

SIGMA-ALDRICH

FLAG[®] 96-well Immunoprecipitation System

User Manual

SIGMA-ALDRICH FLAG[®] 96-well Immunoprecipitation System User Manual

FLAG and ANTI-FLAG are 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.

All other trademarks are the property of their respective owners.

© 2010, Sigma-Aldrich Biotechnology LP. All rights reserved.

1.0 TABLE OF CONTENTS

1.0 Table of Contents	3
2.0 System Components	4
Immunoprecipitation Vector Kit (Product # COIP-P)	4
FLAG 96-well Immunoprecipitation Kit (Product # HT-COIP-1)	4
Immunoprecipitation Detection Kit (Product # COIP-D)	4
3.0 Introduction	5
4.0 FLAG 96-well Immunoprecipitation System Overview	6
5.0 Immunoprecipitation Vector Kit (Product # COIP-P)	7
5.1 Reagents Provided (sufficient for 96 samples).....	7
5.2 Additional Reagents and Equipment (Not Supplied).....	7
5.3 Expression Vectors.....	8
5.4 Control Vectors	10
5.5 Verification Primers	11
6.0 FLAG 96-well Immunoprecipitation Kit (Product # HT-COIP-1)	12
6.1 Reagents Provided (sufficient for 96 samples).....	12
6.2 Additional Reagents and Equipment (Not Supplied).....	12
7.0 Immunoprecipitation Detection Kit (Product # COIP-D)	13
7.1 Reagents Provided (sufficient for 96 samples).....	13
7.2 Additional Reagents and Equipment (Not Supplied).....	13
8.0 Protocols for Studying Protein-Protein Interactions with the ANTI-FLAG M2 Plate	14
8.1 Overview of the Analysis of Protein-Protein Interaction with the ANTI-FLAG M2 Plate by ELISA or Western blotting	14
8.2 General Notes	15
8.2.1 Plasmid DNA for transfection	15
8.2.2 Transfection reagent	15
8.2.3 Cell lysates used for co-immunoprecipitation assay	15
8.2.4 Detection of endogenous binding partners during Western blotting or ELISA	16
8.3 Preparation of Cell Extract.....	16
8.4 Immunoprecipitation with ANTI-FLAG M2 plate	16
8.5 Detection of the Bound Protein by ELISA or Western Blotting.....	17
8.5.1 ELISA	17
8.5.2 Western Blotting.....	18
9.0 Troubleshooting Guide	19
10.0 References	20
Appendix 1: Construction of fusion genes using standard cloning techniques	21
Appendix 2: Screening of recombinant plasmids by colony PCR	22
Procedure for Screening Recombinant Plasmids.....	22
Appendix 3: Capturing the Amino-terminal FLAG-BAP Fusion Protein with the ANTI-FLAG M2 plate and evaluating the pNPP substrate integrity	24

2.0 SYSTEM COMPONENTS

Immunoprecipitation Vector Kit (Product # COIP-P)

- pFLAG-CMV™-2 Expression Vector (10 µg)
- pc-Myc-CMV-2 Expression Vector (10 µg)
- pFLAG-CMV-2-p53 control plasmid (20 µg)
- pc-Myc-CMV-2-Large T antigen control plasmid (10 µg)
- pc-Myc-CMV-2-BAP control plasmid (10 µg)
- Verification Primer-MF (2 nmol)
- Verification Primer-MR (2 nmol)
- Technical Bulletins

Storage Conditions: Store all components at -20°C

FLAG 96-well Immunoprecipitation Kit (Product # HT-COIP-1)

- ANTI-FLAG M2 High Sensitivity Capture Plate (1 plate)
- Amino-terminal FLAG-BAP Fusion Protein (2.5 µg)
- 10× Wash Buffer (50 ml)
- Lysis Buffer (50 ml)
- 2× Sample Buffer (6 ml)
- SealPlate Film (6)
- User Manual

Storage Conditions: Store all components at 4°C

Immunoprecipitation Detection Kit (Product # COIP-D)

- Monoclonal Anti-c-Myc, Clone 9E10, Alkaline Phosphatase Conjugate (0.25 ml)
- Tris-Buffered Saline, pH 8.0, with 3% Nonfat Milk (1 packet)
- SIGMA FAST™ p-Nitrophenyl Phosphate Tablets (5 tablet sets)
- Technical Bulletins

Storage Conditions: Store all components at 4°C

3.0 INTRODUCTION

Researchers studying protein-protein interactions commonly use one of several different two-hybrid systems. While these systems are of some use in determining protein binding partners, all of the current two-hybrid systems generate false positives. Sigma's FLAG-96-well Immunoprecipitation Kit provides a unique method to validate protein-protein interactions in a convenient 96-well format. This system enables potential binding partners for the protein of interest to be tagged with FLAG and c-Myc epitopes, followed by capture and detection using the ANTI-FLAG M2 High Sensitivity Capture Plate. The system offers the following advantages compared to currently available techniques:

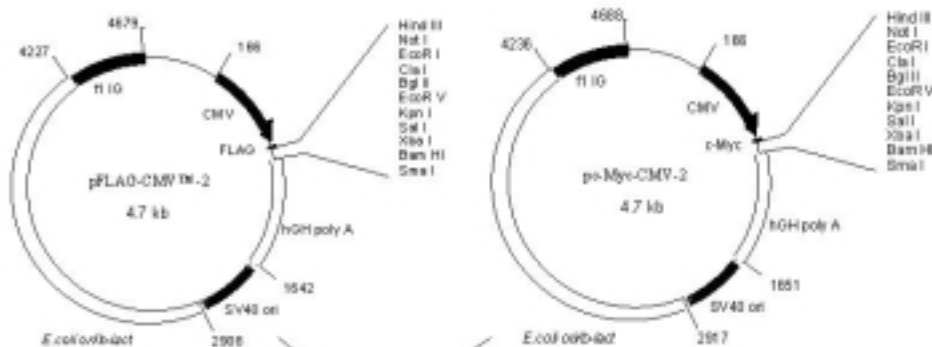
1. **Rapid 96-well format**—The 96-well plate configuration is very effective at handling multiple samples and greatly increases the throughput of the assay. Additionally, this is an ideal system for optimizing *in vitro* binding conditions.
2. **No leaching of ANTI-FLAG M2 antibody**—The ANTI-FLAG M2 antibody is covalently coupled to the surface of the well and the elution conditions used are relatively mild, so very little antibody leaches from the plate. This results in little to no interference from denatured antibody heavy and light chain bands during Western blot analysis.
3. **Rapid protocol with quantitative results**—Because an ELISA format can be used with this system, the protein-protein interaction assay is rapid (less than 4 hours), sensitive, and quantitative.

4.0 FLAG 96-WELL IMMUNOPRECIPITATION SYSTEM OVERVIEW

Figure 1. Outline of the procedure

Clone two potentially interacting proteins into the two mammalian epitope-tagged expression vectors provided in the Immunoprecipitation Vector Kit. One vector generates an N-terminally tagged FLAG fusion protein while the other vector produces an N-terminally tagged c-Myc fusion protein.

- ↓
- Co-express the proteins in appropriate mammalian host cells
- ↓
- Prepare cell extracts with Lysis Buffer
- ↓
- Dilute the cell lysate with 1× Wash Buffer (if necessary)
- ↓
- Incubate cell lysate in the 96-well ANTI-FLAG M2 High Sensitivity Capture Plate
- ↓
- Wash the ANTI-FLAG M2 High Sensitivity Capture Plate with 1× Wash Buffer
- ↓
- Analyze captured proteins by ELISA or Western blotting



1) Clone gene X in pFLAG-CMV™-2 Expression Vector
Clone gene Y in pc-Myc-CMV-2 Expression Vector

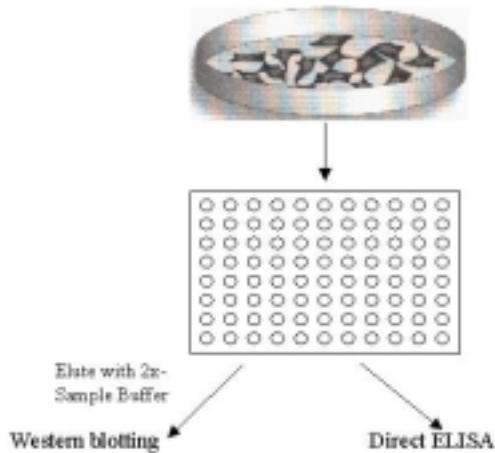
2) Co-express plasmids in mammalian cells

3) Lyse cells

4) Incubate with Anti-FLAG® M2 High Sensitivity Capture Plate

5) Wash

6) Analysis



5.0 IMMUNOPRECIPITATION VECTOR KIT (PRODUCT # COIP-P)

This kit includes the transient mammalian expression vectors pFLAG-CMV[™]-2 and pc-Myc-CMV-2, which are designed for the construction and expression of N-terminally tagged fusion proteins with a FLAG epitope (DYKDDDDK)¹ or a c-Myc epitope (EQKLISEEDL)², respectively. The control plasmids pFLAG-CMV-2-p53, pc-Myc-CMV-2-Large T antigen and pc-Myc-CMV-2-BAP are also supplied. The control plasmids can be used as positive- and negative-binding partners in the immunoprecipitation analysis (see **Section 5.4**). **Figures 2** through **4** and **Table I** illustrate the features of the supplied vectors.

5.1 Reagents Provided (sufficient for 96 samples)

Part	Product #	Amount
pFLAG-CMV [™] -2 Expression Vector	E7398	1 × 10 µg
pc-Myc-CMV-2 Expression Vector	P9236	1 × 10 µg
pFLAG-CMV [™] -2-p53 Control Plasmid	P9986	1 × 20 µg
pc-Myc-CMV-2-Large T antigen Control Plasmid	P9861	1 × 10 µg
pc-Myc-CMV-2-BAP Control Plasmid	P9736	1 × 2.5 µg
Verification Primer-MF	P2987	1 × 2 nmol
Verification Primer-MR	P3112	1 × 2 nmol

5.2 Additional Reagents and Equipment (Not Supplied)

- Pipets (single- or multi-channel)
- Pipette tips
- Microcentrifuge tubes
- Microcentrifuge
- Reagent reservoir
- Vortexer
- MilliQ[®] or sterile water
- DNA plasmid purification kit
- Transfection reagent
- Agarose gel electrophoresis system
- Thermal cycler
- PCR tubes or multi-well PCR plate
- *Taq* DNA polymerase
- Restriction enzymes
- Phosphatase suitable for DNA
- Competent *E. coli* cells
- LB-ampicillin agar plates
- LB medium with 100 µg/ml Ampicillin
- 15 ml culture tubes
- 37°C incubator

5.3 Expression Vectors

Figure 2. Circular maps of the expression vectors

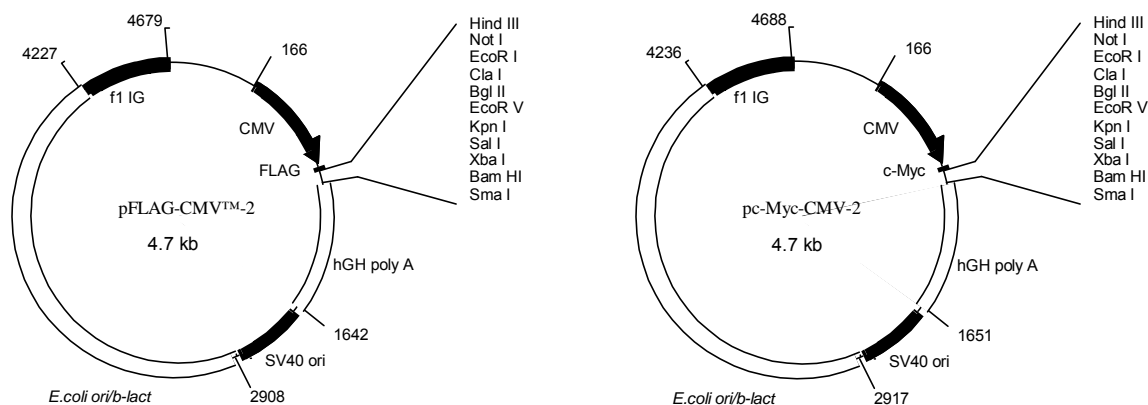
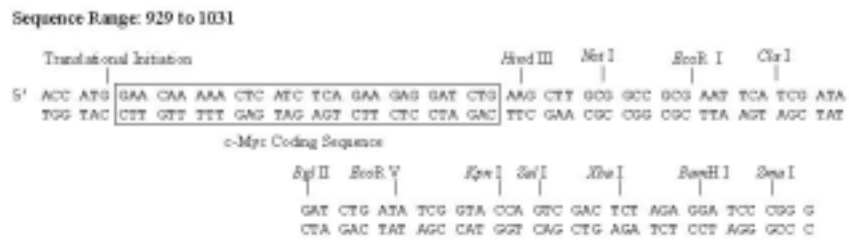
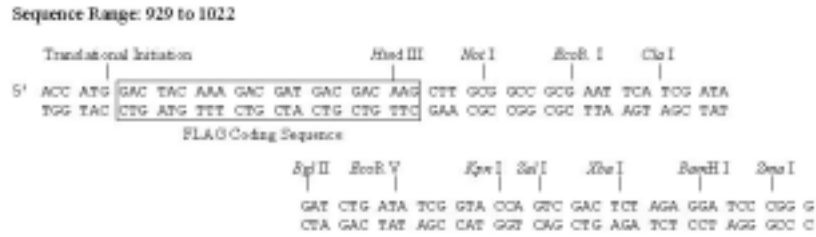


Table I. Expression vector features

Feature	pFLAG-CMV™-2 (4679 bp)	pc-Myc-CMV-2 (4688 bp)
CMV promoter	166-916	166-916
Verification Primer-MF	825-854	825-854
Translational initiation	932-934	932-934
FLAG and c-Myc sequences	935-958	935-967
Multiple cloning region	956-1022	965-1031
hGH poly A	1023-1642	1032-1651
Verification Primer-MR	1080-1103	1089-1112
SV40 ori	1661-2005	1670-2014
pBR322 ori/beta-lactamase	2908-4092	2917-4101
f1 intragenic region	4227-4679	4236-4688

Note: The only difference between the two expression vectors is the epitope tags.

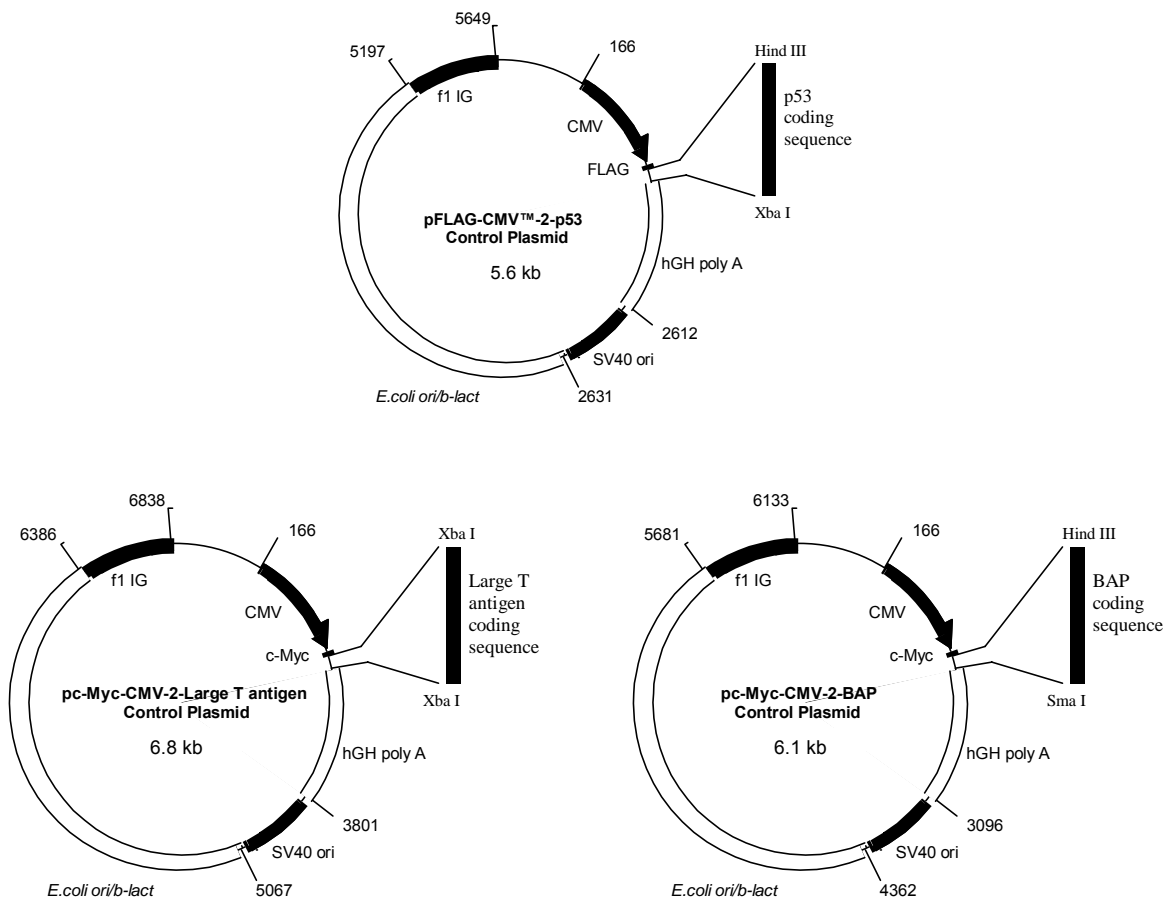
Figure 3. Nucleotide sequences of the epitope tags and multiple cloning region of the pFLAG-CMV™-2 (*upper*) and pc-Myc-CMV-2 (*lower*) expression vectors



5.4 Control Vectors

The Immunoprecipitation Vector Kit provides positive and negative control vectors for co-immunoprecipitation (**Figure 4**). pFLAG-CMV[™]-2-p53 Control Plasmid expresses the N-terminally-tagged FLAG fusion protein containing the wild-type C-terminal domain of mouse p53 (residues 73-390).³ pc-Myc-CMV-2-Large T antigen Control Plasmid expresses the N-terminally-tagged c-Myc fusion protein containing the C-terminal domain of Large T-antigen (residues 86-708).⁴ pc-Myc-CMV-2-BAP Control Plasmid expresses the N-terminally-tagged c-Myc fusion protein containing bacterial alkaline phosphatase (BAP). Mammalian cells co-transfected with control plasmids pFLAG-CMV[™]-2-p53 and pc-Myc-CMV-2-Large T antigen will express tagged fusion proteins that interact, providing a positive control for the co-immunoprecipitation analysis.⁵ Cells co-transfected with pFLAG-CMV[™]-2-p53 and pc-Myc-CMV-2-BAP express tagged fusion proteins that do not interact, demonstrating a negative result in the co-immunoprecipitation analysis.

Figure 4. Circular maps of control vectors



5.5 Verification Primers

Verification primer-MF and Verification primer-MR are supplied with this kit and allow the researcher to screen for the presence, orientation, and reading frame of the DNA insert by colony PCR or DNA sequencing (**Table II**).

Table II. Vector primers and sequences		
Vector primers	Product #.	Sequence
Verification Primer-MF	P2987	5' AAT GTC GTA ATA ACC CCG CCC CGT TGA CGC 3'
Verification Primer-MR	P3112	5' TAT TAG GAC AAG GCT GGT GGG CAC 3'

Notes:

1. Please see **Appendix 1** for procedures involved in cloning recombinant plasmids.
2. Please see **Appendix 2** for procedures involved in screening recombinant plasmids by Colony PCR.
3. If the genes of interest are already cloned into expression vectors that contain FLAG or c-Myc tags, it is not necessary to subclone them into the expression vectors provided in this kit.

6.0 FLAG 96-WELL IMMUNOPRECIPITATION KIT (PRODUCT # HT-COIP-1)

The FLAG-96-well Immunoprecipitation Kit provides all of the unique reagents required to study protein-protein interactions using a 96-well ANTI-FLAG M2 High Sensitivity Capture Plate (ANTI-FLAG M2 plate). The ANTI-FLAG M2 plate is designed to capture recombinant FLAG fusion proteins obtained from cell lysates of transfected cells.

6.1 Reagents Provided (sufficient for 96 samples)

Part	Composition	Product #	Amount
ANTI-FLAG M2 High Sensitivity Capture Plate	—	P2983	1 plate
Lysis Buffer	20 mM Tris-HCl (pH 7.4), 1 M NaCl, 1 mM DTT, 1.0% Triton [®] X-100	L1413	50 ml
10× Wash Buffer	500 mM Tris-HCl (pH 7.4), 1.5 M NaCl	W4511	50 ml
2× Sample Buffer	125 mM Tris-HCl (pH 6.8), 4% SDS, 20% (v/v) glycerol, 0.004% bromphenol blue	B2556	6 ml
Amino-terminal FLAG-BAP Fusion Protein	—	A1974	2.5 µg
SealPlate Film	—	Z36, 965-9	5

6.2 Additional Reagents and Equipment (Not Supplied)

- Pipets (single- or multi-channel)
- Pipette tips
- Microcentrifuge tubes
- Reagent reservoir
- Microcentrifuge at 4°C
- Vortexer
- MilliQ or sterile water
- Tissue culture plates
- Tissue culture medium
- Tissue culture cells, such as COS-7
- Tissue culture incubator
- Mammalian protease inhibitor cocktail
- Ice-cold PBS
- Cell scrapers
- Kit for measuring total protein concentration
- 4°C incubator
- Plate washer
- Rotary shaker

7.0 IMMUNOPRECIPITATION DETECTION KIT (PRODUCT # COIP-D)

The Immunoprecipitation Detection Kit provides all of the unique reagents required to detect protein-protein interactions using a plate-based colorimetric assay. The substrate p-Nitrophenyl Phosphate Disodium (pNPP) is hydrolyzed by Anti-c-Myc Alkaline Phosphatase-conjugated antibody to produce a yellow-colored solution that can be read spectrophotometrically at 405 nm. Tris Buffered Saline (pH 8.0) with 3% Nonfat Milk is used to dilute the Anti-c-Myc Alkaline Phosphatase Conjugate. The pNPP reaction is terminated with 3 N NaOH and then read at 405 nm. The SIGMA *FAST*[™] pNPP tablet set provided in the kit produces a ready-to-use solution after the addition of 5 ml of distilled or deionized water.

7.1 Reagents Provided (sufficient for 96 samples)

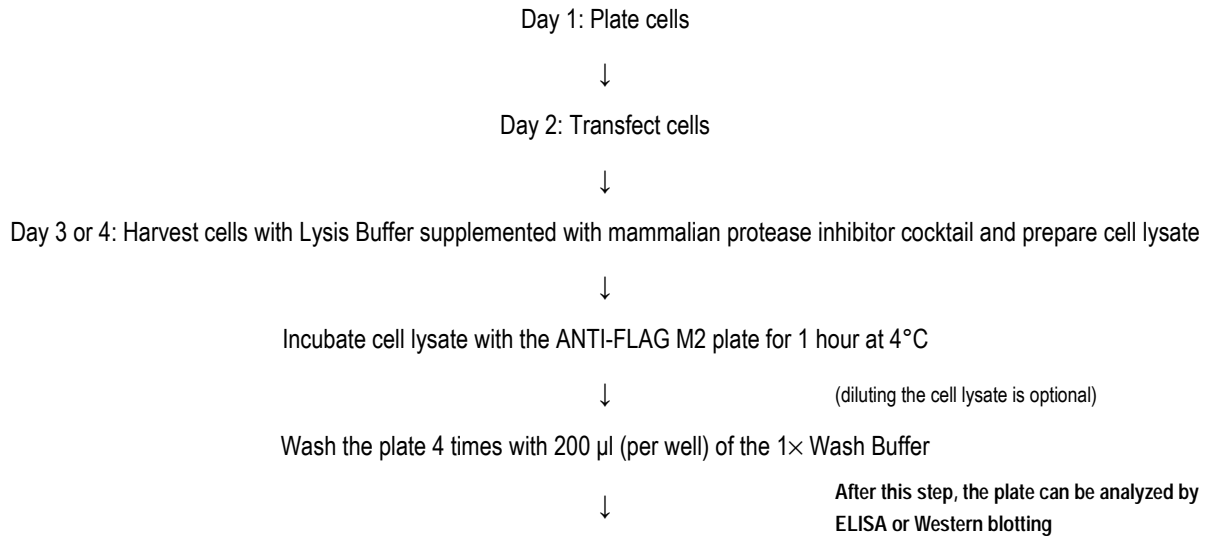
Kit Components	Product #	Quantity
Monoclonal Anti-c-Myc, Clone 9E10, Alkaline Phosphatase Conjugate	A5963	1 × 0.25 ml
Tris Buffered Saline, pH 8.0, with 3% Nonfat Milk	T8793	1 packet
SIGMA <i>FAST</i> [™] p-Nitrophenyl Phosphate Tablets	N1891	5 tablet sets

7.2 Additional Reagents and Equipment (Not Supplied)

- Pipets (single- or multi-channel)
- Pipette tips
- Microcentrifuge tubes
- Reagent reservoir
- Microcentrifuge at 4°C
- Vortexer
- Plate washer
- Plate mixer
- Rotary shaker
- MilliQ or sterile water
- 3 N NaOH
- 50 ml conical tubes
- Spectrophotometer capable of measuring absorbance at 405 nm
- β-Mercaptoethanol
- Bulldog clamps
- Boiling water bath
- SDS-PAGE gels and apparatus
- Protein transfer apparatus
- Protein transfer membrane
- Tris-glycine buffer
- TBST
- TBS
- Anti-mouse IgG peroxidase conjugate
- ANTI-FLAG M2 monoclonal antibody peroxidase (HRP) conjugate
- ECL Plus[™] Reagent

8.0 PROTOCOLS FOR STUDYING PROTEIN-PROTEIN INTERACTIONS WITH THE ANTI-FLAG M2 PLATE

8.1 Overview of the Analysis of Protein-Protein Interaction with the ANTI-FLAG M2 Plate by ELISA or Western blotting



Analysis with ELISA:

Add 200 µl of Anti-c-Myc Alkaline Phosphatase Conjugate (diluted 1:100 in Blocking Buffer) to each well and incubate for 1 hour at room temperature on a rotary shaker

↓

Wash the plate 4 times with 200 µl (per well) 1× Wash Buffer

↓

Incubate each well with 200 µl pNPP substrate

↓

Incubate at room temperature for 15 to 30 min or until the yellow color develops

↓

Stop the reaction with 3 N NaOH

↓

Measure A₄₀₅ with a plate reader

Analysis with Western Blotting:

Add 60 µl of 2× Sample Buffer (supplemented with 10% β-Mercaptoethanol) to each well and incubate at room temperature for 30 minutes (with shaking on a vortexer)

↓

Transfer the contents of each well to microcentrifuge tube and boil for 5 minutes

↓

Western Blotting

8.2 General Notes

Please read this section carefully before proceeding with the experiment.

8.2.1 Plasmid DNA for transfection

To obtain maximum transfection efficiency, the DNA used for transfections must be of high quality (free of protein, RNA, and chemical contamination, with an $A_{260}:A_{280}$ ratio of 1.8-1.9). To produce high-quality plasmid DNA, we recommend Sigma's GeneElute Endo-Free Plasmid Purification Kit (Product # PLEX 15).

8.2.2 Transfection reagent

The quality of a transfection reagent dictates the sensitivity and specificity of an assay using the FLAG 96-well Immunoprecipitation Kit. Generally, high transfection efficiency yields a more sensitive assay, requiring a smaller sample volume. The Immunoprecipitation Vector Kit includes three control plasmids for optimizing and monitoring the efficiency of transfection. To obtain high-efficiency transfections, we recommend Sigma's ESCORTII Transfection Reagent (Product # L6037) for commonly used cell lines.

8.2.3 Cell lysates used for co-immunoprecipitation assay

- 1. Expression level of the fusion proteins in the cell lysate.** The concentration of the expressed fusion protein is critical to the success of the co-immunoprecipitation analysis. The expression of the FLAG and c-Myc fusion proteins should be verified by Western blotting with ANTI-FLAG M2 antibody and anti-c-Myc antibody, prior to the co-immunoprecipitation analysis. In general, a high concentration of fusion protein in the cell lysate will yield more sensitive results. For each tag, the control plasmids supplied in the Immunoprecipitation Vector Kit can be used as positive controls in transfection and Western blot analysis. FLAG-p53 can be detected with ANTI-FLAG M2 antibody, while the c-Myc-Large T antigen and c-Myc-BAP can be detected with monoclonal anti-c-Myc antibody 9E10.
- 2. Amount of cell lysate used in the co-immunoprecipitation assay.** The optimum amount of cell lysate added to the wells of the ANTI-FLAG M2 plate may vary with each experimental system and should be determined empirically. The ANTI-FLAG M2 plate-based immunoprecipitation is very sensitive, especially in ELISA. When analyzing the interaction of FLAG-p53 and c-Myc-Large T antigen, the c-Myc-Large T antigen can be detected in as little as 5 μ g of the cell lysate from co-transfected COS-7 cells (as compared to the negative control cell lysate).
Note: The volume of the cell lysate used should not exceed 200 μ l since the wells of the ANTI-FLAG M2 plate are coated up to the 200 μ l level with the ANTI-FLAG M2 antibody.
- 3. Dilution of the cell lysate.** The relative binding affinity for different protein complexes may vary. Since high salt and detergent levels are present in the Lysis Buffer, it may be necessary to dilute the cell lysate with 1 \times Wash Buffer to minimize disruption of the protein-protein interactions.
Note: For the positive control interaction between FLAG-p53 and c-Myc-Large T antigen, dilution of the cell lysate is not necessary.
- 4. Negative-binding protein partner.** Inclusion of a negative-binding partner along with the FLAG-tagged protein construct is critical for verifying the specific protein-protein interaction observed. Over-expression of a fusion protein may cause nonspecific protein-protein interactions between the proteins being tested. The c-Myc-BAP vector provided in the Immunoprecipitation Vector Kit can be used as a negative-binding protein partner.

8.2.4 Detection of endogenous binding partners during Western blotting or ELISA

Although this kit and protocol are designed for detecting the c-Myc-tagged binding partner from co-transfected cell lysates using anti-c-Myc antibody, you may use the ANTI-FLAG M2 plate to capture endogenous binding partners that interact with the FLAG-fusion protein from cells transfected only with the FLAG-fusion protein. In this situation, optimum dilution and incubation times should be determined for the specific antibodies used to detect the endogenous binding partners during Western blotting or ELISA. Since the ANTI-FLAG M2 plate is coated with monoclonal ANTI-FLAG M2 antibody, either polyclonal antibody or antibody conjugates should be used for ELISA-based detection to avoid problems with cross-reactivity.

Note: This protocol has been optimized for analyzing the interaction of FLAG-p53 and c-Myc-Large T antigen expressed in COS-7 cells. To ensure that the system works in your hands, we recommend including the control plasmids provided in the Immunoprecipitation Vector Kit in your experiment.

8.3 Preparation of Cell Extract

1. Aspirate growth medium from the transfected cells to be assayed. Rinse the cells twice with ice-cold PBS.
2. Add 1 ml of ice-cold Lysis Buffer supplemented with mammalian protease inhibitors (10 μ l per 1 ml Lysis Buffer, Product # P8340) to a 10-cm plate (you can adjust the volume of the Lysis Buffer according to the size of the culture plate used).
3. Scrape the cells off the plate with cell scrapers (Product # C2802) and transfer them to microcentrifuge tubes.
4. Incubate for 1 hour on ice, occasionally mixing with a vortexer.
5. Microcentrifuge the cells for 15 minutes at 4°C and transfer the supernatant to a new tube. If the supernatant is not to be used immediately, store the cell lysate at -70°C.

Note: The supernatant may be viscous during transfer, due to the genomic DNA released from the nuclei of lysed cells. Genomic DNA can be removed from the supernatant with a 1 ml pipet tip prior to the transfer.

6. Determine the total protein concentration of the supernatant using standard techniques. We recommend using Sigma's BCA Protein Assay. (Bicinchoninic Acid Kit, Cat. No BCA-1).

8.4 Immunoprecipitation with ANTI-FLAG M2 plate

Note: For protein complexes where the interaction occurs at salt and detergent concentrations different from the 1 \times Lysis Buffer, use the 1 \times Wash Buffer to dilute the cell lysate. If dilution is required, first add the required amount of the 1 \times Wash Buffer to the plate and then add the cell lysate. The ANTI-FLAG M2 plate is resistant to the following detergents: 5% Tween[®]-20, 5% Triton[®] X-100, 0.1% Igepal CA-630, 0.1% CHAPS, and 0.2% digitonin. The ANTI-FLAG M2 plate can also be used with 1.0 M NaCl.

1. Rehydrate the plate by adding 200 μ l of 1 \times Wash Buffer to each well.
2. Incubate the plate for 2 minutes at room temperature.

Note: If the entire plate is not used at once, cover the unused wells firmly with SealFilm to prevent contamination from the used wells.

3. Remove the liquid from each well manually or with a plate washer. Tap the plate firmly onto a paper towel until most of the solution is removed.

4. Incubate various amounts of cell lysate on the plate using a rotary shaker (with gentle rocking) for 1 hour at 4°C. Cover the plate with the lid.
Note: The ANTI-FLAG M2 plate is coated with 200 µl of ANTI-FLAG M2 antibody. We recommend that the volume of cell lysate added to each well should not exceed 200 µl.
5. This kit includes Amino-terminal FLAG-BAP Fusion Protein as a positive control for immunoprecipitation on the plate. For a positive control, add 0.5 µg of FLAG-BAP fusion protein (Product # A1974) diluted with 150 µl Lysis Buffer (Product # L1413) to a single well. This control sample should be treated identically to all other samples.
6. Wash each well on the plate 4 times with 200 µl of 1× Wash Buffer. The plate can be washed either manually or with a plate washer. After washing, remove excess liquid by firmly tapping the plate onto a paper towel.
Note: After this step, the captured proteins can be analyzed by two methods: ELISA (Section 8.5.1) or Western blotting (Section 8.5.2).

8.5 Detection of the Bound Protein by ELISA or Western Blotting

After the FLAG and c-Myc-tagged protein complexes have been captured (Section 8.4), they can be detected using either ELISA or Western blotting.

8.5.1 ELISA

1. Dilute Monoclonal Anti-c-Myc, Clone 9E10, Alkaline Phosphatase Conjugate (Product # A5963) 1:100 with Blocking Buffer (Tris Buffered Saline [pH 8.0] with 3% Nonfat Milk). Add 200 µl of the diluted antibody to each well. Cover the plate with the lid to minimize evaporation.
2. Incubate the plate for 1 hour at room temperature on a rotary shaker with gentle agitation.
3. During the incubation, prepare the pNPP substrate solution (provided in the Detection Kit) and a 3 N NaOH Stopping Solution:
 - a. Prepare enough pNPP solution for each well being used in the assay. To dissolve 1 tablet set, add 5 ml of MilliQ water to a sterile 50 ml conical tube and add one SIGMA FAST™ p-Nitrophenyl Phosphate Tablet set (one buffer tablet and one pNPP tablet; Product # N1891) to the water.
 - b. Vortex until the tablets dissolve. The pNPP substrate solution should be used within one hour.
4. At the end of incubation, remove the liquid from each well and wash the wells 4 times with 200 µl each of 1× Wash Buffer.
5. Add 200 µl of pNPP substrate to each well.
6. Incubate the plate at room temperature for 15 to 30 minutes or until a yellow color develops.
Note: The time required to develop the yellow color will vary for different cell lysates. The higher the concentration of fusion proteins in the cell lysate, the quicker the yellow color will develop.
7. Stop each reaction with 50 µl of the 3 N NaOH Stopping Solution. Mix gently by pipetting.
Note: Avoid introducing bubbles into the wells. If bubbles are present, remove them by gently tapping the plate prior to reading the absorbance.
8. Using a spectrophotometer, measure the absorbance of each well at 405 nm.

8.5.2 Western Blotting

1. Add 60 µl of 2× Sample Buffer (supplemented with 10% β-mercaptoethanol) to each sample well on the plate. Seal the plate with SealPlate Film (Product # Z36, 965-9) and clamp the plate to a Vortex Genie 2 Mixer (or an equivalent device) with two bulldog clamps.

Notes:

1. **Warm the 2× Sample Buffer (Product # B2566) at 37°C for a few minutes before use.**
2. **Do not cover the plate with a plastic lid, since this combination may be too thick for the bulldog clamps.**
2. Shake the plate for 30 minutes at room temperature on the vortexer at a setting of 3.
3. Transfer the entire contents of each well to separate microcentrifuge tubes.
4. Boil each sample for 5 minutes.
5. Load approximately 50 µl of each sample onto a standard SDS-PAGE gel.
6. Transfer the proteins from the gel to a membrane such as PVDF using Tris-Glycine Buffer (25 mM Tris, 192 mM glycine, 15% v/v methanol [pH 8.3]; Product # T4904) and the settings recommended by the manufacturer of your transfer apparatus.
7. Block the membrane with Blocking Buffer at room temperature for 30 to 60 minutes.
8. Transfer the membrane to a tray containing a 1:2,000 dilution of ANTI-FLAG M2 Monoclonal Antibody Peroxidase (HRP) Conjugate (Product # A8592) or a 1:100 dilution (in Blocking Buffer) of Monoclonal anti-c-Myc Unconjugated, Clone No. 9E10 (Product # M5546).
9. Incubate at room temperature for 1 hour with gentle agitation.
10. Wash the membrane three times (5 minutes per wash) with TBST at room temperature with gentle agitation.
11. For membranes probed with ANTI-FLAG M2 Monoclonal Antibody Peroxidase (HRP) Conjugate, develop the membrane with ECL Plus™ Reagent according to the manufacturer's instructions.
12. For membranes probed with Monoclonal anti-c-Myc Unconjugated, Clone No. 9E10, incubate the membrane with anti-mouse IgG peroxidase conjugate (Product # A9044) as the secondary antibody, diluted 1:80,000 in Blocking Buffer.
13. Incubate the membrane at room temperature for 30 minutes with gentle agitation.
14. Wash the membrane three times (10 minutes per wash) with TBST at room temperature with gentle agitation.
15. Rinse the membrane once with TBS (Product # T6664) to remove residual Tween-20.
16. Develop the membrane with ECL Plus™ Reagent according to the manufacturer's instructions.

9.0 TROUBLESHOOTING GUIDE

Problem	Possible Cause	Solution
Precipitation of the cell lysate after mixing with 2× Sample Buffer for Western blotting.	You may experience a precipitation problem when you mix undiluted cell lysate from certain mammalian cell lines with 2× Sample Buffer. The Lysis Buffer is very efficient at solubilizing proteins from cells, which sometimes results in too much protein in the lysate and causes aggregation.	Add 6 -8 M urea and 10% β-mercaptoethanol to the 2× Sample Buffer and then mix with the cell lysate. Heat at 100°C for 10 minutes before loading on the SDS-PAGE gel.
Samples exhibit no signal or low signal in both Western blotting and ELISA after capturing on the ANTI-FLAG M2 plate.	The concentration of the fusion protein in the cell lysates may be low.	Check the expression level of the fusion proteins in the co-transfected cell lysate by Western blotting using ANTI-FLAG M2 antibody to detect the FLAG-tagged fusion protein or anti-c-Myc antibody to detect the c-Myc-tagged fusion protein. If expression level is low, optimize the conditions for transfection with the control plasmids.
	High salt and detergent present in the Lysis Buffer may interrupt protein-protein interactions.	Dilute the cell lysate 1- to 5-fold with 1× Wash Buffer (e.g., 1 volume of cell lysate mixed with 1 to 4 volumes of 1× Wash Buffer) before immunoprecipitation with ANTI-FLAG M2 plate.
	In the ELISA analysis, the pNPP substrate may have degraded.	Prepare a fresh stock of pNPP and use it within 1 hour. To determine the integrity of the substrate, see Appendix 3 .
Samples exhibit high signal in ELISA.	Too much cell lysate added to the ANTI-FLAG M2 plate.	Dilute the cell lysate. You can do a titration series of the cell lysate to be analyzed and find the optimum amount of cell lysate to generate an A ₄₀₅ reading between 0.5 and 1.0.
High background during ELISA.	Incubation time with pNPP substrate was too long.	Decrease the incubation time with pNPP.
	It is unlikely that the background is contributed by nonspecific interactions between the plate and cellular proteins, since the ANTI-FLAG M2 plate is highly specific. It is likely that the pNPP substrate incubation time was too long or the substrate was previously degraded.	Decrease the incubation time with the pNPP substrate or prepare fresh substrate. To examine the integrity of the substrate, please see Appendix 3 .
Low yield of eluted protein for Western blotting analysis after capturing proteins on the ANTI-FLAG M2 plate.	Loss of the 2× Sample Buffer (with 10% β-mercaptoethanol) during the elution step	Do not vortex the plate too vigorously, which may spill the 2× Sample Buffer Sample Buffer (with 10% β-mercaptoethanol) into adjacent wells. Set the vortexer at a speed that provides adequate motion for covering the upper portions of each well.
	No β-mercaptoethanol in the 2× Sample Buffer	10% β-mercaptoethanol is critical for efficient elution of bound proteins. Add this to your 2× Sample Buffer.

10.0 REFERENCES

1. Hopp, T.V., Prickett, K.S., Price, R.T., Libby R.T., March, C.J., Cerretti, D. Urdal, D.L. and Conlon, P.J. *Bio/Technology* **6**, 1204-1210 (1988).
2. Evan, G .I., Lewis, G.K., Ramsay, G., and Bishop, V.M. *Mol. Cell. Biol.* **5**, 3610-3616 (1985).
3. Iwabuchi, K.M., Li, B., Bartel, P. and Fields, S. *Oncogene* **8**, 1693-1696 (1993).
4. Li, B., and Fields, S., *FASEB J.* **7**, 957-963 (1993).
5. Sambrook, J., Fritsch, E.F., and Maniatis, T. *Molecular Cloning: A Laboratory Manual* (Cold Spring Harbor Laboratory, Cold Spring Harbor, NY) (1989).

APPENDIX 1: CONSTRUCTION OF FUSION GENES USING STANDARD CLONING TECHNIQUES

A brief outline of the recommended cloning technique is supplied below (for more information, please refer to Sambrook *et al.*, 1989).⁵ To test the interaction between protein X and Y, the gene for protein X is cloned into the pFLAG-CMV[™]-2 vector and the gene for protein Y is cloned into the pc-Myc-CMV-2 vector.

1. DNA encoding protein X and protein Y can be generated either by PCR amplification or by cutting them out of another plasmid.
2. Purify the DNA fragments with a DNA purification kit designed for PCR analysis. We recommend using Sigma's GenElute[™] PCR DNA Purification Kit (Product # GEN-PCR).
3. Digest both vectors with the appropriate restriction enzyme(s), dephosphorylate, and purify the DNA.
4. Ligate the vector and insert.
5. Transform the ligation mixtures into competent *E. coli* cells.
6. Plate the transformed *E. coli* on LB-Ampicillin agar plates.
7. Screen clones using colony PCR or restriction analysis to identify those containing the correct insert.

APPENDIX 2: SCREENING OF RECOMBINANT PLASMIDS BY COLONY PCR

This procedure describes colony PCR using the vector primers provided in the kit. We recommend that you use one gene-specific primer and one of the vector primers to perform colony PCR (e.g., a forward gene-specific primer and the Verification Primer-MR, or Verification Primer-MF and a reverse gene-specific primer). This method allows the identification of recombinant clones with the right orientation. In addition to your experimental sample, we recommend that you perform a positive control PCR using one of the control plasmids as the template in conjunction with the vector primers. This control reaction will confirm that the PCR was successful. If control plasmids and the vector primers are used for the positive control PCR, the expected PCR fragments are shown in the table below:

Plasmid	Product #	Sizes of Expected PCR fragment (kb)
pFLAG-CMV™-2-p53 Control Plasmid	P9986	1.3 kb
pc-Myc-CMV-2-Large T antigen Control Plasmid	P9861	2.4 kb
pc-Myc-CMV-2-BAP Control Plasmid	P9736	1.7 kb

Procedure for Screening Recombinant Plasmids

1. Pick single colonies from the LB-Ampicillin agar plate using a sterile pipet tip.
2. Deposit the bacteria at the bottom of individual PCR tubes or the wells of a multiwell PCR plates (Product # Z37, 490-3) containing 20 µl of water (Product # W1754) and vortex.
3. Aliquot 5 µl of the colony mixture into separate PCR tubes or a multiwell PCR plate and set it aside. The rest of the mixture can be left at room temperature until the PCR is completed.
4. Prepare a PCR master mix with Sigma's ReadyMix™ Taq (Product # P4600). Prepare enough PCR master mix for all of the PCR tubes. For each PCR (30 µl total volume), mix the following reagents in a sterile microcentrifuge tube:
 - a. 15 µl of 2× PCR ReadyMix™ Taq
 - b. 1 µl of Forward primer (20 µM Verification Primer-MF)
 - c. 1 µl of Reverse primer (20 µM Verification Primer-MR)
 - d. 8 µl of Water

Note: If gene-specific primers are used in the PCR reaction, we recommend that you make a stock concentration of 20 µM and use 1 µl of this stock per PCR.
5. Mix gently and briefly centrifuge each sample to collect all of the components at the bottom of the tube.
6. Add 25 µl of the PCR master mix (prepared in **Step 4**) into the PCR tubes or the multiwell PCR plate containing 5 µl of the colony mix.

7. Start the PCR using a thermal cycler with a heated lid. The amplification parameters should be optimized for each unique combination of primers, template, and thermal cycler. If the vector primers are used, the cycling parameters should be:
 - a. Incubate at 94°C for 5 minutes
 - b. 25 cycles of: 94°C for 40 seconds, 55°C for 50 seconds, and 72°C for 2 minutes.
 - c. Extend at 72°C for 10 minutes

Note: The proposed extension time is suitable for the amplification of PCR products 0.2 to 2 kb in length. You can alter the extension time in proportion to the length of the expected PCR fragments (generally, 1 kb/minute).
8. When the PCR is complete, analyze 15 µl of the amplified DNA from each PCR by electrophoresis in an agarose/EtBr gel.
9. When positive PCR products are identified, match them to the original PCR tubes or multiwell PCR plates containing the colony mix.
10. Take the positive colony mix and use it to inoculate 3 ml of fresh LB medium containing 100 µg/ml ampicillin in a 15 ml culture tube.
11. Incubate the culture at 37°C overnight.
12. Isolate plasmid DNA. We recommend using Sigma's GenElute™ plasmid Purification Kit (Product # PLN-70).

Note: The vector primers can also be used for DNA sequencing.

APPENDIX 3: CAPTURING THE AMINO-TERMINAL FLAG-BAP FUSION PROTEIN WITH THE ANTI-FLAG M2 PLATE AND EVALUATING THE pNPP SUBSTRATE INTEGRITY

1. Dilute the Amino-Terminal FLAG-BAP Fusion Protein (Product # A1974) or nontransfected control cell lysate to a concentration of 2.5 ng/μl with 1× Wash Buffer.
2. To rehydrate the wells, add 200 μl of 1× Wash Buffer to each of 6 wells and incubate for 2 minutes.
3. Remove the solution from each well and firmly tap the plate on a paper towel until most of the 1× Wash Buffer has been removed.
4. Add 200 μl of Amino-Terminal FLAG-BAP Fusion Protein solution (prepared in **Step 1**) into each well (500 ng/well).
5. Cover the plate with a lid.
6. Incubate the plate at 4°C for 1 hour on an orbital shaker operating at 100 rpm.
Note: During the incubation, prepare the pNPP substrate solution.
 - a. Do not prepare the pNPP solution until the plate is incubating, as it must be used within 60 minutes.
 - b. Add 5 ml of MilliQ water to a sterile 50 ml conical tube.
 - c. Add one SIGMA FAST™ p-Nitrophenyl Phosphate Tablet set (one buffer tablet and one pNPP tablet; Product # N1891) to the water.
 - d. Vortex until the tablets dissolve.
7. Remove the solution from each well and firmly tap the plate on a paper towel until most of the solution has been removed.
8. Wash each well 4 times with 200 μl of 1× Wash Buffer.
9. Remove the solution and tap the plate dry onto a paper towel.
10. Add 200 μl of the pNPP solution to each well.
11. Incubate at room temperature for 15 minutes.
12. Immediately at the end of the incubation, add 50 μl of 3 N NaOH to each well (keep the pipette tips submerged in the wells).
13. Mix the solution several times by pipetting. This step stops the color development and distributes the yellow color evenly in each well.
Note: Avoid the formation of bubbles. If bubbles form, disrupt them with a clean pipette tip.
14. Using a spectrophotometer, measure the A_{405} with a plate reader. If the control immunoprecipitation worked correctly and the pNPP solution is fresh (and prepared properly), the A_{405} reading should be between 0.5 and 0.7.