# Fas polyclonal antibody

Fas/APO-1/CD95 is a 36 kDa cell surface type I-membrane glycoprotein with an apparent molecular weight of 44 kDa on SDS-PAGE. Fas belongs to the tumor necrosis factor receptor (TNFR) family, which includes TNFR1, TNFR2, CD27, CD30 and CD40. Binding of Fas-ligand (FasL) to Fas or crosslinking of Fas by anti-Fas monoclonal antibodies leads to apoptosis of Fas expressing cells. Both Fas and TNFR contain a 70 amino acid cytoplasmic "death domain" responsible for transmitting signals for apoptosis. Apoptosis induced by Fas-Fas ligand interaction is inhibited by a number of cytosolic factors such as the ICE/CED3 family inhibitor CrmA, and the short and long forms of FLIP. Insertion of an early transposable element in intron 2 of the Fas gene, as in the mouse lpr mutations, causes lymphoproliferation and autoimmunity in mutant mice, suggesting the involvement of Fas in clonal deletion of autoreactive T cells.

This antibody is covered by our Worry-Free Guarantee.

Citations: 2

View Online »

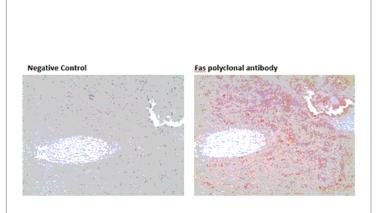
# **Ordering Information**

**Order Online** »

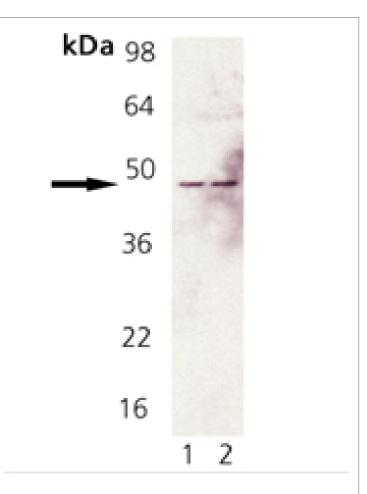
ADI-AAP-221-D	50µg
ADI-AAP-221-F	200μg

Manuals, SDS & CofA

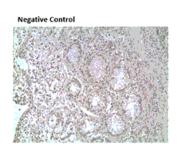
**View Online »** 

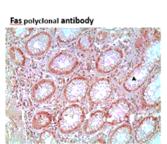


Immunohistochemistry analysis of formalin-fixed, paraffin-embedded human spleen tissue stained with Fas polyclonal antibody, (Prod. No. ADI-AAP-221) at  $10\mu g/ml$ .



Western Blot Analysis of Fas pAb : Lane 1: Rat Brain Tissue Extract, Lane 2: Mouse Brain Tissue Extract.





Immunohistochemistry analysis of formalin-fixed, paraffin-embedded human colon tumor tissue stained with Fas polyclonal antibody, (Prod. No. ADI-AAP-221) 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 CD95, APO-1, TNFRSF 6

**Application** IHC, WB

**Application Notes**Detects a band of ~45kDa by Western blot.

**Formulation** Liquid. In PBS, pH 7.2, containing 0.09% sodium azide.

GenBank ID M83649

**Host** Rabbit

**Immunogen** Synthetic peptide corresponding to a portion of mouse

Fas.

Purity Detail Peptide affinity purified.

**Recommendation Dilutions/Conditions** Immunohistochemistry (1:50)Western Blot (1:1,000,

ECL)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 Human, Mouse, Rat

UniProt ID P25446

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

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com