[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.

lsotype lgG1

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions Western Blot (5µg/ml, ECL)Suggested dilutions/conditions

may not be available for all 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

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com