

Figure 5. MCM6 IF results

MCM6 antibody detects MCM6 protein at nucleus by immunofluorescent analysis.

Sample: U2OS cells were fixed in 4% paraformaldehyde at RT for 15 min.

Green: MCM6 protein stained by MCM6 antibody diluted 1:500.

Blue: Hoechst 33342 staining.

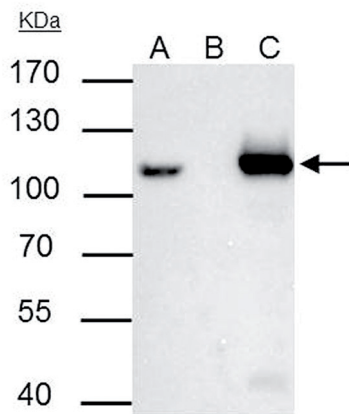


Figure 6. MCM6 IP results

MCM6 antibody immunoprecipitates MCM6 protein in IP experiments. IP Sample: 293T whole cell extract A : 30 µg whole cell extract of MCM6 protein expressing 293T cells B : Control with 3 µg of pre-immune rabbit IgG C : Immunoprecipitation of MCM6 by 3 µg of MCM6 antibody. The immunoprecipitated MCM6 protein was detected by western blot using the MCM6 antibody diluted 1:1,000.