

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- κ 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](#).

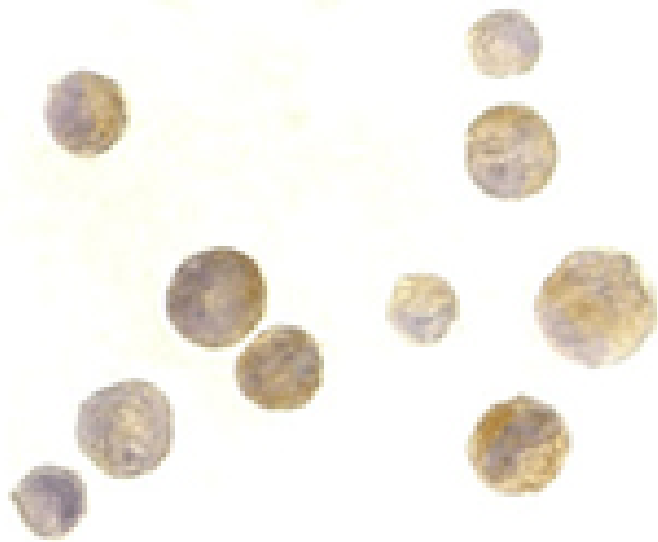
Ordering Information

[Order Online »](#)

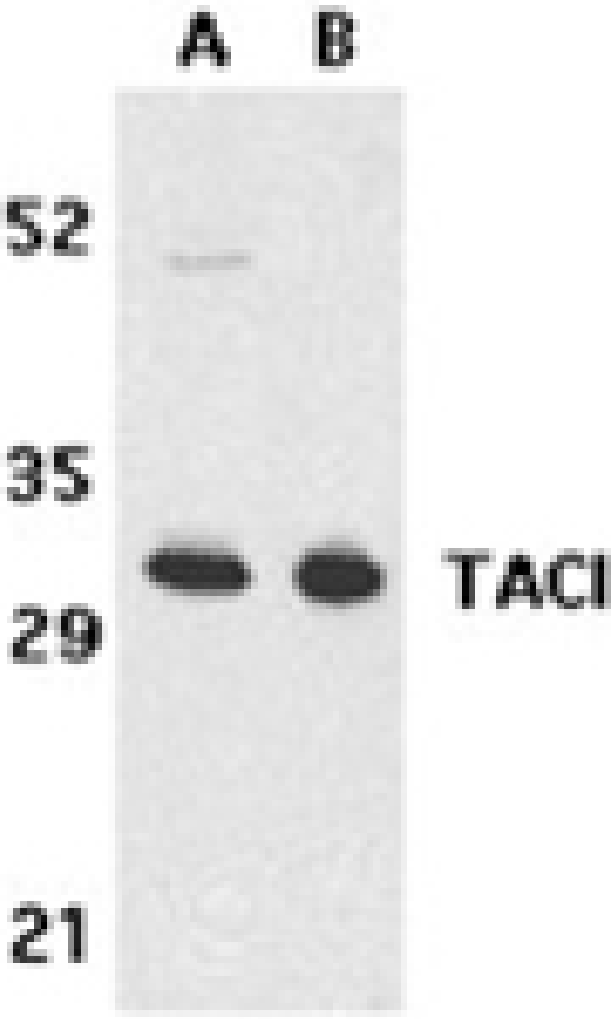
ADI-905-312-100	100 μ g
-----------------	-------------

Manuals, SDS & CofA

[View Online »](#)



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



Western blot analysis of TACI in K562 (A) and U937 (B) cell lysates with TACI antibody at 5 μ g/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 Dilutions/Conditions	Immunocytochemistry (2µg/ml)Western Blot (2-5µg/ml)Suggested 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

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-
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)