

# 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 »](#)

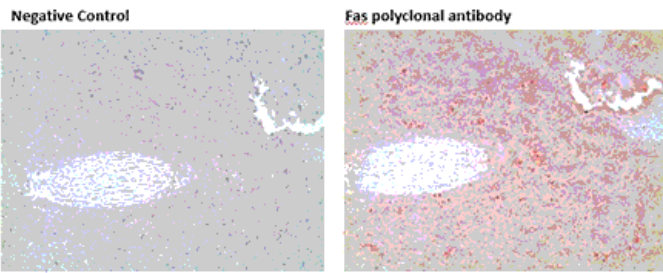
Ordering Information

[Order Online »](#)

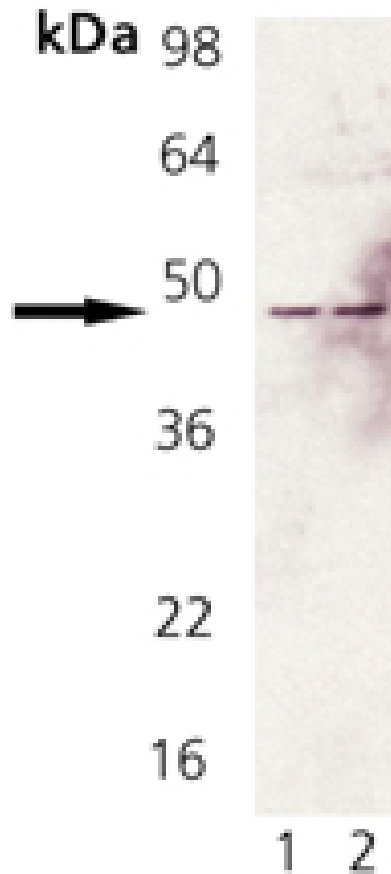
ADI-AAP-221-F	200µg
---------------	-------

Manuals, SDS & CofA

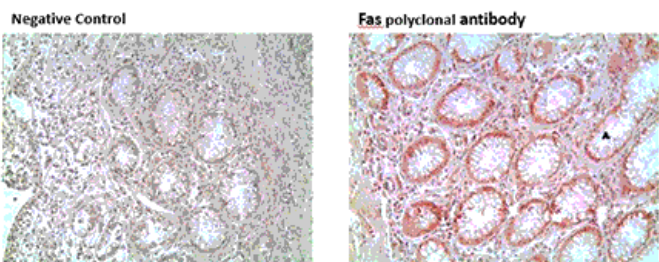
[View Online »](#)



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



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)