

Product Information

FLAG[®] Tagged Protein Immunoprecipitation Kit

Product Code **FLAGIPT-1**

TECHNICAL BULLETIN

Product Description

Immunoprecipitation is a powerful technique for the isolation of proteins or protein complexes. Immunoprecipitation consists of multiple ordered steps: cell lysis, binding of specific antigen to an antibody, antibody-antigen complex precipitation, precipitant wash and antigen dissociation from the immune complex.

Epitope-tagged proteins can be affinity purified and immunoprecipitated, using highly specific antibodies raised against their epitope. The use of such antibodies facilitates subsequent biochemical and immunological analysis. The FLAG[®] epitope system relies on the small FLAG octapeptide (N-Asp-Tyr-Lys-Asp-Asp-Asp-Asp-Lys-C), which allows fusion proteins to retain their original conformation and function. The hydrophilic character of FLAG increases the likelihood that it will be located on the surface of the fusion protein where it is accessible to antibodies.

The FLAG Tagged Protein Immunoprecipitation Kit allows a rapid and efficient immunoprecipitation and elution of an active FLAG-tagged protein. The immunoprecipitation is performed with ANTI-FLAG[®]-M2 affinity gel, which is a highly specific monoclonal antibody covalently attached to agarose resin. The use of affinity resin allows an efficient binding of FLAG-tagged proteins without the need for preliminary steps and calibrations. The immunoprecipitated FLAG-tagged proteins can be efficiently eluted from the resin with acidic conditions or by competition with the FLAG peptide. The immunoprecipitated proteins can be detected for their size, post-translational modifications and interactions on gel electrophoresis, and by activity assays.

Reagents Provided

Sufficient for 50 immunoprecipitation reactions

- Lysis Buffer, 50 ml
Product Code L 3412
50 mM Tris HCl, pH 7.4, with 150 mM NaCl,
1 mM EDTA, and 1% TRITON[™] X-100
- 10x Wash Buffer, 30 ml
Product Code W 0390
0.5 M Tris HCl, pH 7.4, with 1.5 M NaCl
- Elution Buffer, 30 ml
Product Code E 6150
0.1 M Glycine, pH 3.5
- 2x Sample Buffer, 1.5 ml
Product Code S 8684
125 mM Tris HCl, pH 6.8, with 4% SDS,
20% (v/v) glycerol, and
0.004 % bromphenol blue
- 3X FLAG Peptide, 1 mg
Product Code F 4799
- ANTI-FLAG M2-Agarose Affinity Gel, 1 ml
Product Code A 2220
- Amino-terminal FLAG-BAP Fusion Protein, 50 µg
Product Code P 7582

Reagents and Equipment Required but Not Provided

(Product Codes are given where appropriate)

- Test tubes
- Hamilton[®] syringe, Product Code Z10,928-2 or a Pasteur pipette, Product Code S 6268
- Shaker
- Microcentrifuge
- Dulbecco's phosphate buffered saline (PBS), Product Code D 8537
- Cells scrapers, Product Code C 2802
- Protease inhibitor cocktail, Product Code P 8340
- SIGMAFAST[™] pNPP substrate tablets set Product Code N 1891, or pNPP liquid substrate system, Product Code N 7653

Precautions and Disclaimer

This product 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

The kit is shipped on wet ice and stored at $-20\text{ }^{\circ}\text{C}$.

Preparation Instructions

Preparation of working solutions

- Thaw all kit solutions and mix to ensure homogeneity.
- Before use, equilibrate the 2x sample buffer to room temperature. Be sure that the solution is homogenous.
- Prepare 5 $\mu\text{g}/\mu\text{l}$ 3X FLAG peptide solution. The 3X FLAG peptide (N-Met-Asp-Tyr-Lys-Asp-His-Asp-Gly-Asp-Tyr-Lys-Asp-His-Asp-Ile-Asp-Tyr-Lys-Asp-Asp-Asp-Lys-C) is acidic. In order to dissolve it properly add 40 μl of 10x Wash Buffer to 1 mg of 3X FLAG peptide. When the peptide is completely dissolved, add 160 μl of distilled water to the sample. Mix well and store aliquots at $-20\text{ }^{\circ}\text{C}$.
- Dilute a portion ($\sim 2\text{ }\mu\text{l}$) of the FLAG-BAP fusion protein stock solution to a concentration of 50 $\text{ng}/\mu\text{l}$ with 1x wash buffer. The diluted solution is stable at $-20\text{ }^{\circ}\text{C}$ for about two months.
- Prepare 5 ml of 1x Wash Buffer for each immunoprecipitation sample by adding 0.5 ml of 10x Wash Buffer to 4.5 ml sterile deionized water and mixing well.

Procedure

Perform all steps at $2\text{-}8\text{ }^{\circ}\text{C}$, unless the procedure specifies otherwise. Use precooled lysis and wash buffers and equipment. **Do not precool** the sample and elution buffers. All centrifugations are done at $2\text{-}8\text{ }^{\circ}\text{C}$ with precooled rotors.

For antigens and protein:protein complexes requiring a special lysis buffer composed of a different percentage of a detergent, it is recommended to use the 10x Wash Buffer as a core buffer. The ANTI-FLAG M2 affinity gel is resistant to the following detergents: 5.0% TWEEN[®] 20, 5.0% TRITON X-100, 0.1% IGEPAL[®] CA-630, 0.1% CHAPS, and 0.2% digitonin. It can also be used with 1.0 M NaCl or 1.0 M urea. Do not use the ANTI-FLAG M2 affinity gel in the presence of SDS, 2-mercaptoethanol, dithiothreitol (DTT), deoxycholate (DOC), or guanidine HCl. This is not a comprehensive list of interferences.

A. Cell lysis

For 70-90% confluent 100 mm dish ($10^6\text{-}10^7$ cells), use 1 ml lysis buffer. If the expression level of the FLAG-tagged protein is relatively low, lyse the cells with a reduced volume of lysis buffer. It is highly recommended to add protease inhibitor cocktail (Product Code P 8340) to the lysis buffer (10 μl per 1 ml of lysis buffer), especially if the lysate is to be stored for further use.

1. Wash cells.
 - a. For adherent cells.

Remove the growth medium from the cells to be assayed. Rinse the cells twice with PBS buffer, being careful not to dislodge any of the cells. Discard PBS. Add lysis buffer ($10^6\text{-}10^7$ cells/ml).
 - b. For cells in suspension.

Collect the cells into an appropriate centrifuge conical test tube. Centrifuge for 5 minutes at $420\text{ } \times g$. Decant the supernatant and discard. Wash the cells twice by re-suspending the cell pellet with PBS and centrifuge for 5 minutes at $420\text{ } \times g$. Decant the supernatant and discard. Resuspend the cell pellet in lysis buffer ($10^6\text{-}10^7$ cells/ml).
2. Incubate the cells 15-30 minutes on a shaker.
3. For adherent cells only, scrape and collect cells. For cells in suspension, proceed to step 4.
4. Centrifuge the cell lysate for 10 minutes at $12,000\text{ } \times g$.
5. Transfer the supernatant to a chilled test tube. For immediate use, keep on ice. If the supernatant is not to be used immediately, store it at $-70\text{ }^{\circ}\text{C}$.

B. FLAG Fusion Protein Immunoprecipitation

The procedure described is for a single immunoprecipitation reaction. 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, it is recommended to use 40 μ l of gel suspension per reaction (~20 μ l of packed gel volume). Smaller amounts of resin (~10 μ l of packed gel volume, which binds >1 μ g FLAG-tagged protein) can be used.

Two control reactions are recommended for the procedure. The first control is immunoprecipitation with FLAG-BAP fusion protein (positive control) and the second is a reagent blank with no protein (negative control).

The immunoprecipitation procedure includes several steps of precipitation and supernatant removal. To avoid resin removal during the wash and elution steps, it is possible to use a column for the immunoprecipitation reaction. Place an empty chromatography spin column (with a frit) in a centrifuge tube for support.

Load the resin on the column and wash the resin using centrifugation to remove the wash solutions. If the column has an outlet plug the binding and elution steps may be performed on the column. An efficient separation of the eluted protein from the resin is performed by a short centrifugation. This procedure is suitable for either small (20 μ l bed volume) or large resin volumes.

1. Thoroughly suspend the ANTI-FLAG M2 affinity gel in the vial, in order to make a uniform suspension of the resin. The ratio of suspension to packed gel volume should be 2:1. Immediately transfer 40 μ l of the resin in its suspension buffer to a fresh test tube to allow a homogenous dispensation of the resin. For resin transfer use a plastic pipette tip with the end enlarged to allow the resin to be transferred.
2. Centrifuge the resin for 30 seconds at 5,000-8,200 x g. In order to let the resin settle in the tube, wait for 1-2 minutes before handling the samples. Remove the supernatant with a narrow-end pipette tip or a Hamilton syringe, being careful not to transfer any resin. Narrow-end pipette tips can be made using forceps to pinch the opening of a plastic pipette tip until it is partially closed.

3. Wash the packed gel twice with 0.5 ml of 1x Wash Buffer. Be sure that most of the wash buffer is removed and no resin is discarded.

Optional Wash Step:

In order to remove any traces of an unbound ANTI-FLAG antibody from the resin suspension, wash the resin with 0.5 ml elution buffer before continuing with the binding step. **Do not leave the resin in elution buffer more than 20 minutes.** Discard the supernatant immediately, being careful to remove all supernatant from the resin, and follow with three washes consisting of 0.5 ml 1x wash buffer each.

If the Optional Wash Step is not performed: Wash the packed gel an additional 2-3 times using 0.5 ml of 1x Wash Buffer for each wash. This will ensure all of the glycerol is removed before the protein is bound to the gel. In case of numerous immunoprecipitation samples, wash the resin needed for all samples together. After washing, divide the resin according to the number of samples tested. Each wash should be performed with a 1x wash buffer at a volume equal to 20 times the total packed gel volume.

4. Add 200-1000 μ l of cell lysate 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-tagged protein in the transfected cells.

For the positive control, add 1 ml 1x Wash Buffer and 4 μ l of 50 ng/ μ l FLAG-BAP fusion protein (~200 ng) to the washed resin. For the negative control, add only 1 ml of Lysis Buffer with **no** protein.

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.

5. Agitate or shake all samples and controls gently (a roller shaker is recommended) for 2 hours. In order to increase the binding efficiency, the binding step may be extended overnight.

6. Centrifuge the resin for 30 seconds at 5,000-8,200 x *g*. Remove the supernatants with a narrow-end pipette tip.
7. Wash the resin three times with 0.5 ml of 1x Wash Buffer. Make sure all the supernatant is removed by using a Hamilton syringe or equivalent device.

C. Elution of the FLAG-fusion protein

Three elution methods are recommended according to protein characteristics or further usage:

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

1. Elution with 3X FLAG peptide
 - a. Prepare 3X FLAG Elution Buffer. Add 3 μ l of 5 μ g/ μ l 3X FLAG peptide solution to 100 μ l of 1x Wash Buffer (150 ng/ μ l final concentration).
 - b. Add 100 μ l of 3X FLAG Elution Buffer to the resin in the test tube.
 - c. Incubate the samples and controls with gentle shaking for 30 minutes at 4 °C.
 - d. Centrifuge the resin for 30 seconds at 5,000-8,200 x *g*. Transfer the supernatants to fresh test tubes using a Hamilton syringe or equivalent device. 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.

2. Elution with Elution Buffer (Product Code E 6150)

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

 - a. Add 100 μ l of Elution Buffer (Product Code E 6150) to each sample and control resin.
 - b. Incubate the samples and controls with gentle shaking for 5 minutes at room temperature.

- c. Centrifuge the resin for 30 seconds at 5,000-8,200 x *g*. Transfer the supernatants to fresh test tubes containing 10 μ l of 10x Wash Buffer, using a Hamilton syringe or equivalent device. 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.

3. Elution with Sample Buffer

The procedure should be performed at room temperature. Sample buffer should be at room temperature before use.

In order to minimize the denaturation of the antibody, no reducing agent (i.e. 2-mercaptoethanol or DTT) is included in the sample buffer in this kit. If reducing conditions are absolutely necessary, a reducing agent may be added. The final concentration of 2-mercaptoethanol or DTT in the 1x sample buffer should be 5% or 50 mM, respectively.

- a. Add 20 μ l of 2x sample buffer to each sample and control.
- b. Boil the samples and controls for 3 minutes.
- c. Centrifuge the samples and controls at 5,000-8,200 x *g* for 30 seconds to pellet any undissolved agarose. Transfer the supernatants to fresh test tubes with a Hamilton syringe or a narrow-end Pasteur pipette. The samples and controls are ready for loading on SDS-PAGE and immunoblotting using ANTI-FLAG or specific antibodies against the tagged protein.

D. Detection of the positive control
(FLAG-BAP fusion protein):

1. Run the samples and controls on an SDS-PAGE. The protein has a molecular weight of 49.3 kDa. It migrates as a 45-55 kDa band by SDS-PAGE depending on the electrophoresis conditions. Immunoblot the gel with ANTI-FLAG or anti-Bacterial Alkaline Phosphatase (BAP) antibodies, or stain the gel with one of the staining methods (silver staining or Coomassie blue).
2. Detect the BAP presence by enzymatic activity. SIGMAFAST™ pNPP substrate tablets set (Product Code N 1891) or pNPP liquid substrate system (Product Code N 7653) are recommended for the detection of BAP activity.

Troubleshooting Guide

Problem	Possible reason	Solution
Low yield of eluted protein	Low binding efficiency	Increase the amount of lysate/protein in the binding step and/or extend the binding duration to overnight.
	Low elution efficiency	Improve the elution in one of two ways: <ul style="list-style-type: none"> • Increase the concentration of 3X FLAG in the elution solution. • Add salt to the Elution Buffer (0.1 M glycine, pH 3.5).
Appearance of multiple bands in SDS-PAGE in samples eluted with sample buffer	ANTI-FLAG-M2 antibody subunits are removed from the resin by SDS or reducing agents.	Use sample buffer with no SDS or reducing agents, and elute the protein by boiling. After the separation of the supernatant from the resin, SDS can be added. Then samples should be boiled and loaded on an SDS-PAGE.

References

1. Brizzard, B.L., et al., *BioTechniques*, **16**, 730 (1994).
2. Knappik, A., and Pluckthun, A., *BioTechniques*, **17**, 754 (1994).
3. Chiang, C.M., and Roeder, R.G., *Pept. Res.*, **6**, 62 (1993).
4. *Current Protocols in Molecular Biology*, Ausubel F.M., et al. (John Wiley and sons Inc., N.Y.1998), pp. 10.15.1 - 10.16.29
5. *Antibodies, A Laboratory Manual*, Harlow E. and Lane D. (Cold Spring Harbor Laboratory Press, NY, 1988), pp. 514-517, 541-542, 547-549.

TRITON is a trademark of the Union Carbide Corp.
 TWEEN is a registered trademark of Uniqema, a business unit of ICI Americas, Inc.
 IGEPAL is a registered trademark of Rhone-Poulenc AG Co.
 Hamilton is a registered trademark of Hamilton Co.
 Coomassie is a registered trademark of Imperial Chemical Industries Ltd.

DJ/MAM 09/04

These products and/or their use are covered by one or more of the following patents:US 5,011,912, US 4,703,004, US 4,782,137, US 4,851,341, EP 150126,EP 335899, JP 1983150,JP 2665359, CA 1307752. Use of these products are subject to the terms of a license provided in the product packaging, a copy of which will be provided upon request. FLAG® and ANTI-FLAG® registered trademarks of Sigma-Aldrich Biotechnology LP. The product designations of pFLAG™, p3XFLAG™, pFLAG-1™,pFLAG-2™,pFLAGSHIFT™,pFLAG-CTS™,pFLAG-ATS™pFLAG-MAC™,pFLAG-CMV™,YEpFLAG™, and FLAG-BAP™ are trademarks of Sigma-Aldrich Biotechnology LP.

LICENSE AGREEMENT

The enclosed DNA expression vector and/or antibody are specifically adapted for a method of producing selected protein molecules covered by one or more of the following patents owned by Sigma-Aldrich Co.: U.S. Patent Nos. 5,011,912, 4,703,004, 4,782,137 and 4,851,341; EP Patent No. 150,126 (Austria, Belgium, Switzerland, France, United Kingdom, Italy, Netherlands and Sweden); EP Patent No. 335,899 (Belgium, Switzerland, Germany, France, United Kingdom, Italy, Luxembourg and Sweden); German Patent No. P3584260.1; Canadian Patent No. 1,307,752; and Japanese Patent Nos. 1,983,150 and 2,665,359. Your payment includes a limited license under these patents to make only the following uses of these products:

A. Vector License: You may use the enclosed vector to transform cells to produce proteins containing the amino acid sequence DYKDDDDK for research purposes provided, however, such research purposes do not include binding an unlicensed antibody to any portion of this amino acid sequence nor using such proteins for the preparation of antibodies having an affinity for any portion of this amino acid sequence.

B. Antibody License: You may only use the enclosed antibody for research purposes to perform a method of producing a protein in which the protein is expressed in a host cell and purified by use of the antibody in accordance with a claim in one of the above patents in force in a country where the use actually occurs so long as: (1) you perform such method with a DNA expression vector licensed from Sigma-Aldrich Co.; and (2) you do not bind (or allow others to bind) an unlicensed antibody to any DYKDDDDK epitope of any fusion protein that is produced by use of the method.

This license does not include any rights under any other patents. You are not licensed to use the vector and/or antibody in any manner or for any purpose not recited above. As used above, the term "unlicensed antibody" means any antibody which Sigma-Aldrich Co. has not expressly licensed pursuant to Paragraph B, above. Sigma-Aldrich Co. hereby expressly retains all rights in the above listed patents not expressly licensed hereunder.

If the terms and conditions of this License Agreement are acceptable to you, then you may open the vessel(s) containing the vector and/or antibody and, through such act of opening a vessel, will have shown your acceptance to these terms and conditions.

If the terms and conditions of this License Agreement are not acceptable to you, then please return the vessel(s) unopened to Sigma-Aldrich Co. for a complete refund of your payment.

For additional licensing information or to receive a copy of any of the above patents, please contact the Sigma-Aldrich Co. licensing department at telephone number 314-771-5765.

Sigma brand products are sold through Sigma-Aldrich, Inc.

Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.