# Wizard<sup>®</sup> Plus Megapreps DNA Purification System

INSTRUCTIONS FOR USE OF PRODUCTS A7300, A7361 AND A7421.



## Megapreps Procedure (500-1,000ml culture volumes)

#### Preparation of Cleared Lysate

- 1. Pellet cells at  $1,500 \times g$  for 20 minutes at room temperature.\*
- 2. Suspend pellet in 30ml Cell Resuspension Solution.
- 3. Add 30ml Cell Lysis Solution. Invert to mix.
- 4. Add 30ml Neutralization Solution. Invert to mix.<sup>†</sup>
- 5. Centrifuge lysate at  $14,000 \times g$  for 15 minutes.\*
- 6. Filter supernatant into a graduated cylinder. Measure volume, then transfer liquid to centrifuge bottle.
- 7. Add 0.5 volume isopropanol. Invert to mix.\*
- 8. Centrifuge at  $14,000 \times g$  for 15 minutes.\*
- 9. Resuspend pellet in 4ml TE buffer.

### Plasmid DNA Purification

- 10. Resuspend resin. Add 20ml resin to DNA from Step 9. Swirl to mix.
- Attach Megacolumn to vacuum manifold. Transfer resin/DNA mixture to Megacolumn. Apply vacuum and release when all liquid has passed through Megacolumn.<sup>†</sup>

### Washing

- 12. Add 25ml Column Wash Solution containing ethanol. Apply vacuum, pulling liquid through column. Release vacuum and repeat.
- 13. Add 5ml of 80% ethanol. Apply vacuum; continue for 1 minute after liquid has passed through column.
- 14. Transfer Megacolumn to 50ml centrifuge tube. Centrifuge at  $1,300 \times g$  for 5 minutes, using a swinging bucket rotor.\*
- 15. Place Megacolumn on manifold. Apply vacuum and continue for 5 minutes.

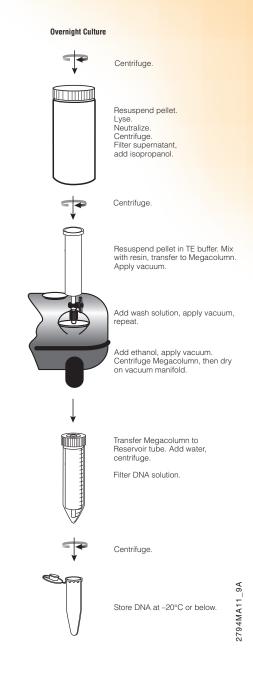
### Elution

- Place Megacolumn in Reservoir tube provided. Add 3.0ml preheated water. Wait 1 minute, then centrifuge at 1,300 × g for 5 minutes to elute DNA. For plasmids ≥10kb, use water preheated to 70°C; for plasmids ≥20kb, use water preheated to 80°C.
- 17. Filter eluted DNA. Centrifuge filtrate at  $14,000 \times g$  for 1 minute.
- 18. Transfer supernatant containing DNA to a new centrifuge tube and store at  $-20^{\circ}$ C or below.

\*Rotor, solutions and centrifugation must be at room temperature. <sup>†</sup>For EndA+ strains and other modifications, additional protocol information in Technical Bulletin #TB140, available online at www.promega.com

#### ORDERING/TECHNICAL INFORMATION:

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





Part #9FB020 © 1999, 2009 Promega Corporation. All Rights Reserved.