

# TRAF5 polyclonal antibody (1847)

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

Citations: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-210-027-R050	50µl
------------------	------

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored at -80°C.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-80°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	Tumor necrosis factor (TNF) receptor-associated factor 5
Application	IHC (PS), WB
Application Notes	Detects a band of ~64kDa by Western blot.
Crossreactivity	Does not cross-react with other members of the TRAF family.
Formulation	Liquid. In PBS containing 0.05% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to aa 1-20 (M1AHSEEQAAPCAFIRQNSG20) of mouse TRAF5 (tumor necrosis factor (TNF) receptor-associated factor 5).
Recommendation Dilutions/Conditions	Immunohistochemistry (paraffin sections, 1:500)Western blot (1:500-1:2,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse
UniProt ID	P70191
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-  
fr@enzolifesciences.com](mailto:info-fr@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)