

# 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 »](#)

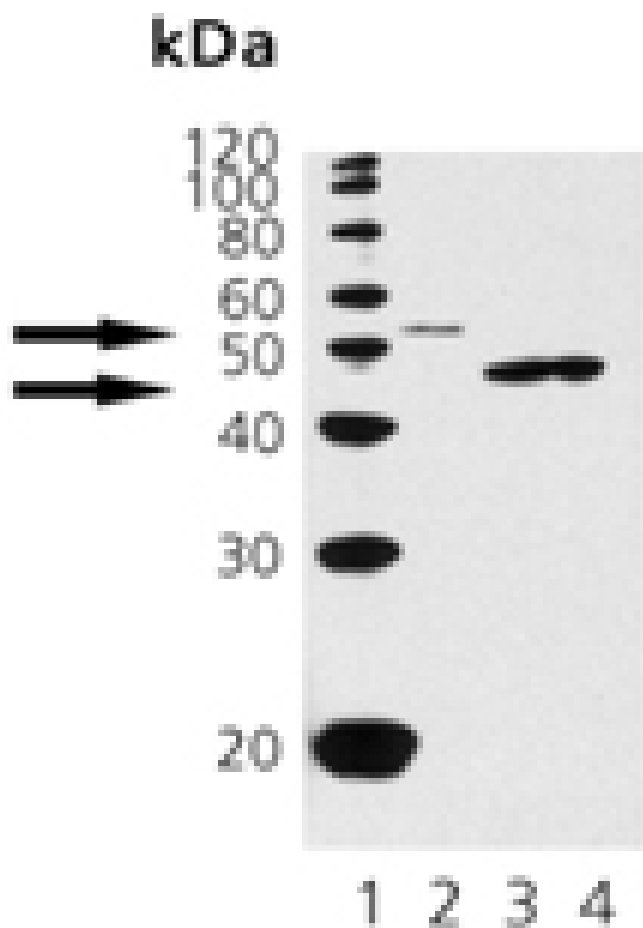
## Ordering Information

[Order Online »](#)

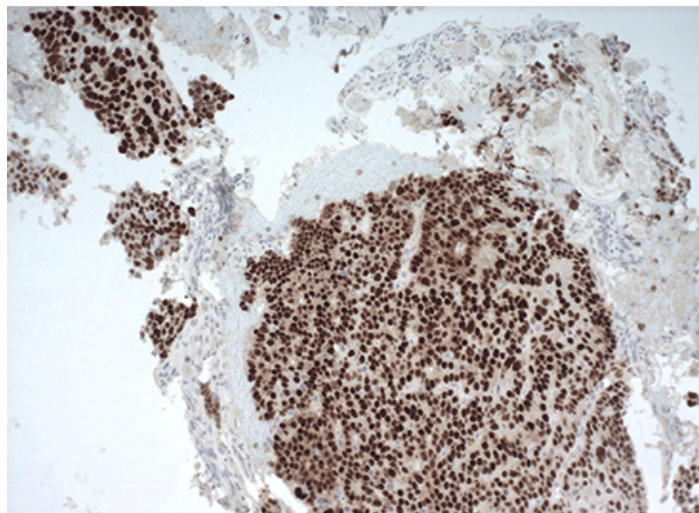
|                 |       |
|-----------------|-------|
| ADI-KAM-CC002-J | 1mg   |
| ADI-KAM-CC002-D | 50µg  |
| ADI-KAM-CC002-F | 200µg |

## Manuals, SDS & CofA

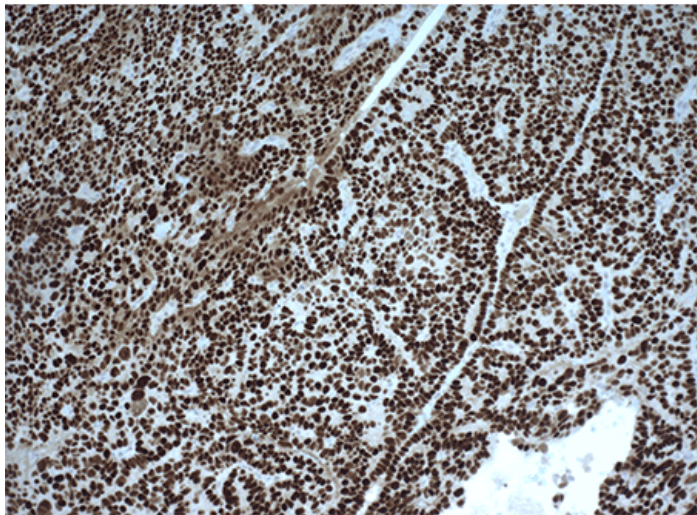
[View Online »](#)



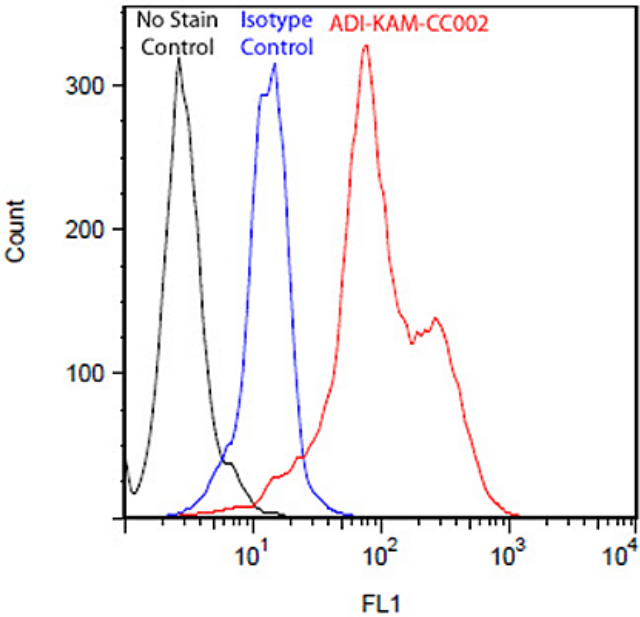
Western blot analysis of p53: Lane 1: MW marker, Lane 2: A431, Lane 3: rat brain, Lane 4: mouse brain.



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  $\mu\text{g/mL}$ .

## 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** Western Blot (1:250 colorimetric, 1:250 ECL) Immunohistochemistry (1:20 –  
**Dilutions/Conditions** 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



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-  
fr@enzolifesciences.com](mailto:info-fr@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)

**Worry-free Guarantee**

**This antibody is covered by our Worry-Free Guarantee.**

For local distributors and detailed product information visit us online: [enzo.com](http://enzo.com) All product names, logos, and brands are the property of their respective owners.