

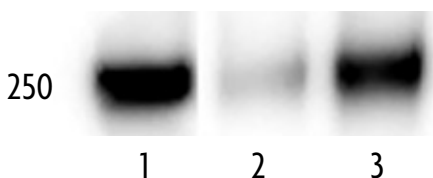
# Nuclear Complex Co-IP Kit

optimized co-immunoprecipitation of nuclear protein complexes

Co-Immunoprecipitation (Co-IP) is a powerful method used to study protein/protein interactions. However, traditional Co-IP methods are not optimal for studying DNA-binding protein complexes as these complexes are often disrupted during the extraction process. Active Motif's Nuclear Complex Co-IP Kit simplifies Co-IP of nuclear protein complexes by providing high-quality reagents for both nuclear extract preparation and immunoprecipitation. The kit's extraction process offers a simple and effective method to obtain and maintain protein complexes contained in nuclear compartments of the cell, especially complexes that had been bound to DNA. And, the versatile Co-IP reagents offer the flexibility to vary the stringency of the buffer compositions, enabling the study of tightly bound or weak protein complexes.

## The Nuclear Complex Co-IP Kit advantage

- Simple and efficient
- Optimized extraction procedure preserves nuclear protein complexes
- Easily alter IP stringency to detect interactions of varying strengths



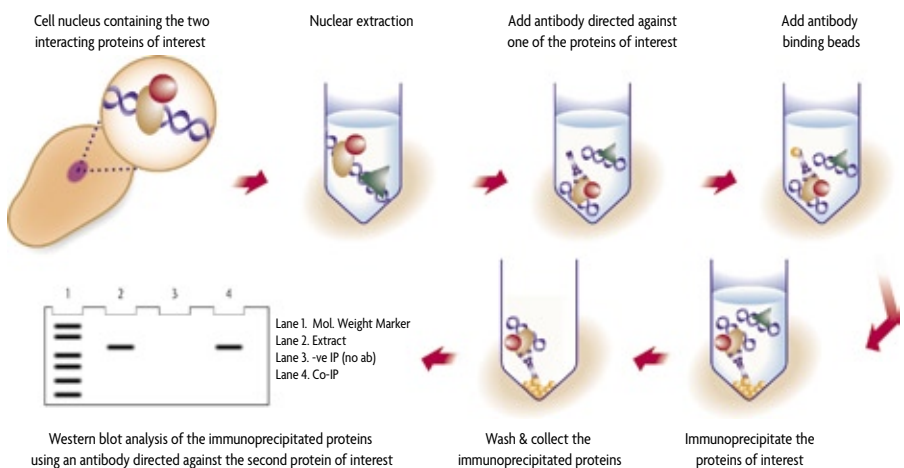
**Figure 2: Western blot analysis of the IP'd p33 subunit of RNA pol II.**

HeLa nuclear extracts were prepared using the kit's extraction reagents. IP with the p33 protein was performed with 100 µg of nuclear extract, followed by Western blot analysis with an RNA pol II mouse mAb. Detection of the p33/RNA pol II complex by the RNA pol II antibody (lane 3) demonstrates that the Co-IP was successful in maintaining the protein complex.

Lane 1: Western blot control

Lane 2: Negative Control (no antibody used in IP)

Lane 3: Co-IP: IP using p33/WB using RNA pol II



**Figure 1: Schematic of Co-Immunoprecipitation procedure.**

Product	Format	Catalog No.
Nuclear Complex Co-IP Kit	50 rxns	54001

### CONTENTS

Each Nuclear Complex Co-IP Kit supplies sufficient reagents to perform 50 co-immunoprecipitation experiments and provides Hypotonic Buffer, PBS, Phosphatase Inhibitors, Protease Inhibitor Cocktail, PMSF, EDTA, Digestion Buffer, Enzymatic Shearing Cocktail, Detergent, IP High and IP Low Buffers, 5M NaCl, BSA and DTT.