

I. Introduction

This protocol is provided for the **Capturem Protein G 24-Well Plate - HC** (Cat. Nos. 635744 & 635745), a single-use, disposable 24-well plate for the simple, rapid purification of antibodies from animal sera, ascites fluid, cell culture media, and other sources. Each well can hold up to 4 ml of sample and requires a minimum elution volume of 1 ml. Depending upon the sample type, species, and antibody isotype being purified, the binding capacity of the Capturem membranes will vary. More information about bed volume and capacities can be found on [our website](#).

II. Materials and Reagents

A. Components

Cat. No. 635744:

- 4 x Capturem Protein G 24-Well Plate - HC

Cat. No. 635745:

- 1 x Capturem Protein G 24-Well Plate - HC

B. Additional Materials Required

1. Purification Buffers

This kit is compatible with all standard buffers typically used for antibody purification, such as phosphate- and acetate-based buffers. Loading serum samples without prior dilution is not recommended. Depending on the sample, transiently transfected cell supernatants (e.g., Expi293 HEK cells) may be loaded without prior dilution in binding buffer.

- **Equilibration/Binding:** The following buffers have been tested and shown to perform equally well:
 - **Recommended:** Pierce Protein G IgG Binding Buffer (Thermo Fisher Scientific, Cat. No. 21019)
 - 20 mM sodium phosphate (pH 6.8), 500 mM sodium phosphate, 500 mM NaCl (pH 6.0)
 - PBS: 10 mM Na₂HPO₄, 1.8 mM KH₂PO₄, 2.7 mM KCl, 0.137 M NaCl (pH 7.4)
 - PBS: 10 mM Na₂HPO₄, 1.8 mM KH₂PO₄, 2.7 mM KCl, 0.137 M NaCl (pH 5.0)
 - TBS: 50 mM Tris, 2.7 mM KCl, 0.137 M NaCl (pH 8.0)
 - 100 mM sodium acetate, 2 M NaCl (pH 5.0)
- **Wash Buffer:** Dulbecco's PBS Buffer w/o Ca²⁺, Mg²⁺ (pH 7.5)
- **Elution Buffer:** 0.1 M glycine (pH 2.5–3.0) or 0.1 M citric acid (pH 2.5–3.0)
- **Neutralization Buffer:** 1 M Tris (pH 8.5)

2. Collection Plates

- **24-well collection plates:** Each purification will require four standard 24-well collection plates (e.g., Thomas Scientific, Cat. No. 1149Q53, 24-Well Polypropylene Storage/Reaction Microplate, 10mL/square well, round bottoms, 44mm height). These plates should be used throughout the protocol to collect flowthrough, wash, and elution samples that will be saved for SDS-PAGE analysis and/or colorimetric protein assays (e.g., Bradford assays). Alternatively, if only the eluates are to be analyzed, the flowthrough from the equilibration and wash steps can be discarded. In this case, only two collection plates are required, one for collecting the flowthrough and washes and a second for the actual eluates.

- **Recommended filtration device:** The Capturem Protein G 24-Well Plate - HC is compatible with standard 96-well plate filtration devices, such as the NucleoVac 96 Vacuum Manifold (Cat. No. 740681).
- **Recommended plate centrifuge:** The Capturem Protein G 24-Well Plate - HC is compatible with standard 96-well plate centrifuges (e.g., Eppendorf Centrifuge 5804 R with deepwell-plate rotor, A-2-DWP).
- **Recommended positive pressure device:** The Capturem Protein G 24-Well Plate - HC is compatible with standard 96-well positive pressure devices (e.g., Waters Positive Pressure-96 Processor, Cat. No. 186006961).
- **Single-channel pipettes** (various).

III. Sample Preparation

1. Follow the standard protocols for preparing antibody samples. We recommend diluting the antibody sample in the range of 1:2–1:50 sample to Equilibration/Binding Buffer (preferably Pierce Protein G IgG Binding Buffer; see Section II.B.1 for other buffer options). We recommend 1:2–1:4 dilution for hybridoma samples; and 1:15 dilution for serum samples, which may differ depending on animal species and may require further optimization for obtaining high yield and purity. Loading the sample (e.g., serum matrix) without prior dilution is not recommended. Loading high-density cell culture supernatants (e.g., Expi293 HEK cells) directly without dilution may work from case to case but could lead to higher well-to-well variability.
2. Samples with visible precipitates must be clarified by centrifugation or filtered through a 0.5-micron filter before loading into the plate.

IV. Antibody Purification

Purification of antibodies from the samples prepared in Section III may be performed using vacuum filtration (Protocol A), centrifugation in a standard 96-well plate centrifuge (Protocol B), or using positive pressure (Protocol C).

A. Protocol: Antibody Purification Using Vacuum Filtration

1. Assemble the Capturem Protein G 24-Well Plate - HC (“Capturem plate” thereafter) in a filtration device together with a 24-well collection plate (supplied by the user—see Section II.B).
2. Add 2 ml of Equilibration/Binding Buffer to each well of the Capturem plate to equilibrate the wells. We recommend Pierce Protein G IgG Binding Buffer (Thermo Fisher Scientific). Use a vacuum (~0.3–0.4 bar) to run the buffer through the Capturem plate wells into the collection plate. Remove the flowthrough and discard it along with the collection plate. Reassemble the Capturem plate in the filtration device together with a new collection plate.

NOTE: The required pressure may vary between 0.2–0.8 bar, depending on the sample. When performing vacuum filtration, use the following formula to convert between different vacuum pressure units:

Pressure Conversion Formula:

$$1 \text{ mbar} = 100 \text{ Pa} = 0.750 \text{ mm Hg} = 14.504 \times 10^{-3} \text{ psi} = 0.987 \times 10^{-3} \text{ atm}$$

3. Load 0.5–4.3 ml of diluted antibody (from Section III) into each well of the equilibrated Capturem plate. Use a vacuum (~0.3–0.4 bar) to run the buffer through the Capturem plate wells into the collection plate. Save the collection plate containing the flowthrough for antibody analysis. Reassemble the Capturem plate in the filtration device together with a new collection plate.
4. Add 4 ml of Wash Buffer (PBS Buffer w/o Ca²⁺, Mg²⁺, pH 7.5) to each well of the Capturem plate. Use a vacuum (~0.3–0.4 bar) to run the buffer through the Capturem plate wells into the collection plate. Save the collection plate containing the flowthrough for antibody analysis.
5. Repeat Step 4 and perform a second wash step with PBS buffer. Use a vacuum (~0.3–0.4 bar) to run the buffer through the Capturem plate wells into the collection plate. Save the collection plate containing the flowthrough for antibody analysis.
6. Elute your antibody from the Capturem plate as follows:
 - If using citric acid:**
 - a. Add 500 µl of Neutralization Buffer to each well of a new collection plate and place the Capturem plate securely on top of the collection plate for the first elution.
 - b. Add 900 µl of 0.1 M citric acid to each well of the Capturem plate. Use a vacuum (~0.3–0.4 bar) to run the buffer through the Capturem plate wells into the collection plate.
 - c. Perform a second elution by adding 900 µl of 0.1 M citric acid to each well of the Capturem plate. Use a vacuum (~0.3–0.4 bar) to run the buffer through the Capturem plate wells into the collection plate. The wells of the collection plate now contain your eluted antibodies.
 - If using glycine:**
 - d. Add 180 µl of Neutralization Buffer to each well of a new collection plate and place the Capturem plate securely on top of the collection plate for the first elution.
 - e. Add 900 µl of 0.1 M glycine to each well of the Capturem plate. Use a vacuum (~0.3–0.4 bar) to run the buffer through the Capturem plate wells into the collection plate.
 - f. Perform a second elution by adding 900 µl of 0.1 M glycine to each well of the Capturem plate. Use a vacuum (~0.3–0.4 bar) to run the buffer through the Capturem plate wells into the collection plate.
7. Measure the amount of antibody in the eluate using UV absorbance at 280 nm or another colorimetric protein analysis method. One OD₂₈₀ unit typically corresponds to 0.73 mg/ml of IgG. The purity of the eluted antibodies can be analyzed by SDS-PAGE, size exclusion chromatography, etc.

B. Protocol: Antibody Purification Using Centrifugation

1. Place the Capturem Protein G 24-Well Plate - HC (“Capturem plate” thereafter) securely on top of a 24-well collection plate (supplied by the user—see Section II.B).
2. Add 2 ml of Equilibration/Binding Buffer, preferably Pierce Protein G IgG Binding Buffer (Thermo Fisher Scientific), to each well of the Capturem plate to equilibrate the wells. Centrifuge at 2,000g for 4 min at room temperature. Make sure wells are drained, remove the flowthrough and discard it along with the collection plate. Place the Capturem plate securely on top of a new 24-well collection plate.

NOTES:

- If wells did not drain fully, centrifuge again at 2,000g for 4 min.
- When performing centrifugation, use the following formula to convert rpm to centrifugal force units (RCF or g) if the centrifuge does not automatically provide this information:

Centrifugal Force Conversion Formula: $RCF \text{ or } g = (1.12) \times (R) \times (\text{rpm}/1,000)^2$

where R is the distance (in millimeters) from the center of the rotor to the end of the spin bucket when held horizontally and away from the center.

3. Load 0.5–4.3 ml of diluted antibody (from Section III) into each well of the equilibrated Capturem plate. Centrifuge at 2,000g for 4 min at room temperature. Save the collection plate containing the flowthrough for antibody analysis. Place the Capturem plate securely on top of a new 24-well collection plate.
4. Add 4 ml of Wash Buffer (PBS Buffer w/o Ca^{2+} , Mg^{2+} , pH 7.5) to each well of the Capturem plate. Centrifuge at 2,000g for 4 min at room temperature. Save the collection plate containing the flowthrough for antibody analysis.
5. Wash a second time by adding 4 ml of Wash Buffer to each well of the Capturem plate. Centrifuge at 2,000g for 4 min at room temperature. Save the collection plate containing the flowthrough for antibody analysis.
6. Elute your antibody from the Capturem plate as follows:

If using citric acid:

- a. Add 500 μl of Neutralization Buffer to each well of a new collection plate and place the Capturem plate securely on top of the collection plate for the first elution.
- b. Add 900 μl of 0.1 M citric acid to each well of the Capturem plate. Centrifuge at 2,000g for 4 min at room temperature.
- c. Perform a second elution by adding 900 μl of 0.1 M citric acid to each well of the Capturem plate. Centrifuge at 2,000g for 4 min at room temperature. The wells of the collection plate now contain your eluted antibodies.

If using glycine:

- d. Add 180 μl of Neutralization Buffer to each well of a new collection plate and place the Capturem plate securely on top of the collection plate for the first elution.
- e. Add 900 μl of 0.1 M glycine to each well of the Capturem plate. Centrifuge at 2,000g for 4 min at room temperature.

- f. Perform a second elution by adding 900 µl of 0.1 M glycine to each well of the Capturem plate. Centrifuge at 2,000g for 4 min at room temperature. The wells of the collection plate now contain your eluted antibodies.
7. Measure the amount of antibody in the eluate using UV absorbance at 280 nm or another colorimetric protein analysis method. One OD₂₈₀ unit typically corresponds to 0.73 mg/ml of IgG. The purity of the eluted antibodies can be analyzed by SDS-PAGE, size exclusion chromatography, etc.

C. Protocol: Antibody Purification Using Positive Pressure

1. Place the Capturem Protein G 24-Well Plate - HC (“Capturem plate” thereafter) securely on top of a 24-well collection plate (supplied by the user—see Section II.B).
2. Add 2 ml of Equilibration/Binding Buffer, to each well of the Capturem plate to equilibrate the wells. Apply positive pressure of 10 psi at room temperature until all liquid has passed through the wells. Make sure wells are drained completely, remove the flowthrough and discard it along with the collection plate. Place the Capturem plate securely on top of a new 24-well collection plate.

NOTE: If wells did not drain fully, apply positive pressure again at 10 psi.

3. Load 0.5–4.3 ml of diluted antibody (from Section III) into each well of the equilibrated Capturem plate and apply 10 psi at room temperature. Save the collection plate containing the flowthrough for antibody analysis. Place the Capturem plate securely on top of a new 24-well collection plate.
4. Add 4 ml of Wash Buffer (PBS Buffer w/o Ca²⁺, Mg²⁺, pH 7.5) to each well of the Capturem plate and apply 10 psi at room temperature. Save the collection plate containing the flowthrough for antibody analysis.
5. Wash a second time by adding 4 ml of Wash Buffer to each well of the Capturem plate and apply 10 psi at room temperature. Save the collection plate containing the flowthrough for antibody analysis.
6. Elute your antibody from the Capturem plate as follows:

If using citric acid:

- a. Add 500 µl of Neutralization Buffer to each well of a new collection plate and place the Capturem plate securely on top of the collection plate for the first elution.
- b. Add 900 µl of 0.1 M citric acid to each well of the Capturem plate. Apply 10 psi at room temperature to elute.
- c. Add 900 µl of 0.1 M citric acid to each well of the Capturem plate. Apply 10 psi at room temperature to elute. The wells of the collection plate now contain your eluted antibodies.

If using glycine:

- d. Add 180 µl of Neutralization Buffer to each well of a new collection plate and place the Capturem plate securely on top of the collection plate for the first elution.
- e. Add 900 µl of 0.1 M glycine to each well of the Capturem plate. Apply 10 psi at room temperature to elute.
- f. Add 900 µl of 0.1 M glycine to each well of the Capturem plate. Apply 10 psi at room temperature to elute. The wells of the collection plate now contain your eluted antibodies.

Capturem™ Protein G 24-Well Plate - HC Protocol-At-A-Glance

7. Measure the amount of antibody in the eluate using UV absorbance at 280 nm or another colorimetric protein analysis method. One OD₂₈₀ unit typically corresponds to 0.73 mg/ml of IgG. The purity of the eluted antibodies can be analyzed by SDS-PAGE, size exclusion chromatography, etc.

Appendix A. Troubleshooting Guide

Table 1. Troubleshooting Guide

Problem	Possible Explanation	Solution
Background bands/poor purity	Nonspecific binding of proteins to the membrane	Add an additional wash step with Wash Buffer (PBS) or TBS
Minimal recovery	The sample contains more antibody than the Protein G 24-Well Plate - HC has capacity for.	Reduce the amount of sample added to each well. If you need to purify more antibody, consider splitting your sample into two separate 24 wells or use a Capturem Protein G Maxiprep column which has a higher capacity (1 mg).
Low yield of purified antibody	Poor binding affinity of the IgG subtype used. Different subtype classes and species bind differently to Protein G. Please, refer to Roque, A. C. A., Silva, C. S. O. & Taipa, M. A. Affinity-based methodologies and ligands for antibody purification: advances and perspectives. <i>J. Chromatogr. A</i> 1160 , 44–55 (2007).	Verify that Protein G-based purification is suitable for the IgG subtype class and species used. Alternatively, consider using Protein A-based purification instead.
Low yield of purified antibody	pH of the binding buffer was altered by the hybridoma, serum, or cell culture medium used.	Check and make sure the pH of the buffer is in an optimal range for your IgG sample.
Low yield of purified antibody	Suboptimal dilution	Make sure to use the optimal dilution of your sample. For hybridoma medium, 1:2 to 1:4 were found to work well, but it is sample dependent.
The antibody does not elute	Elution conditions too mild	<ul style="list-style-type: none"> Follow the instructions use recommended elution buffer (0.1 M glycine, pH 2.5–3.0) Make sure the pH of the elution buffer is 2.5-3.0 if you are using a different buffer than the one recommended.
Well does not fully drain	Clogging due to particles or very viscous sample	<ul style="list-style-type: none"> Preclarify the solution by centrifugation or using a 0.5-micron filter Repeat spin at 2,000g for 4 min Reapply positive pressure at 10 psi.

Capturem™ Protein G 24-Well Plate - HC Protocol-At-A-Glance

Contact Us	
Customer Service/Ordering	Technical Support
tel: 800.662.2566 (toll-free)	tel: 800.662.2566 (toll-free)
fax: 800.424.1350 (toll-free)	fax: 800.424.1350 (toll-free)
web: takarabio.com	web: takarabio.com
e-mail: ordersUS@takarabio.com	e-mail: techUS@takarabio.com

Notice to Purchaser

Our products are to be used for research purposes only. They may not be used for any other purpose, including, but not limited to, use in drugs, *in vitro* diagnostic purposes, therapeutics, or in humans. Our products may not be transferred to third parties, resold, modified for resale, or used to manufacture commercial products or to provide a service to third parties without prior written approval of Takara Bio USA, Inc.

Your use of this product is also subject to compliance with any applicable licensing requirements described on the product's web page at takarabio.com. It is your responsibility to review, understand and adhere to any restrictions imposed by such statements.

© 2019 Takara Bio Inc. All Rights Reserved.

All trademarks are the property of Takara Bio Inc. or its affiliate(s) in the U.S. and/or other countries or their respective owners. Certain trademarks may not be registered in all jurisdictions. Additional product, intellectual property, and restricted use information is available at takarabio.com.

This document has been reviewed and approved by the Quality Department.