

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 ubiquitinated protein is then selectively degraded by the 26S complex, a multicatalytic cytosolic and nuclear protease. The Ub-proteasome 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: 37

[View Online »](#)

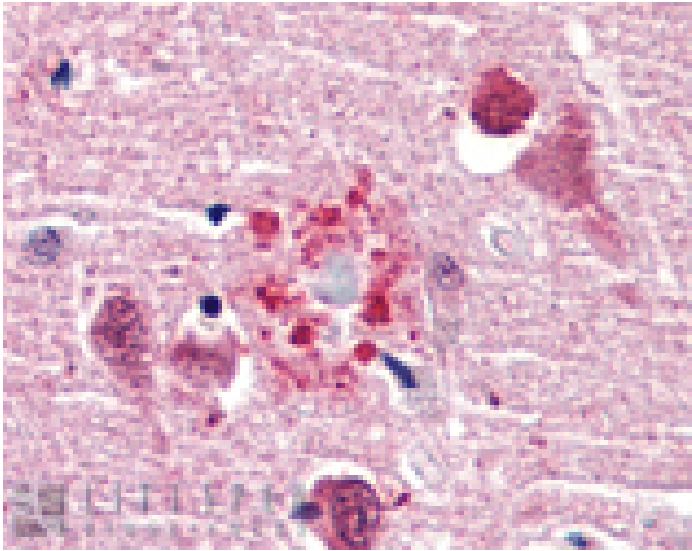
Ordering Information

[Order Online »](#)

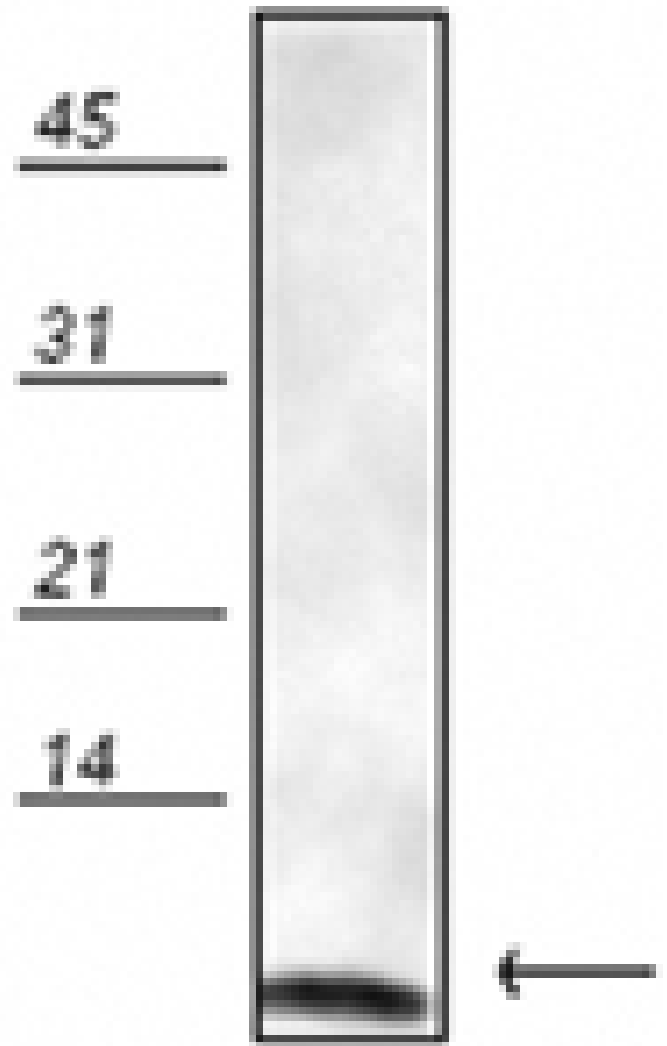
ADI-SPA-200-D	50µg
ADI-SPA-200-F	200µg

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 Notes	Detects 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 .



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)