

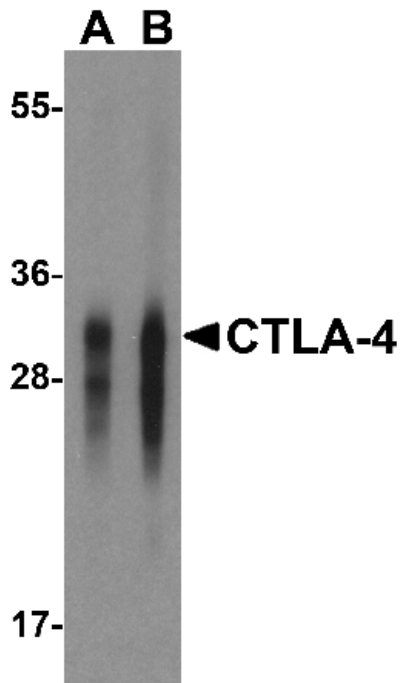
# CTLA4 monoclonal antibody (1E6)

The cytotoxic T-lymphocyte-associated protein 4 (CTLA-4), also known as CD152, is a member of the immunoglobulin superfamily that is expressed by activated T cells and transmits an inhibitory signal to T cells. CTLA-4 also functions as an immune checkpoint protein, and anti-CTLA-4 antibodies have been successfully used in the treatment of cancer. Mutations in the CTLA-4 gene have been implicated in multiple autoimmune diseases.

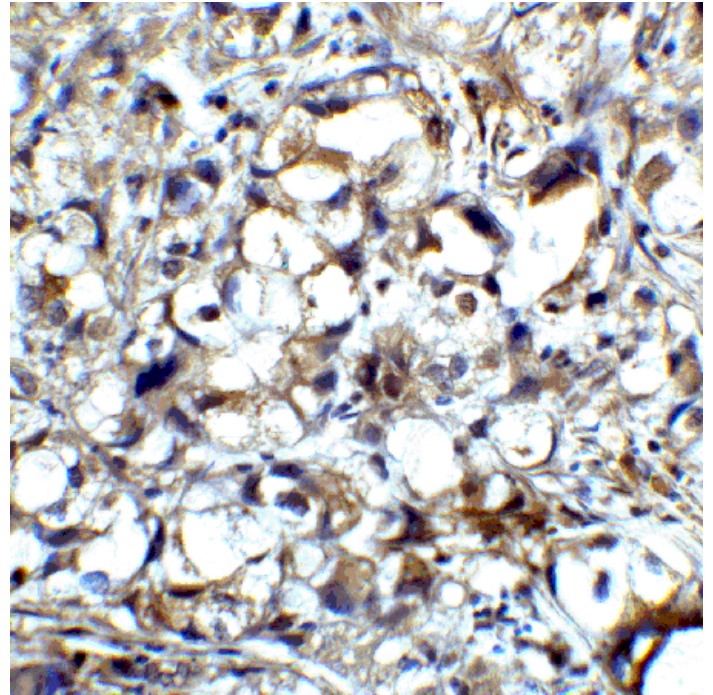
This antibody is covered by our [Worry-Free Guarantee](#).

**Manuals, SDS & CofA**

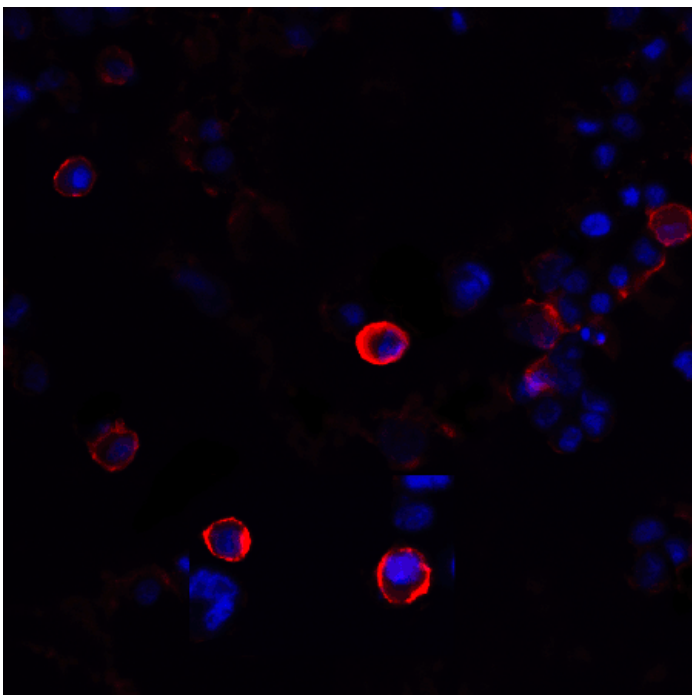
[View Online »](#)



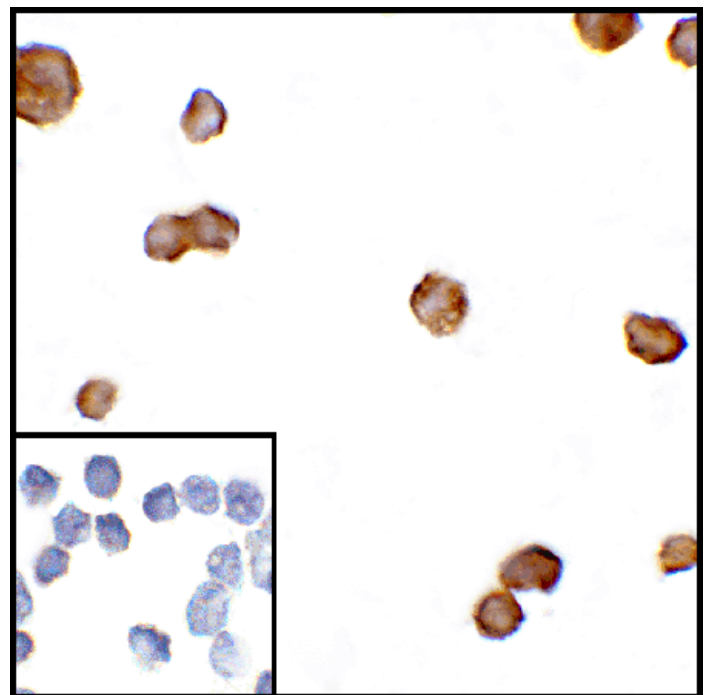
Western blot analysis of CTLA-4 in overexpressing HEK293 cells. CTLA-4 antibody at 0.125 and 0.25  $\mu\text{g}/\text{mL}$ .



Immunohistochemistry of CTLA-4 in human lymph node tissue with CTLA-4 antibody at 5  $\mu\text{g}/\text{mL}$ .



Immunofluorescence of CTLA-4 in transfected HEK293 cells with CTLA-4 antibody at 20  $\mu\text{g}/\text{mL}$ . Red: CTLA4 Antibody [1E6] (RF16011) Blue: DAPI staining



Immunocytochemistry of CTLA-4 in transfected HEK293 cells with CTLA-4 antibody at 5  $\mu\text{g}/\text{mL}$ . Lower left: Immunocytochemistry in transfected HEK293 cells with control mouse IgG antibody at 5  $\mu\text{g}/\text{mL}$ .

## Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored as recommended.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	Cytotoxic T-lymphocyte associated protein 4, CD152
Application	ICC, IF, IHC (PS), WB
Clone	1E6
Formulation	Liquid. In PBS containing 0.02% sodium azide and 50% glycerol.
Host	Mouse
Immunogen	Extracellular domain of human CTLA-4.
Isotype	IgG1
Purity Detail	Protein A affinity purified.
Species Reactivity	Human
UniProt ID	P16410
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>

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)