

# TAK1 polyclonal antibody

TAK1 (TGF-beta-activated kinase-1) is a 65 kDa serine/threonine kinase that is a member of the MAPKKK family. TAK1 is involved in the regulation of transcription by the TGF-beta super family, as its kinase activity is stimulated by TGF-beta and bone morphogenetic protein (BMP).

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

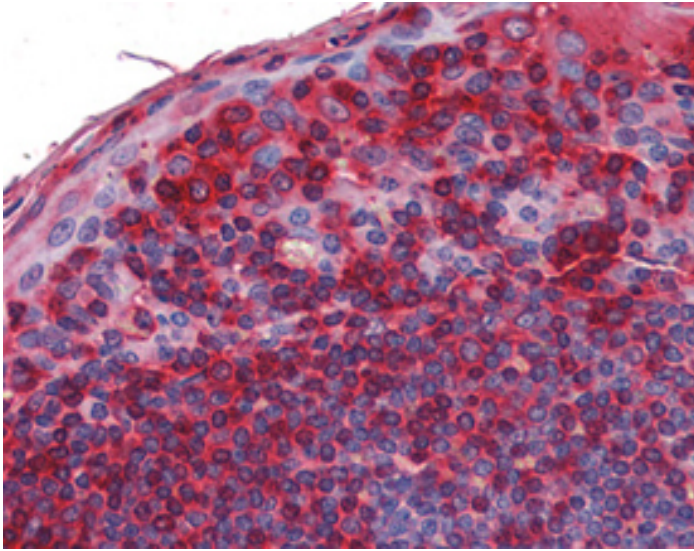
## Ordering Information

[Order Online »](#)

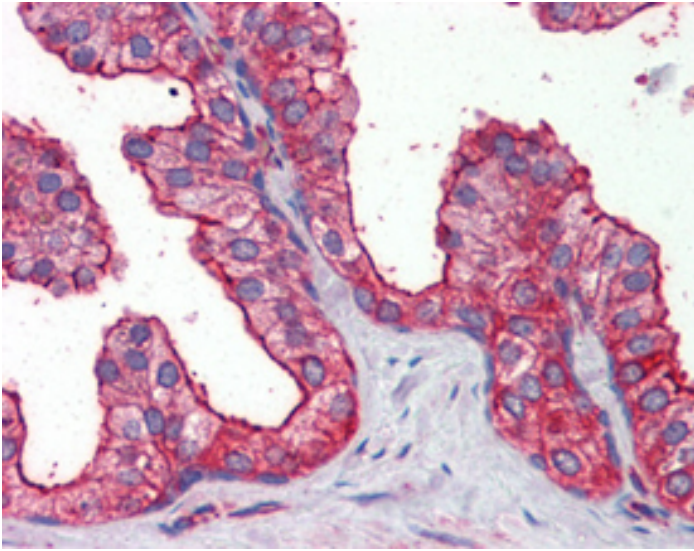
ADI-KAS-ST009-E	100µg
-----------------	-------

## Manuals, SDS & CofA

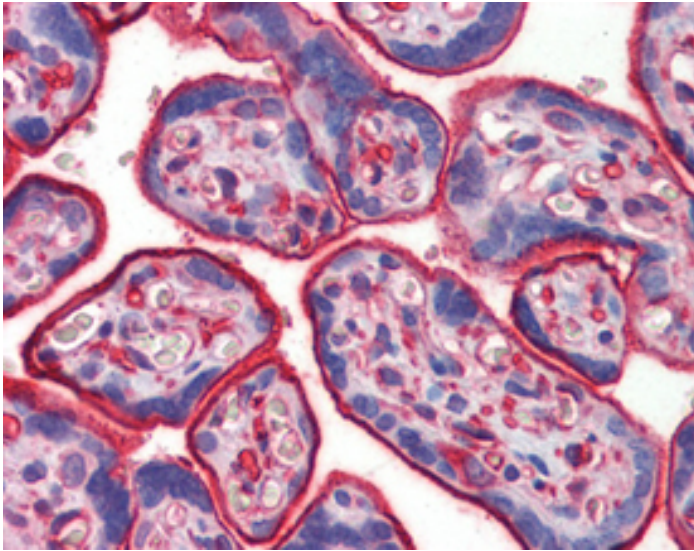
[View Online »](#)



IHC Analysis: Human tonsil tissue stained with TAK1, pAb at 10µg/ml.



IHC Analysis: Human prostate tissue stained with TAK1, pAb at 10µg/ml.



IHC Analysis: Human placenta tissue stained with TAK1, pAb at 10µg/ml.

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

## Product Details

Alternative Name	TGF-β-activated kinase-1
Application	IHC (PS), WB
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.
GenBank ID	D76466
Host	Rabbit
Immunogen	Synthetic peptide corresponding to a portion of mouse TAK1.
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1µg/ml, colorimetric)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from rabbit serum.
Species Reactivity	Bovine, Chicken, Dog, Guinea pig, Hamster, Human, Monkey, Mouse, Porcine, Rabbit, Rat, Sheep, Xenopus
UniProt ID	Q62073
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)