

# FOXP3 (mouse) polyclonal antibody

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

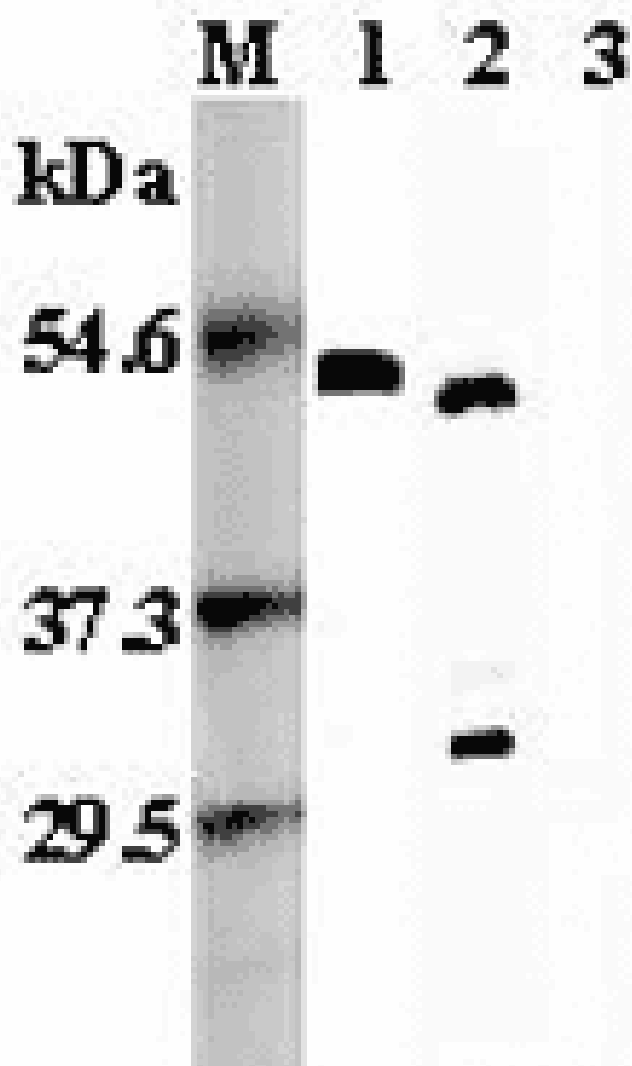
## Ordering Information

[Order Online »](#)

ALX-210-436-C100	100µg
------------------	-------

## Manuals, SDS & CofA

[View Online »](#)



**Figure:** Western blot analysis of mouse FOXP3 with PAb to FOXP3 (mouse) (Prod. No. ALX-210-436) at 1:5'000 dilution.

1. Recombinant mouse FOXP3 protein (His tagged)
2. Transfected mouse FOXP3 full length cell lysate (HEK 293)
3. Transfected vector only cell lysate

# Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

# Product Details

Alternative Name	Forkhead Box protein P3, Scurfin
Application	ELISA, WB
Application Notes	Detects a band of ~46-49kDa by Western blot.
Formulation	Liquid. In PBS, pH 7.4. Contains no preservatives.
Host	Rabbit
Immunogen	Recombinant mouse FOXP3 (forkhead box protein P3).
Purity Detail	Purified IgG.
Recommendation Dilutions/Conditions	ELISA (direct and indirect: 1:5,000-1:10,000)Western Blot (1:2,000-1:5,000 using ECL; suggested blocking and dilution buffer is TBST containing 0.05% Tween 20 and 5% skim milk; suggested incubation time is 1 hour at room temperature)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Mouse
Specificity	Recognizes both natural and recombinant FOXP3.
UniProt ID	Q99JB6
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)