

Wizard® DNA Clean-Up System

INSTRUCTIONS FOR USE OF PRODUCT A7280.

Quick
PROTOCOL

Purification With a Vacuum Manifold

Before You Begin

Thoroughly mix the Wizard® DNA Clean-Up Resin before removing an aliquot. If crystals or aggregates are present, dissolve by warming the resin to 37°C for 10 minutes. The resin itself is insoluble. Cool to 25–30°C before use.

The sample volume must be between 50 and 500µl. If the sample volume is less than 50µl, bring the volume up to at least 50µl with sterile water.

Binding of DNA

1. Use one Wizard® Minicolumn for each sample.
2. Attach the provided Syringe Barrel to the Luer-Lok® extension of each Minicolumn. Insert the tip of the Minicolumn/Syringe Barrel assembly into the vacuum manifold.
3. Add 1ml of Wizard® DNA Clean-Up Resin to a 1.5ml microcentrifuge tube. Add the sample to the Clean-Up Resin and mix by inversion.
4. Pipet the resin/DNA mix into the Syringe Barrel. Apply a vacuum to draw the solution through the Minicolumn. Break the vacuum to the Minicolumn.

Washing

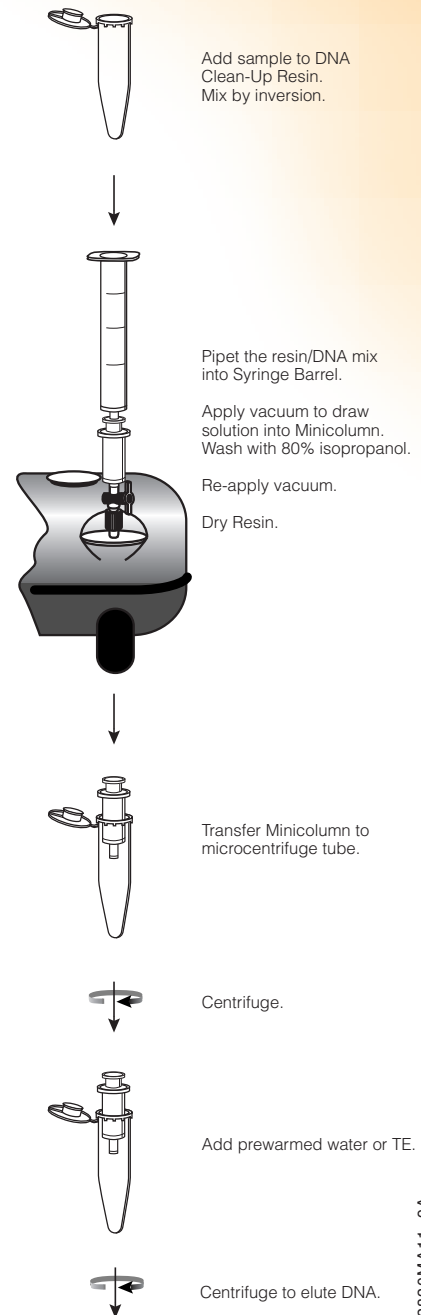
5. Add 2ml of 80% isopropanol to the Syringe Barrel, and re-apply a vacuum to draw the solution through the Minicolumn.
6. Dry the resin by continuing to draw a vacuum for 30 seconds after the solution has been pulled through the column. Do not dry the resin for more than 30 seconds. Remove the Syringe Barrel and transfer the Minicolumn to a 1.5ml microcentrifuge tube.

Centrifuge the Minicolumn at maximum speed in a microcentrifuge for 2 minutes to remove any residual isopropanol.

Elution

7. Transfer the Minicolumn to a new microcentrifuge tube. Apply 50µl of pre-warmed (65–70°C) water or TE buffer to the Minicolumn and wait 1 minute. Centrifuge the Minicolumn for 20 seconds at maximum speed to elute the bound DNA.
8. Remove and discard the Minicolumn. The purified DNA may be stored in the microcentrifuge tube at 4°C or –20°C.

Additional protocol information is available in Technical Bulletin #TB141, available online at: www.promega.com



2820MA11_9A

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 #9FB021

Wizard® DNA Clean-Up System

INSTRUCTIONS FOR USE OF PRODUCT A7280.

Quick
PROTOCOL

Purification Without a Vacuum Manifold (Using 3ml Luer-Lok® Syringes)

Before You Begin

Thoroughly mix the Wizard® DNA Clean-Up Resin before removing an aliquot. If crystals or aggregates are present, dissolve by warming the resin to 37°C for 10 minutes. The resin itself is insoluble. Cool to 25–30°C before use.

The sample volume must be between 50 and 500µl. If the sample volume is less than 50µl, bring the volume up to at least 50µl with sterile water.

Binding of DNA

1. Use one Wizard® Minicolumn for each sample. Remove and set aside the plunger from a 3ml disposable syringe. Attach the Syringe Barrel to the Luer-Lok® extension of each Minicolumn.
2. Add 1ml of Wizard® DNA Clean-Up Resin to a 1.5ml microcentrifuge tube. Add the sample to the Clean-Up Resin and mix by inversion.
3. Pipet the Wizard® DNA Clean-Up Resin containing the bound DNA into the Syringe Barrel. Insert the syringe plunger slowly and gently push the slurry into the Minicolumn with the syringe plunger.

Washing

4. Detach the syringe from the Minicolumn and remove the plunger from the syringe. Reattach the Syringe Barrel to the Minicolumn. Pipet 2ml of 80% isopropanol into the syringe. Reinsert the plunger and push the solution through the Minicolumn.
5. Remove the Syringe Barrel and transfer the Minicolumn to a 1.5ml microcentrifuge tube. Centrifuge the Minicolumn at maximum speed in a microcentrifuge for 2 minutes to dry the resin.

Elution

6. Transfer the Minicolumn to a new microcentrifuge tube. Apply 50µl of pre-warmed (65–70°C) water or TE buffer to the Minicolumn and wait 1 minute. Centrifuge the Minicolumn for 20 seconds at maximum speed to elute the bound DNA.
7. Remove and discard the Minicolumn. The purified DNA may be stored in the microcentrifuge tube at 4°C or –20°C.

Additional protocol information is available in Technical Bulletin #TB141, available online at: www.promega.com



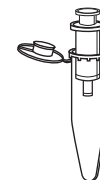
Add sample to DNA Clean-Up Resin. Mix by inversion.



Attach Syringe Barrel to Minicolumn.

Pipet the resin/DNA into Syringe Barrel.

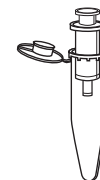
Insert syringe plunger and push slurry into Minicolumn.



Transfer Minicolumn to microcentrifuge tube.



Centrifuge.



Transfer Minicolumn to new microcentrifuge tube.

Add prewarmed water or TE.



Centrifuge to elute DNA.

2824MA11_9A

ORDERING/TECHNICAL INFORMATION:

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



Promega