

# Cas9 monoclonal antibody (7A9)

Cas9 (CRISPR associated protein 9) is an RNA-guided DNA endonuclease enzyme associated with the CRISPR (Clustered Regularly Interspersed Palindromic Repeats) adaptive immunity system in *Streptococcus pyogenes* and other bacteria. Cas9 is used to interrogate and cleave foreign DNA, such as invading bacteriophage DNA or plasmid DNA. This is accomplished by unwinding foreign DNA and checking for if it is complementary to the 20 base pair spacer region of the guide RNA. If the DNA substrate is complementary to the guide RNA, Cas9 cleaves the invading DNA.

Cas9 is becoming a prominent tool in the field of genome editing because it can cleave nearly any sequence complementary to the guide RNA. Because the target specificity of Cas9 stems from the guide RNA, engineering Cas9 to target new DNA is straightforward, and versions of Cas9 that bind but do not cleave cognate DNA can be used to control transcriptional activation and repression.

This antibody is covered by our [Worry-Free Guarantee](#).

## Ordering Information [Order Online »](#)

ENZ-ABS294-0100	100µg
-----------------	-------

## Manuals, SDS & CofA [View Online »](#)

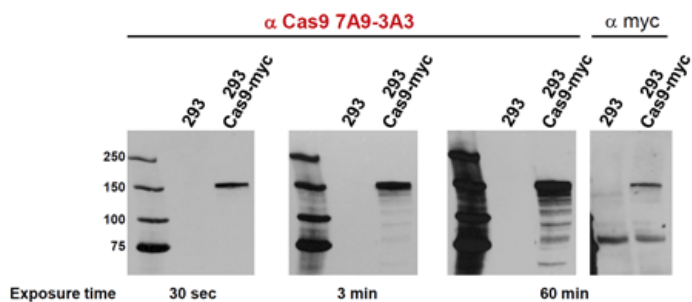


Figure 1. Western blot. HEK293 cells were transiently transfected with pMLM3613myc and lysed in hot Laemmli buffer 40 hours post transfection. Proteins were separated by 7.5% SDS-PAGE and transferred to nitrocellulose. Membranes were blocked in 3% dry milk (NFDM) in PBS-Tween for 1 h at RT and incubated with the indicated antibodies diluted in 0.5% NFDM in PBS-T for 2 hours at RT, followed by incubation with HRP-coupled anti-mouse antibody together with antiBLUE protein marker antibody for 1 hour at RT. Proteins were visualized by ECL.

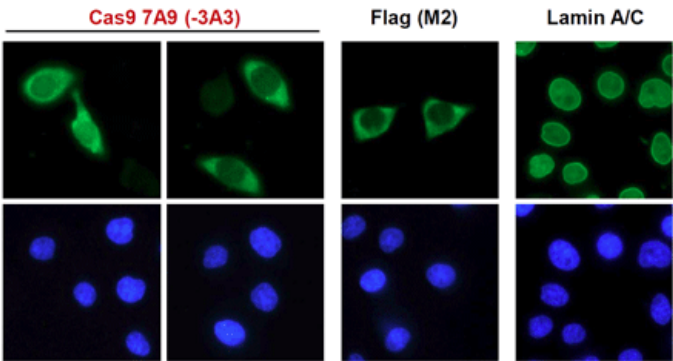


Figure 2. Immunofluorescence analysis. HeLa cells were transiently transfected with an N-terminally Flag-tagged *S. pyogenes* Cas9 expression vector. 48 hours post transfection cells were fixed in 3.7% formaldehyde, permeabilized in 0.5% Triton-X and blocked in 2% BSA in PBS for 2 hours at RT. Cells were stained with the indicated antibodies at 4C o/n, followed by incubation with anti mouse-AF488 coupled secondary antibody for 1 h at RT. Nuclei were counter-stained with Hoechst 33342. Images were taken with an Axiovert 135 TV fluorescence microscope. For presentation purposes, images were adjusted for brightness and contrast.

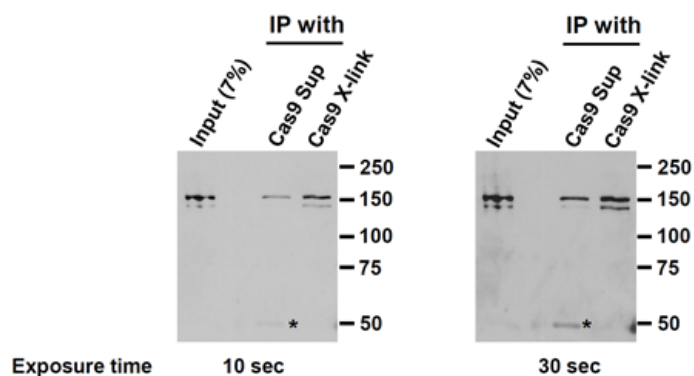


Figure 3. Immunoprecipitation. HEK293T cells expressing N-terminally Flag-tagged *S. pyogenes* Cas9 were lysed 72 hours post transfection by resuspending the cells in Hunt buffer and subjecting them to three freeze-thaw cycles in liquid nitrogen/on ice. Proteins were immunoprecipitated from 100 µg of whole cell lysate either for 1 hour at 4°C with crude Cas9 clone 7A9 cell culture supernatant (IgG concentration n.d., lane 3) followed by incubation for 1 hour at 4°C with a 1:1 mixture of protein A/G sepharose beads, or for 2 hours at 4°C with Cas9 clone 7A9 monoclonal antibody crosslinked to a 1:1 mixture of protein A/G sepharose beads. Beads were washed two times with Hunt buffer and once with TBS. Bound proteins were eluted by boiling in Laemmli buffer, separated by 7.5% SDS-PAGE and transferred to nitrocellulose. Membranes were blocked in 3% dry milk (NFDM) in PBS-Tween for 1 h at RT and incubated with Cas9 7A9 monoclonal antibody diluted in 0.5% NFDM in PBS-T for 1 hour at RT, followed by incubation with HRP-coupled anti-mouse antibody. Proteins were visualized by ECL. Asterisk indicates IgG heavy chain.

# Handling & Storage

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

# Product Details

Alternative Name	CRISPR-associated endonuclease Cas9/Csn1, CRISPR associated protein 9, SpyCas9
Application	IF, IP, WB
Clone	7A9
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Mouse
Immunogen	N-terminal fragment of Streptococcus pyogenes Cas9.
Isotype	IgG1κ
Purity Detail	Protein G affinity purified.
Recommendation Dilutions/Conditions	Immunofluorescence (1:500-1:2,000)Immunoprecipitation (1:20)Western Blot (1:1,000)Optimal conditions must be determined individually for each application.
Source	Purified from mouse ascites.
Specificity	Recognizes both Cas9 and dCas9 (nuclease deficient Cas9).
UniProt ID	Q99ZW2
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)