

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 »](#)

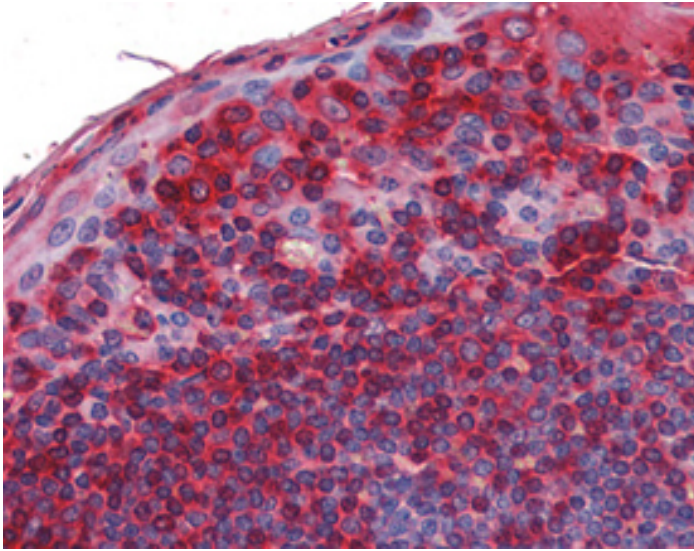
Ordering Information

[Order Online »](#)

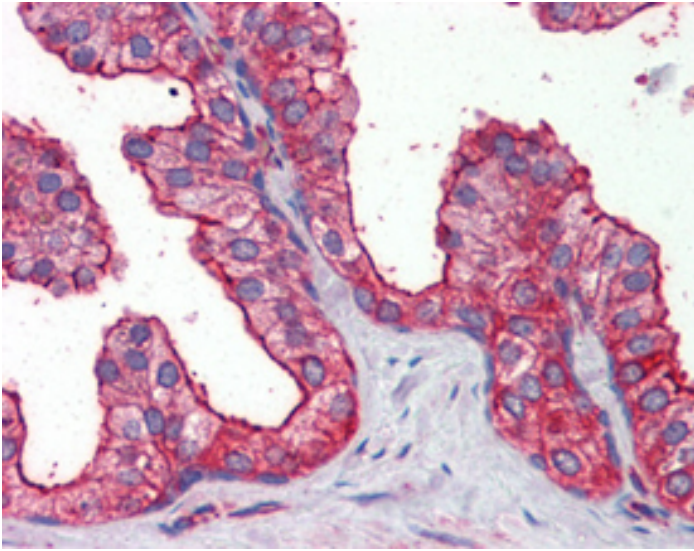
| | |
|-----------------|-------|
| ADI-KAS-ST009-E | 100µg |
|-----------------|-------|

Manuals, SDS & CofA

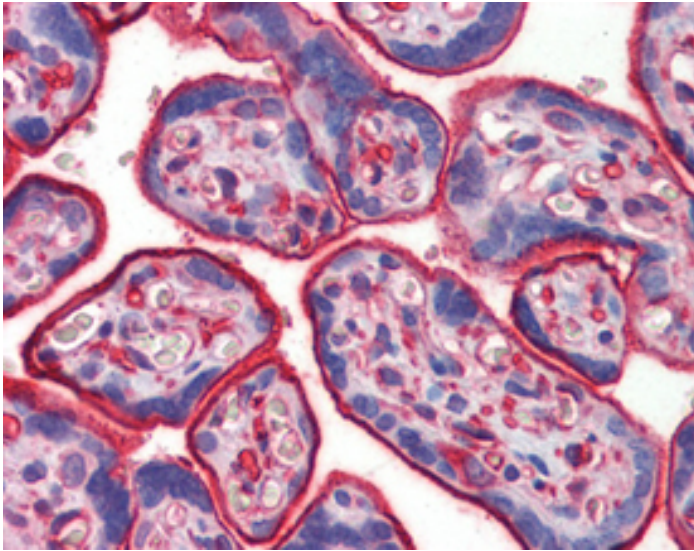
[View Online »](#)



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.

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-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 Worry-Free Guarantee . |



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)