

# [pSer315]p53 monoclonal antibody (FPS315)

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](#).

## Ordering Information

[Order Online »](#)

ADI-KAM-CC239-E	100µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage      -20°C

Shipping      Blue Ice

**Regulatory Status**      RUO - Research Use Only

## Product Details

Alternative Name      TRP53, Tumor protein p53

Application      WB

Application Notes      Western Blot (5µg/ml, ECL)  
Optimal conditions must be determined individually for each application.

Clone      FPS315

Formulation      Liquid. In PBS, pH 7.2, containing 50% glycerol.

Host      Mouse

Immunogen      Synthetic peptide corresponding to a portion of human p53 phosphorylated at Ser315.

Isotype      IgG1

Purity Detail      Protein A affinity purified.

Recommendation      Western Blot (5µg/ml, ECL) Suggested dilutions/conditions may not be available for all  
Dilutions/Conditions      applications. Optimal conditions must be determined individually for each application.

Species Reactivity      Human

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)