

# TACI polyclonal antibody

TACI (transmembrane activator and CAML interactor), also known as TNF receptor superfamily member 13B, is a lymphocyte-specific member of the TNF-R family. TACI and its ligands (BAFF, APRIL) regulate humoral immune responses, activate NF- $\kappa$ B and c-jun N-terminal kinase, and play a role in B cell associated autoimmune diseases. Mutations in TACI are associated with a subset of common variable immunodeficiency (CVID).

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

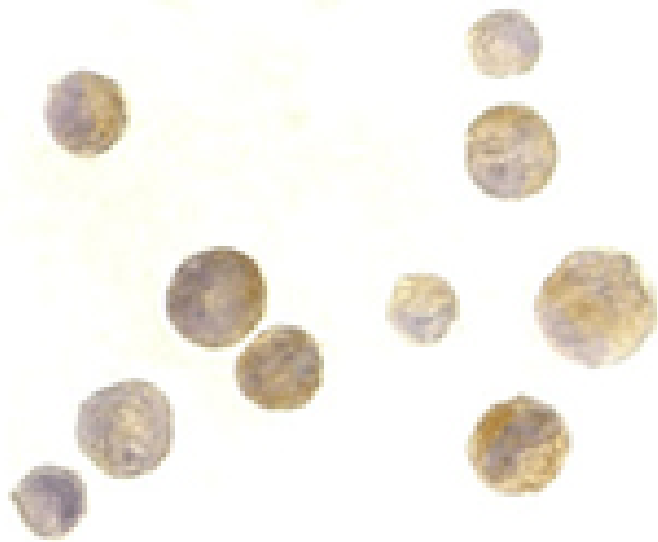
## Ordering Information

[Order Online »](#)

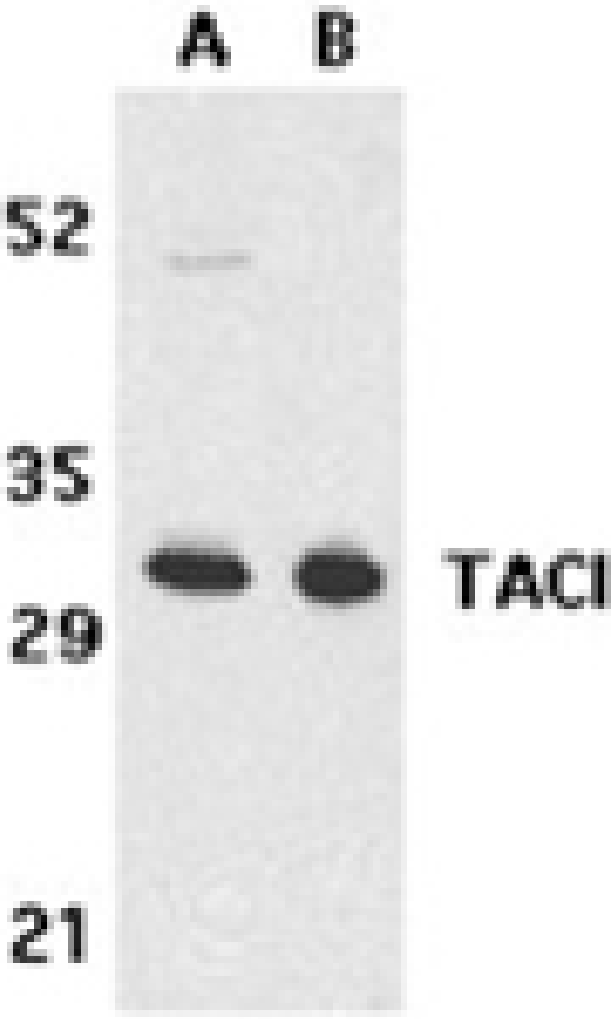
ADI-905-312-100	100 $\mu$ g
-----------------	-------------

## Manuals, SDS & CofA

[View Online »](#)



Immunocytochemistry analysis of TACI in K562 cells with TACI antibody at 2µg/ml.



Western blot analysis of TACI in K562 (A) and U937 (B) cell lysates with TACI antibody at 5µg/ml.

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** TNFRSF 13B, Transmembrane activator and CAML interactor, CD267

**Application** ICC, WB

**Application Notes** Detects a band of ~30kDa by Western blot.

**Formulation** Liquid. In PBS containing 0.02% sodium azide.

**Host** Rabbit

**Immunogen** Synthetic peptide corresponding to the sequence near the N-terminus of human TACI.

**Purity Detail** Peptide affinity purified.

**Recommendation** Immunocytochemistry (2µg/ml)Western Blot (2-5µg/ml)Suggested dilutions/conditions  
**Dilutions/Conditions** may not be available for all applications.Optimal conditions must be determined individually for each application.

**Species Reactivity** Human

**UniProt ID** O14836

**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)