

Ubiquitin monoclonal antibody (P4G7-H11)

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 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: 14

[View Online »](#)

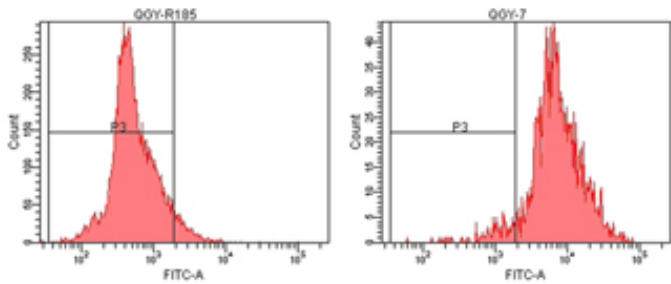
Ordering Information

[Order Online »](#)

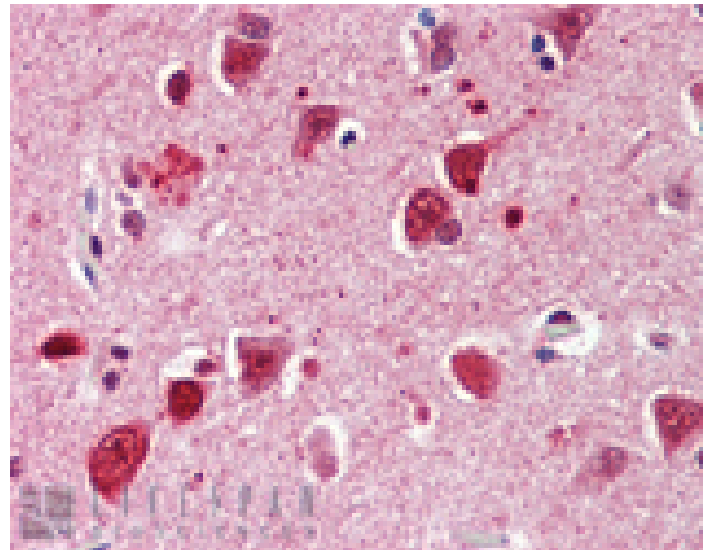
ADI-SPA-203-D	50µg
ADI-SPA-203-F	200µg

Manuals, SDS & CofA

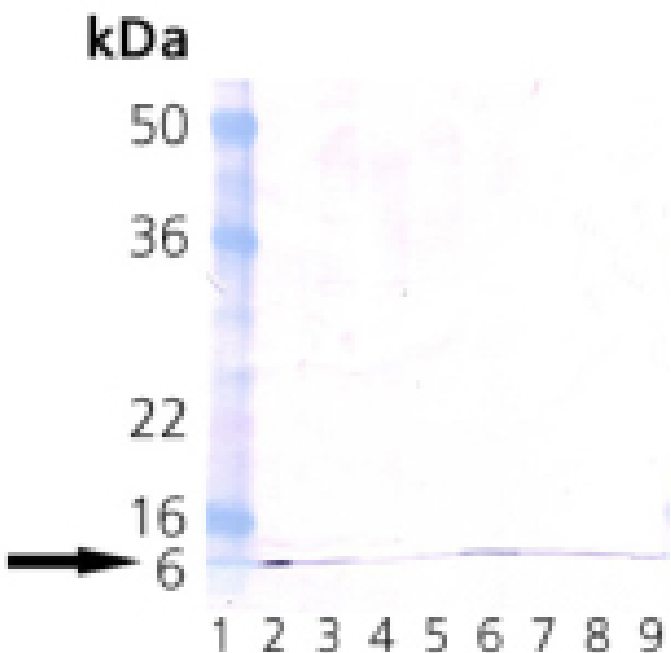
[View Online »](#)



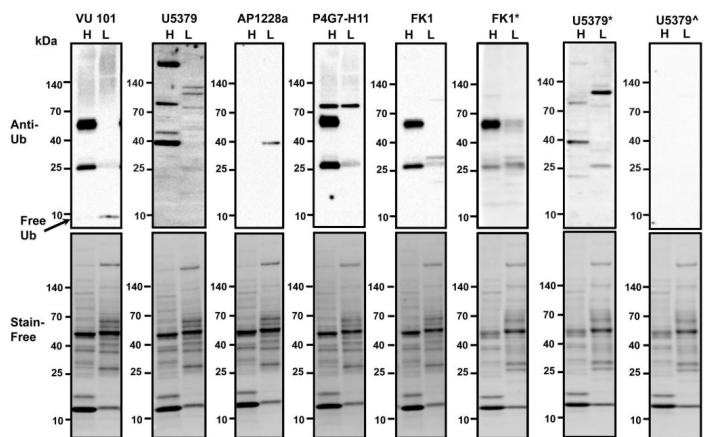
Flow cytometry analysis of human hepatoma QGY cells analyzed by flow cytometry using isotype control antibody (left) or Ubiquitin Monoclonal Antibody (P4G7-H11) (right) at 10 μ g/mL



Immunohistochemistry analysis of human brain, cortex tissue stained with Ubiquitin, mAb (P4G7-H11) at 5-10 μ g/ml.



Western blot analysis of Ubiquitin: Lane 1: MW marker, Lane 2: Ubiquitin, Lane 3: HeLa, Lane 4: RK-13, Lane 5: 3T3, Lane 6: PC-12, Lane 7: Mouse Brain, Lane 8: Rat Brain, Lane 9: MDBK



Comparison of anti-ubiquitin antibodies. Heart and liver lysates (20 μ g each) were investigated by Western blotting using five commercially available anti-ubiquitin antibodies (VU101, U5379, AP1228a, P4G7-H11, FK1). Arrow shows location of free unbound ubiquitin. Stain-free staining of total proteins loaded was used as the normalization control. H, heart; L, liver. BSA was used as the blocking reagent for the blot labeled FK1* while non-fat milk was used as the blocking reagent in all the other blots shown. All antibodies were used at a dilution of 1:1000 except for blots labeled U5379* and U5379[^] which were used at dilutions of 1:100 and 1:2000 respectively.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Western Blotting Inaccuracies with Unverified Antibodies: Need for a Western Blotting Minimal Reporting Standard (WBMRS). *PLoS One* (2015)

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Flow Cytometry, IHC (PS), WB

Application Notes Detects a band of ~6.5kDa by Western blot.

Clone P4G7-H11

Formulation Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.

Gene/Protein Identifier NP_776558 (RefSeq)

Host Mouse

Immunogen Native bovine Ubiquitin.

Isotype IgG1

Purity Detail Protein G affinity purified.

Recommendation Dilutions/Conditions Flow Cytometry (10µg/ml)Western Blot (1:1,000, colorimetric)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Source Purified from ascites.

Species Reactivity Bovine, Hamster, Human, Mouse, Rabbit, Rat

UniProt ID P0CG53

Last modified: May 29, 2024



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)