Ubiquitin polyclonal antibody

Ubiquitin (Ub) plays a very important role in regulated non-lysosomal ATP dependent protein degradation. The protein to be degraded is conjugated to Ub and the ubiquinated protein is then selectively degraded by the 26S complex, a multicatalytic cytosolic and nuclear protease. The Ubproteasome proteolytic pathway, which is a complex process, is implicated to be of great importance for regulating numerous cellular processes.

This antibody is covered by our Worry-Free Guarantee.

Citations: 36

View Online »

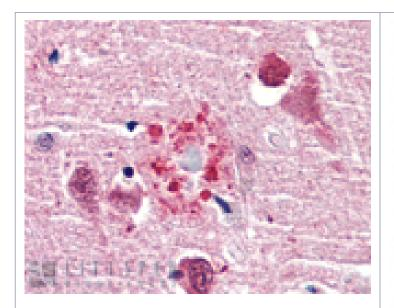
Ordering Information

Order Online »

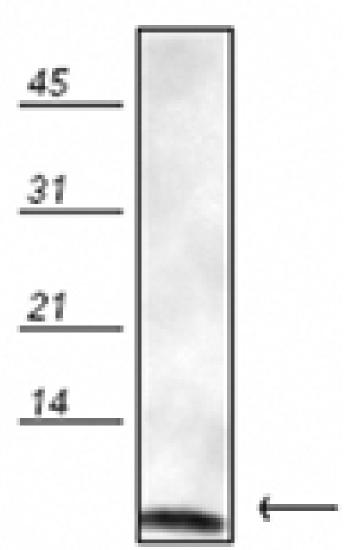
ADI-SPA-200-D	50µg
ADI-SPA-200-F	200μg

Manuals, SDS & CofA

View Online »



Immunohistochemistry analysis of human brain, cortex, senile plaque tissue stained with Ubiquitin, pAb at 10µg/ml.



Western blot analysis: Ubiquitin protein probed with Ubiquitin pAb.

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ICC, IHC (PS), WB

Application NotesDetects a band of ~8kDa by Western blot.

Formulation Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09%

sodium azide.

Gene/Protein Identifier NP_776558 (RefSeq)

Host Rabbit

Immunogen Native bovine Ubiquitin.

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions Western Blot (1:5,000, ECL)Suggested dilutions/conditions

may not be available for all applications. Optimal conditions

must be determined individually for each application.

Source Purified from rabbit serum.

Species Reactivity Bovine, Chicken, Dog, Drosophila, Fish, Guinea pig,

Hamster, Human, Monkey, Mouse, Porcine, Rabbit, Rat,

Sheep, Xenopus, Yeast

UniProt ID P0CG53

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024



info-

eu@enzolifesciences.com