

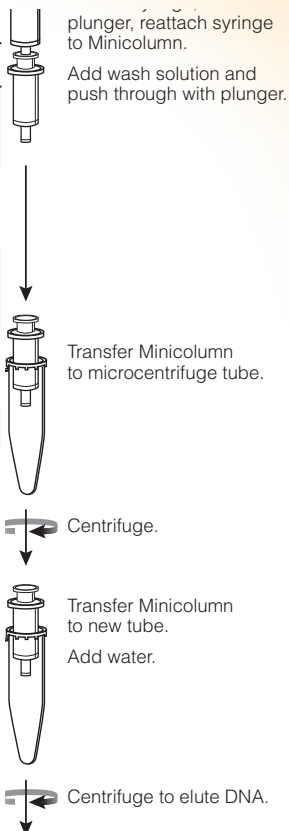
# Wizard® Plus Minipreps DNA Purification System

INSTRUCTIONS FOR USE OF PRODUCTS A7100, A7500, A7510, A7141 AND A7211.

**Quick**  
PROTOCOL

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

Prepare Cleared Lysate	Culture Volume		
	1–3ml	3–5ml	5–10ml
1. Pellet cells.	1–2 minutes 10,000 × g <sup>†</sup>	10 minutes 10,000 × g <sup>†</sup>	10 minutes 1,400 × g
2. Suspend pellet in Cell Resuspension Solution.	200µl	300µl	400µl
3. Add Cell Lysis Solution. Invert 4 times to mix.	200µl	300µl	400µl
4. Add Neutralization Solution. Invert 4 times to mix.*	200µl	300µl	400µl
5. Centrifuge lysate for 5 minutes at 10,000 × g.			



### Plasmid DNA Purification

- Remove plunger from 3ml Luer-Lok® syringe (Becton-Dickinson Cat.# 9585). Attach syringe barrel to Luer-Lok® extension of Minicolumn.
- Resuspend resin. Add 1ml resin to each Minicolumn/syringe assembly. Carefully transfer cleared lysate (from Step 5) to resin in each assembly.
- Insert plunger and push resin and lysate into Minicolumn.\*

### Washing

- Detach syringe from Minicolumn; remove plunger from syringe barrel. Reattach barrel to Minicolumn.
- Add 2ml Column Wash Solution containing ethanol. Insert the plunger and push the Column Wash Solution through the Minicolumn.
- Remove syringe and transfer the Minicolumn to a 1.5ml microcentrifuge tube. Centrifuge at 10,000 × g for 2 minutes.

### Elution

- Transfer Minicolumn to a new microcentrifuge tube.
- Add 50µl of Nuclease-Free Water to the Minicolumn and wait 1 minute.  
For plasmids ≥10kb, use water preheated to 70°C; for plasmids ≥20kb, use water preheated to 80°C.
- Centrifuge at 10,000 × g for 20 seconds at room temperature.
- Remove and discard Minicolumn. Store DNA at –20°C or below.

<sup>†</sup>Maximum speed on a microcentrifuge.

\*For EndA<sup>+</sup> strains and other modifications, additional protocol information is available in Technical Bulletin #TB117, available from Promega or online at [www.promega.com](http://www.promega.com)

### ORDERING/TECHNICAL INFORMATION:

[www.promega.com](http://www.promega.com) • Phone 608-274-4330 or 800-356-9526 • Fax 608-277-2601



**Promega**

Printed in USA Revised 4/09. 8196BI-PC-CP  
Part #9FB016

# Wizard® Plus Minipreps DNA Purification System

INSTRUCTIONS FOR USE OF PRODUCTS A7100, A7500, A7510, A7141 AND A7211.

**Quick**  
PROTOCOL

## Purification With a Vacuum Manifold

Prepare Cleared Lysate	Culture Volume		
	1–3ml	3–5ml	5–10ml
1. Pellet cells.	1–2 minutes 10,000 × <i>g</i> <sup>†</sup>	10 minutes 10,000 × <i>g</i> <sup>†</sup>	10 minutes 1,400 × <i>g</i>
2. Suspend pellet in Cell Resuspension Solution.	200µl	300µl	400µl
3. Add Cell Lysis Solution. Invert 4 times to mix.	200µl	300µl	400µl
4. Add Neutralization Solution. Invert 4 times to mix.*	200µl	300µl	400µl
5. Centrifuge lysate for 5 minutes at 10,000 × <i>g</i> .			

### Plasmid DNA Purification

- Resuspend resin. Add 1ml resin to each Minicolumn/syringe assembly. Carefully transfer cleared lysate (from Step 5) to resin in each assembly.
- Open stopcocks. Apply vacuum to pull liquid through column. Release vacuum when all liquid has passed through column.\*

### Washing

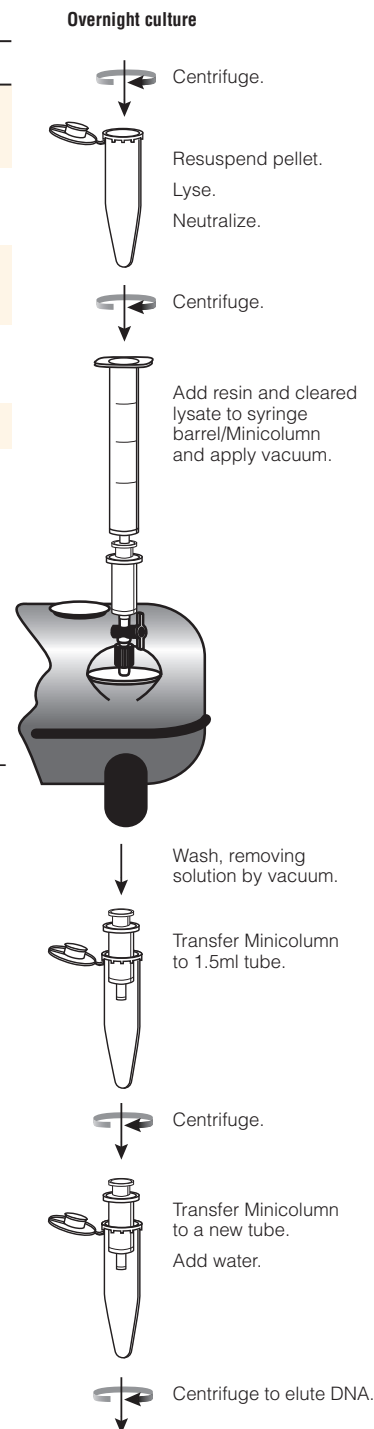
- Add 2ml Column Wash Solution containing ethanol. Apply vacuum to pull liquid through column. Continue vacuum for additional 30 seconds to dry resin.
- Remove syringe barrel; transfer Minicolumn to a 1.5ml microcentrifuge tube. Centrifuge at 10,000 × *g* for 2 minutes.

### Elution

- Transfer Minicolumn to a new microcentrifuge tube.
- Add 50µl of Nuclease-Free Water to the Minicolumn and wait 1 minute.  
For plasmids ≥10kb, use water preheated to 70°C; for plasmids ≥20kb, use water preheated to 80°C.
- Centrifuge at 10,000 × *g* for 20 seconds at room temperature.
- Remove and discard Minicolumn. Store DNA at –20°C or below.

<sup>†</sup>Maximum speed on a microcentrifuge.

\*For EndA<sup>+</sup> strains and other modifications, additional protocol information is available in Technical Bulletin #TB117, available from Promega or online at [www.promega.com](http://www.promega.com)



2776MA10\_9A

2776MA10\_9A



**Promega**

### ORDERING/TECHNICAL INFORMATION:

[www.promega.com](http://www.promega.com) • Phone 608-274-4330 or 800-356-9526 • Fax 608-277-2601

Printed in USA Revised 4/09. 8196BI-PC-CP  
Part #9FB016

© 1999, 2009 Promega Corporation. All Rights Reserved.