Wizard® SV 96 Plasmid DNA Purification System

INSTRUCTIONS FOR USE OF PRODUCTS A2250, A2255 AND A2291.



Production of Cleared Lysate and DNA Binding

- 1. Pellet cells at $1,500 \times g$ for 15 minutes.
- 2. Suspend each pellet in 250µl of Cell Resuspension Solution.
- 3. Add 250 μ l of Cell Lysis Solution per well. Tap the plate to mix.

Optional: Add 10µl of Alkaline Protease Solution per well. Mix and incubate 3 minutes.

- 4. Add 350µl of Neutralization Solution per well and mix.
- 5. Transfer samples to Lysate Clearing Plate on manifold. Apply a vacuum. Release vacuum when liquid has cleared the Lysate Clearing Plate. Remove the Clearing Plate.
- 6. Add 500µl of Neutralization Solution per well to the Binding Plate.
- 7. Apply vacuum for 1 minute and release.

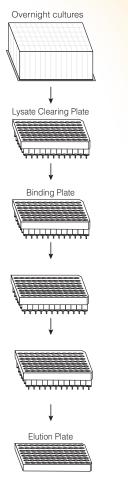
Washing

- 8. Add 1.0ml of Column Wash Solution (CWA) containing ethanol to each well of the Binding Plate. Apply vacuum for 1 minute.
- 9. Release vacuum and repeat Step 8. After second wash continue vacuum for 10 minutes.
- 10. Release vacuum. Connect the vacuum line to the Manifold Collar.

Elution

- 11. Place Elution Plate on Manifold Bed. Add the Manifold Collar.
- 12. Place the Binding Plate on the collar. Add 100µl of water to each well. Wait 1 minute, then apply vacuum for 1 minute.
- 13. Release vacuum, remove Binding Plate and collar. Store DNA at 4°C or -20°C after covering the Elution Plate with a plate sealer.

Additional protocol information is available in Technical Bulletin #TB272, 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 #9FB023

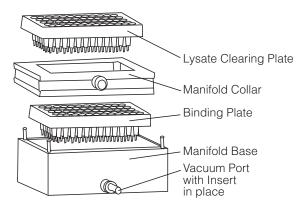
Wizard® SV 96 Plasmid DNA Purification System

INSTRUCTIONS FOR USE OF PRODUCTS A2250, A2255 AND A2291.

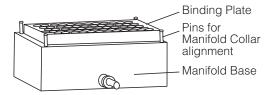


Setup for Vac-Man® 96 Vacuum Manifold

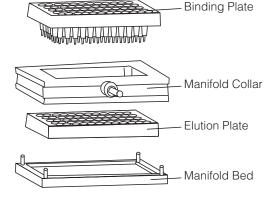
A. Lysate Clearing and DNA Binding Apparatus



B. Washing Apparatus



C. Elution Apparatus



SESEMANE/OA

ORDERING/TECHNICAL INFORMATION:

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

