

# **Anti-*myc* Antibody** **Anti-*myc*-HRP Antibody**

**Catalog nos. R950-25, R951-25**

**Version D**

080101  
28-0138



[www.invitrogen.com](http://www.invitrogen.com)  
[tech\\_service@invitrogen.com](mailto:tech_service@invitrogen.com)



# Table of Contents

|  |            |
|--|------------|
| <b>Table of Contents .....</b>                         | <b>iii</b> |
| <b>Overview .....</b>                                  | <b>1</b>   |
| <b>Western and Dot Blot.....</b>                       | <b>2</b>   |
| <b>Enzyme-Linked Immunosorbent Assay (ELISA) .....</b> | <b>4</b>   |
| <b>Immunoprecipitation .....</b>                       | <b>7</b>   |
| <b>Technical Service .....</b>                         | <b>8</b>   |
| <b>Product Qualification .....</b>                     | <b>10</b>  |
| <b>References .....</b>                                | <b>11</b>  |



# Overview

## Introduction

The Anti-*myc* and Anti-*myc*-HRP Antibody allows detection of recombinant proteins containing the *c-myc* epitope (Evan, G. I., *et al.*, 1985). The Anti-*myc* Antibody recognizes the sequence:

**-Glu-Gln-Lys-Leu-Ile-Ser-Glu-Glu-Asp-Leu-**

## Contents

The table below supplies information on the concentration of antibody, buffer, and amount supplied.

The Anti-*myc* antibody is a mouse monoclonal IgG<sub>1</sub> antibody. Anti-*myc*-HRP antibody was prepared by crosslinking the Anti-*myc* antibody with horseradish peroxidase using glutaraldehyde.

For MSDS information, please refer to page 8.

| Antibody              | Concentration | Buffer           | Amount              |
|-----------------------|---------------|------------------|---------------------|
| Anti- <i>myc</i>      | ~1 mg/ml*     | PBS, 0.01% azide | 50 µl (25 westerns) |
| Anti- <i>myc</i> -HRP | *             | PBS              | 50 µl (25 westerns) |

\*For exact concentration please refer to the label on the tube.

## Shipping/Storage

The Anti-*myc* and Anti-*myc*-HRP antibody are shipped and stored at +4°C.

For long-term storage, aliquot the antibody and store at -20°C or -80°C. Repeated freezing and thawing is not recommended as it may result in loss of antibody activity.

## Specificity of the Antibody

The Anti-*myc* and Anti-*myc*-HRP antibody have been tested in both immunoblotting and ELISA procedures. In western blot experiments with purified protein, 500 ng of recombinant protein gave a strong signal using chemiluminescent or alkaline phosphatase detection reagents.

## Recommended Dilutions

We recommend the following dilutions of the supplied solution of antibody for these applications:

- For western blots, dilute 1:5000 into Phosphate-Buffered Saline (PBS) containing 0.05% Tween-20 and 5% nonfat, dry milk (PBSTM).
- For ELISA assays, serially dilute into PBSTM.

If you use a different buffer for washing and blocking your blots, then dilute as described above with that buffer. You may use other blocking agents such as bovine serum albumin (BSA) or gelatin.



If you are using the unconjugated antibody, please review the following information.

- If you use alkaline phosphatase-conjugated secondary antibody, do not use PBS. Phosphate inhibits alkaline phosphatase. Use Tris-Buffered Saline (TBS) instead.
- If you use horseradish peroxidase-conjugated secondary antibody, be sure to wash the western blot or microtiter wells thoroughly before adding the secondary antibody. Azide will inhibit horseradish peroxidase activity.

# Western and Dot Blot

---

## Introduction

This procedure can be used for detection of fusion protein expression. We have included a general protocol for your convenience. The table below outlines the basic steps of a western blot.

| Step | Description   |
|------|---|
| 1    | Run an SDS polyacrylamide gel of the purified or partially purified protein or cell lysate with appropriate controls  |
| 2    | Transfer the proteins electrophoretically to a nylon or nitrocellulose membrane                                       |
| 3    | Probe the blot with Anti- <i>myc</i> and Anti- <i>myc</i> -HRP Antibody   |
| 4    | Incubate the blot with appropriate secondary antibodies for chemiluminescent, alkaline phosphatase, or HRP detection. |
| 5    | Generate signal using the appropriate detection reagents.   |

---

## Suggested Solutions

We use chemiluminescence to detect binding of the Anti-*myc* and Anti-*myc*-HRP antibody to the recombinant protein. Other detection methods can be used to detect your protein. The following materials and solutions are needed for immunoblotting:

- Phosphate-Buffered Saline (137 mM NaCl, 2.7 mM KCl, 10 mM Na<sub>2</sub>HPO<sub>4</sub>, 1.8 mM KH<sub>2</sub>PO<sub>4</sub>)
  - Phosphate-Buffered Saline + Tween 20 (PBST: PBS plus 0.05% Tween-20, v/v)
  - Blocking buffer (PBST + 5% nonfat, dry milk, w/v)
  - Secondary Antibody: Anti-Mouse IgG (whole molecule) HRP
- 

*continued on next page*

## Western and Dot Blot, continued

---

### Immunoblotting Protocol

Prepare an SDS polyacrylamide gel (either Tris/Glycine or Tris/Tricine) designed to resolve your recombinant protein or you may purchase Novex<sup>®</sup> pre-cast gels (please call Technical Service for more information). Prepare your samples for electrophoresis. For information about SDS-polyacrylamide gel electrophoresis, please see Ausubel *et al.*, 1994. Load at least 50 to 500 ng of your recombinant protein onto the gel in order to get a good signal.

1. Load your samples and electrophorese your SDS polyacrylamide gel.
2. Transfer proteins to nitrocellulose membrane electrophoretically. We use 25 mM Tris (pH 8.3), 192 mM glycine, 20% v/v methanol as a transfer buffer.
3. Run at 100V, 150 mA (100V, 240 mA at the finish) for 1 hour. Be sure to have a cooling system in place and operational with these electrophoretic settings. You may also transfer overnight at 30V, 40 mA (30V, 90 mA at the finish).
4. Remove the nitrocellulose membrane and incubate it in 10 ml blocking buffer. Gently agitate using a rocker platform for 1 hour at room temperature.
5. Wash the nitrocellulose membrane in 20 ml PBST 2X for 5 minutes each with gentle agitation.
6. Transfer membrane to a tray containing the Anti-*myc* or the Anti-*myc*-HRP antibody diluted 1:5000 in 10 ml blocking buffer (2  $\mu$ l of antibody diluted into 10 ml blocking buffer). Incubate with gentle agitation for 1-2 hours.  
**Note:** Overnight incubation may be preferred, since longer incubations may increase the sensitivity of detection.
7. Transfer membrane to a tray containing 20 ml PBST and wash for 2 x 5 minutes with gentle agitation. If you are using the Anti-*myc*-HRP antibody, proceed to detection.
8. If you are using the Anti-*myc* antibody, transfer the membrane to a tray containing the secondary antibody. Dilute the secondary antibody according to the manufacturer's recommendation into blocking buffer. Incubate with gentle agitation for 1 hour.
9. Wash for 2 x 5 minutes in PBST as described in Step 7.

### Detection Reaction

We use chemiluminescence to detect the fusion proteins. Please follow the manufacturer's instructions. Other detection methods are suitable.

---

### Dot Blot Protocol

This protocol is used to quickly detect the presence of recombinant protein. Be sure to spot equivalent amounts of protein for each sample.

1. Make serial dilutions of samples (purified or partially purified protein or cell lysates) in 10 mM Tris-HCl, 25 mM EDTA, pH 8.0. The lowest dilution should have at least 30 ng of protein present. Nitrocellulose membrane can bind approximately 100  $\mu$ g protein per cm<sup>2</sup>.
  2. Spot 1  $\mu$ l of each sample onto nitrocellulose paper, or alternatively, use a slot blot apparatus.
  3. Allow membrane to air-dry.
  4. Proceed with **Immunoblotting Protocol**, steps 4-9, then perform the detection reaction of choice.
-

# Enzyme-Linked Immunosorbent Assay (ELISA)

---

## Introduction

In addition to western blots, the Anti-*myc* and Anti-*myc*-HRP antibody can also be used in immunoassays. A sample protocol is provided for your convenience. Other protocols are suitable. For more information, please refer to *Antibodies* (Harlow and Lane, 1988) and *Current Protocols in Molecular Biology* (Ausubel *et al.*, 1994), unit 11.2.

---

## Guidelines

Please consider the items below when setting up your ELISA.

- Include controls for cross-reactivity and nonspecific binding to host cell proteins
  - Be sure that all experimental conditions are kept constant to ensure reproducibility
  - Determine optimal dilution of the Anti-*myc* or Anti-*myc*-HRP antibody for use with your antigen
  - Always include a standard curve with each plate
  - Analyze samples in duplicate
  - Be sure that the concentration of antigen falls within the dynamic range of the standard curve.
- 

## Detection of Antigen

ELISAs can be used to detect 1 ng/ml to 1 µg/ml antigen in a bacterial cell lysate. Sandwich ELISAs are more sensitive than direct ELISAs.

---

## Buffers

We have used the following buffers with these antibodies.

- PBS: 137 mM NaCl, 2.7 mM KCl, 10 mM NaHPO<sub>4</sub>, 1.8 mM KH<sub>2</sub>PO<sub>4</sub>
  - PBST: PBS, 0.05% Tween-20
  - PBSTM: PBST, 5% nonfat, dry milk
- 

*continued on next page*

## ELISA, continued

---

### General Procedure

The procedure below is an example of a direct ELISA to detect the amount of antigen in a bacterial cell lysate. For details, please refer to Unit 11.2 of *Current Protocols in Molecular Biology*. Please note that other protocols are suitable.

1. Prepare a 10 µg/ml solution of antigen in PBS. This will be used for your standard curve so it should be as pure as possible.
  2. Take a polystyrene 96-well plate and serially dilute the antigen solution across the columns so that each row has a different dilution of antigen. Use PBSTM as the dilution buffer and 50 µl as the final volume in each well. Leave column 12 blank as a background control. Repeat with a second plate. **Note:** We recommend that you include controls such as a lysate that does not contain the antigen of interest.
  3. Cover the plates and incubate overnight at 4°C or 2 hours at room temperature to allow antigen to bind to the plate.
  4. Remove antigen solution and wash with wells with PBS three times.
  5. Add 0.2 ml PBSTM to each well and incubate for 1 hour at room temperature.
  6. Remove PBSTM and wash with PBST three times.
  7. Serially dilute the Anti-*myc* or Anti-*myc*-HRP antibody across the rows. Use PBSTM as the diluting buffer and 50 µl as the final volume in each well. Start with a 1:500 or 1:1000 dilution in row A.
  8. Cover and incubate plates at room temperature for 2 hours.
  9. Remove antibody and wash wells three times with PBST. If you used Anti-*myc*-HRP antibody, proceed to **ELISA Detection Reaction**, below.
  10. For unconjugated Anti-*myc* antibody, add 50 µl of diluted anti-mouse HRP-conjugated secondary antibody to each well. Use PBSTM as the dilution buffer. For the appropriate dilution of secondary antibody, please see the manufacturer's instructions.
  11. Incubate at room temperature for 30 to 60 minutes.
  12. Remove antibody and wash wells four times with PBS.
  13. Proceed to **ELISA Detection Reaction**, below.
- 

### ELISA Detection Reaction

The protocol uses a sensitive chromogenic substrate for detection of HRP-labeled reagents. Other detection reagents are suitable, please refer to *Antibodies* (Harlow and Lane, 1988). This protocol makes enough substrate solution for up to two 96-well microtiter plates. Pre-made substrate solution is available from other vendors (Sigma, Catalog no. T8665).

1. Dissolve 0.1 mg of TMB (3, 3', 5, 5'-tetramethylbenzidine, Sigma Catalog no. T2885) in 0.1 ml of dimethylsulfoxide (DMSO).
  2. Add 9.9 ml of 0.1 M sodium acetate, pH 6.0.
  3. Filter through Whatman No. 1 paper or equivalent.
  4. Add hydrogen peroxide to a final concentration of 0.01%.
  5. Add 50 µl of the substrate solution to each well.
  6. Incubate 10-30 minutes at room temperature. Positives appear pale blue.
  7. Add 50 µl of 1 M H<sub>2</sub>SO<sub>4</sub> to each well. Positives now appear bright yellow.
  8. Read the results at 450 nm using a spectrophotometer.
- 

*continued on next page*

## ELISA, continued

---

### Analysis of Experiment

Plot absorbance versus known antigen concentration on semilog paper to analyze each antibody dilution. For a working dilution of antibody, choose the dilution that provides the maximum sensitivity over a linear range of antigen concentrations and a minimum binding ( $< 0.05$  absorbance units) for background.

---

### Analyzing Lysates

Once you have identified the optimal working dilution of the Anti-*myc* or Anti-*myc*-HRP antibody, you are ready to analyze your lysates.

1. Prepare an  $\sim 10$   $\mu\text{g/ml}$  solution of your lysate in PBS.
  2. Take a polystyrene 96-well plate and serially dilute the lysates across the columns so that each row has a different dilution of antigen. Use PBSTM as the diluting buffer and 50  $\mu\text{l}$  as the final volume in each well. Leave column 12 blank as a background control. Repeat with a second plate.
  3. Cover the plates and incubate overnight at 4°C or 2 hours at room temperature.
  4. Remove antigen solution and wash with wells with PBS three times.
  5. Add 0.2 ml PBSTM to each well and incubate for 1 hour at room temperature.
  6. Remove PBSTM and wash with PBST three times
  7. Add 50  $\mu\text{l}$  of the appropriate dilution of Anti-*myc* or Anti-*myc*-HRP antibody in PBSTM.
  8. Cover and incubate plates at room temperature for 2 hours.
  9. Remove antibody and wash wells three times with PBST.
  10. Add 50  $\mu\text{l}$  of diluted anti-mouse HRP-conjugated secondary antibody to each well. Use PBSTM as the dilution buffer. For the appropriate dilution of secondary antibody, please see the manufacturer's instructions.
  11. Incubate the blot for 30-60 minutes at room temperature.
  12. Remove antibody and wash wells four times with PBS.
  13. Proceed to **ELISA Detection Reaction**, previous page.
-

# Immunoprecipitation

---

## Immuno- precipitation Protocol

The Anti-*myc* antibody can be used to immunoprecipitate fusion proteins that contain the *c-myc* epitope. The procedure below describes a general immunoprecipitation protocol using the Anti-*myc* antibody. Other protocols are suitable. For more details, please refer to *Antibodies* (Harlow and Lane, 1988).

1. Before beginning, you will need to prepare the appropriate lysis buffer. Many lysis buffers are suitable. We recommend using either:
  - RIPA buffer (150 mM NaCl, 1% NP-40, 0.5% DOC, 0.1% SDS, 50 mM Tris, pH 7.5) or
  - NP-40 lysis buffer (150 mM NaCl, 1% NP-40, 50 mM Tris, pH 8.0)When preparing the lysis buffer, you include a cocktail of protease inhibitors such as 10-100  $\mu$ M leupeptin, 1-10 mM EDTA, 1  $\mu$ M pepstatin, and 0.2-1 mM PMSF (from 100 mM stock in 2-propanol).
2. Lyse cells using one of the lysis buffers above or one of your choice. For mammalian and insect cells, use approximately  $5 \times 10^6$  to  $1 \times 10^7$  cells/ml of lysis buffer. For *E. coli* and yeast cells, use approximately  $1 \times 10^9$  cells/ml of lysis buffer.
3. Centrifuge the lysate for 20 minutes at 10,000 x g at +4°C.
4. Carefully transfer supernatant to a sterile microcentrifuge tube and place on ice.
5. Add 50  $\mu$ l of Protein-G Sepharose<sup>®</sup> resin slurry (50% slurry in lysis buffer) per 1 ml of supernatant to pre-clear the lysate.
6. Rock at +4°C for 1 hour.
7. Centrifuge for 1 minute at 10,000 x g at +4°C.
8. Transfer supernatant to a sterile microcentrifuge tube and place on ice.
9. Add 1-2  $\mu$ g (typically, 1-2  $\mu$ l) of the Anti-*myc* antibody and 50  $\mu$ l of the Protein-G Sepharose<sup>®</sup> resin slurry to the supernatant. Rock for 2-24 hours at +4°C.
10. Centrifuge for 1 minute at 10,000 x g at +4°C.
11. Remove supernatant.
12. Wash the resin 2X with 500  $\mu$ l lysis buffer.
13. The fusion protein immune complexes may now be used in the appropriate assay.
14. For SDS polyacrylamide gels, add 50  $\mu$ l of SDS-PAGE sample buffer (2% SDS, 10% glycerol, 100 mM DTT, 60 mM Tris pH 6.8, and 0.001% bromophenol blue) to the resin. Boil for 5 minutes. Spin down the resin and load supernatant onto gel. Samples can be used immediately or stored at -20°C. Always boil samples before loading.

---

Sepharose<sup>®</sup> is a registered trademark of Amersham Pharmacia Biotech.

## Technical Service

---

### World Wide Web



Visit the Invitrogen Web Resource using your World Wide Web browser. At the site, you can:

- Get the scoop on our hot new products and special product offers
- View and download vector maps and sequences
- Download manuals in Adobe® Acrobat® (PDF) format
- Explore our catalog with full color graphics
- Obtain citations for Invitrogen products
- Request catalog and product literature

Once connected to the Internet, launch your Web browser (Internet Explorer 5.0 or newer or Netscape 4.0 or newer), then enter the following location (or URL):

**<http://www.invitrogen.com>**

...and the program will connect directly. Click on underlined text or outlined graphics to explore. Don't forget to put a bookmark at our site for easy reference!

---

### Contact us

For more information or technical assistance, please call, write, fax, or email. Additional international offices are listed on our Web page ([www.invitrogen.com](http://www.invitrogen.com)).

#### United States Headquarters:

Invitrogen Corporation  
1600 Faraday Avenue  
Carlsbad, CA 92008 USA  
Tel: 1 760 603 7200  
Tel (Toll Free): 1 800 955 6288  
Fax: 1 760 602 6500  
E-mail:  
[tech\\_service@invitrogen.com](mailto:tech_service@invitrogen.com)

#### Japanese Headquarters

Invitrogen Japan K.K.  
Nihonbashi Hama-Cho Park Bldg. 4F  
2-35-4, Hama-Cho, Nihonbashi  
Tel: 81 3 3663 7972  
Fax: 81 3 3663 8242  
E-mail: [jpinfo@invitrogen.com](mailto:jpinfo@invitrogen.com)

#### European Headquarters:

Invitrogen Ltd  
3 Fountain Drive  
Inchinnan Business Park  
Paisley PA4 9RF, UK  
Tel (Free Phone Orders): 0800 269 210  
Tel (General Enquiries): 0800 5345 5345  
Fax: +44 (0) 141 814 6287  
E-mail: [eurotech@invitrogen.com](mailto:eurotech@invitrogen.com)

---

### MSDS Requests

To request an MSDS, please visit our Web site ([www.invitrogen.com](http://www.invitrogen.com)) and follow the instructions below.

1. On the home page, go to the left-hand column under 'Technical Resources' and select 'MSDS Requests'.
  2. Follow instructions on the page and fill out all the required fields.
  3. To request additional MSDSs, click the 'Add Another' button.
  4. All requests will be faxed unless another method is selected.
  5. When you are finished entering information, click the 'Submit' button. Your MSDS will be sent within 24 hours.
- 

*continued on next page*

## Technical Service, continued

---

### Emergency Information

In the event of an emergency, customers of Invitrogen can call the 3E Company, 24 hours a day, 7 days a week for disposal or spill information. The 3E Company can also connect the customer with poison control or with the University of California at San Diego Medical Center doctors.

3E Company  
Voice: 1-760-602-8700

---

### Limited Warranty

Invitrogen is committed to providing our customers with high-quality goods and services. Our goal is to ensure that every customer is 100% satisfied with our products and our service. If you should have any questions or concerns about an Invitrogen product or service, please contact our Technical Service Representatives.

Invitrogen warrants that all of its products will perform according to the specifications stated on the certificate of analysis. The company will replace, free of charge, any product that does not meet those specifications. This warranty limits Invitrogen Corporation's liability only to the cost of the product. No warranty is granted for products beyond their listed expiration date. No warranty is applicable unless all product components are stored in accordance with instructions. Invitrogen reserves the right to select the method(s) used to analyze a product unless Invitrogen agrees to a specified method in writing prior to acceptance of the order.

Invitrogen makes every effort to ensure the accuracy of its publications, but realizes that the occasional typographical or other error is inevitable. Therefore Invitrogen makes no warranty of any kind regarding the contents of any publications or documentation. If you discover an error in any of our publications, please report it to our Technical Service Representatives.

**Invitrogen assumes no responsibility or liability for any special, incidental, indirect or consequential loss or damage whatsoever. The above limited warranty is sole and exclusive. No other warranty is made, whether expressed or implied, including any warranty of merchantability or fitness for a particular purpose.**

---

## Product Qualification

---

### Functional QC

Both antibodies are functionally tested by western blot using the protocols on page 2.

#### **Anti-*myc* Antibody**

The antibody must react specifically with 100 ng of an *E. coli* expressed fusion protein containing a *myc* epitope. Western blots must reveal a strong signal, with no non-specific background, after 10 minutes of color development.

#### **Anti-*myc*-HRP Antibody**

The antibody must react specifically with 50 ng of an *E. coli* expressed fusion protein containing a *myc* epitope. Western blots must reveal a strong signal, with no non-specific background, after development with a chemiluminescent substrate followed by a 1 minute exposure to an x-ray film

---

## References

---

- Ausubel, F. M., Brent, R., Kingston, R. E., Moore, D. D., Seidman, J. G., Smith, J. A. and Struhl, K. (1990) *Current Protocols in Molecular Biology*. Greene Publishing Associates and Wiley-Interscience, New York.
- Evan, G. I., Lewis, G. K., Ramsay, G., and Bishop, V. M. (1985) Isolation of Monoclonal Antibodies Specific for *c-myc* Proto-oncogene Product. *Mol. Cell. Biol.* **5**: 3610-3616.
- Harlow, E. and Lane, D. (1988) *Antibodies: A Laboratory Manual*. Cold Spring Harbor Laboratory. Cold Spring Harbor, NY.
- For Immunoprecipitation, see Chapter 11, pp. 421-470.
- For Immunoblotting (westerns), see Chapter 12, pp. 471-510.
- For Immunoassays (ELISA), see Chapter 14, pp. 553-612.
- Sambrook, J., Fritsch, E. F. and Maniatis, T. (1989) *Molecular Cloning: A Laboratory Manual*. Cold Spring Harbor Laboratory Press. Plainview, New York.

---

©1999-2001 Invitrogen Corporation. All rights reserved.