

# Aurum™ Ion Exchange Mini Kits and Columns

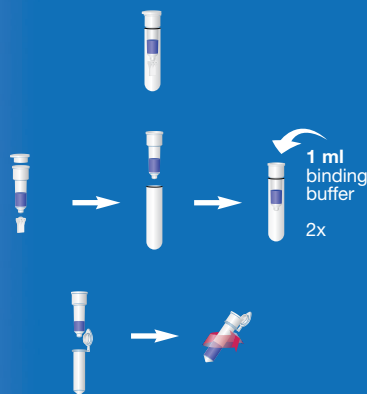


## Protocol Overview

For complete protocol, see instruction manual.

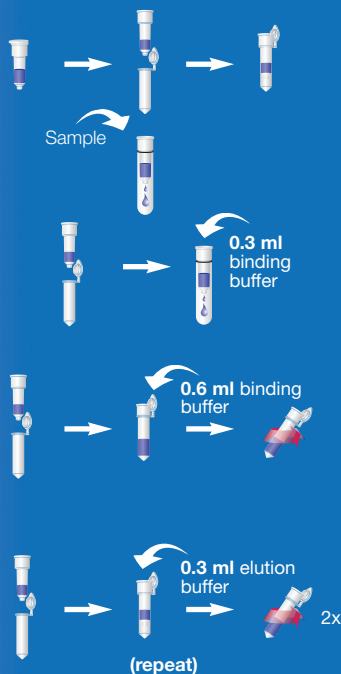
### Column Setup

1. Place AEX or CEX column in a test tube for 5 min to allow resin to settle.
2. Remove cap and break tip from column. Return column to test tube to start gravity flow.
3. Wash column with 1 ml AEX or CEX binding buffer. Repeat.
4. Place column in empty 2.0 ml collection tube and centrifuge for 10 sec at 1,000 x g to dry resin bed. Discard collection tube.



### Sample Binding

5. Place column in a clean 2.0 ml collection tube labeled "unbound".
6. Load sample onto the column and allow it to filter through the column.
7. Place column in another clean 2.0 ml collection tube and wash column with 0.3 ml AEX or CEX binding buffer. Unbound fractions can be discarded or combined for further analysis.
8. Place column in a clean 2.0 ml collection tube. Wash with additional 0.6 ml AEX or CEX binding buffer by centrifuging for 20 sec at 1,000 x g. Discard collection tube.



### Sample Collection

9. Place column in a clean 2.0 ml collection tube labeled "bound #1".
10. Add 0.3 ml AEX or CEX elution buffer to column and centrifuge for 10 sec at 1,000 x g. Repeat.
11. Repeat steps 9 and 10 with a new 2.0 ml collection tube labeled "bound #2".



Aurum Ion Exchange Mini Kits and Columns  
Catalog #732-6705 (AEX) and 732-6702 (CEX)

For more information, call Technical Support at  
1-800-4BIORAD (1-800-424-6723). Visit us on the  
Web at [www.bio-rad.com](http://www.bio-rad.com)