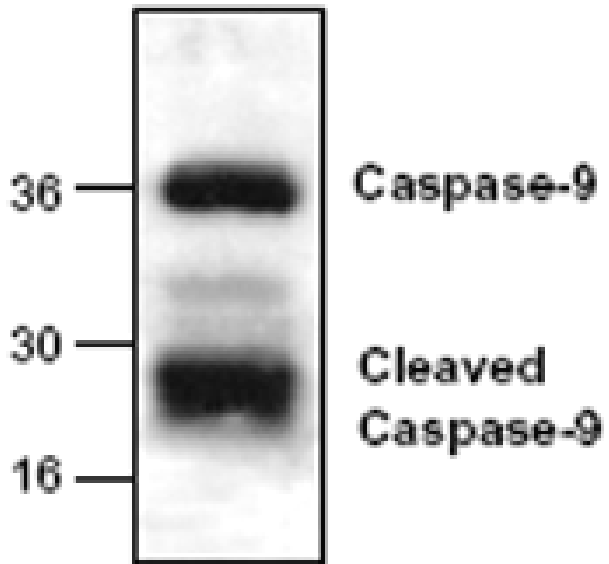


# Caspase-9 polyclonal antibody

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

Manuals, SDS & CofA

[View Online »](#)



**Figure 1:** Western blot analysis of caspase 9 with rat kidney tissue lysate using Caspase-9, pAb (prod. No. ALX-210-815).

## Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Application	WB
Application Notes	Detects procaspase-9 and a caspase-9 cleavage product by Western blot.
Formulation	Liquid. In PBS, pH 7.2, containing 30% glycerol, 0.5% BSA, 5mM EDTA, and 0.03% proclin.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to aa 287-306 of human caspase-9.
Purity Detail	Epitope-affinity purified.
Recommendation Dilutions/Conditions	Western Blot (0.5-4µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Bovine, Dog, Hamster, Human, Monkey, Mouse, Porcine, Rabbit, Rat
UniProt ID	P55211
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>

Last modified: May 29, 2024  
Enzo Life Sciences, Inc.  
Phone: 800.942.0430  
info-  
usa@enzolifesciences.com

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

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

France  
Phone: +33 472 440 655  
info-  
fr@enzolifesciences.com

Germany  
Phone: +49 7621 5500 526  
info-  
de@enzolifesciences.com

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