

Pierce Magnetic c-Myc-Tag IP/Co-IP Kit

Pub. No. MAN0016845
Pub. Part No. 2162548.3
Rev. A.0

88844

Number	Description
88844	<p>Pierce Magnetic c-Myc-Tag IP/Co-IP Kit, contains sufficient reagents to perform 40 reactions using 25µL of magnetic beads</p> <p>Kit Contents:</p> <p>c-Myc-tagged Positive Control (23633), 500µL, 1 mg/mL, <i>Escherichia coli</i> extract containing c-Myc-tagged GST fusion protein</p> <p>Application Set (88844X):</p> <p>Pierce Anti-c-Myc Magnetic Beads, 1mL, supplied at 10mg/mL in PBS containing 0.05% Tween™-20 Detergent and 0.05% sodium azide</p> <p>Mag c-Myc IP/Co-IP Buffer-1, 50mL, 25mM Tris, 0.15M NaCl, 1mM EDTA, 1% NP40, 5% glycerol; pH 7.4</p> <p>Mag c-Myc IP/Co-IP Buffer-2 (20X), 20mL, TBS Tween buffer (500mM Tris, 3M NaCl, 1% Tween-20 Detergent; pH 7.5)</p> <p>Lane Marker Sample Buffer, Non-reducing (5X), 5mL, 0.3M Tris·HCl, 5% SDS, 50% glycerol, lane marker tracking dye; pH 6.8</p> <p>Elution Buffer, 5mL, pH 2.0</p> <p>Storage: c-Myc-tagged Positive Control ships separately on dry ice; upon receipt store at -70°C. Application Set shipped at ambient temperature; upon receipt store at 4°C.</p> <p>For research use only. Not for use in diagnostic procedures.</p>

Table of Contents

Introduction	2
Important Product Information.....	2
Additional Materials Required.....	3
Procedure for Lysis of Mammalian Cells	3
Procedure for IP of c-Myc-Tagged Proteins	4
A. Manual IP/Co-IP.....	4
B. Elution of c-Myc-Tagged Protein.....	4
C. Automated IP/Co-IP.....	5
Troubleshooting	6
Additional Information Available on Our Website.....	6
Frequently Asked Questions for the KingFisher Instrument.....	7
Related Thermo Scientific Products	7

Introduction

The Thermo Scientific™ Pierce™ Magnetic c-Myc-Tag IP/Co-IP Kit provides a simple and fast method to study c-Myc-tagged proteins with advantages over the traditional immunoprecipitation (IP) procedure using Protein A/G magnetic beads. The high-affinity, anti-c-Myc antibody-coupled beads enable IP of c-Myc-tagged proteins or co-immunoprecipitation (co-IP) of their interacting partners without antibody contamination. The Thermo Scientific™ Pierce™ Anti-c-Myc Magnetic Beads are used for IP of specific c-Myc-tagged proteins expressed in human *in vitro* expression systems and mammalian and bacterial cell lysates. The anti-c-Myc antibody coupled to the blocked magnetic beads is a high-affinity mouse IgG₁ monoclonal antibody that recognizes the c-Myc-epitope tag (EEQK LIS EEDLLRKRREQLKHKLEQLRNSCA) derived from the human c-Myc protein. For immunoprecipitation, the beads are added to sample containing c-Myc-tagged proteins. The bound c-Myc-tagged proteins are dissociated from the beads using a low-pH 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.

Table 1. Characteristics of the Thermo Scientific Pierce Anti-c-Myc Magnetic Beads.

Composition:	High-affinity mouse IgG ₁ monoclonal antibody 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:	≥ 10 μg GST-c-Myc (26kDa fusion protein)/mg of beads <i>or</i> ≥ 10 μg GST-c-Myc (26kDa fusion protein)/100 μL of bead suspension

Important Product Information

- Do not centrifuge, dry or freeze the magnetic beads, as this can cause the beads to aggregate and lose binding activity.
- For best results, determine optimal conditions for expression of c-Myc-tagged fusion protein before attempting IP.
- Include a non-transfected lysate as a negative control to identify nonspecific binding of proteins to the anti-c-Myc magnetic beads. The c-Myc-tagged Positive Control assists in verifying whether the anti-c-Myc magnetic beads can successfully capture the c-Myc-tagged protein.
- Cell lysate with expressed c-Myc-tagged protein may be prepared from mammalian cells using the procedure provided or from bacterial cells using a bacterial lysis protocol (e.g., Thermo Scientific™ B-PER™ II Bacterial Protein Extraction Reagent, Product No. 78260). The buffers provided in this kit are compatible with samples prepared from bacterial lysate.
- To minimize protein degradation, include protease inhibitors (e.g., Thermo Scientific™ Halt™ Protease Inhibitor Single-Use Cocktail, EDTA-free, Product No. 78425) when preparing cell lysates.
- Binding capacity and elution recovery will vary depending on the c-Myc fusion protein and the elution method used.
- 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. The c-Myc antibody will not leach from the beads when eluting with IgG Elution Buffer, pH 2.0.
- Pierce Anti-c-Myc Magnetic Beads are compatible with IP and Western blot analysis.
- Do not use cell lysate containing dithiothreitol (DTT). DTT may cause the c-Myc antibody to leach from the beads.

Additional Materials Required

- Phosphate-buffered saline (PBS, 100mM sodium phosphate, 100mM NaCl; pH 7.2, Product No. 28372)
- DTT (Product No. 20290)
- 1M Tris·HCl, pH 8.5
- Sample containing c-Myc-tagged protein

For Automated IP:

- KingFisher Flex System with 96 deep well head (Product No. 5400630)
- 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)

For Manual IP:

- 1.5mL microcentrifuge tubes
- Magnetic stand (e.g., Thermo Scientific™ MagnaBind™ Magnet for 6 × 1.5mL microcentrifuge tubes, Product No. 21359)
- End-over-end rocker or rotator

Procedure for Lysis of Mammalian Cells

Note: For optimal results, use a protease inhibitor cocktail such as Halt Protease Inhibitor Single-Use Cocktail, EDTA-free, (Product No. 78425) when preparing cell lysate.

Protocol I: Lysis of Cell Monolayer (Adherent) Mammalian Cells

1. Carefully remove culture medium from confluent cells.
2. Wash cells once with ice-cold PBS.
3. Add the appropriate volume of Mag c-Myc IP/Co-IP Buffer-1 (Table 2) to the cells. Incubate on ice for 5 minutes with periodic mixing.
4. Transfer the lysate to a microcentrifuge tube and centrifuge at $\sim 13,000 \times g$ for 10 minutes to pellet cell debris.
5. Transfer the supernatant to a new tube for protein concentration determination (e.g., Thermo Scientific™ BCA Protein Assay Kit, Product No. 23225) and further analysis.

Table 2. Recommended Mag c-Myc IP/Co-IP Buffer-1 volume for different standard culture plates.

<u>Plate Size/Surface Area</u>	<u>Volume of Mag c-Myc IP/Co-IP Buffer-1</u>
100 × 100mm	500-1000µL
100 × 60mm	250-500µL
6-well plate	200-400µL per well
24-well plate	100-200µL per well

Protocol II: Lysis of Cell Suspension Cultures

1. Centrifuge the cell suspension at $1000 \times g$ for 5 minutes to pellet the cells. Discard the supernatant.
2. Wash cells once by suspending the cell pellet in PBS. Pellet cells by centrifugation at $1000 \times g$ for 5 minutes.
3. Add ice-cold Mag c-Myc IP/Co-IP Buffer-1 to the cell pellet. Use 500µL of Mag c-Myc IP/Co-IP Buffer-1 per 50mg of wet cell pellet (i.e., 10:1 v/w). If using a large amount of cells, first add 10% of the final volume of Mag c-Myc IP/Co-IP Buffer-1 to the pellet and mix by pipetting up and down. Then add the remaining volume of Mag c-Myc IP/Co-IP Buffer-1 to the cell suspension.

4. Incubate on ice for 5 minutes with periodic mixing. Remove cell debris by centrifugation at $\sim 13,000 \times g$ for 10 minutes.
5. Transfer the supernatant to a new tube for protein concentration determination (e.g., BCA Protein Assay Kit, Product No. 23225) and further analysis.

Procedure for IP of c-Myc-Tagged Proteins

A. Manual IP/Co-IP

Note: The amount of lysate needed and incubation time required are dependent on the expression level of the c-Myc-tagged protein and require optimization for each specific system. For co-IP experiments, the buffer system must be optimized to maintain the protein:protein interaction.

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

1. Place 25 μ L of Pierce Anti-c-Myc Magnetic Beads into a 1.5mL microcentrifuge tube.
2. Add 175 μ L of Mag c-Myc IP/Co-IP Buffer-1 to the beads and gently vortex to mix.
3. Place the tube into a magnetic stand to collect the beads against the side of the tube. Remove and discard the supernatant.
4. Add 1mL of Mag c-Myc IP/Co-IP Buffer-1 to the beads. Invert the tube several times or gently vortex to mix for 1 minute. Collect beads with a magnetic stand. Remove and discard the supernatant.
5. Add the sample containing the c-Myc-tagged protein to the pre-washed magnetic beads and incubate at room temperature for 30 minutes with mixing.

Note: To prepare the positive control, dilute 30 μ L of positive control into 270 μ L of Mag c-Myc IP/Co-IP Buffer-1.

6. Collect the beads with a magnetic stand, remove the unbound sample and save for analysis.
7. Add 300 μ L of Mag c-Myc IP/Co-IP Buffer-2, diluted 1:20 with water, to the tube and gently mix. Collect the beads and discard the supernatant. Repeat this wash twice.
8. Add 300 μ L of ultrapure water to the tube and gently mix. Collect the beads on a magnetic stand and discard the supernatant.

B. Elution of c-Myc-Tagged Protein

Note: If the eluted c-Myc-tagged protein will be used for functional applications, use Elution Protocol 1. If the protein is sensitive to the low pH, use Thermo Scientific™ Pierce™ c-Myc Peptide (Product No. 20170) to competitively elute the c-Myc-tagged protein. Elute in 100 μ L of Pierce c-Myc Peptide (0.5mg/mL peptide in Tris-buffered saline containing 0.05% Tween-20 Detergent) for 5-10 minutes at 37°C. For electrophoretic analysis, use Elution Protocol 2.

• Elution Protocol 1

1. Add 100 μ L of Elution Buffer to the beads. Incubate the beads at room temperature with mixing for 5-10 minutes.
2. Separate the beads on a magnetic stand and save the supernatant containing the target antigen. To neutralize the low pH, add 15 μ L of 1M Tris·HCl, pH 8.5 for each 100 μ L of eluate.
3. For reducing-gel analysis, prepare Reducing Sample Buffer by mixing 2.5 μ L of 2M DTT with 20 μ L of Lane Marker Non-reducing Sample Buffer (5X).
4. Add 77.5 μ L of elution sample to the 22.5 μ L of prepared Reducing Sample Buffer. Heat sample at 95-100°C for 5-10 minutes in a heat block.

• Elution Protocol 2

1. Prepare Non-reducing Sample Buffer by mixing 200 μ L of Lane Marker Non-reducing Sample Buffer (5X) with 800 μ L of ultrapure water. The 1X solution may be stored at 4°C for up to one year.
2. Add 100 μ L of 1X Non-reducing Sample Buffer to the beads. Vortex briefly to resuspend the beads and then incubate at 95-100°C for 5-10 minutes in a heat block.

Note: Using Non-reducing Sample Buffer can minimize interference from co-eluting antibody fragments.

3. For reducing-gel analysis, add 2.5 μ L of 2M DTT to the 100 μ L sample.

C. Automated IP/Co-IP

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

1. Download the “c-Myc-Tag-Immunoprecipitation” protocol from the Thermo Scientific website (<http://www.thermoscientific.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 plates according to Table 3.

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

Plate #	Plate Name	Content	Volume	Time/Speed
1	Beads	Anti-c-Myc Magnetic Beads	25µL	5 seconds
		Mag c-Myc IP/Co-IP Buffer-1	175µL	
2	Bead Wash	Mag c-Myc IP/Co-IP Buffer-1	1000µL	1 minute/slow
3	Bind	Sample containing c-Myc-tagged Protein	300µL	30 minutes/slow
4	Wash 1	Mag c-Myc IP/Co-IP Buffer-2, diluted 1:4 with ultrapure water	300µL	30 seconds/slow
5	Wash 2	Mag c-Myc IP/Co-IP Buffer-2, diluted 1:4 with ultrapure water	300µL	30 seconds/slow
6	Wash 3	Ultrapure Water	300µL	30 seconds/slow
7	Elution	Elution Buffer	100µL	10 minutes/slow
8	Tip Plate	KingFisher 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.
 - If low-pH Elution Buffer is selected for elution, neutralize the pH by adding 15µL of 1M Tris·HCl, pH 8.5 for each 100µL of eluate upon run completion.
 - If using SDS-PAGE Sample Buffer in a heated elution, install the KingFisher Flex Heating Block (see manual for proper installation) to heat samples at 95-100°C for 5-10 minutes.
4. Select the protocol using the arrow keys on the instrument keypad and press Start. See the KingFisher Flex Instrument User Manual for detailed information.
 5. Slide open the door of the instrument’s protective cover.
 6. Load plates into the instrument according to the protocol request, placing each plate in the same orientation. Confirm each action by pressing Start.
 7. 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
Little or no c-Myc-tagged protein is detected	Tagged protein degraded	Include protease inhibitors (e.g., Halt Protease Inhibitor Single-Use Cocktail, EDTA-free, Product No. 78425) in the lysis buffer Use new lysate or lysate stored at -80°C
	No or minimal tagged protein was expressed	Verify protein expression by SDS-PAGE or Western blot analysis of the lysate using c-Myc-tagged Positive Control as a reference
		Increase the amount of lysate used for IP/co-IP
		Use a more sensitive detection system such as Thermo Scientific™ SuperSignal™ West Femto Chemiluminescent Substrate (Product No. 34095)
Magnetic beads aggregated	Magnetic beads were frozen or centrifuged	Handle the beads as directed in the instructions
Failure to co-IP interacting protein	Wash conditions were too stringent for the weak or transient interaction	Reduce the number of washes and/or lower the ionic strength of wash buffer
	Interacting protein was expressed at a low level	Apply additional protein sample
		Use a more sensitive detection system
	Buffer system was not optimal for the protein:protein interaction	Optimize the co-IP buffer
	Insufficient amount of magnetic beads	Increase magnetic bead volume to 50-100µL per reaction
Insufficient sample loaded on the gel to detect by Western blot analysis	Elute sample in 30% acetonitrile/0.5% formic acid, then dry down using a Thermo Scientific™ Speedvac™ Vacuum Concentrator. Resuspend the sample in SDS-PAGE Sample Buffer and load entire elution fraction on to gel	

Additional Information Available on Our Website

- Frequently Asked Questions
- Tech Tip #43: Protein stability and storage
- Visit www.thermoscientific.com/kingfisher for information on the KingFisher Products.
- In the U.S.A., purchase KingFisher Supplies from Fisher Scientific. Contact your local Thermo Fisher Scientific office to purchase KingFisher Supplies outside the U.S.A.

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 plates and KingFisher 96 Plates may 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 and the eluted proteins stick to the wells?	Eluted proteins and proteins conjugated to beads can nonspecifically bind to plastics. The addition of detergent (e.g., 0.05%-0.1% Tween-20 Detergent) to the Binding/Wash Buffer prevents the protein conjugated to the bead from sticking. Also including a small amount of detergent in the elution buffer (e.g., 0.05% Tween-20 Detergent) or silanizing the elution plate can minimize sticking
Are the reagent volumes in each well critical?	For best results, keep the specified volumes within defined limits to avoid spillover

Related Thermo Scientific Products

88842-3	Pierce Anti-c-Myc Magnetic Beads
88831-2	His Pur™ Ni-NTA Magnetic Beads
88821-2	Pierce Glutathione Magnetic Beads
88836-7	Pierce Anti-HA Magnetic Beads
23620	Pierce c-Myc-Tag IP/Co-IP Kit
20168-9	Pierce Anti-c-Myc Agarose
MA1980	Anti-c-Myc Antibody
20170	Pierce c-Myc Peptide
20290	DTT (Dithiothreitol)
78260	B-PER II Bacterial Protein Extraction Reagent
82033	Pierce Agarose qIP Protein Interaction Kit, Tluc and Myc tags
82036	Pierce Magnetic qIP Protein Interaction Kit, Tluc and Myc tags

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 thermo.fisher.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.