

# 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).

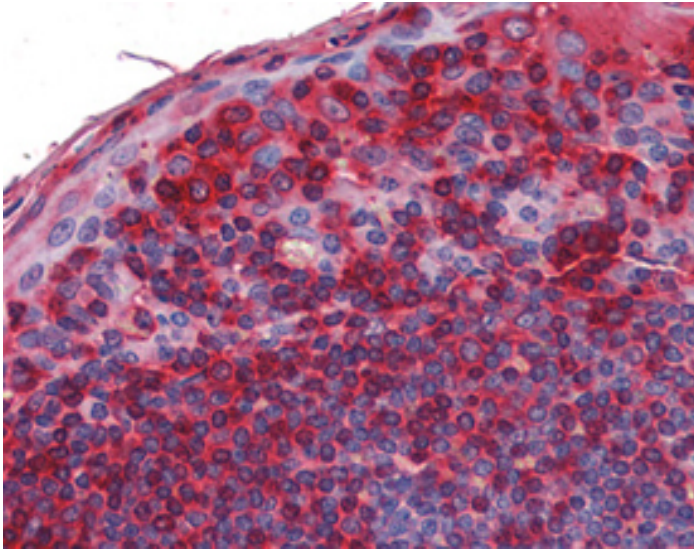
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

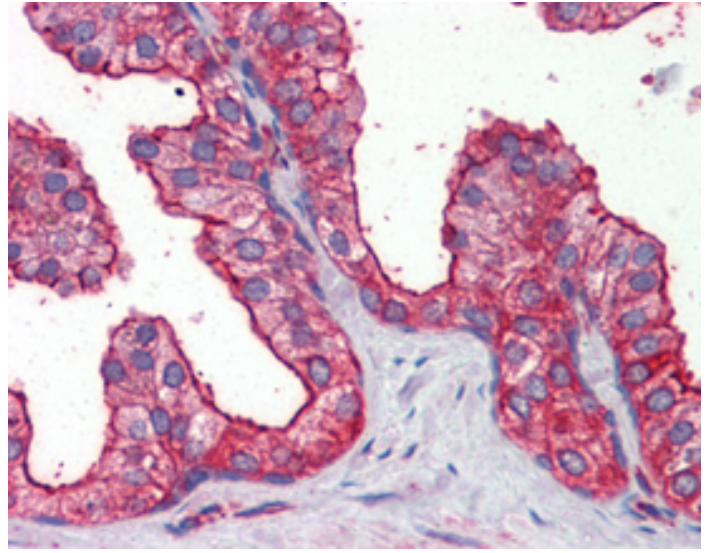
[View Online »](#)

Manuals, SDS & CofA

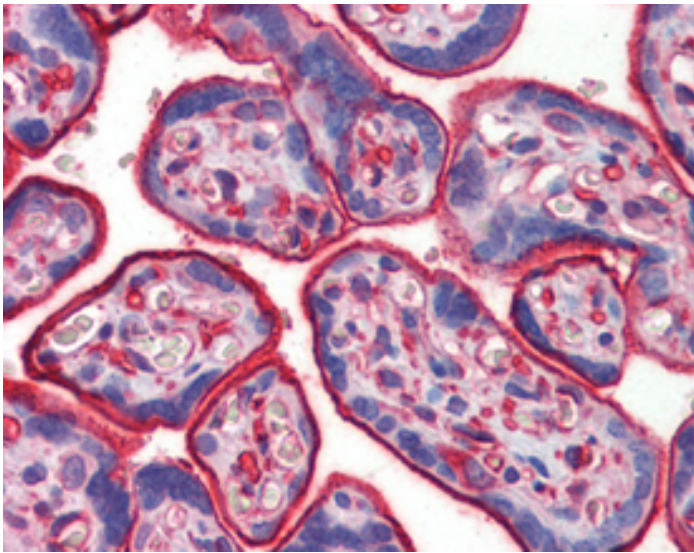
[View Online »](#)



IHC Analysis: Human tonsil tissue stained with TAK1, pAb at 10 $\mu$ g/ml.



IHC Analysis: Human prostate tissue stained with TAK1, pAb at 10 $\mu$ g/ml.



IHC Analysis: Human placenta tissue stained with TAK1, pAb at 10 $\mu$ g/ml.

## 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** TGF- $\beta$ -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 (1 $\mu$ g/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 [Worry-Free Guarantee](#)

Last modified: May 29, 2024



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)