Wizard® Plus Maxipreps DNA Purification System

INSTRUCTIONS FOR USE OF PRODUCTS A7270, A7401 AND A7421.



Maxipreps Procedure (100–500ml culture volumes)

Preparation of Cleared Lysate

- 1. Pellet cells at 5,000 \times g for 10 minutes at room temperature.*
- 2. Suspend pellet in 15ml Cell Resuspension Solution.
- 3. Add 15ml Cell Lysis Solution. Invert to mix.
- 4. Add 15ml Neutralization Solution. Invert to mix.⁺
- 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 2ml TE buffer.

Plasmid DNA Purification

- 10. Resuspend resin. Add 10ml resin to DNA from Step 9. Swirl to mix.
- 11. Attach Maxicolumn to vacuum manifold. Transfer resin/DNA mixture to Maxicolumn. Apply vacuum, releasing when all liquid has passed through the column.[†]

Washing

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

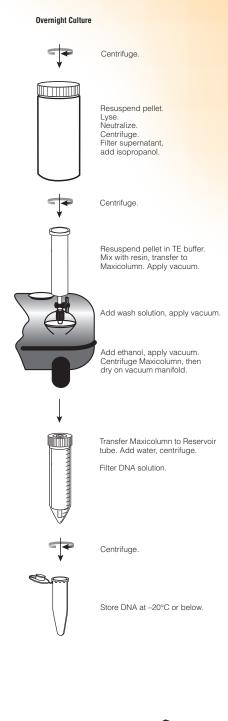
Elution

- 16. Place Maxicolumn in Reservoir tube provided. Add 1.5ml preheated water. Wait 1 minute, then centrifuge at 1,300 × g for 5 minutes to elute DNA. For plasmids \geq 10kb, use water preheated to 70°C; for plasmids \geq 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° C or below.

*Rotor, solutions and centrifugation must be at room temperature. [†]For EndA+ strains and other modifications, additional protocol information in Technical Bulletin #TB139, 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 #9FB018 © 1999, 2009 Promega Corporation. All Rights Reserved.