# 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.

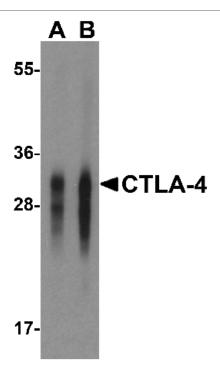
## **Ordering Information**

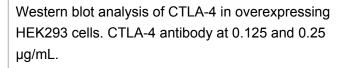
Order Online »

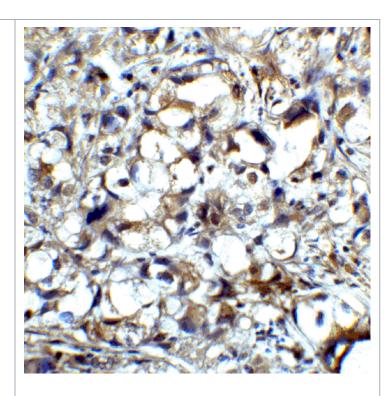
**ENZ-ABS696-0100** 100μg

Manuals, SDS & CofA

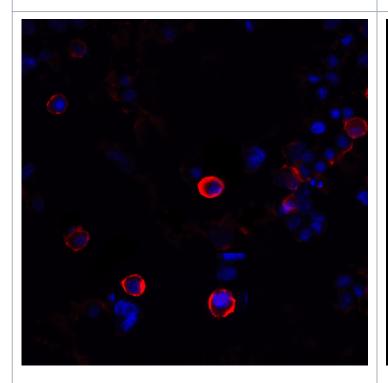
**View Online** »



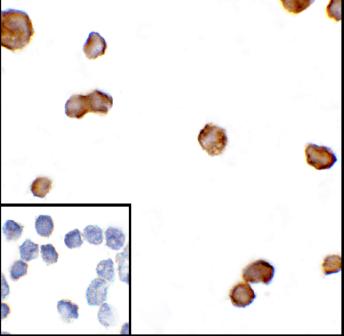




Immunohistochemistry of CTLA-4 in human lymph node tissue with CTLA-4 antibody at 5  $\mu$ g/mL.



Immunofluorescence of CTLA-4 in transfected HEK293 cells with CTLA-4 antibody at 20 µg/mL. Red: CTLA4 Antibody [1E6] (RF16011) Blue: DAPI staining



Immunocytochemistry of CTLA-4 in transfected HEK293 cells with CTLA-4 antibody at 5 ug/mL. Lower left: Immunocytochemistry in transfected HEK293 cells with control mouse IgG antibody at 5 ug/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.

lsotype lgG1

Purity Detail Protein A affinity purified.

Species Reactivity Human

UniProt ID P16410

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•

