

p53 monoclonal antibody (PAb122)

p53 is a tumor suppressor gene expressed in a wide variety of tissue types and is involved in regulating cell growth, replication, and apoptosis. It binds to Mdm2, SV40 T antigen and human Papilloma virus E6 protein. p53 senses DNA damage and facilitates its repair. Mutation involving p53 is found in a wide variety of malignant tumors, including breast, ovary, bladder, colon, lung, and melanoma.

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

Citations: 4

[View Online »](#)

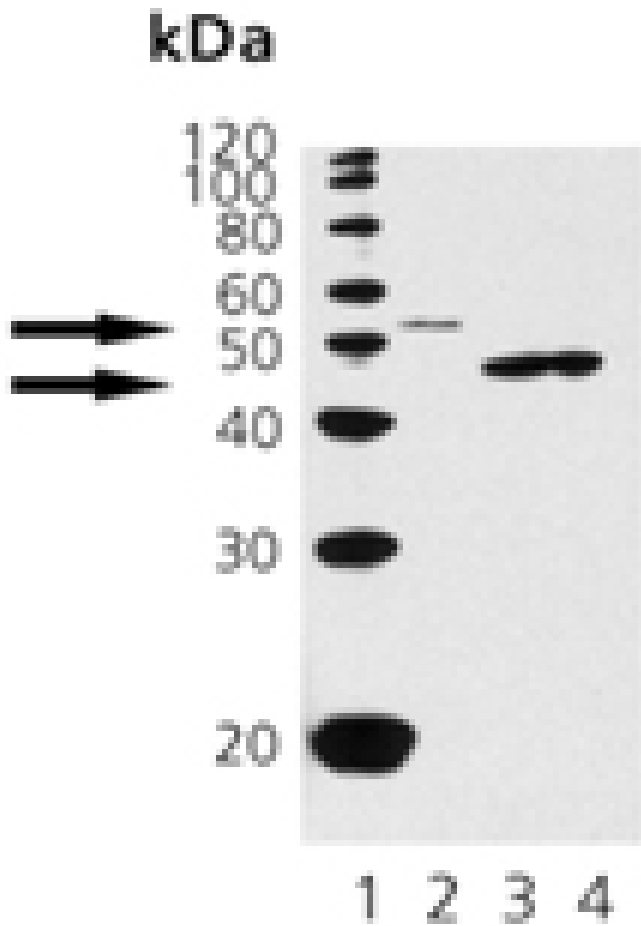
Ordering Information

[Order Online »](#)

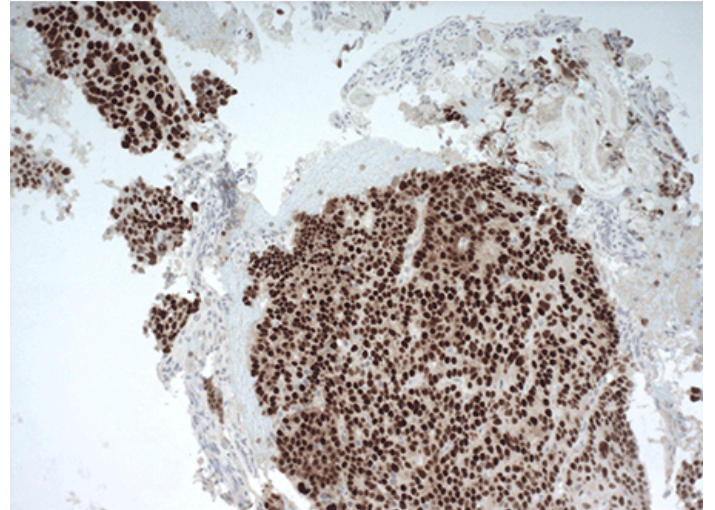
ADI-KAM-CC002-D	50µg
ADI-KAM-CC002-F	200µg

Manuals, SDS & CofA

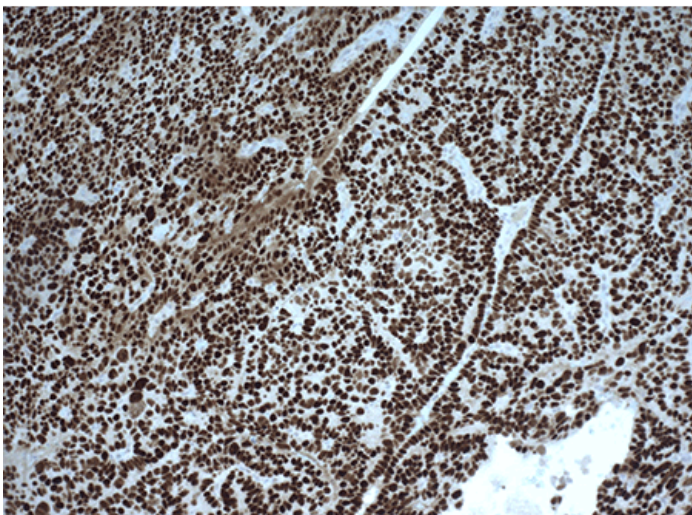
[View Online »](#)



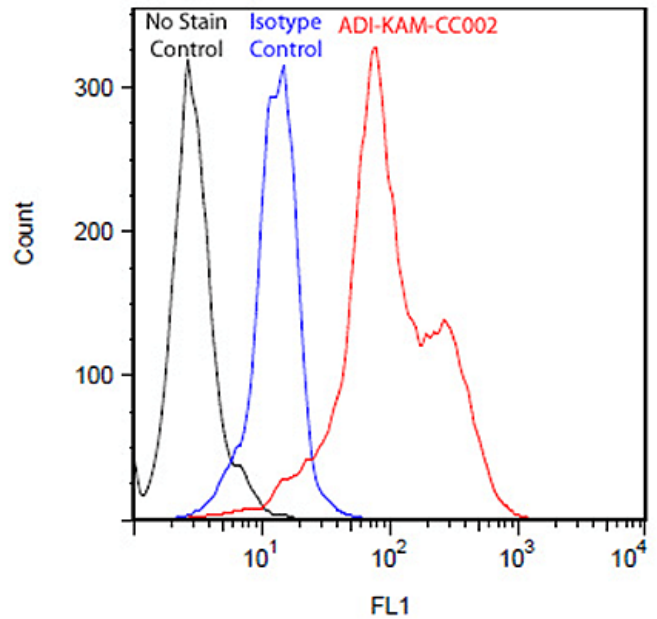
Western blot analysis of p53: Lane 1: MW marker, Lane 2: A431, Lane 3: rat brain, Lane 4: mouse brain.



Endometrial carcinoma stained by immunohistochemistry with p53 monoclonal antibody (PAb122) (ENZ-ABS849, 1:20 Dilution).



Endometrial carcinoma stained by immunohistochemistry with p53 monoclonal antibody (PAb122) (ENZ-ABS849, 1:20 Dilution).



Flow cytometry analysis of 10^6 Jurkat cells stained using p53 monoclonal antibody (PAb122), (Prod. No. ENZ-ABS849), at a concentration of 50 μ g/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 TRP53, Tumor protein p53

Application Flow Cytometry, ICC, IHC, IP, WB

Application Notes Detects a band of ~48/53kDa by Western blot.

Clone pAb122

Crossreactivity Weakly cross-reacts with dog p53.

Formulation Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.

Host Mouse

Immunogen SV40 transformed 3T3 cells.

Isotype IgG2b

Purity Detail Protein G affinity purified.

Recommendation Dilutions/Conditions Western Blot (1:250 colorimetric, 1:250 ECL) Immunohistochemistry (1:20 – 1:100) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Human, Mouse, Rat

UniProt ID

P02340

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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com