

Wizard® SV 96 Plasmid DNA Purification System

INSTRUCTIONS FOR USE OF PRODUCTS A2250, A2255 AND A2291.

Quick
PROTOCOL

Production of Cleared Lysate and DNA Binding

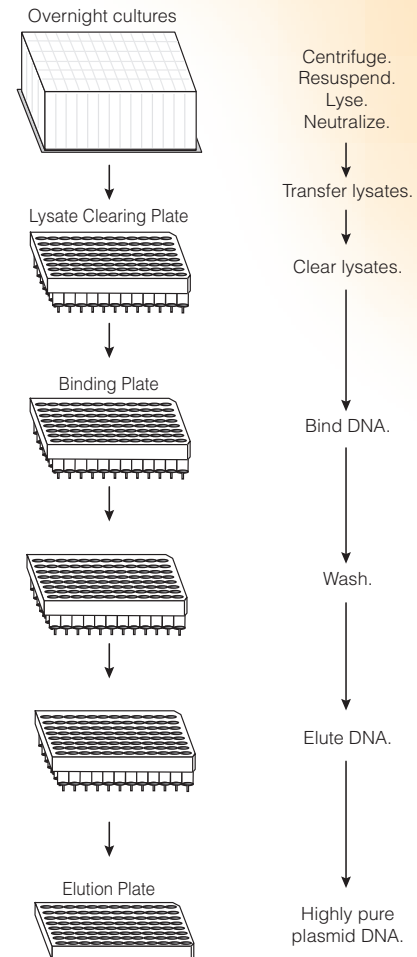
1. Pellet cells at $1,500 \times g$ for 15 minutes.
2. Suspend each pellet in 250 μ l Cell Resuspension Solution.
3. Add 250 μ l Cell Lysis Solution per well. Tap the plate to mix.
Optional: Add 10 μ l Alkaline Protease Solution per well. Mix and incubate 3 minutes.
4. Add 350 μ l Neutralization Solution 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 500 μ l Neutralization Solution per well to the Binding Plate.
7. Apply vacuum for 1 minute and release.

Washing

8. Add 1.0ml Wash Solution 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 100 μ l 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



Promega

Printed in USA. Revised 3/09
Part #9FB023

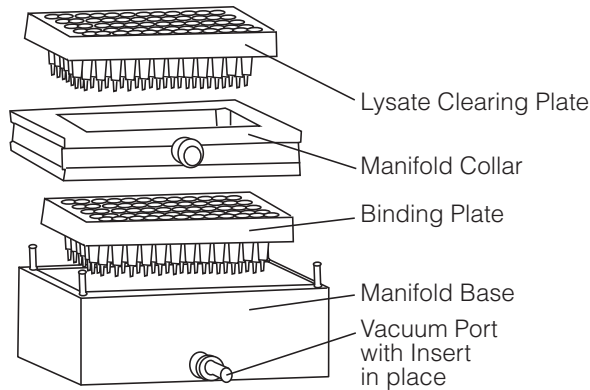
Wizard® SV 96 Plasmid DNA Purification System

INSTRUCTIONS FOR USE OF PRODUCTS A2250, A2255 AND A2291.

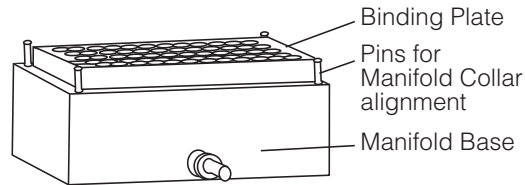
Quick
PROTOCOL

Set-up for Vac-Man® 96 Vacuum Manifold

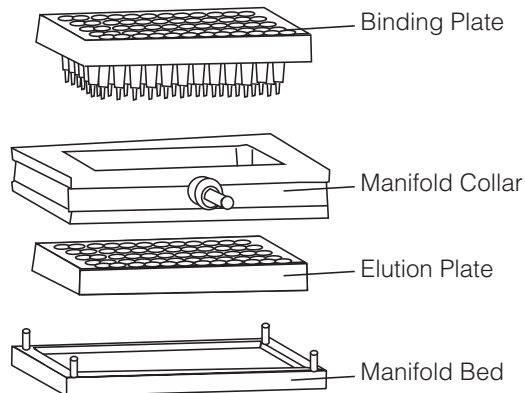
A. Lysate Clearing and DNA Binding Apparatus



B. Washing Apparatus



C. Elution Apparatus



2626MA05/9A

ORDERING/TECHNICAL INFORMATION:

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

© 1999–2009 Promega Corporation. All Rights Reserved.



Promega

Printed in USA. Revised 3/09
Part #9FB023