

Wizard® SV 96 PCR Clean-Up System

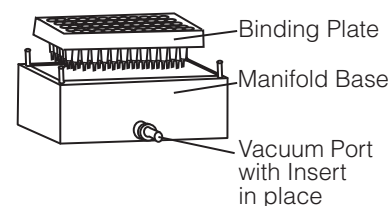
INSTRUCTIONS FOR USE OF PRODUCTS A9340, A9341, A9342 AND A9345.

Quick
PROTOCOL

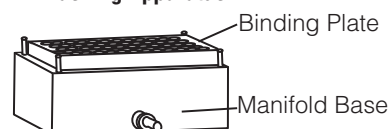
PCR Clean-Up Protocol

1. Prepare the vacuum manifold as shown in the figure. Place the Binding Plate in the vacuum Manifold Base. Orient the Binding Plate in the vacuum manifold with the numerical column headers toward the vacuum port. Attach the vacuum line to the vacuum port on the Manifold Base.
2. Add an equal volume of Membrane Binding Solution to each PCR sample in a 96-well plate (e.g., 100µl of Membrane Binding Solution to each 100µl PCR sample).
3. Mix by pipetting and transfer entire sample volume to the wells of the Binding Plate sitting on the vacuum manifold. Incubate for 1 minute at room temperature.
4. Apply vacuum at 15 inches Hg until the sample passes through the Binding Plate, approximately 30 seconds. Release vacuum.
5. Add 200µl of freshly made 80% ethanol to each well of the Binding Plate. Incubate for 1 minute at room temperature. Apply the vacuum until the 80% ethanol passes through the plate, approximately 30 seconds. Release vacuum.
6. Repeat Step 5 for a total of three 200µl 80% ethanol washes.
7. After the wells of the Binding Plate have emptied from the final wash, continue to apply the vacuum for an additional 4 minutes to allow the binding matrix to dry.
8. Turn off the vacuum. Release the vacuum line from the Manifold Base and snap it into the vacuum port in the Vacuum Manifold Collar. Remove the Binding Plate from the Manifold Base and blot by gently tapping onto a clean paper towel to remove residual ethanol.
9. Place a 96-well, U-bottom Collection Plate in the Manifold Bed and position the Vacuum Manifold Collar on top. Orient the U-bottom Collection Plate with the numerical column headers toward the vacuum port.
10. Position the Binding Plate on top of the Manifold Collar and the Collection Plate as shown in the figure. The Binding Plate tips must be centered on the Collection Plate wells, and both plates must be in the same orientation. Add 100µl Nuclease-Free Water to each well of the Binding Plate and incubate for 1 minute at room temperature. Apply vacuum until the solution passes through the plate, approximately 1 minute.
11. Release the vacuum and remove the Binding Plate. Carefully remove the Manifold Collar. If droplets are present on the walls of the Collection Plate, briefly centrifuge the plate to collect the droplets on the bottom of the wells. Eluate volumes may vary but are approximately 75µl.

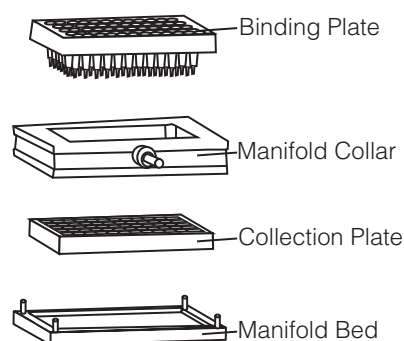
A. PCR Product Binding Apparatus



B. Washing Apparatus



C. Elution Apparatus



See additional protocol information in Technical Bulletin #TB311, available online at

www.promega.com

ORDERING/TECHNICAL INFORMATION:

www.promega.com • Phone 608-274-4330 or 800-356-9526 • Fax 608-277-2601



Promega

Printed in USA. Revised 3/09

Part #9FB071

© 2002–2009 Promega Corporation. All Rights Reserved.

3744MA06_2A