

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 of Membrane Binding Solution per 10mg of gel slice. Vortex and incubate at 50–65°C until gel slice is completely dissolved.



Prepare gel slice or PCR product.

B. Processing PCR Amplifications

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

Binding of DNA

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



Add dissolved gel mixture or prepared PCR product to SV Minicolumn assembly.



Centrifuge.

Washing

5. Add 700µl of Membrane Wash Solution (ethanol added). Centrifuge at 16,000 × g for 1 minute. Discard flowthrough and reinsert Minicolumn into Collection Tube.
6. Repeat Step 4 with 500µl of Membrane Wash Solution. Centrifuge at 16,000 × 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.



Wash, removing solution by centrifugation.

Elution

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



Elute DNA.

3760MA07_2A

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 10µl of 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

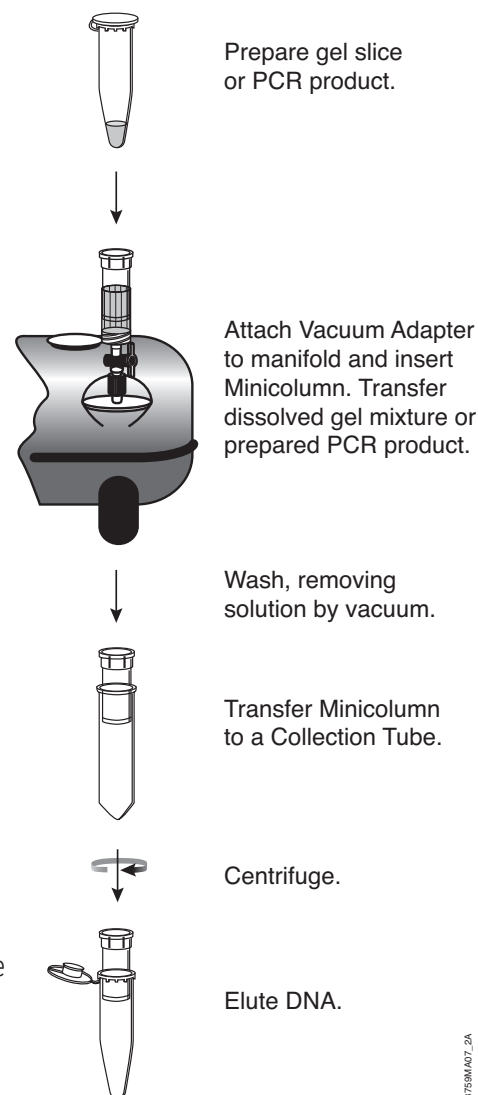
2. Attach Vacuum Adapter to manifold port and insert SV Micolumn into Adapter.
3. Transfer dissolved gel mixture or prepared PCR product to the Micolumn. Incubate at room temperature for 1 minute.
4. Apply vacuum to pull liquid through Micolumn. Release vacuum when all liquid has passed through Micolumn.ollection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

Washing

5. Add 700µl of Membrane Wash Solution (ethanol added). Apply a vacuum to pull solution through Micolumn.
6. Turn off vacuum and repeat Step 4 with 500µl of Membrane Wash Solution. Apply a vacuum to pull solution through Micolumn.
7. Transfer Micolumn to a Collection Tube. Centrifuge at 16,000 × g for 5 minutes.
8. 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

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



3759MA07_2A

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