

Protocol for Protein G Agarose Columns

Catalog Nos.: 53037, 53039

Name: Protein G Agarose Columns

Format: 5 columns, 30 columns

Store at: -20°C

The following protocol is designed to be used with Active Motif's Protein G Agarose Columns for use in chromatin immunoprecipitation (ChIP). Each column is pre-packed with 30 μ l Protein G agarose beads and is ready to use. Simply add your ChIP reaction to the column and proceed with your ChIP incubation and wash steps within the column. The Protein G Agarose Columns can be adapted to work with any available ChIP protocol. The Protein G Agarose is supplied in each column has a binding capacity of 10 μ g IgG/ μ l bead. Columns can be stored at -20°C until ready to use. Do not reuse columns.

Protein G Agarose Columns for ChIP

1. Remove the desired number of columns from the freezer and thaw at room temperature.
2. Briefly spin the column at 1,250 x g for 10 seconds to collect any beads that may have shifted within the column.
3. ChIP reactions containing the chromatin and antibody of interest can be prepared using your desired protocol.
4. Add the ChIP reaction to the Protein G Agarose Column. Cap the column tightly to ensure it is sealed. If desired, use parafilm to seal the cap. (Do not remove the tab from the bottom of the column)
5. Incubate the ChIP reaction in the column according to your protocol.
6. Following the incubation, spin the column at 1,250 x g for 1 minute to collect liquid from inside of the cap.
7. The columns are designed to fit within an empty 1 ml pipet tip box which will serve as a stand for wash steps (see image below). Remove the cap and the tab from the bottom of the column and place the columns in the empty tip box. Add the appropriate wash buffer to the column and allow flow-through to occur by gravity.
8. Proceed with your ChIP protocol.

