

Aurum™ Plasmid Mini Kit: Spin Format



Protocol Overview

For complete protocol, consult instruction manual.

Growth and Isolation

1. Grow 1–5 ml bacterial culture overnight or ≥ 16 hr.
2. Measure A_{600} (if higher yield required).
3. Transfer an appropriate volume of culture to a capped 2 ml tube. Centrifuge and decant supernatant.
4. Add 250 μ l resuspension solution; vortex.
5. Add 250 μ l lysis solution; invert 6–8x.
6. Add 350 μ l neutralization solution; invert 6–8x.
7. Centrifuge 5 min to pellet cell debris.

250 μ l
resuspension
solution



Vortex

250 μ l
lysis solution



Invert 6–8x

350 μ l
neutralization
solution



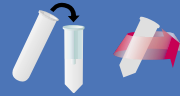
Invert 6–8x



Purification

8. Transfer cleared lysate (supernatant) to mini spin/vac column.
9. Centrifuge 1 min to bind plasmid DNA. Decant flow-through.
10. Add 750 μ l wash solution and centrifuge 1 min. Decant flow-through.
11. Centrifuge additional **1 min** to remove residual wash solution.

Transfer cleared lysate



750 μ l
wash
solution



Collection of Purified Samples

12. Transfer mini spin/vac column to a clean 1.5–2.0 ml capped tube.
13. Add 50 μ l elution solution. Let stand 1 min and then centrifuge 1 min to elute.
14. Purified DNA is ready to use or can be stored at 4°C.

50 μ l
elution
solution



BIO-RAD

Aurum Plasmid Mini Kit: Cat. #732-6400

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

Aurum™ Plasmid Mini Kit: Vacuum Format



Protocol Overview

For complete protocol, consult instruction manual.

Growth and Isolation

1. Grow 1–5 ml bacterial culture overnight or ≥ 16 hr.
2. Measure A_{600} (if higher yield required).
3. Transfer an appropriate volume of culture to a capped 2 ml tube. Centrifuge and decant supernatant.
4. Add 250 μ l resuspension solution; vortex.
5. Add 250 μ l lysis solution; invert 6–8x.
6. Add 350 μ l neutralization solution; invert 6–8x.
7. Centrifuge 5 min to pellet cell debris.

250 μ l
resuspension
solution



Vortex

250 μ l
lysis solution



Invert 6–8x

350 μ l
neutralization
solution



Invert 6–8x



Transfer
cleared lysate

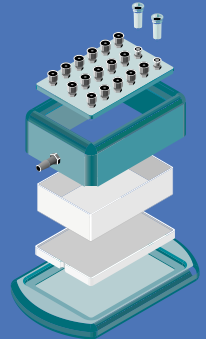


Purification on Aurum or Comparable Manifold

(See exploded view for proper setup of manifold.)

8. Transfer cleared lysate (supernatant) to mini spin/vac column.
9. Apply vacuum at -20 to -23" Hg to bind plasmid DNA. Turn vacuum off.
10. Add 750 μ l wash solution and reapply vacuum until all liquid has passed through column.

750 μ l
wash solution



11. Transfer mini spin/vac column to a 2 ml wash tube. Spin **1 min** to remove residual wash.



Collection of Purified Samples

12. Transfer mini spin/vac column to a clean 1.5–2.0 ml capped tube.
13. Add 50 μ l elution solution. Let stand 1 min and then centrifuge 1 min to elute.
14. Purified DNA is ready to use or can be stored at 4°C.

50 μ l
elution
solution



BIO-RAD

Aurum Plasmid Mini Kit: Cat. #732-6400

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