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 »

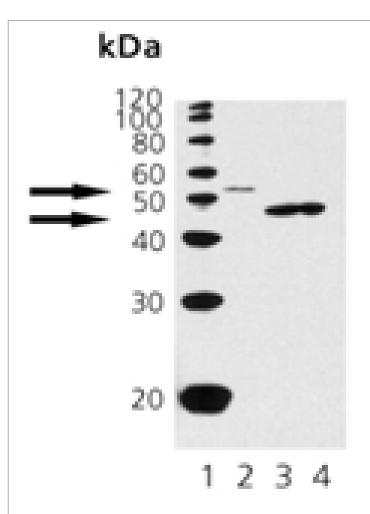
Ordering Information

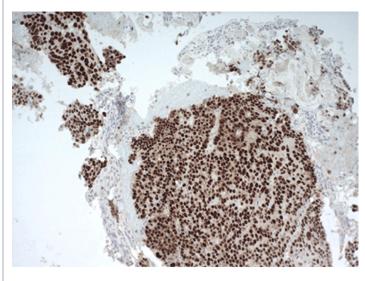
Order Online »

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

Manuals, SDS & CofA

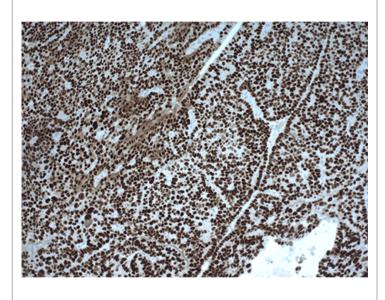
View Online »



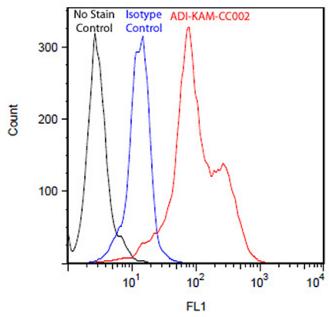


Endometrial carcinoma stained by immunohistochemistry with p53 monoclonal antibody (PAb122) (ENZ-ABS849, 1:20 Dilution).

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).



Flow cytometry analysis of 10⁶ Jurkat cells stained using p53 monoclonal antibody (PAb122), (Prod. No. ENZ-ABS849), at a concentration of 50 µg/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.

lgG2b

Purity Detail Protein G affinity purified.

Recommendation Dilutions/ConditionsWestern 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

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024

