

Wizard® SV Gel and PCR Clean-Up System

INSTRUCTIONS FOR USE OF PRODUCTS A9280, A9281, A9282, AND A9285.

Quick
PROTOCOL

DNA Purification by Centrifugation

Gel Slice and PCR Product Preparation

A. Dissolving the Gel Slice

1. Following electrophoresis, excise DNA band from gel and place gel slice in a 1.5ml microcentrifuge tube.
2. Add 10µl Membrane Binding Solution per 10mg of gel slice. Vortex and incubate at 50–65°C until gel slice is completely dissolved.

B. Processing PCR Amplifications

1. Add an equal volume of Membrane Binding Solution to the PCR amplification.

Binding of DNA

1. Insert SV Minicolumn into Collection Tube.
2. Transfer dissolved gel mixture or prepared PCR product to the Minicolumn assembly. Incubate at room temperature for 1 minute.
3. Centrifuge at $16,000 \times g$ for 1 minute. Discard flowthrough and reinsert Minicolumn into Collection Tube.

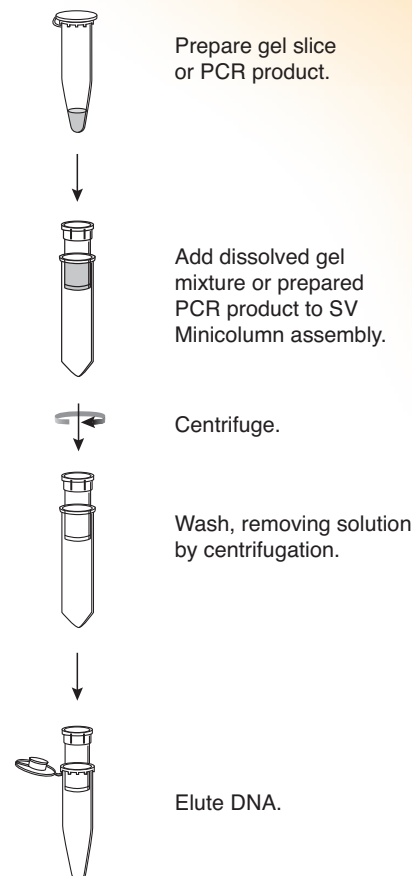
Washing

4. Add 700µl Membrane Wash Solution (ethanol added). Centrifuge at $16,000 \times g$ for 1 minute. Discard flowthrough and reinsert Minicolumn into Collection Tube.
5. Repeat Step 4 with 500µl Membrane Wash Solution. Centrifuge at $16,000 \times g$ for 5 minutes.
6. Empty the Collection Tube and recentrifuge the column assembly for 1 minute with the microcentrifuge lid open (or off) to allow evaporation of any residual ethanol.

Elution

7. Carefully transfer Minicolumn to a clean 1.5ml microcentrifuge tube.
8. Add 50µl of Nuclease-Free Water to the Minicolumn. Incubate at room temperature for 1 minute. Centrifuge at $16,000 \times g$ for 1 minute.
9. Discard Minicolumn and store DNA at 4°C or –20°C.

Additional protocol information is available in Technical Bulletin #TB308, available online at: www.promega.com



3760MA07_2A

ORDERING/TECHNICAL INFORMATION:

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

© 2002, 2004, 2005 and 2009 Promega Corporation. All Rights Reserved.



Promega

Printed in USA. Revised 11/09
Part #9FB072

Wizard® SV Gel and PCR Clean-Up System

INSTRUCTIONS FOR USE OF PRODUCTS A9280, A9281, A9282 AND A9285.

Quick
PROTOCOL

DNA Purification by Vacuum

Gel Slice and PCR Product Preparation

A. Dissolving the Gel Slice

1. Following electrophoresis, excise DNA band from gel and place gel slice in a 1.5ml microcentrifuge tube.
2. Add 10 μ l Membrane Binding Solution per 10mg of gel slice. Vortex and incubate at 50–65°C until gel slice is completely dissolved.

B. Processing PCR Amplifications

1. Add an equal volume of Membrane Binding Solution to the PCR amplification.

Binding of DNA

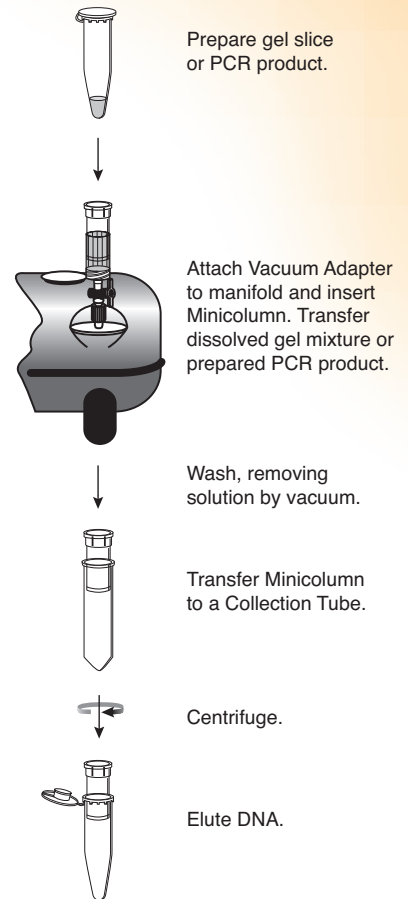
1. Attach Vacuum Adapter to manifold port and insert SV Minicolumn into Adapter.
2. Transfer dissolved gel mixture or prepared PCR product to the Minicolumn. Incubate at room temperature for 1 minute.
3. Apply vacuum to pull liquid through Minicolumn. Release vacuum when all liquid has passed through Minicolumn.

Washing

4. Add 700 μ l Membrane Wash Solution (ethanol added). Apply a vacuum to pull solution through Minicolumn.
5. Turn off vacuum and repeat Step 4 with 500 μ l Membrane Wash Solution. Apply a vacuum to pull solution through Minicolumn.
6. Transfer Minicolumn to a Collection Tube. Centrifuge at 16,000 $\times g$ for 5 minutes.
7. Empty the Collection Tube and recentrifuge the column assembly for 1 minute with the microcentrifuge lid open (or off) to allow evaporation of any residual ethanol.

Elution

8. Carefully transfer Minicolumn to a clean 1.5ml microcentrifuge tube.
9. Add 50 μ l of Nuclease-Free Water to the Minicolumn. Incubate at room temperature for 1 minute. Centrifuge at 16,000 $\times g$ for 1 minute.
10. Discard Minicolumn and store DNA at 4°C or –20°C.



Additional protocol information is available in Technical Bulletin #TB308, 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 11/09
Part #9FB072

© 2002, 2004, 2005 and 2009 Promega Corporation. All Rights Reserved.

3759MA07_2A