



G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ [technical@GBiosciences.com](mailto:technical@GBiosciences.com)

A Geno Technology, Inc. (USA) brand name

# Calmodulin Resin

(Cat. # 786-282, 786-1305)



think proteins! think G-Biosciences [www.GBiosciences.com](http://www.GBiosciences.com)

## INTRODUCTION

Calmodulin is immobilized on 4% agarose using the cyanogen bromide method to make Calmodulin Resin. This resin is ideal for the purification of calmodulin binding proteins that are involved in many biological pathways, including glycogen metabolism, neurotransmission and cytoskeletal control. In addition, a growing use is the isolation of recombinant proteins that are fused to the calmodulin-binding peptide (CBP).

## ITEMS SUPPLIED (Cat. # 786-282)

Cat. #	Description	Size
786-282	Calmodulin Resin*	10ml resin
786-1305	Calmodulin Resin*	50ml resin

\*Calmodulin Resin is supplied as 50% slurry in 20% ethanol

## STORAGE CONDITIONS

It is shipped at ambient temperature. Upon arrival, store it refrigerated at 4°C. **DO NOT FREEZE**. This product is stable for 1 year at 4°C.

## SPECIFICATIONS

Ligand Density: 0.9-1.2mg calmodulin/ml resin  
Binding capacity: 1-3mg/ml (approx.)  
Bead Structure: 4% agarose  
Bead Size: 50-160µm

## ADDITIONAL ITEMS REQUIRED

- Disposable columns
- Binding Buffer: 50mM Tris, pH 7.5, 0.05-0.2M NaCl, 2mM CaCl<sub>2</sub>
- Elution Buffer: 50mM Tris, pH 7.5, 0.05-0.2M NaCl, 2mM EGTA  
**NOTE:** For the binding buffer, the hydrophobic sites are exposed in the presence of Ca<sup>2+</sup>, however in some cases there is an increase in non-specific binding, which can be eliminated by the presence of low salt concentrations (0.05-2.0M NaCl). EDTA may be used in place of EGTA in the elution buffer, however it is less efficient.
- Regeneration reagents: 0.1M NaHCO<sub>3</sub>, pH 8.6, 2mM CaCl<sub>2</sub>; 1M NaCl, 2mM CaCl<sub>2</sub>; 0.1M acetate buffer, pH 4.4, 2mM CaCl<sub>2</sub>; 20% ethanol

## PREPARATION BEFORE USE

*Sample preparation:* Common lysis buffers are compatible with the resin, but must contain 2mM CaCl<sub>2</sub>. The following list is the maximum compatible levels of some common reagents: 50-300mM NaCl/ KCl/ NH<sub>3</sub>SO<sub>4</sub>, 5mM DTT, 10mM β-mercaptoethanol, 0.1% TX-100/ NP-40. Avoid EDTA and EGTA.

## PROTOCOL

1. Add an appropriate amount of calmodulin resin to a suitable tube (suitable to hold 7 columns volumes (CV)). Allow resin to settle and carefully decant the storage buffer.
2. Equilibration Step: Re-suspend the resin in 5CV of binding buffer and allow resin to settle. Decant the supernatant and repeat step 2 once. Finally add an equal volume of binding buffer to the resin.

### Column Method

**NOTE:** Reaction can be performed at 4°C or room temperature. Ensure all reagents and components are at the same temperature.

1. Add 10% volume of binding buffer to the column and pipette in the desired amount of calmodulin resin. Allow the column to drain.
2. Gently load an appropriate volume of sample. Allow column to drain under gravity.
3. Wash Step: Wash the column with 10CV of binding buffer to remove unbound material.
4. Elution Step: Elute the bound proteins in a stepwise manner with 0.5-2ml aliquots of elution buffer.
5. Identify the CBP-tagged protein fractions using a suitable protein assay.

### Batch Binding Method

1. Add the equilibrated Calmodulin Resin directly to the sample lysate and allow binding for several hours to overnight with mechanical rotation at 4°C.
2. After binding, pour the resin into a column and wash with at least 10CV of binding buffer, or until there is no protein in the flow-through (measure absorbance at 280nm or use a protein assay (NI-Protein Assay Cat. # 786-005))
3. Elute the protein with 10CV of elution buffer in a stepwise manner with 0.5-2ml aliquots of elution buffer.
4. Identify the CBP-tagged protein fractions using a suitable protein assay (NI-Protein Assay Cat# 786-005).

## COLUMN REGENERATION

1. Wash the resin with 3CV of 0.1M NaHCO<sub>3</sub>, pH 8.6 containing 2mM EGTA.
2. Wash with 3CV of 1M NaCl containing 2mM CaCl<sub>2</sub>.
3. Wash with 3CV of 0.1M acetate buffer, pH 4.4 containing 2mM CaCl<sub>2</sub>.
4. Wash with binding buffer and store washed resin at 4°C in 20% ethanol.

**NOTE:** Do not regenerate resin more than 3 times.

## RELATED PRODUCTS

Download our Protein Purification Handbook.

<http://info.gbiosciences.com/complete-protein-purification-handbook>

For other related products, visit our website at [www.GBiosciences.com](http://www.GBiosciences.com) or contact us.



[www.GBiosciences.com](http://www.GBiosciences.com)