

Technical Manual

| | |
|---|---|
| I. Description | 1 |
| II. Equipments and Reagents Required but not Supplied | 1 |
| III. Instructions for Use | 2 |
| IV. Reagents Compatibility Table | 5 |
| V. Application Examples | 6 |
| VI. Troubleshooting | 8 |
| VII. Related Products..... | 9 |

I. Description

DYKDDDDK or FLAG® tag is an octapeptide that can be added to a protein using recombinant DNA technology. The tag sequence can be fused to the coding sequence of a target protein to facilitate detection and purification. GenScript’s **MonoRab™ Anti-DYKDDDDK Affinity Resin (Cat. No. L00766)** is designed for optimal purification of DYKDDDDK-tagged proteins. Based on our new rabbit monoclonal antibody technology (MonoRab™), this resin offers high binding capacity, high purity, and resistance to high concentrations of salt or acid. The binding capacity of MonoRab™ Anti-DYKDDDDK Affinity Resin is 1.5 mg DYKDDDDK-tagged protein per mL of settled resin. The recombinant proteins expressed in bacterial or mammalian cells cultures specifically bind to the anti-DYKDDDDK monoclonal antibodies coupled to the resin, resulting in highly pure eluates. The resin can withstand stringent washing steps, especially with buffers containing high concentration of salt. The resin is resistant to up to 2 M NaCl, and to buffers with pH as low as 2.5. Table 1 lists the main characteristics of the MonoRab™ Anti-DYKDDDDK Affinity Resin.

Table 1. Characteristics of MonoRab™ Anti-DYKDDDDK Affinity Resin

| | |
|------------------------|---|
| Product content | 50% settled resin in TBS with 50% glycerol and 0.02% sodium azide |
| Matrix | 4% cross-linked agarose |
| Average bead size | 90 μm |
| Ligand | MonoRab™ monoclonal antibody against DYKDDDDK tag |
| Binding capacity | Approximately 1.5 mg DYKDDDDK-tagged protein (Size: 50KDa) per ml settled resin |
| Storage and stability | Store at -20 °C for up to 12 months. |
| Resin reuse | The resin can be recycled for at least 4 times. If maintained as per manufacturer’s instructions, the resin can be reused 10 times with minimum loss of binding capability. |
| Elution method | Two methods can be implied for elution: 1. Alkaline, Elution under native conditions by competition with peptide 2. Elution with SDS-PAGE loading buffer*. |
| Reagents compatibility | Compatible with commonly used reagents at certain concentrations (Please refer to reagent compatibility table on page 5) |

*When eluting with SDS-PAGE loading buffer, the heavy chain of the MonoRab™ antibody will be denatured and released from the resin.

II. Equipments and Reagents Required but not Supplied

- Distilled water
- Micropipettors
- Microcentrifuge tubes
- Vortex mixer
- Reagent reservoirs
- Empty Columns
- Serological pipettes (5 ml, 10 ml)
- Cell lysis buffer
- Protease inhibitor reagents
- Buffers (Table 2 lists the necessary buffers for use with Anti-DYKDDDDK MonoRab™ Affinity Resin.)

Table 2. Buffers necessary for the use of MonoRab™ Anti-DYKDDDDK Affinity Resin

| Purpose | Buffer | Formulation |
|----------------------------|---------------------------------|--|
| Equilibration and Washing | Tris-buffered saline, TBS | 50 mM Tris-HCl, 150 mM NaCl, pH 7.4 |
| | Alkaline elution buffer pH 12.0 | 0.1 M Tris, 0.5 M NaCl, pH 12.0 |
| Elution | Competitive elution buffer | DYKDDDDK or FLAG® peptide in TBS with concentration of 100-500 µg/ml |
| | PAGE gel sample buffer | 10% glycerol, 1% lithium dodecyl sulfate (LDS), 0.2 M triethanolamine-Cl pH 7.6, 1% Ficol®-400, 0.00625% phenol red, 0.00625% Coomassie G250, 0.5 mM EDTA disodium |
| Regeneration of used resin | Regeneration buffer | 0.1 M Tris HCl, 0.5 M NaCl, pH 12.0 |

III. Instruction for Use

1. Sample Preparation

For optimal results, follow the recommendations below for sample preparation.

- 1) Prepare the sample according to the protein's biophysical characteristics. Optimize lysis conditions to minimize factors interfering with protein binding (See *Reagent Compatibility Table*).
- 2) To prevent protein degradation during the purification process, perform sample preparation on ice and/or add protease inhibitors to the sample during cell lysis.
- 3) During cell lysis, add endonucleases to reduce sample viscosity caused by the release of chromosomal DNA or RNA.
- 4) Sample should not contain any insoluble particles. Filter the sample or centrifuge with high speed for 10-15 min at 4 °C to remove the insoluble materials before binding procedure.
- 5) Avoid frequent freeze-thaw cycles. Make lysate/sample aliquots and store at -80°C.

2. Resin Preparation

- 1) Place an empty column on a firm support; rinse the column once with 1x TBS.
- 2) Thoroughly resuspend the resin by gentle inversion and immediately load appropriate volume of the slurry into the column. Wide bore pipette tips are recommended for easy resin slurry transfer.
- 3) Equilibrate the resin by washing with 3 bed volumes of TBS, repeat for a total of 3 times. Allow the buffer to drain from the column by gravity; do not let the resin run dry.

3. Binding Procedures

The MonoRab™ Anti-DYKDDDDK Affinity Resin can purify DYKDDDDK-tagged proteins by different methods including column chromatography, batch binding, or by immunoprecipitation (IP). If the sample volume is around 50 mL with less than 4 mg DYKDDDDK-tagged protein (Size: 50KDa), column binding format with 2-4 ml of settled resin is recommended. For every additional 1.5 mg protein, use an additional 1 ml of settled resin. For a larger sample volume, batch binding format is recommended for quick and efficient purification. For a small volume of sample with low protein expression level, apply the immunoprecipitation method (See *Section III.7 Immunoprecipitation of DYKDDDDK-tagged Protein*)

3.1 Column Chromatography

- 1) Load the prepared sample onto the column under gravity flow at room temperature. Attach a buffer reservoir to the top of the column for loading large volume. Collect and reload flowthrough several times for maximal binding. Lower flow rate may facilitate better target protein binding.
- 2) Wash the column with 10 - 20 bed volumes of TBS to reduce non-specific bindings.
- 3) Allow the column to drain completely and proceed to elution procedure. (See *Section III.4 Elution of DYKDDDDK-tagged protein*)

3.2 Batch Binding

- 1) Resuspend prepared resin (as described in *Section III.2*) in the column completely with prepared sample containing target protein. For every 50 mL sample with less than 4 mg target protein, use 4 mL of settled resin.
- 2) Transfer the resuspended resin to the remaining sample. Incubate the sample with the resin at room temperature for at least 30 min with rotation.
- 3) After incubation, load the sample with resin onto another column to collect the resin. Preserve the flow through of the sample for further use.
- 4) Wash the resin with 10-20 bed volumes of TBS to remove any non-specific binding.
- 5) Allow the column to drain completely and proceed to elution procedure. (See *Section III.4 Elution of DYKDDDDK-tagged protein*)

Notes:

1. The binding time can be extended empirically. The incubation can be done overnight at 4°C with end-to-end rotation. To prevent protein degradation, protease inhibitor should be added to the sample.
2. Preserve the flow through after binding for further SDS-PAGE analysis or Western Blot detection to determine the binding efficiency of the DYKDDDDK-tagged protein.

4. Elution of DYKDDDDK-tagged Protein

Two elution methods are described below (sections 4.1 and 4.2), are recommended. However, according to protein characteristics and downstream applications, one or more elution buffers may be used to optimize recovery. The selected elution method should preserve protein structure integrity and biological function, and should be compatible with downstream applications. All the methods can be carried out at room temperature; lower temperature may be necessary depending on special requirements. After the final wash step, immediately load elution buffer onto the column. See *Section II, Equipment and Reagents Required but not Supplied*, for recipes of recommended elution buffers.

4.1 Elution with alkaline elution buffer , pH 12.0

Elute the bound DYKDDDDK-tagged protein with 6 bed volumes of Alkaline elution buffer (pH 12.0), into vials containing 1/20 bed volume of 1 M HCl to neutralize the eluate. For example, for 1 ml settled resin, use 6 × 1 ml elution buffer to elute sequentially and collect the eluate with six vials containing 50 µl 1M HCl, respectively. Do not leave the column in the alkaline elution buffer for more than 15 minutes; re-equilibrate the column immediately after elution. (See *Section III.5*)

4.2 Elution with competitive elution buffer

Competitive binding with added DYKDDDDK or FLAG[®] peptide can elute bound target protein. The concentration of DYKDDDDK or FLAG[®] peptide in competitive elution buffer may vary from 100-500 µg/ml. Load 2-3 bed volumes of competitive elution buffer into the column by gravity flow; when there is about 1 bed volume of elution buffer left on top of the resin, cap the bottom of column and incubate at room temperature for 30-60 minutes; open the end of the column and collect the eluate.

5. Re-equilibration

After elution, the resin should be washed with TBS buffer to remove residual elution buffer that may denature the antibodies immobilized on the resin. It is recommended to wash the resin 3 times each with 5-10 bed volumes of TBS buffer. After the final wash, allow TBS buffer to drain completely before adding 2 bed volumes of TBS to the resin for next use.

6. Regeneration of MonoRab[™] Anti-DYKDDDDK Affinity Resin

The MonoRab[™] Anti-DYKDDDDK Affinity Resin can be reused multiple times to purify the same protein without regeneration. If the target DYKDDDDK-tagged protein to be purified is different, the resin must be regenerated using the following protocol:

- 6.1 Wash the resin with 2 bed volumes of 0.1 M Tris HCl, 0.5 M NaCl, pH 12.0.
- 6.2 Re-equilibrate the resin with 3-5 bed volumes of TBS.
- 6.3 For short-term storage, the resin can be stored in TBS containing 0.02% sodium azide at 2-8 °C, or stored in TBS containing 50% glycerol and 0.02% sodium azide at -20 °C. For most cases, the resin can be stored for a month in these conditions.

The resin can be recycled for at least 4 times. If maintained properly, the resin can be reused 10 times with minimum loss of binding capability.

7. Immunoprecipitation of DYKDDDDK-tagged Protein

For target proteins to be purified from a small volume of starting sample (e.g., 1–2 ml of cell lysate), the immunoprecipitation procedure can be applied.

- 7.1 Resuspend the resin to form a uniform slurry and transfer 20-100 µl of the slurry into a 1.5 ml vial.

Note: Wide bore pipette tips are recommended for transferring resin slurry.

- 7.2 Add 500 µl TBS into the vial and gently mix the resin. Centrifuge at 6,000 × g for 30 seconds, remove supernatant carefully. Repeat this step two more times. Carefully remove as much supernatant as possible after each wash without disturbing the settled resin.
- 7.3 Add sample to the washed resin. If the sample volume is less than 1 ml, add lysis buffer or TBS to make the total sample volume to 1 ml.
- 7.4 Mix by end-to-end rotation on a tube rotator for at least 1 hour at room temperature.

Note: For optimal elution, the incubation time can be extended. In cases where the protein is highly unstable, lowering the incubating temperature during elution is recommended.

- 7.5 Centrifuge at 6,000×g for 30 seconds. Carefully remove supernatant and add TBS to wash the resin three times. Remove as much supernatant as possible without disturbing the resin, proceed to elution procedure.
- 7.6 Extreme pH condition or DYKDDDDK peptide can be applied to elute the DYKDDDDK-tagged protein. The resin bound with target protein also can be applied directly for SDS-PAGE analysis. Choose one of the following elution methods according to the characteristics of target protein and downstream application.
- 1) Elution with alkaline elution buffer, pH 12.0
 Add 60-300 μ l (3 column volumes) of alkaline elution buffer into the washed resin and use a wide bore pipette tip to gently resuspend the resin. Incubate at room temperature for 5 minutes, mix gently by tapping the tube once or twice during the incubation period. After incubation, centrifuge at 6,000 \times g for 30 seconds. Carefully transfer supernatant into a new vial containing 3-15 μ l 1M HCl for further application.
 - 2) Elution with competitive elution buffer
 Add 60-300 μ l (3 column volumes) of 100-500 μ g/ml DYKDDDDK peptide elution buffer into the washed resin and use a wide bore pipette tip to gently resuspend the resin. Incubate at room temperature for 5 minutes, mix gently by tapping the tube once or twice during the incubation period. After incubation, centrifuge at 6,000×g for 30 seconds. Carefully transfer supernatant into a new vial for further application.
 - 3) Elution with PAGE gel sample buffer
 In order to minimize the denaturation and elution of the MonoRab™ antibody immobilized on the resin, no reducing reagents (β -mercaptoethanol or DTT) should be included in the sample buffer. Reducing reagents will dissociate the heavy and light chains of the antibody. If reducing condition are absolutely necessary, a reducing agent may be added. Please refer to *Section IV Reagents Compatibility, Table 3*.

Add 40-100 μ l (1 column volume) of PAGE gel sample buffer to the washed resin and mix well. Boil for 3 minutes and centrifuge at 8,000×g for 30 seconds. Carefully transfer supernatant into a new vial.

IV. Reagents Compatibility Table

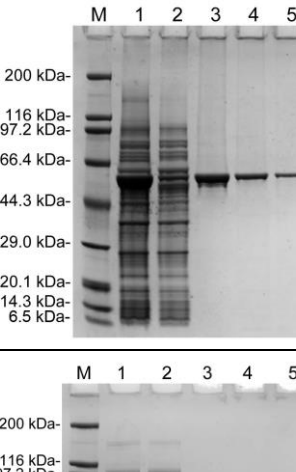
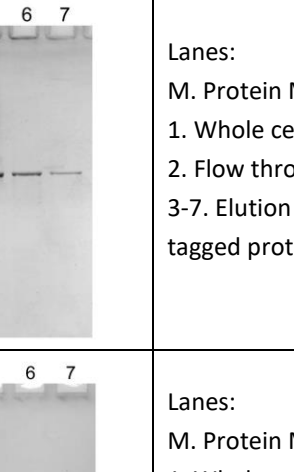
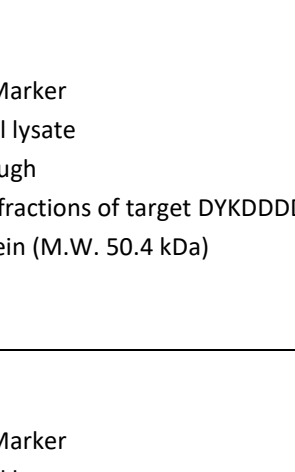
The tolerable concentration of listed reagents are tested by addition of these reagents at indicated concentrations.

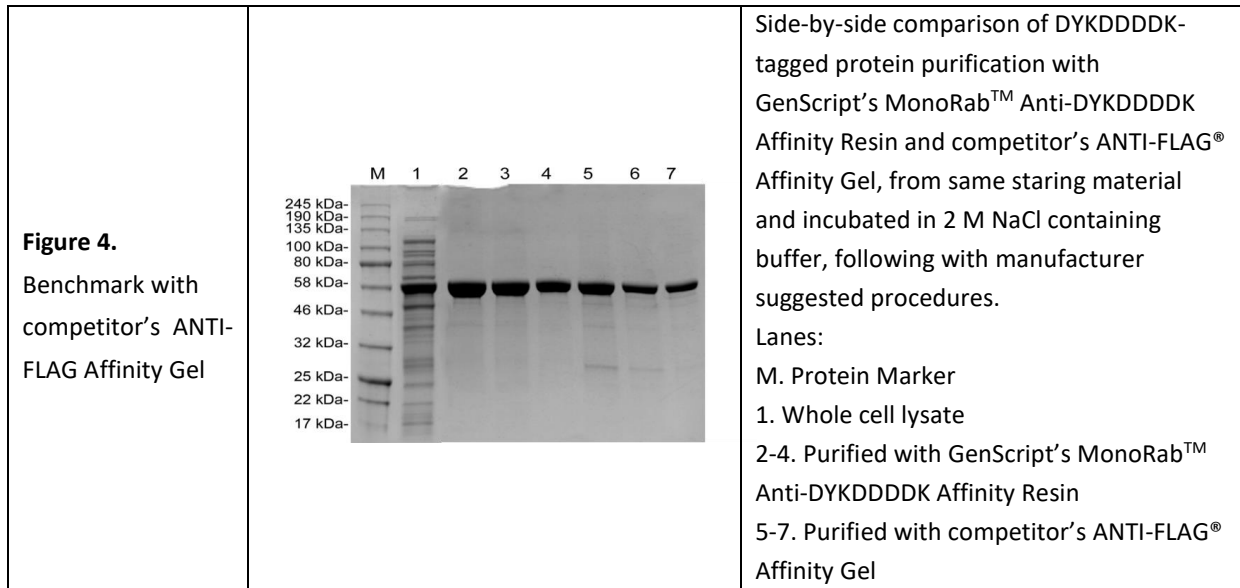
Table 3. Reagents Compatibility

| Reagent | Maximum Tolerable concentration | Note |
|--------------|---------------------------------|--|
| EDTA | 50 mM | Higher concentration of chelating agent will reduce purification efficiency with less target protein recovery. |
| β -ME | 5 mM | Reducing agents will reduce disulfide bonds in the antibody on the resin. Avoid reducing agents or keep at low concentration (<5 mM) during purification process. When the reducing agents reach the maximum tolerable concentration mentioned here, there will be a band of antibody heavy chain (~50 kDa) in SDS-PAGE analysis. The resin cannot be reused if samples containing higher concentration of reducing reagents are applied to the resin. |
| DTT | 80 mM | |
| Tween 20 | 10% | The concentration of detergents should not exceed 10%. |
| Triton X-100 | 10% | |

| | | |
|----------|---------------|---|
| SDS | Not suggested | This detergent will denature the MonoRab™ Anti-DYKDDDDK antibody on the resin. |
| NP-40 | 10% | Higher concentration will reduce purification efficiency with less target protein recovery. |
| HCl | pH 2.5 | More acid will reduce the binding capacity of the resin by destroying the MonoRab™ Anti-DYKDDDDK antibody on the resin. |
| Glycerol | 20% | Higher concentration will interfere with the binding of DYKDDDDK-tagged protein. |
| NaCl | 2 M | Higher concentration will reduce purification efficiency with less target protein recovery. |

V. Application Examples

| Description | Figure | Figure legend |
|---|--|---|
| <p>Figure 1. Purification of DYKDDDDK-tagged protein from <i>E. coli</i> lysate and elution with alkaline elution buffer , pH 12.0</p> |  | <p>Lanes: M. Protein Marker 1. Whole cell lysate 2. Flow through 3-7. Elution fractions of target DYKDDDDK-tagged protein (M.W. 50.4 kDa)</p> |
| <p>Figure 2. Purification of DYKDDDDK-tagged protein from <i>E. coli</i> lysate and Elution with 300 µg/ml DYKDDDDK peptide</p> |  | <p>Lanes: M. Protein Marker 1. Whole cell lysate 2. Flow through 3-7. Elution fractions of target DYKDDDDK-tagged protein (M.W. 50.4 kDa)</p> |
| <p>Figure 3. Purification of DYKDDDDK-tagged protein from mammalian cell lysate and elution with alkaline elution buffer , pH 12.0</p> |  | <p>Lanes: M. Protein Marker 1. Whole cell lysate 2. Flow through 3-7. Elution fractions of target DYKDDDDK-tagged protein (M.W. 50.4 kDa)</p> |



VI. Troubleshooting

| Problem | Possible Cause | Solution |
|---|---|---|
| Large amount of DYKDDDDK-tagged protein found in the flow through | Binding time is not enough | If using batch method, increase the binding time experimentally; If using column method, use a lower flow rate when loading samples. |
| | Column is overloaded | Reduce the amount of the sample added to the resin or increase the amount of resin. |
| | DYKDDDDK-tag is not accessible to resin. | Expose the epitope tag by adding low amount of denaturant to the protein extract (dialysis may be needed before applying onto resin), or fuse DYKDDDDK tag to the other terminus of the target protein. |
| | Resin has not been regenerated since last purification. | Perform resin regeneration procedure prior to binding. |
| | Reagent compatibility problem | Dialyze the sample against TBS before purification procedure. |
| Very few or no DYKDDDDK-tagged protein exists in the eluate. | The target protein has been degraded. | <ol style="list-style-type: none"> 1. Use freshly prepared sample 2. Perform purification at lower temperature, such as 4 °C 3. Include protease inhibitors to the sample during cell lysis and binding steps. |
| | Protein is not completely eluted | Change elution methods according to the instructions in <i>Section III.4.</i> |
| | No target protein expressed | Confirm the presence of target DYKDDDDK-tagged protein in cell lysate with Western blot before purification. |
| | Very low protein expression level | <ol style="list-style-type: none"> 1. Use larger volume of cell lysate. 2. Optimize expression conditions to raise the protein expression level. |
| Multiple protein bands found in the eluate. | The protein is not stable at room temperature. | Purify the target protein at lower temperature, such as 4 °C. |
| | Protein degradation due to proteases activity during purification process | Add protease inhibitors to the cell lysate. |
| | Non-specific binding | <ol style="list-style-type: none"> 1. Prepare cell lysate again. 2. Add additional wash steps. |

VII. DYKDDDDK-tag Related Products

- A00187 THE™ DYKDDDDK Tag Antibody, mAb, Mouse
- A01428 THE™ DYKDDDDK Tag Antibody [HRP], mAb, Mouse
- A01429 THE™ DYKDDDDK Tag Antibody [Biotin], mAb, Mouse
- A01632 THE™ DYKDDDDK Tag Antibody [FITC], mAb, Mouse
- A01868 MonoRab™ DYKDDDDK Tag Antibody, mAb, Rabbit
- A01869 MonoRab™ DYKDDDDK Tag Antibody [HRP], mAb, Rabbit
- A01870 MonoRab™ DYKDDDDK Tag Antibody [Biotin], mAb, Rabbit
- A01871 MonoRab™ DYKDDDDK Tag Antibody [FITC], mAb, Rabbit
- A00170 DYKDDDDK-tag Antibody, pAb, Rabbit
- RP10586 DYKDDDDK Peptide
- M00676 4x LDS Sample Buffer
- L00455W DYDDDK Tag Antibody Plate (White, 96-well)

Limited Use Label license: This product may be the subject of one or more patents filed by GenScript Corporation. The purchase of this product conveys to the buyer the non-transferable right to use the purchased amount of the product and components of the product in research conducted by the buyer (whether the buyer is an academic or for-profit entity). The buyer may not sell or otherwise transfer (a) this product (b) its components or (c) materials made using this product or its components to a third party or otherwise use this product or its components or materials made using this product or its components for any commercial purposes. For commercial use, please contact GenScript at product@genscript.com.

For laboratory or further manufacturing use only.

GenScript USA Inc.

860 Centennial Ave.,
Piscataway, NJ 08854

Tel: 1-877-436-7274, 1-732-885-9188

Fax: 732-210-0262, 732-885-5878

Email: product@genscript.com

Web: www.genscript.com