

Aurum™ Vacuum Manifold: 96-Well Purification



Protocol Overview

For complete protocol, consult instruction manual.

Tubing provided is 4 ft long and must be cut into appropriate pieces before proceeding.

Vacuum Setup Conditions

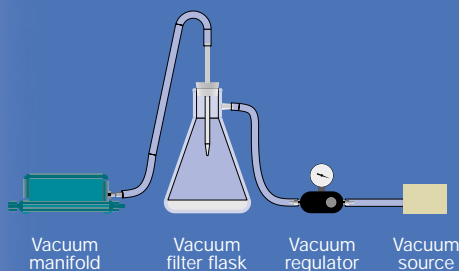
Vacuum source to manifold

1. Use tubing to connect vacuum source to **right side** of vacuum regulator unit.
2. Use tubing to connect **left side** of vacuum regulator to filter flask.

Rubber stopper of vacuum filter flask must contain a hole to accommodate serological (or comparable) pipet.

Vacuum trap must contain additional sidearm for connection to regulator.

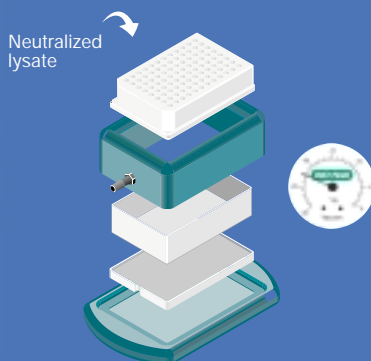
3. Use tubing to connect filter flask to nozzle of manifold top.



Vacuum Purification Conditions

When ready to purify

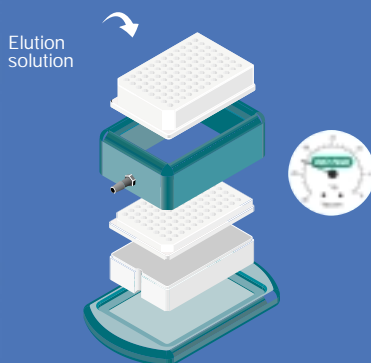
4. Place A stage in center of manifold base.
5. Place waste collection tray on top of A stage.
6. Place manifold top on base to ensure complete contact.
7. Place desired Aurum 96-well plate on manifold top and proceed with appropriate protocol.



Vacuum Elution Conditions

When ready to elute

8. Remove A stage and waste collection tray from manifold.
9. Place B stage in center of manifold base.
10. Place clean 96-well collection plate on top of B stage.
11. Replace manifold top on base to ensure complete contact.
12. Proceed with appropriate elution step as recommended by protocol.



* The Aurum vacuum manifold and stages should be rinsed with water after each use.

BIO-RAD

Aurum Vacuum Manifold: Cat. #732-6470

For more information, call Technical Service at 1-800-4BIORAD (1-800-424-6723). Visit us on the Web at www.bio-rad.com/aurum/

Aurum™ Vacuum Manifold: Column Purification



Protocol Overview

For complete protocol, consult instruction manual.

Tubing provided is 4 ft long and must be cut into appropriate pieces before proceeding.

Vacuum Setup Conditions

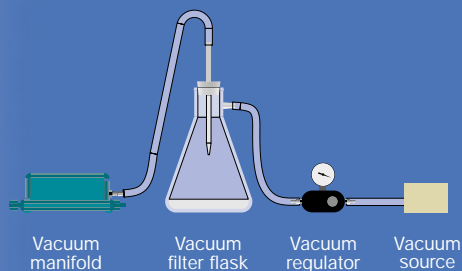
Vacuum source to manifold

1. Use tubing to connect vacuum source to **right side** of vacuum regulator unit.
2. Use tubing to connect **left side** of vacuum regulator to filter flask.

Rubber stopper of vacuum filter flask must contain a hole to accommodate serological (or comparable) pipet.

Vacuum trap must contain additional sidearm for connection to regulator.

3. Use tubing to connect filter flask to nozzle of manifold top.



Vacuum Purification Conditions

When ready to purify

4. Place A stage in center of manifold base.
5. Place waste collection tray on top of A stage.
6. Place manifold top on base to ensure complete contact.
7. Place the column adapter plate (CAP) on manifold top.

Ensure luer is end up.

Ensure unused luer fittings are tightly closed.

8. Place desired Aurum vacuum column(s) into open luer ends.
9. Proceed with appropriate purification protocol for column processing.



Spin Elution Conditions

When ready to elute

10. Proceed with appropriate spin elution step as recommended by protocol.



* The Aurum vacuum manifold and stages should be rinsed with water after each use.

BIO-RAD

Aurum Vacuum Manifold: Cat. #732-6470

For more information, call Technical Service at 1-800-4BIORAD (1-800-424-6723). Visit us on the Web at www.bio-rad.com/aurum/