

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 »](#)

Ordering Information

[Order Online »](#)

BML-SA293-0050	50µg
----------------	------

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

Last modified: May 29, 2024