

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

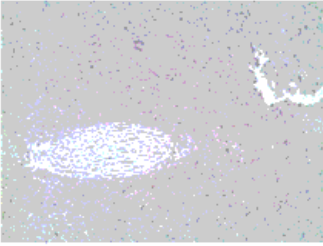
[Order Online »](#)

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

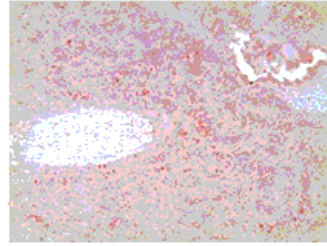
Manuals, SDS & CofA

[View Online »](#)

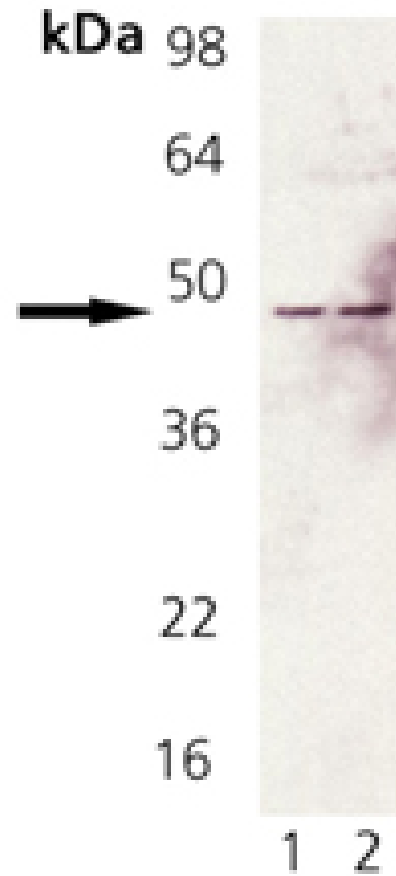
Negative Control



Fas polyclonal antibody

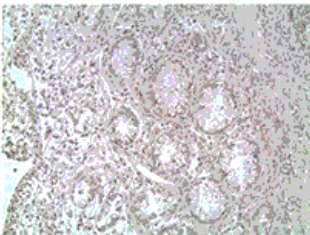


Immunohistochemistry analysis of formalin-fixed, paraffin-embedded human spleen tissue stained with Fas polyclonal antibody, (Prod. No. ADI-AAP-221) at 10µg/ml.

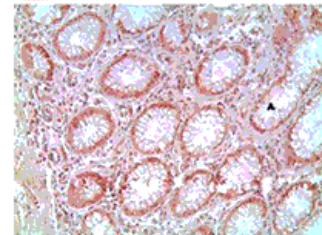


Western Blot Analysis of Fas pAb : Lane 1: Rat Brain Tissue Extract, Lane 2: Mouse Brain Tissue Extract.

Negative Control



Fas polyclonal antibody



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 Immunohistochemistry (1:50)Western Blot (1:1,000, ECL)Suggested dilutions/conditions
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](#).



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)