

Product Information

EZview™ Red ANTI-FLAG® M2 Affinity Gel

Catalog Number **F2426**

Storage Temperature $-20\text{ }^{\circ}\text{C}$

TECHNICAL BULLETIN

Product Description

EZview™ Red ANTI-FLAG® M2 Affinity Gel is a highly visible, red colored ANTI-FLAG M2 agarose affinity gel, designed for use in immunoprecipitation experiments. The affinity resin contains ANTI-FLAG M2 monoclonal antibody covalently attached to crosslinked 4% agarose beads.

EZview Red ANTI-FLAG M2 Affinity Gel functions to bind FLAG® fusion proteins from cell lysates and other biological samples in the same manner as the standard non-colored ANTI-FLAG M2 Monoclonal Antibody Agarose Affinity Gel (Catalog Number A2220). FLAG fusion proteins, bound to the ANTI-FLAG Affinity Gel, are recovered by centrifugation. The red color gives the affinity gel enhanced visibility in order to ease manipulations, such as repetitive washing, recovery of the affinity resin beads, and recovery of the bound FLAG fusion proteins. The enhanced visibility results in less tedious sample processing and greater experimental reproducibility for more accurate quantitation of the proteins of interest.

EZview Red ANTI-FLAG M2 Affinity Gel is supplied as an ~50% slurry suspension in phosphate buffered saline (PBS), containing 50% glycerol and 0.0015% (15 ppm) Kathon® CG/IPCII, as an antimicrobial preservative.

Binding Specificity: FLAG octapeptide (N-Asp-Tyr-Lys-Asp-Asp-Asp-Lys-C), at N-terminal, C-terminal, or Met-N-terminal locations of a fusion protein.

Binding capacity for the FLAG fusion protein, bacterial alkaline phosphatase, is ≥ 0.6 mg per ml of packed resin.

Equipment and Reagents to be Supplied by User

- Cells to be used for preparation of lysate
- Lysis buffer (50 mM Tris HCl, pH 7.4, 150 mM NaCl, 1 mM EDTA, 1% TRITON® X-100), CelLytic™ M (C2978) or CelLytic B (B7435, B7310, or C8740)
- Tris Buffered Saline (TBS): 50 mM Tris HCl, 150 mM NaCl, pH 7.4
- Vortex Mixer
- Protease Inhibitor Cocktail (P8340 or P2714)
- Pipette tips (200 μ l) (P5161)
- Pipette tips, wide orifice (200 μ l) (P1678)
- Pipette tips (1000 μ l) (P1665)
- Pipette (200 μ l) (Z368113)
- Pipette (1000 μ l) (Z368121)
- Microcentrifuge Tubes (T9661)
- 2 \times Laemmli Sample Buffer (S3401)
- 3X FLAG Peptide (F4799) or Glycine HCl (G2879)

Precautions and Disclaimer

EZview Red ANTI-FLAG M2 Affinity Gel is for R&D use only, not for drug, household, or other uses. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

Storage/Stability

EZview Red ANTI-FLAG M2 Affinity Gel should be stored in 50% glycerol at $-20\text{ }^{\circ}\text{C}$ for maximum stability. The unopened product is stable for one year when stored as indicated. Do not freeze in the absence of glycerol.

Procedure

Note: It is recommended the entire technical bulletin be read before use, especially the Reagent Compatibility Table at the end of this bulletin. There are many different procedures and variations for performing immunoprecipitation experiments. The specific method should be chosen by the investigator to suit the particular experiment. See Reference 1 for additional information and procedures.

The generic procedure below is intended to be an example or starting point and may not be appropriate for all situations. The procedure below is written for one sample and is appropriate for most mammalian tissue culture cell lines. For multiple immunoprecipitation reactions, calculate the volume of reagents needed according to the number of samples to be processed. For easy performance of immunoprecipitation reactions, use 40 μ l of gel suspension per reaction (~20 μ l packed gel volume). Smaller amounts of resin (~10 μ l packed gel volume, which bind >1 μ g FLAG fusion protein) can be used.

Manipulations should be carried out on ice or at 2–8 °C, unless otherwise specified.

Immunoprecipitation of FLAG fusion proteins

1. Carefully mix EZview Red ANTI-FLAG M2 Affinity Gel beads until completely and uniformly suspended. Immediately aliquot 40 μ l of the 50% slurry into a clean 1.5 ml microcentrifuge tube on ice. To dispense beads, use a wide orifice pipette tip or cut ~1 mm off the tip of a regular pipette tip, to enlarge the opening and allow unrestricted flow of the bead suspension.
2. Wash/equilibrate beads: Add 500 μ l of TBS (50 mM Tris HCl, 150 mM NaCl, pH 7.4) to the tube, vortex, and centrifuge in a microcentrifuge for 30 seconds at 8,200 \times g (10,000 rpm in an Eppendorf® 5415C microcentrifuge). Carefully remove the supernatant with pipette (or carefully aspirate supernatant) and set tube with the bead pellet on ice.
3. Repeat wash as above. After removing supernatant, set washed bead pellet on ice.
Note: In case of multiple immunoprecipitation samples, the amount of resin needed for all samples can be washed together. Each wash should be performed with TBS at a volume of at least 20 times the total packed gel volume. The washed resin can then be aliquoted for the desired number of samples.
4. Clarify lysate of denatured protein and cell debris by centrifugation for 10 minutes 8,200 \times g at 2–8 °C. Add 200–1,000 μ l cell lysate supernatant to the washed resin. If necessary, bring the final volume to 1 ml by adding lysis buffer. The volume of cell lysate to be used depends on the expression level of FLAG fusion protein in the transfected cells.
5. Gently agitate or shake all samples and controls (a roller shaker is recommended) for 1–2 hours at 2–8 °C. The binding step may be extended overnight to ensure maximum binding.
6. Centrifuge in microcentrifuge for 30 seconds at 8,200 \times g. Place on ice. Aspirate supernatant carefully or remove with a pipette. The supernatant may be saved for analysis, if desired. Place the tube with the bead pellet on ice.
7. Wash the bead pellet by adding 500 μ l of TBS. Vortex briefly and incubate with thorough, gentle mixing at 2–8 °C for 5 minutes. Centrifuge in microcentrifuge for 30 seconds at 8,200 \times g. Aspirate supernatant carefully (or remove with a pipette) and place the tube with the bead pellet on ice.
8. Repeat washes two more times as in step 7. After removing the final wash supernatant, the bound antigen can be eluted and analyzed as desired (see Analysis of Results, below).

Analysis of Results

Elution of the FLAG fusion protein

Three elution methods are suggested:

- A. Protein elution under native conditions by competition with 3X FLAG peptide. The elution efficiency is very high using this method.
- B. Elution under acidic conditions with 0.1 M glycine HCl, pH 3.5. This is a fast and efficient elution method. Immediate neutralization of the eluted protein with wash buffer may help preserve its activity.
- C. Elution with sample buffer for gel electrophoresis and immunoblotting.

Choice of method will depend on protein characteristics and downstream applications.

A. Elution with 3X FLAG peptide:

1. Prepare 3X FLAG elution concentrate by dissolving 3X FLAG peptide (Catalog Number F4799) in 0.5 M Tris HCl, pH 7.5, 1 M NaCl at a concentration of 25 μ g/ μ l. Dilute 5-fold with water to prepare a 3X FLAG stock solution containing 5 μ g/ μ l 3X FLAG peptide. For 3X FLAG elution solution, add 3 μ l of 5 μ g/ μ l 3X FLAG peptide stock solution to 100 μ l of TBS (150 ng/ μ l final concentration).
2. Add 100 μ l of 3X FLAG elution solution (150 ng/ μ l) to each sample and control resin.
3. Incubate the samples and controls with gentle shaking for 30 minutes at 2–8 °C.

4. Centrifuge the resin for 30 seconds at $8,200 \times g$. Transfer the supernatants to fresh test tubes. Be careful not to transfer any resin.

For immediate use, store the supernatants at 2–8 °C. Store at –20 °C for long-term storage.

B. Elution with 0.1 M glycine HCl at pH 3.5:

Note: The procedure should be performed at room temperature. **Do not leave the resin in this buffer more than 20 minutes.**

1. Add 100 μ l of 0.1 M glycine HCl, pH 3.5, buffer to each sample and control resin.
2. Incubate the samples and controls with gentle shaking for 5 minutes at room temperature.
3. Centrifuge the resin for 5 seconds at $8,200 \times g$. Immediately transfer the supernatants to fresh test tubes containing 10 μ l of 0.5 M Tris HCl, pH 7.4, 1.5 M NaCl, to neutralize samples. Be careful not to transfer any resin.

For immediate use, store the supernatant at 2–8 °C. Store at –20 °C for long-term storage.

C. Elution with SDS-PAGE Sample Buffer:

Notes: In order to minimize the denaturation and elution of the antibody, no reducing agent (such as 2-mercaptoethanol or DTT) should be included in the sample buffer. The addition of reducing agents will result in the dissociation of the heavy and light chains of the immobilized M2 antibody (25 and 50 kDa bands). However, if reducing conditions are absolutely necessary, a reducing agent may be added. The final concentration of 2-mercaptoethanol or DTT in the 1 \times sample buffer (62.5 mM Tris HCl, pH 6.8, 2% SDS, 10% (v/v) glycerol, and 0.002% bromophenol blue) should be 5% or 50 mM, respectively.

SDS in the sample buffer will denature the M2 antibody, and, consequently, the EZview Red ANTI-FLAG M2 Affinity Gel beads cannot be reused after treatment with the SDS-PAGE sample buffer.

1. Add 20 μ l of 2 \times Laemmli Sample Buffer (Catalog Number S3401) (125 mM Tris HCl, pH 6.8, 4% SDS, 20% (v/v) glycerol, and 0.004% bromophenol blue) to each sample and control.
2. Boil the samples and controls for 5 minutes.
3. Vortex briefly and centrifuge the samples and controls at $8,200 \times g$ for 30 seconds. The supernatants may be transferred to fresh tubes or may be directly loaded onto SDS-PAGE for subsequent analysis by staining, autoradiography or immunoblotting, using ANTI-FLAG antibodies or specific antibodies against the fusion protein.

Enzyme assays

Enzyme assays, such as kinase assays, can be performed by adding assay mixture and substrate directly into the sample tube. The bead pellet first should be equilibrated in assay buffer by pre-washing in assay buffer before the assay.

Troubleshooting

See troubleshooting guide table on next page.

Because of the enhanced visibility of the affinity resin beads, it is easy to see if beads are accidentally removed during wash steps. If this happens, simply put the wash supernatant back into the tube and repeat the centrifugation step to pellet the resin again.

Controls: For a positive control, add 1 ml TBS and 4 μ l of 50 ng/ μ l FLAG-BAP™ fusion protein (~200 ng) to a comparable tube of washed resin. The amount of FLAG-BAP fusion protein to be precipitated depends on the detection method. 200 ng of protein is sufficient for an activity assay or for an immunoblot analysis. For SDS-PAGE analysis with Coomassie® blue or silver staining, use 1 μ g of FLAG-BAP fusion protein.

For a negative control to monitor non-specific background, use a comparable lysate volume from a lysate of cells not expressing the FLAG fusion protein. Also, FLAG peptide can be used as a competitor in the immunoprecipitation to show specificity for the FLAG fusion protein.

Troubleshooting Guide

Problem	Possible Cause	Solution
No signal is observed.	FLAG fusion protein is not present in the sample.	<ul style="list-style-type: none"> • Make sure the protein of interest contains the FLAG tag by immunoblot or dot blot analyses. • Prepare fresh lysates. Avoid using frozen lysates. • Use appropriate protease inhibitors in the lysate or increase their concentrations to prevent degradation of the FLAG fusion protein.
	Washes are too stringent.	<ul style="list-style-type: none"> • Reduce the number of washes. • Avoid adding high concentrations of NaCl to the mixture. • Use solutions that contain less or no detergent.
	Incubation times are inadequate.	<ul style="list-style-type: none"> • Increase the incubation times with the affinity resin (from several hours to overnight).
	Interfering substance is present in sample.	<ul style="list-style-type: none"> • Lysates containing high concentrations of dithiothreitol (DTT), 2-mercaptoethanol, or other reducing agents may destroy antibody function, and must be avoided. • Excessive detergent concentrations may interfere with the antibody-antigen interaction. Detergent levels in buffers may be reduced by dilution.
	Detection system is inadequate.	<p>If Western blotting detection is used:</p> <ul style="list-style-type: none"> • Check primary and secondary antibodies using proper controls to confirm binding and reactivity. • Verify that the transfer was adequate by staining the membrane with Ponceau S. • Use fresh detection substrate or try a different detection system.
Background is too high.	Proteins bind non-specifically to the ANTI-FLAG monoclonal antibody, the resin beads, or the microcentrifuge tubes.	<ul style="list-style-type: none"> • Pre-clear lysate with Mouse IgG-Agarose (A0919) and/or EZview Red Protein A Affinity Gel (P6486) to remove non-specific binding proteins. • After suspending beads for the final wash, transfer entire sample to a clean microcentrifuge tube before centrifugation.
	Washes are insufficient.	<ul style="list-style-type: none"> • Increase the number of washes. • Prolong duration of the washes, incubating each wash for at least 15 minutes. • Increase the salt and/or detergent concentrations in the wash solutions. • Centrifuge at lower speed to avoid non-specific trapping of denatured proteins from the lysate during the initial centrifugation of the affinity resin complexes.

Reagent Compatibility Table

Reagent	Effect	Comments
Chaotropic agents (such as urea, guanidine HCl)	Denatures the immobilized M2 antibody	Do not use any reagent that contains these types of components since they will denature the M2 antibody on the resin and destroy its ability to bind FLAG fusion proteins. Low concentrations of urea (1 M or less) can be used.
Reducing agents (such as DTT, DTE, 2-mercaptoethanol)	Reduces the disulfide bridges holding the M2 antibody chains together	Do not use any reagent that contains these types of components since they will reduce the disulfide linkages in the M2 antibody on the resin and destroy its ability to bind the FLAG fusion proteins.
Deoxycholate	Interferes with M2 binding to FLAG fusion proteins	Do not use any reagent that contains this detergent since it will inhibit the M2 antibody from binding to FLAG fusion proteins.
Sodium dodecyl sulfate (SDS)	Denatures the immobilized M2 antibody	Do not use any reagent that contains this detergent in the lysis and washing buffers since it will denature the M2 antibody on the resin and destroy its ability to bind FLAG fusion proteins. It is included in the sample buffer for removal of protein from the affinity resin after immunoprecipitation, but the resin cannot be re-used.
TWEEN® 20	Reduces non-specific protein binding to the resin	May be used up to a concentration of 5%. Do not exceed.
TRITON® X-100	Reduces non-specific protein binding to the resin	May be used up to a concentration of 5%. Do not exceed.
IGEPAL® CA-630 (NP-40)	Reduces non-specific protein binding to the resin	May be used up to a concentration of 0.1%. Do not exceed.
CHAPS	Reduces non-specific protein binding to the resin	May be used up to a concentration of 0.1%. Do not exceed.
Digitonin	Reduces non-specific protein binding to the resin	May be used up to a concentration of 0.2%. Do not exceed.
Sodium chloride	Reduces non-specific protein binding to the resin by reducing ionic interactions	May be used up to a concentration of 1.0 M. Do not exceed.
0.1 M glycine HCl, pH 3.5	Elutes FLAG fusion protein from the resin	Do not leave the affinity resin in glycine-HCl for longer than 20 minutes. Longer incubation times will begin to denature the M2 antibody.

Related Products

- Mammalian FLAG expression kits (Catalog Numbers FLMAS and FLMC)
- Mammalian expression vectors (Catalog Numbers E7398, E8770, E1775, E3762, E2275, and E2400)
- Mammalian sequencing primers (Catalog Numbers P5350 and P5475)
- ANTI-FLAG M1 Monoclonal Antibody (Catalog Number F3040)
- ANTI-FLAG M1 Agarose Affinity Gel (Catalog Number A4596)
- ANTI-FLAG M2 Monoclonal Antibody (Catalog Number F3165)
- ANTI-FLAG M2 Affinity Gel (Catalog Number A2220)
- ANTI-FLAG M5 Monoclonal Antibody (Catalog Number F4042)
- FLAG Immunoprecipitation Kit (Catalog Number FLAGIPT1)
- Bicinchoninic Acid (BCA) Kit for Protein Determination (Catalog Number BCA1)
- QuantiPro™ BCA Assay Kit (Catalog Number QPBCA)
- FLAG-BAP control fusion proteins (Catalog Numbers P7582, P7457, and P5975)
- Protease inhibitor cocktails (for general, bacterial, mammalian, fungal & yeast, plant, tissue culture, and histidine-tagged) (Catalog Numbers P2714, P8465, P8340, P8215, P9599, P1860, and P8849)
- Phosphatase inhibitor cocktails (Catalog Numbers P2850 and P5726)
- FLAG peptide (Catalog Number F3290)
- 3X FLAG peptide (Catalog Number F4799)
- ANTI-FLAG M2-Peroxidase (HRP) Conjugate (Catalog Number A8592)
- ANTI-FLAG M2-Alkaline Phosphatase Conjugate (Catalog Number A9469)
- Dithiothreitol (DTT) (Catalog Number D9779)
- 2-Mercaptoethanol (Catalog Number M7154)
- Glycine hydrochloride (Catalog Numbers G2879)
- EZview Red Protein A Affinity Gel (Catalog Numbers P6486)

References

1. Harlow, E., and Lane, D. Antibodies, A Laboratory Manual, pp. 423-470, 513-517 (Cold Spring Harbor Laboratory Press, NY, 1988).
2. Brizzard, B.L., *et al.*, *BioTechniques*, **16**, 730-735 (1994).
3. Knappik, A., and Pluckthun, A., *BioTechniques*, **17**, 754-761 (1994).
4. Chiang, C.M., and Roeder, R.G., *Pept. Res.*, **6**, 62-64 (1993).
5. Ausubel F.M., *et al.*, Current Protocols in Molecular Biology, John Wiley and Sons Inc., NY (1998), pp. 10.15.1-10.16.29

ANTI-FLAG and FLAG are registered trademarks of Sigma-Aldrich® Biotechnology LP and Sigma-Aldrich Co.

EZview, CellLytic, QuantiPro, and FLAG-BAP are trademarks of Sigma-Aldrich Biotechnology LP and Sigma-Aldrich Co.

Kathon is a registered trademark of Rohm and Haas Company.

Eppendorf is a registered trademark of Eppendorf-Netheler-Hinz GmbH.

TRITON is a registered trademark of Union Carbide Corporation.

TWEEN is a registered trademark of Croda International PLC.

IGEPAL is a registered trademark of Rhodia Operations.

Coomassie is a registered trademark of Imperial Industries PLC.

TD,DWF,KMR,CMH,KTA,MAM 08/10-1