

Pierce™ Protein A Magnetic Beads

Pub. No. MAN0011856

Rev. B.0

Pub Part No. 2162512.3

88845 88846

Number	Description
88845	Pierce Protein A Magnetic Beads , 1mL, supplied at 10mg/mL in PBS containing 0.05% Tween™-20 Detergent and 0.05% NaN ₃
88846	Pierce Protein A Magnetic Beads , 5mL, supplied at 10mg/mL in PBS containing 0.05% Tween-20 Detergent and 0.05% NaN ₃

Storage: Upon receipt store at 4°C. Product shipped with an ice pack.

Table of Contents

Introduction	1
Important Product Information	2
Procedure for Manual Antibody Purification.....	2
Procedure for Automated Antibody Purification.....	3
Procedure for Manual Immunoprecipitation.....	4
Procedure for Automated Immunoprecipitation	5
Troubleshooting.....	6
Additional Information Available on Our Website.....	7
Frequently Asked Questions for the KingFisher Instrument	7
Related Thermo Scientific Products	7

Introduction

The Thermo Scientific™ Pierce™ Magnetic Beads provide a fast and convenient method for both manual and automated magnetic isolation of proteins using affinity binding. Pierce Protein A Magnetic Beads (Table 1) are typically used for isolating antibodies from serum, cell culture supernatant or ascites and for immunoprecipitation and co-immunoprecipitation of antigens from cell or tissue extracts. For antibody purification, the beads are incubated with the antibody solution and then magnetically separated from the supernatant. For immunoprecipitation, the beads are added to an antigen-containing sample to which antibody has been added and allowed to incubate to form the antibody-antigen complex. The bound antibodies or antigens are dissociated from the beads using an elution buffer and removed from the solution manually using a magnetic stand or by automation using an instrument such as the Thermo Scientific™ KingFisher™ Flex or KingFisher Duo Instrument. Automated instruments are especially useful for large-scale screening of multiple samples.

The Pierce Protein A Magnetic Beads contain a recombinant Protein A (molecular weight 44.7kDa; apparent molecular weight by SDS-PAGE ~47kDa) with five Fc-binding domains per protein. These domains can bind to antibodies from many different species, including mouse, human, rabbit, pig, dog and cat. The proprietary blocking agent on the magnetic beads minimizes or eliminates nonspecific binding on the bead surface when working with complex biological samples. Detailed instructions and optimized buffer components are provided for best results.

Table 1. Characteristics of the Thermo Scientific Pierce Protein A Magnetic Beads.

Composition:	Recombinant Protein A monolayer covalently coupled to a blocked magnetic bead surface
Magnetization:	Superparamagnetic (no magnetic memory)
Mean Diameter:	1 μm (nominal)
Density:	2.0g/cm ³
Bead Concentration:	10mg/mL
Binding Capacity:	≥ 40μg rabbit IgG/mg of bead

Important Product Information

- Do not centrifuge, dry or freeze the Pierce Magnetic Beads. Centrifuging, drying or freezing will cause the beads to aggregate and lose binding activity. To ensure good dispersal of beads for optimal antibody binding, it is important to include 0.025% to 0.1% non-ionic (e.g., Tween-20 Detergent) or zwitterionic (e.g., CHAPS) detergent in the binding buffer and mix the beads during incubation.
- To minimize protein degradation, include protease inhibitors (e.g., Thermo Scientific Halt Protease Inhibitor Single-Use Cocktail, EDTA-free, Product No. 78425) in preparation of cell lysates.
- A low-pH elution may be used for single-use applications. Optimal time for low-pH elution is 10 minutes; exceeding 10 minutes may result in nonspecific binding and yield reduction.
- When using rabbit antibodies (primary or secondary) in downstream Western blot applications, perform elution in SDS-PAGE sample buffer at room temperature. For all other antibody species, boiling the beads in SDS-PAGE sample buffer is acceptable for single-use applications.

Note: Boiling the magnetic beads can cause bead aggregation and loss of binding activity.

- Pierce Protein A Magnetic Beads are compatible with small-scale antibody purification and immunoprecipitation and analyses by Western blot.
- Protein A is preferred over Protein G for the isolation of polyclonal IgG from rabbit, pig, dog and cat serum. Protein A is not recommended for the isolation of mouse IgG₁; human IgG₃; or rat, goat and cow antibodies because of weak binding. For more information, see Tech Tip #34: Binding characteristics of Protein A, Protein G, Protein A/G and Protein L from our website.

Procedure for Manual Antibody Purification

A. Additional Materials Required

- 1.5mL microcentrifuge tubes
 - Sample: Serum, concentrated cell culture supernatant or concentrated ascites
- Note:** Samples can be concentrated using the Pierce Concentrators 20mL/30K, Product No. 88529 or 88531.
- Binding/Wash Buffer: Tris-buffered saline (TBS, Product No. 28360) containing 0.05% Tween-20 Detergent
 - Elution Buffer: IgG Elution Buffer (Product No. 21028) or 0.1M glycine, pH 2.0
 - Neutralization Buffer: 1M Tris, pH 8.5
 - Magnetic stand (e.g., Thermo Scientific™ DynaMag™-2 Magnet, Product No. 12321D)

B. Antibody Purification from Serum, Cell Culture Supernatant or Ascites

Note: To ensure homogeneity, mix the beads thoroughly before use by repeated inversion, gentle vortexing or using a rotating platform.

1. Place 50µL (0.50mg) of Pierce Protein A Magnetic Beads into a 1.5mL microcentrifuge tube. Add 150µL of Binding/Wash Buffer to the beads and gently vortex to mix.
2. Place the tube into a magnetic stand to collect the beads against the side of the tube. Remove and discard the supernatant.
3. Add 1mL of Binding/Wash Buffer to the tube. Invert the tube several times or gently vortex to mix for 1 minute. Collect beads with magnetic stand, then remove and discard the supernatant.
4. Dilute 10µL of sample with 490µL of Binding/Wash Buffer.

Note: Sample volume can be modified according to user preference. If the sample volume is < 500µL, dilute it to a final volume of 500µL with Binding/Wash Buffer.

5. Add the diluted sample to the tube containing pre-washed magnetic beads and gently vortex or invert to mix.
6. Incubate the samples at room temperature with mixing for 1 hour.
7. Collect the beads with a magnetic stand, then remove and discard the supernatant.

8. Add 500µL of Binding/Wash Buffer to the tube, mix well, collect the beads with a magnetic stand and discard the supernatant. Repeat this wash twice.
9. Add 100µL of Elution Buffer to the tube, mix well and incubate 10 minutes at room temperature with occasional mixing.
10. Collect the beads with a magnetic stand and then remove and save the supernatant that contains the eluted antibody. To neutralize the low pH, add 15µL of Neutralization Buffer for each 100µL of eluate.

Note: The minimum volume of beads recommended for antibody purification is 50µL.

Procedure for Automated Antibody Purification

A. Additional Materials Required

- KingFisher Flex System with 96 deep well head (Product No. 5400630) or KingFisher 96 System (Product No. 5400500)
- Thermo Scientific™ Microtiter Deep Well 96 Plate, V-bottom, polypropylene (100-1000µL; Product No. 95040450)
- KingFisher Flex 96 Tip Comb for Deep Well Magnets (Product No. 97002534)
- Binding/Wash Buffer: Tris-buffered saline (TBS, Product No. 28360) containing 0.05% Tween-20 Detergent
- Elution Buffer: IgG Elution Buffer (Product No. 21028) or 0.1M glycine, pH 2.0
- Neutralization Buffer: 1M Tris, pH 8.5

B. Preparation of Instrument and Plate Set-Up

Note: The following protocol is designed for use with the KingFisher Flex Instrument. The protocol can be modified according to customer needs using the Thermo Scientific™ BindIt™ Software provided with the instrument.

1. Download the “Antibody Purification” protocol from the Thermo Scientific website (thermofisher.com/bindit-protocols) into the BindIt Software on an external computer.
2. Transfer the protocol to the KingFisher Flex Instrument from an external computer. See the BindIt Software User Manual for detailed instructions on importing protocols.
3. Set up the plates according to Table 2.

Table 2. Pipetting instructions for the antibody purification protocol using the Thermo Scientific Microtiter Deep Well 96 Plates.

Plate #	Plate Name	Content	Volume
1	Beads	Protein A beads	50µL
		Binding/Wash Buffer	150µL
2	Bead Wash	Binding/Wash Buffer	1000µL
3	Bind	Sample	10µL
		Binding/Wash Buffer	490µL
4	Wash 1	Binding/Wash Buffer	500µL
5	Wash 2	Binding/Wash Buffer	500µL
6	Wash 3	Water	500µL
7	Elution	Elution Buffer	100µL
8	Tip Plate	KingFisher Flex 96 Tip Comb for Deep Well Magnets	-

Notes:

- If using less than 96 wells, fill the same wells in each plate. For example, if using wells A1 through A12, use these same wells in all plates.
- To ensure bead homogeneity, mix the vial thoroughly by repeated inversion, gentle vortexing or rotating platform before adding the beads to Plate 1.
- Combine the Tip Comb with a Deep Well 96 Plate. See the KingFisher Instrument user manual for detailed instructions.

-
- Sample volume can be modified according to user preference. If the sample volume is < 500 μ L, dilute to a final volume of 500 μ L with Binding/Wash Buffer.

C. Executing the Antibody Purification Protocol on the KingFisher Flex Instrument

1. Select the protocol using the arrows on the instrument keypad and press Start. See the KingFisher Flex Instrument User Manual for detailed information.
2. Slide open the door of the instrument's protective cover.
3. Load the plates into the KingFisher Flex Instrument according to the protocol request, placing each plate in the same orientation. Confirm each action by pressing Start.
4. After sample processing, remove plates as instructed by the instrument's display. Press Start after removing each plate. Press Stop after all plates are removed. Upon completion, if desired, neutralize the low pH by adding 15 μ L of Neutralization Buffer for each 100 μ L of eluate.

Procedure for Manual Immunoprecipitation

A. Additional Materials Required

- 1.5mL microcentrifuge tubes
- Binding Buffer: Buffer used to prepare antigen sample (e.g., Thermo Scientific Pierce IP Lysis Buffer, Product No. 87788)
- Wash Buffer: Tris-buffered saline (TBS, Product No. 28360) containing 0.05% Tween-20 Detergent and 0.5M NaCl
- Elution Buffer: IgG Elution Buffer (Product No. 21028) or 0.1M glycine, pH 2.0
- Neutralization Buffer: 1M Tris, pH 8.5
- Cell Lysis Buffer (used to adjust IP reaction volume)
- Alternative Elution Buffer: SDS-PAGE reducing sample buffer
- Antibody for immunoprecipitation
- Antigen sample
- Magnetic stand (e.g., Thermo Scientific™ DynaMag™-2 Magnet; Product No. 12321D)

B. Immunoprecipitation

Note: This protocol is a general guideline for immunoprecipitation and will require optimization for each application.

1. Combine the antigen sample with 5-10 μ g of antibody. Adjust the reaction volume to 500 μ L with the Cell Lysis Buffer. Incubate the reaction for 1-2 hours at room temperature or overnight at 4°C with mixing.
2. Place 25 μ L (0.25mg) of Pierce Protein A Magnetic Beads into a 1.5mL microcentrifuge tube.
3. Add 175 μ L of Wash Buffer to the beads and gently vortex to mix.
4. Place the tube into a magnetic stand to collect the beads against the side of the tube. Remove and discard the supernatant.
5. Add 1mL of Wash Buffer to the tube. Invert the tube several times or gently vortex to mix for 1 minute. Collect beads with magnetic stand. Remove and discard the supernatant.
6. Add the antigen sample/antibody mixture to a 1.5mL microcentrifuge tube containing pre-washed magnetic beads and incubate at room temperature for 1 hour with mixing.
7. Collect the beads with a magnetic stand and then remove the flow-through and save for analysis.
8. Add 500 μ L of Wash Buffer to the tube and gently mix. Collect the beads and discard the supernatant. Repeat wash twice.
9. Add 500 μ L of ultrapure water to the tube and gently mix. Collect the beads on a magnetic stand and discard the supernatant.

10. Elution: Add 100µL of Elution Buffer to the tube. Incubate the tube at room temperature with mixing for 10 minutes. Magnetically separate the beads and save the supernatant containing the target antigen. To neutralize the low pH, add 15µL of Neutralization Buffer for each 100µL of eluate.

Alternative Elution: Add 100µL of SDS-PAGE reducing sample buffer to the tube and heat samples at 96-100°C in a heating block for 10 minutes. Magnetically separate the beads and save the supernatant containing the target antigen.

Note: If you will be performing a Western blot using rabbit antibodies (primary or secondary) do not heat the samples. Incubate at room temperature for 10 minutes with mixing.

Procedure for Automated Immunoprecipitation

A. Additional Materials Required

- KingFisher Flex System with 96 deep well head (Product No. 5400630)
- Microtiter Deep Well 96 Plate, V-bottom, polypropylene (100-1000µL; Product No. 95040450)
- KingFisher Flex 96 Tip Comb for Deep Well Magnets (Product No. 97002534)
- 1.5mL microcentrifuge tubes
- Binding Buffer: Buffer used to prepare antigen sample (e.g., Pierce IP Lysis Buffer, Product No. 87788)
- Wash Buffer: Tris-buffered saline (TBS, Product No. 28360) containing 0.05% Tween-20 Detergent and 0.5M NaCl
- Elution Buffer: IgG Elution Buffer (Product No. 21028) or 0.1M glycine, pH 2.0
- Alternative Elution Buffer: SDS-PAGE reducing sample buffer
- Neutralization Buffer: 1M Tris, pH 8.5
- Antibody for immunoprecipitation
- Antigen sample
- Cell Lysis Buffer (used to prepare the antigen sample)

B. Instrument Preparation and Plate Set-Up

Note: The following protocol is designed for general use with the KingFisher Flex Instrument. The protocol can be modified according to your needs using the BindIt Software provided with the instrument.

1. Combine antigen sample with 5-10µg of immunoprecipitation antibody per sample. Incubate 1-2 hours at room temperature or overnight at 4°C with mixing.
2. Download the “Immunoprecipitation” protocol from the Thermo Fisher Scientific website (thermofisher.com/bindit-protocols) into the BindIt™ Software on an external computer.
3. Transfer the protocol to the KingFisher Flex Instrument from an external computer. See BindIt Software User Manual for detailed instructions on importing protocols.
4. Set up plates according to Table 3.

Table 3. Pipetting instructions for the immunoprecipitation protocol using the Thermo Scientific Microtiter Deep Well 96 Plates.

Plate #	Plate Name	Content	Volume	Time/Speed
1	Beads	Protein A Beads	25µL	5 seconds
		Binding Buffer	175µL	
2	Bead Wash	Binding Buffer	1000µL	1 minute/Slow
3	Bind	Antibody/Antigen Sample	500µL	1 hour/Slow
4	Wash 1	Wash Buffer	500µL	30 seconds/Slow
5	Wash 2	Wash Buffer	500µL	30 seconds/Slow
6	Wash 3	Ultrapure Water	500µL	30 seconds/Slow
7	Elution	Elution Buffer	100µL	10 minutes/Medium
8	Tip Plate	KingFisher Flex 96 Tip Comb for Deep Well Magnets	–	10 seconds/Fast

Notes:

- If less than 96 wells are used, fill the same wells in each plate. For example, if using wells A1 through A12, use these same wells in all plates.
- To ensure bead homogeneity, mix the vial thoroughly by repeated inversion, gentle vortexing or rotating platform before adding the beads to Plate 1.
- Combine the Tip Comb with a Deep Well 96 Plate. See the instrument user manual for detailed instructions.
- The beads can be eluted into 100µL of 0.1M glycine, pH 2.0 or 100µL SDS-PAGE reducing sample buffer. If using SDS-PAGE reducing sample buffer in a heated elution, install the KingFisher Flex Heating Block (see manual for proper installation) to heat samples at 96-100°C for 10 minutes.
- If you select SDS-PAGE reducing sample buffer for elution and will be performing a Western blot using rabbit antibodies (primary or secondary) **do not** heat the samples. Incubate at room temperature for 10 minutes.
- If low-pH elution buffer is selected for elution, neutralize the pH using 15µL of Neutralization Buffer for each 100µL of eluate upon run completion.
- To limit evaporation, select “Mix” and “Slow” speed under the subheading “Heating Action.”

C. Executing Automated Immunoprecipitation Protocol

1. Select the protocol using the arrow keys on the instrument keypad and press Start. See the KingFisher Flex Instrument User Manual for detailed information.
2. Slide open the door of the instrument’s protective cover.
3. Load plates into the instrument according to the protocol request, placing each plate in the same orientation. Confirm each action by pressing Start.
4. After the samples are processed, remove the plates as instructed by the instrument’s display. Press Start after removing each plate. Press Stop after all the plates are removed.

Troubleshooting

Problem	Possible Cause	Solution
Low amount of protein was recovered	The protein degraded	Add protease inhibitors
	Not enough magnetic beads were used	Increase the amount of magnetic bead used for capture
	Sample had an insufficient amount of target protein	Increase amount of antigen sample
Protein does not elute	Elution conditions were too mild	Increase incubation time with elution buffer
Bands at ~50kDa appeared on the Western blot	Elution was performed in Lane Marker Sample Buffer at temperatures above room temperature and rabbit antibody was used in the Western blot detection	Perform elution at room temperature when using a rabbit antibody for Western blot detection
Multiple nonspecific bands	Nonspecific protein bound to the magnetic beads	Add 50-350mM NaCl to the Binding/Wash Buffers
Recovered protein was inactive	Elution conditions were too stringent	Use a milder elution buffer (e.g., Pierce Gentle Ag/Ab Elution Buffer, Product No. 21027)
Magnetic beads aggregated	Magnetic beads were frozen or centrifuged	Handle the beads as directed in the instructions
	Buffer was incompatible with magnetic beads	

Additional Information Available on Our Website

- Frequently Asked Questions
- Tech Tip #43: Protein stability and storage
- Tech Tip #34: Binding characteristics of Protein A, Protein G, Protein A/G and Protein L
- Visit thermofisher.com/kingfisher for information on the KingFisher Products

Frequently Asked Questions for the KingFisher Instrument

Question	Answer
Which plates are compatible with the KingFisher Flex Instrument?	The KingFisher Flex Instrument is compatible with the KingFisher 24 Deep Well Plates, Microtiter Deep Well 96 Plates and KingFisher 96 and 96 PCR Plates
Is it possible to concentrate samples during the run?	Both Deep Well 96 Plates and KingFisher 96 Plates can be used during the same run. Therefore, it is possible to start the processing using larger volumes (in a deep well plate) and elute the purified sample to a smaller volume (in a KingFisher 96 Plate)
Is it possible to heat the samples during the run?	The heating block is located inside the instrument and can be used automatically during the sample process. All plates compatible with the KingFisher Flex Instrument can be heated using specially designed, interchangeable heating blocks
Why do the beads stick to the plastic tips and wells or the eluted proteins stick to the wells?	Proteins conjugated to beads and eluted proteins can nonspecifically bind to plastics. Adding detergent to Binding/Wash Buffer prevents the protein conjugated to the bead from sticking (0.05%-0.1% Tween-20 Detergent). Also include a small amount of detergent in the Elution Buffer (e.g., 0.05% Tween-20 Detergent) or silanize the elution plate
Are the reagent volumes in each well critical?	For best results, keep the specified volumes within defined limits to avoid spillover

Related Products

88802-3	Pierce Protein A/G Magnetic Beads
88847-8	Pierce Protein G Magnetic Beads
88849-50	Pierce Protein L Magnetic Beads
88816-7	Pierce Streptavidin Magnetic Beads
88826-7	Pierce NHS-Activated Magnetic Beads
88804	Pierce Classic Magnetic IP/Co-IP Kit
88805	Pierce Crosslink Magnetic IP/Co-IP Kit
88828	Pierce Direct Magnetic IP/Co-IP Kit
24615	Imperial™ Protein Stain
34075	SuperSignal™ West Dura Extended Duration Substrate
XP04200BOX	Novex™ Tris-Glycine protein gels (see thermofisher.com/proteingels for a complete listing)

For research use only. Not for use in diagnostic procedures.

Products are warranted to operate or perform substantially in conformance with published Product specifications in effect at the time of sale, as set forth in the Product documentation, specifications and/or accompanying package inserts (“Documentation”). No claim of suitability for use in applications regulated by FDA is made. The warranty provided herein is valid only when used by properly trained individuals. Unless otherwise stated in the Documentation, this warranty is limited to one year from date of shipment when the Product is subjected to normal, proper and intended usage. This warranty does not extend to anyone other than Buyer. Any model or sample furnished to Buyer is merely illustrative of the general type and quality of goods and does not represent that any Product will conform to such model or sample.

NO OTHER WARRANTIES, EXPRESS OR IMPLIED, ARE GRANTED, INCLUDING WITHOUT LIMITATION, IMPLIED WARRANTIES OF MERCHANTABILITY, FITNESS FOR ANY PARTICULAR PURPOSE, OR NON INFRINGEMENT. BUYER’S EXCLUSIVE REMEDY FOR NON-CONFORMING PRODUCTS DURING THE WARRANTY PERIOD IS LIMITED TO REPAIR, REPLACEMENT OF OR REFUND FOR THE NON-CONFORMING PRODUCT(S) AT SELLER’S SOLE OPTION. THERE IS NO OBLIGATION TO REPAIR, REPLACE OR REFUND FOR PRODUCTS AS THE RESULT OF (I) ACCIDENT, DISASTER OR EVENT OF FORCE MAJEURE, (II) MISUSE, FAULT OR NEGLIGENCE OF OR BY BUYER, (III) USE OF THE PRODUCTS IN A MANNER FOR WHICH THEY WERE NOT DESIGNED, OR (IV) IMPROPER STORAGE AND HANDLING OF THE PRODUCTS.

Unless otherwise expressly stated on the Product or in the documentation accompanying the Product, the Product is intended for research only and is not to be used for any other purpose, including without limitation, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses, or any type of consumption by or application to humans or animals.

Current product instructions are available at thermofisher.com. For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2017 Thermo Fisher Scientific Inc. All rights reserved. Tween is a trademark of Croda International PLC. All (other) trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.