

# p53 monoclonal antibody (PAb421)

PAb421 antibody induces allosteric changes in certain tumor-derived p53 mutants and restores tumor suppressor activity. Microinjection of PAb421 into SW480 colorectal carcinoma cells restored transcriptional activation function to the resident mutant p53 (Arg to His 273, Pro to Ser 309). The epitope for PAb421 is contained within the "p53 activator peptide" 46. Thus the antibody recognizes this peptide and has been used to examine cellular uptake of the p53 activator peptide.

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

Citations: 8

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Handling	Avoid freeze/thaw cycles.
Long Term Storage	+4°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	TRP53, Tumor protein p53
Application	ICC, IHC, IP, WB
Clone	PAb421
Formulation	Liquid. In 50 mM sodium phosphate buffer, pH 7.5, containing 0.2% gelatin and 0.09% sodium azide.
Host	Mouse
Immunogen	Synthetic peptide corresponding to aa 371-380 of human p53.
Isotype	IgG2a
Purity Detail	Purified.
Recommendation Dilutions/Conditions	Western Blot (10µg/ml) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse, Rat
UniProt ID	P04637

## 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)

Last modified: May 29, 2024