

# Active Rap1 Pull-Down and Detection Kit

16120

2273.2

Number	Description
16120	<b>Active Rap1 Pull-Down and Detection Kit</b> , contains sufficient reagents for 30 pull-down reactions

**Kit Contents:****Box 16120X (these items ship together on dry ice; upon receipt store at -70°C):**

**GST-RalGDS-RBD Fusion Protein**, 1 × 600µg, contains 1-2mg/mL in 50mM Tris•HCl, pH 7.2, 150mM NaCl, 0.5% Triton® X-100, 5mM MgCl<sub>2</sub>, 1mM DTT, protease inhibitors and 10% glycerol; ~40kDa; GST-RalGDS-RBD interacts with Rap1 from human, mouse, and rat origin, store at -70°C

**GDP (100mM)**, 50µL, 100X, in sterile water, store at -70°C (or -20°C)

**GTPγS (10mM)**, 50µL, 100X, in sterile water, store at -70°C (or -20°C)

**Box 16120Y (these items ship together with an ice pack; upon receipt store products as directed):**

**Glutathione Resin**, 3.0mL, supplied as a 50% slurry containing 0.05% sodium azide, store at 4°C

**1X Lysis/Binding/Wash Buffer**, 100mL, 25mM Tris•HCl, pH 7.2, 150mM NaCl, 5mM MgCl<sub>2</sub>, 1% NP-40 and 5% glycerol, store at 4°C

**2X SDS Sample Buffer**, 1.5mL, 125mM Tris•HCl, pH 6.8, 2% glycerol, 4% SDS (w/v) and 0.05% bromophenol blue, store at 4°C

**Anti-Rap1 Antibody**, 50µL (5 units), rabbit monoclonal IgG; Anti-Rap1 antibody reacts with Rap1A and Rap1B of human, rat, and mouse, store at -20°C.

**Note:** One unit of Anti-Rap1 antibody is defined as the amount of antibody required to efficiently detect Rap1 in 40µg NIH3T3 whole cell lysate by Western blotting (8.5 × 7.5cm membrane).

**Spin Cups**, 30 each, maximum capacity 850µL, store at room temperature or 4°C

**Collection Tubes**, 90 each, store at room temperature or 4°C

**Table of Contents**

Introduction .....	2
Important Product Information .....	2
Additional Materials Required.....	2
Procedure for Pull-Down and Detection of GTP-Rap1 .....	3
A. Cell Lysis.....	3
B. <i>In vitro</i> GTPγS or GDP Treatment (Optional).....	3
C. Affinity Precipitation of Activated Rap1.....	3
D. Western Blot Analysis.....	4
Troubleshooting.....	5
Additional Information .....	5
Related Thermo Scientific Products .....	6
General References.....	6

---

## Introduction

The Thermo Scientific Active Rap1 Pull-Down and Detection Kit is a simple and fast tool to monitor Rap1 small GTPase activation. The kit provides a GST-fusion protein of the Rap1-binding domain (RBD) from human RalGDS along with glutathione agarose resin to specifically pull down active Rap1 and an anti-Rap1 antibody for Western blot detection. Also included are two control nucleotides, GTP $\gamma$ S and GDP, which can be used to generate positive and negative control lysates, respectively. Each kit is functionally tested to ensure component performance.

Small GTP-binding proteins (or GTPases) serve as molecular switches in signaling transduction pathways. The small GTPase Rap1 (~24kDa) is a member of the Ras superfamily of small GTPases. Rap1 was originally characterized as an antagonist of Ras-induced cell transformation and has been implicated in regulation of cell morphogenesis and cell-cell adhesion. Rap1 is activated by a diverse array of extracellular stimuli and secondary messengers. The resulting signal cascade can result in both the activation and inhibition of ERK (extracellular signal-regulated kinase). Like other small GTPases, Rap1 is active when bound to GTP and inactive when bound to GDP.

## Important Product Information

- Rap1-GTP is quickly hydrolyzed to Rap1-GDP; prepare lysate just before use for each assay, if possible. Otherwise, immediately freeze the lysate at -70°C after preparation.
- For optimal pilot experiments, use 500 $\mu$ g to 1 mg of total lysate per assay.
- Lysis/Binding/Wash buffer is compatible with Thermo Scientific Pierce BCA (Product No. 23227) and Pierce 660nm (Product No. 22660) Protein Assays but not the Coomassie (Bradford) Protein Assay.
- For best results always use protease inhibitors for lysis, and keep lysates on ice between steps.
- For best results when performing the Western blotting procedure, use Pierce™ Goat Anti-Rabbit IgG (H+L), Peroxidase Conjugated (Product No. 31460) and Thermo Scientific SuperSignal West Pico Chemiluminescent Substrate (Product No. 34080) (refer to Additional Information Section, Figure 1). If similar products from other vendors are used, the Western blotting procedure will require optimization.

## Additional Materials Required

- Protease inhibitors (e.g., Thermo Scientific Halt Protease Inhibitor Single-Use Cocktail – EDTA Free, Product No. 78425)
- Pierce BCA Protein Assay Reagent (Product No. 23227) or Pierce 660nm Protein Assay (Product No. 22660)
- $\beta$ -mercaptoethanol (Product No. 35602) or dithiothreitol (DTT) (Product No. 20291)
- Polyacrylamide gel, 12% or 4-20% (Thermo Scientific Precise Protein Gels; see catalog or website)
- Nitrocellulose (Product No. 88014) or PVDF (Product No. 88585) membrane
- Tris-buffered saline (TBS; 25mM Tris•HCl, pH 7.5, 150mM NaCl; Product No. 28379 or 28358)
- Tween™-20 Detergent (Product No. 28320)
- BSA, Fraction V
- Nonfat Dry Milk
- Pierce Goat Anti-Rabbit IgG (H+L), Peroxidase Conjugated (Product No. 31460)
- SuperSignal™ West Pico Chemiluminescent Substrate (Product No. 34080)
- Thermo Scientific CL-XPosure X-ray Film (Product No. 34090 or 34091) or a CCD camera
- 0.5M EDTA, pH 8.0
- 1M MgCl<sub>2</sub>
- Sodium azide (NaN<sub>3</sub>)
- Electrophoresis Apparatus
- Variable-speed Bench-top Microcentrifuge

---

## Procedure for Pull-Down and Detection of GTP-Rap1

### A. Cell Lysis

**Note:** Add protease inhibitors to Lysis/Binding/Wash Buffer before use.

- **For adherent cells:**

1. Carefully remove the culture medium and gently rinse the cells once with ice-cold TBS.
2. Add 0.5-1.0mL Lysis/Binding/Wash Buffer per 75cm<sup>2</sup> flask or 0.3-0.5mL Lysis/Binding/Wash Buffer per 100mm plate with cells at 80-90% confluency.
3. Scrape the cells and transfer to a microcentrifuge tube. Vortex the tube briefly and incubate on ice for 5 minutes.
4. Centrifuge at 16,000 × g at 4°C for 15 minutes.
5. Transfer the supernatant (total lysate) to a new tube.

- **For non-adherent cells:**

1. Pellet cells from one 75cm<sup>2</sup> flask (approx. 1-2 × 10<sup>7</sup> cells) at 100 × g for 5 minutes and then resuspend cells in 10mL ice-cold TBS.
2. Pellet the cells at 100 × g for 5 minutes and carefully remove TBS.
3. Add 0.5-1.0mL Lysis/Binding/Wash Buffer to the cell pellet and resuspend the pellet.
4. Transfer the sample to a microcentrifuge tube and incubate on ice for 5 minutes.
5. Centrifuge at 16,000 × g at 4°C for 15 minutes.
6. Transfer the supernatant (total lysate) to a new tube.

### B. *In vitro* GTPγS or GDP Treatment (Optional)

Perform the following treatments, GTPγS (positive control) and GDP (negative control) to ensure the pull-down procedures are working properly. Use 500μg of cell lysate for each treatment. For best results, aliquot GTPγS and GDP at first use to minimize freeze/thaw cycles.

1. For 500μL lysate, add 10μL 0.5M EDTA pH 8.0 (for a final concentration of 10mM) and vortex the sample.
2. Add 5μL of 10mM GTPγS (for a final concentration of 0.1mM) or 5μL 100mM GDP (for a final concentration of 1mM) and vortex the sample.
3. Incubate the mixture at 30°C for 30 minutes with constant agitation.
4. Terminate the reaction by placing the sample on ice and adding 32μL of 1M MgCl<sub>2</sub> (for a final concentration of 60mM) and vortex the sample.

### C. Affinity Precipitation of Activated Rap1

1. Save a sample of the cell lysate for protein assay using the Pierce BCA or 660nm Protein Assay.
2. Place a spin cup into a collection tube for each sample.
3. Swirl the bottle of Glutathione Resin to thoroughly resuspend the agarose beads. Add 100μL of the 50% resin slurry to the spin cup with collection tube. Centrifuge the tubes at 6000 × g for 10-30 seconds.
4. Discard the flow-through. Add 400μL of Lysis/Binding/Wash Buffer to each tube with resin. Invert the tubes gently several times. Centrifuge the tubes at 6000 × g for 10-30 seconds. Discard the flow-through.
5. Thaw the GST-RalGDS-RBD on ice and immediately make 20μg aliquots. Store aliquots for later use at -70°C.
6. Add 20μg of GST-RalGDS-RBD to the spin cup containing the glutathione resin.
7. Immediately transfer up to 700μL of the cell lysate (containing at least 500μg of total proteins) to the spin cup, close the cap and vortex the sample.

8. Seal cap of the collection tube with laboratory film to prevent leakage, which may result from the presence of detergent in the lysate, and vortex the sample.
9. Incubate the reaction mixture at 4°C for 1 hour with gentle rocking.
10. Centrifuge the spin cup with collection tube at 6000 × g for 10-30 seconds.
11. Remove the laboratory film and transfer the spin cup to a new collection tube.
12. To wash resin, add 400µL of Lysis/Binding/Wash Buffer, invert the tube three times, and centrifuge at 6000 × g for 10-30 seconds. Decant the buffer. Repeat this wash step two additional times.
13. Transfer the spin cup to a new collection tube.
14. Prepare 50µL of reducing sample buffer for each pull-down reaction by mixing 1 part β-mercaptoethanol to 20 parts 2X SDS Sample Buffer (e.g., mix 2.5µL of β-mercaptoethanol to 50µL of 2X SDS Sample Buffer), or by adding dithiothreitol (DTT) to a final concentration of 200mM.
15. Add 50µL 2X reducing sample buffer to the resin. Vortex the sample and incubate at room temperature for 2 minutes.
16. Centrifuge the tube at 6000 × g for 2 minutes. Remove and discard the spin cup containing the resin.
17. Heat the eluted samples for 5 minutes at 95-100°C. Samples may be electrophoresed on a gel or stored at -20°C until use. Apply at least 25µL per lane for a 10 × 10cm mini-gel (12% or 4-20% acrylamide gel provides the best separation).

#### D. Western Blot Analysis

##### Notes:

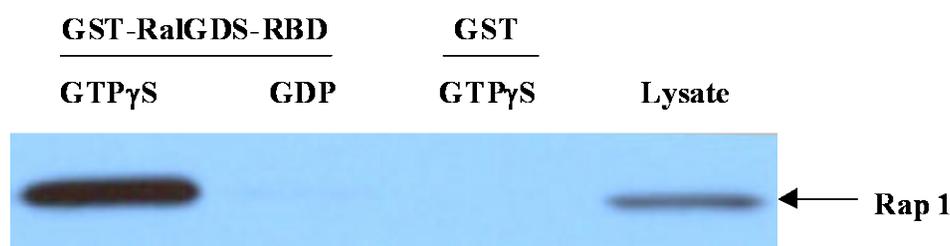
- This procedure has been optimized for use with SuperSignal West Pico Chemiluminescent Substrate (see Important Product Information section).
  - Include unfractionated cell lysate as a control to verify that the Western blot analysis is functioning properly.
  - Perform all blocking, probing and washing incubation steps using constant agitation.
1. Separate the proteins by SDS-PAGE and transfer to nitrocellulose membrane.  
**Note:** Using a PVDF membrane may result in a high background signal.
  2. Block the membrane in TBS containing 3% BSA at room temperature for 2 hours.  
**Note:** Do not block with nonfat dry milk. Nonfat dry milk will significantly reduce the Rap1 signal on the Western blot.
  3. Rinse the membrane with TBS containing 0.05% Tween-20 (TBST).
  4. Prepare a solution containing the Anti-Rap1 Antibody (1:1000 dilution) in 5% BSA and 0.1% NaN<sub>3</sub> in TBST. An example of a 1:1000 dilution is to add 10µL of the stock antibody solution to 10mL of buffer.
  5. Incubate the membrane in the anti-Rap1 antibody solution at 4°C overnight.  
**Note:** If the number of pull-down reactions per blot is low, the diluted anti-Rap1 antibody solution can be re-used up to three times with no performance loss. Store the diluted anti-Rap1 antibody solution at 4°C for up to two months.
  6. Wash the membrane five times for 5 minutes each with TBST.
  7. Dilute the anti-rabbit IgG-HRP conjugate in TBST containing 5% BSA [e.g., if using Pierce Goat Anti-Rabbit IgG (H+L), Peroxidase Conjugated (Product No. 31460), dilute at 1:20,000 to 1:100,000].
  8. Incubate the membrane in the diluted anti-rabbit IgG HRP conjugate at room temperature for 1 hour.
  9. Wash the membrane five times for 5 minutes each with TBST.
  10. Incubate the membrane with chemiluminescent substrate at room temperature (e.g., SuperSignal West Pico Chemiluminescent Substrate).
  11. Immediately expose the membrane to X-ray film or a CCD camera.  
**Note:** The Rap1 band is located at ~24kDa.

## Troubleshooting

Problem	Cause	Solution
No activated Rap1 isolated	No activated Rap1 was present in lysates	Include GTP $\gamma$ S-treated lysate as positive control for pull-down
	Insufficient activated Rap1	Increase the amount of lysate used for detection
	GST-RalGDS-RBD was not added	Add GST-RalGDS-RBD to the reactions
	Degraded GST-RalGDS-RBD	Avoid multiple freeze/thaw cycles of GST-RalGDS-RBD
	Degraded lysate proteins	Add protease inhibitors to the Lysis/Binding/Wash Buffer before lysing the cells
	Incorrect secondary antibody was used for detection	Use goat anti-rabbit IgG
	Detection system was not functioning properly or requires optimization	Consult the instructions for the detection system being used
No signal with GTP $\gamma$ S or strong signal with GDP	GTP $\gamma$ S or GDP were no longer functional	Aliquot GTP $\gamma$ S or GDP after first thaw and store at -70°C; avoid repeated freeze/thaw cycles of the resuspended solution
	Incorrect concentration of EDTA or MgCl <sub>2</sub>	Prepare new solutions with correct concentrations
Western blot resulted in high background	Inadequate blocking and/or washing	Optimize blocking; increase length, number or volume of washes
	Secondary antibody concentration was too high	Use the secondary antibody at a higher dilution

## Additional Information

Rap1 is active when bound to GTP and inactive when bound to GDP. Upon binding to GTP, Rap1 interacts with downstream effectors such as RalGDS. Furthermore, binding of Rap1 to the Rap1-binding domain (RBD) from RalGDS inhibits intrinsic and GAP-enhanced GTPase activity of Rap1. Therefore, the RalGDS-RBD can be used as a probe to specifically isolate active or GTP-Rap1 (Figure 1).



**Figure 1. Western blot of control reactions.** NIH3T3 cell lysates (500 $\mu$ g) were treated *in vitro* with GTP $\gamma$ S or GDP to activate or inactivate Rap1 (refer to optional section B). The lysates (500 $\mu$ g) were then incubated with GST-RalGDS-RBD and a Glutathione Resin. GTP $\gamma$ S-treated lysate was also incubated with GST alone in the presence of a Glutathione Resin (negative control). Fifty percent of the volume of the eluted samples (24 $\mu$ L) and 30 $\mu$ g of cell lysate were separated by 4-20% SDS-PAGE, transferred to a nitrocellulose membrane, and probed with Anti-Rap1 Antibody. Pierce Goat Anti-Rabbit IgG (H+L), Peroxidase Conjugated (Product No. 31460; 1:20,000 dilution) was used as the secondary antibody. The detection was performed with SuperSignal West Pico Chemiluminescent Substrate (Product No. 34080) and followed by exposure to X-ray film.

## Related Thermo Scientific Products

25200-44	Precise™ Protein Gels (see catalog or website for a complete listing)
21065	Pierce Background Eliminator Kit
23236	Pierce Coomassie Plus (Bradford) Protein Assay Reagent
23227	Pierce BCA Protein Assay Reagent Kit
22660	Pierce 660nm Protein Assay Reagent, 750mL
28320	Surfact-Amps™ 20 Detergent Solution, 6 × 10mL
28379	BupH™ Tris Buffered Saline Packs, 10 packs, each makes 500mL
28358	Tris Buffered Saline, 20X, 500mL
78425	Halt™ Protease Inhibitor Single-Use Cocktail, EDTA-free (100X), 24 × 100µL microtubes
34079	SuperSignal West Pico Chemiluminescent Substrate, 500mL
34090	CL-XPosure™ Film (5" × 7" sheets), 100 sheets/pkg
34091	CL-XPosure Film (8" × 10"), 100 sheets/pkg
20291	Dithiothreitol, (DTT) No-Weigh™ Format, 7.7 mg DTT/Tube × 48 tubes
88014	Nitrocellulose Membrane, 0.45µm, 7.9cm × 10.5cm
88585	PVDF Membrane, 0.45µm, 7.9cm × 10.5cm
21059	Restore® Western Blot Stripping Buffer, 500mL

## General References

- Asha, H., *et al.* (1999). The Rap1 GTPase functions as a regulator of morphogenesis *in vivo*. *EMBO J* **18**(3):605-15.
- Knox, A.L. and Brown, N.H. (2002). Rap1 GTPase regulation of adherens junction positioning and cell adhesion. *Science* **295**:1285-8.
- Stork, P.J.S. and Schmitt, J.M. (2002). Crosstalk between cAMP and MAP kinase signaling in the regulation of cell proliferation. *Trends in Cell Biology* **12**(6):258-66.
- Zwartkruis, F.J. and Bos, J.L. (1999). Ras and Rap1: two highly related small GTPases with distinct function. *Exp Cell Res* **253**:157-65.

Parafilm is a trademark of Bemis Company, Inc.

Tween is a trademark of Croda International Plc.

Triton is a trademark of The Dow Chemical Company.

This product ("Product") is 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") and to be free from defects in material and workmanship. Unless otherwise expressly authorized in writing, Products are supplied for research use only. 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 the original purchaser of the Product ("Buyer").

**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 replacement of or refund for the non-conforming Product(s).**

There is no obligation to replace 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.

Current product instructions are available at [www.thermoscientific.com/pierce](http://www.thermoscientific.com/pierce). For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2012 Thermo Fisher Scientific Inc. All rights reserved. Unless otherwise indicated, all trademarks are property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.