

---

## CRISPR/Cas9 Monoclonal Antibody [7A9]

(Catalog # A-9000)

---

### Background

The discovery of CRISPR (Clustered Regularly Interspaced Short Palindromic Repeats) and Cas9 (CRISPR associated system or CRISPR associated protein 9 nuclease) found in bacteria to work as a defense mechanism against foreign DNA has proven to be an invaluable tool to target and modify a genetic sequence in gene editing and genome engineering applications. The system, known as CRISPR/Cas9, allows for sequence-specific cleavage of a targeted genomic locus by delivering the RNA-guided nuclease (Cas9) and appropriate guide RNAs (gRNA) into a cell. In addition, Protospacer Adjacent Motif (PAM) sequence immediately following the specificity sequence is necessary for successful binding of the Cas9 nuclease.

### Concentration

1 mg/ml

### Description

Mouse monoclonal antibody raised against CRISPR/Cas9, clone 7A9, generated with synthesized peptide corresponding to sequence of *Streptococcus pyogenes* (*S. pyogenes*) CRISPR-associated endonuclease Cas9/Csn1. This Anti-Cas9 mAb can detect CRISPR/Cas9 expression in target cells by WB, IF, IP, or ELISA to confirm and verify whether gRNA and Cas9 vectors are successfully transfected.

### Specificity

Recognizes both Cas9 and dCas9 (nuclease deficient Cas9)

### Reactivity

Species Independent

### Isotype

IgG1/Kappa

### Immunogen

Recombinant N-terminal fragment of *S. pyogenes* Cas9 protein expressed in *E. coli*.

### Formulation

PBS, 30% Glycerol

### Storage

Store at 4°C for short-term storage (1-2 weeks). For long-term storage, aliquot and then store at -20°C. Avoid multiple freeze/thaw cycles.

### Purity

Protein G Purified

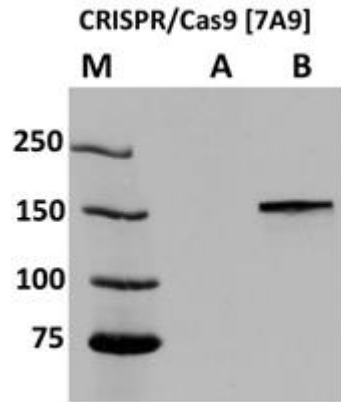
### Alternative Names

Anti-Cas9, Anti-CRISPR, Anti-CRISPR/Cas9, CRISPR antibody, Cas9 antibody, Cas9 7A9, 7A9-3A3

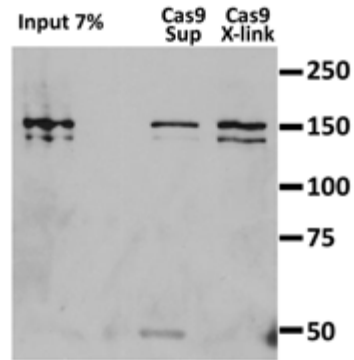
### Application

WB (1:200-1:500), IP (2 µg/10<sup>6</sup> cells), IF (1:200-1:500), ELISA (1:1000-1:2000)

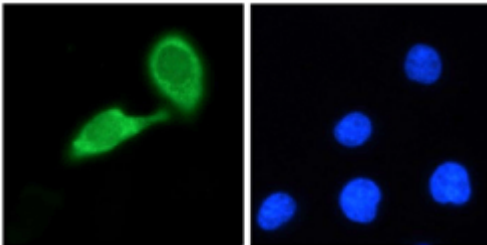
---



Western Blot: Shown are the results of WB on protein extracts from untransfected (A) and transfected (B) HEK293 cells using the Anti-CRISPR-Cas9 mAb.



Immunoprecipitation: HEK293T expressing N-terminally Flag-tagged S.pyogenes Cas9 were lysed 72 hours post transfection. Proteins were immunoprecipitated from 100 µg of whole cell lysate for 1H at 4



Immunofluorescence: HeLa cells were transiently transfected with an N-terminally Flag-tagged S. pyogenes Cas9 expression vector. The cells were stained with the Anti-Cas9 monoclonal antibody followed by anti-mouse-AF488 coupled secondary antibody. Nuclei were counter-stained with Hoechst 33342.