minipreps DNA Purification System + Vacuum Adapters*	50 preps	A1340
	250 preps	A1470
Wizard® <i>Plus</i> Minipreps DNA Purification System*	50 preps	A7100
	100 preps	A7500
	250 preps	A7510
Wizard® Minipreps DNA Purification Resin*	250ml	A7141
Wizard® <i>Plus</i> Midipreps DNA Purification System*	25 preps	A7640
Wizard® <i>Plus</i> Maxipreps DNA Purification System*	10 preps	A7270
Wizard® <i>Plus</i> Megapreps DNA Purification System*	5 preps	A7300
Wizard® DNA Clean-Up System*	100 preps	A7280
Wizard® PCR Preps DNA Purification System*	50 preps	A7170
Wizard® Enviro Start-Up Kit	110V	A3050
	220V	A3060
Maxwell® RSC Enviro Start-Up Kit	110V	A3070
	220V	A3080

*For Laboratory Use.



6. Summary of Changes

The following changes were made to the 5/23 revision of this document:

1. Added Cat. # A3110.
2. Updated Section 5.
3. Moved document to a new template.

© 1996–2023 Promega Corporation. All Rights Reserved.

Maxwell, Vac-Man and Wizard are registered trademarks of Promega Corporation. PureYield is a trademark of Promega Corporation.

Luer-Lok is a registered trademark of Becton, Dickinson and Company

Products may be covered by pending or issued patents or may have certain limitations. Please visit our Web site for more information.

All prices and specifications are subject to change without prior notice.

Product claims are subject to change. Please contact Promega Technical Services or access the Promega online catalog for the most up-to-date information on Promega products.