

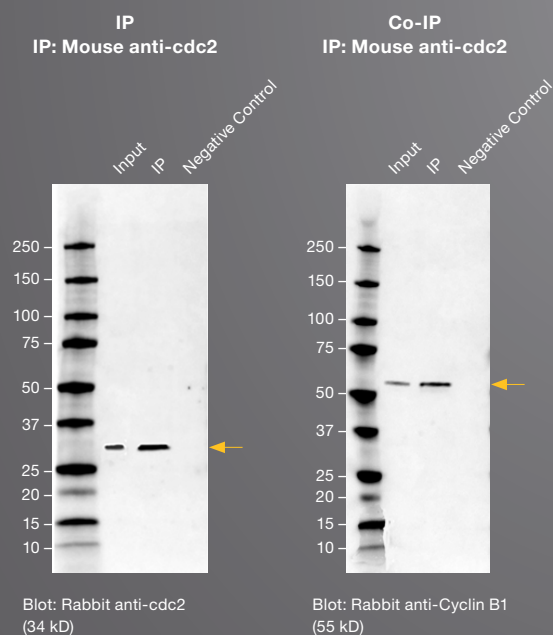


# Six Tips to Improve Your Co-IP Results

## Achieve Successful Co-IP with SureBeads Magnetic Beads

Co-immunoprecipitation (Co-IP) is a good tool for identifying target protein complexes and interacting partners. Though highly informative, Co-IP is also a challenging technique to master. Below are real results from our Co-IP experiments with SureBeads Magnetic Beads.

We've included six useful tips and a protocol to help you see the same success.



**Successful IP and Co-IP results.** HEK293 whole cell lysates were immunoprecipitated with *cdc2* using the SureBeads IP protocol. Cyclin B1 was successfully co-immunoprecipitated with *cdc2*.

## Tips



### 1. Samples

- Select biologically relevant samples that have your target protein complex
- Avoid high concentrations, harsh detergents like SDS, and additives that will cause protein complex dissociation



### 2. Immunoprecipitation

- Maintain protein complexes by using freshly prepared lysates
- Select capture reagent based on the host species of the capture antibody. Protein A has higher affinity for rabbit IgGs. Protein G has higher affinity for mouse IgGs



### 3. Unidirectional Co-IP

- Due to differences in abundance, you may observe Co-IP in one direction, rather than two. When this occurs, immunoprecipitate the component that has the highest percentage of its total population bound to its partner

continues

## Tips



### 4. Other Antibodies

- Try another antibody if you do not observe Co-IP. Only unbound protein will immunoprecipitate if the epitope of the capture antibody overlaps the protein-protein interaction site



### 5. Positive and Negative Controls

- To show that the IP worked, include a blot for just the immunoprecipitated component (positive control)
- To show that the Co-IP worked, include a blot for just the SureBeads Magnetic Beads or IgGs (negative control)



### 6. Analysis

- To avoid antibody detection in the western blot, use a detection antibody from a different host species than the antibody used in the IP



SureBeads Magnetic Rack and SureBeads Magnetic Beads.

## Ordering Information

Catalog #	Description
1614013	<b>SureBeads Protein A Magnetic Beads</b> , 3 ml
1614023	<b>SureBeads Protein G Magnetic Beads</b> , 3 ml
1614916	<b>SureBeads Magnetic Rack</b> , 16-tube holder
1614011	<b>SureBeads Protein A Magnetic Beads</b> , 1 ml
1614021	<b>SureBeads Protein G Magnetic Beads</b> , 1 ml
1614833	<b>SureBeads Starter Kit Protein A and G</b> , Protein A and G beads and magnetic rack
1614813	<b>SureBeads Starter Kit Protein A</b> , two Protein A beads and magnetic rack
1614823	<b>SureBeads Starter Kit Protein G</b> , two Protein G beads and magnetic rack

Visit [bio-rad.com/SureBeadsPubs](http://bio-rad.com/SureBeadsPubs) for publications about SureBeads Magnetic Beads.

Bio-Rad is a trademark of Bio-Rad Laboratories, Inc. in certain jurisdictions. SureBeads are manufactured with magnetic bead technology from JSR Life Sciences. All trademarks used herein are the property of their respective owner.

## Co-IP with SureBeads Magnetic Beads Protocol

### Sample Preparation

- Determine the best cell line for target of interest.
- Grow at least  $3.6 \times 10^7$  cells (two 80–95% confluent T-175 flasks).
- Split cultures 24–48 hours prior to harvest.
- Grow cells to mid-log growth (determine corresponding confluency or density).
- Lyse cells using this lysis buffer:
  - 20 mM Tris (pH 7.5)
  - 150 mM NaCl
  - 1 mM EDTA
  - 1 mM EGTA
  - 1% Triton X-100
  - Add fresh protease inhibitor before use
- IP freshly prepared lysates to avoid complex dissociation.
- IP using SureBeads Magnetic Beads.

### Co-Immunoprecipitation

- Select SureBeads Protein A or Protein G Magnetic Beads appropriate for the antibody used for IP.
- Vortex SureBeads to resuspend them and transfer 100  $\mu$ l (1 mg at 10 mg/ml) of SureBeads to tube.
- Magnetize beads and discard supernatant.
- Wash 3x with 1 ml PBS-T (1x PBS + 0.1% Tween-20).
- Add 7.5  $\mu$ g of antibody. Rotate 20 min at room temperature.
- Wash 3x with 1 ml PBS-T.
- Add 1 mg of antigen-containing lysate. Rotate 1 hr at room temperature.
- Magnetize beads and wash 3x with 1 ml PBS-T.
- Elute with 20  $\mu$ l of 2x Laemmli Buffer. Incubate for 5 min at 90°C.
- Magnetize and transfer supernatant to new vial (IP sample).
- Elute again with 20  $\mu$ l of 2x Laemmli Buffer. Incubate for 5 min at 90°C.
- Magnetize and transfer supernatant to new vial (IP sample).
- Run SDS-PAGE.

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** [bio-rad.com](http://bio-rad.com) **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 01 877 89019 **Belgium** 32 03 710 53 00 **Brazil** 55 11 3065 7550 **Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 36 01 459 6192 **Denmark** 45 04 452 10 00 **Finland** 35 08 980 422 00 **France** 33 01 479 593 00 **Germany** 49 089 3188 4393 **Hong Kong** 852 2789 3300 **Hungary** 36 01 459 6190 **India** 91 124 4029300 **Israel** 972 03 963 6050 **Italy** 39 02 49486600 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 0 318 540 666 **New Zealand** 64 9 415 2280 **Norway** 47 0 233 841 30 **Poland** 36 01 459 6191 **Portugal** 351 21 4727717 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 36 01 459 6193 **Spain** 34 091 49 06 580 **Sweden** 46 08 555 127 00 **Switzerland** 41 0617 17 9555 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 971 4 8187300 **United Kingdom** 44 01923 47 1301

