

# 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 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: 13

[View Online »](#)

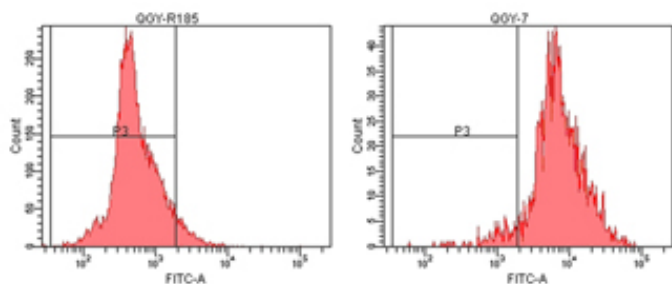
## Ordering Information

[Order Online »](#)

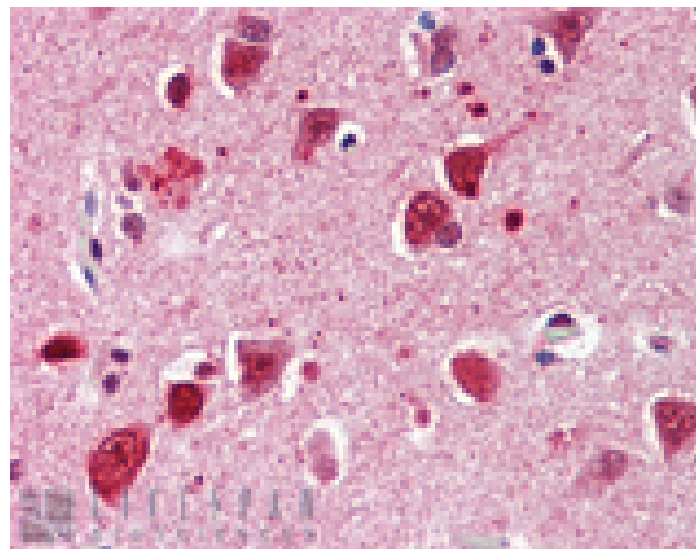
ADI-SPA-203-D	50µg
ADI-SPA-203-F	200µg

## Manuals, SDS & CofA

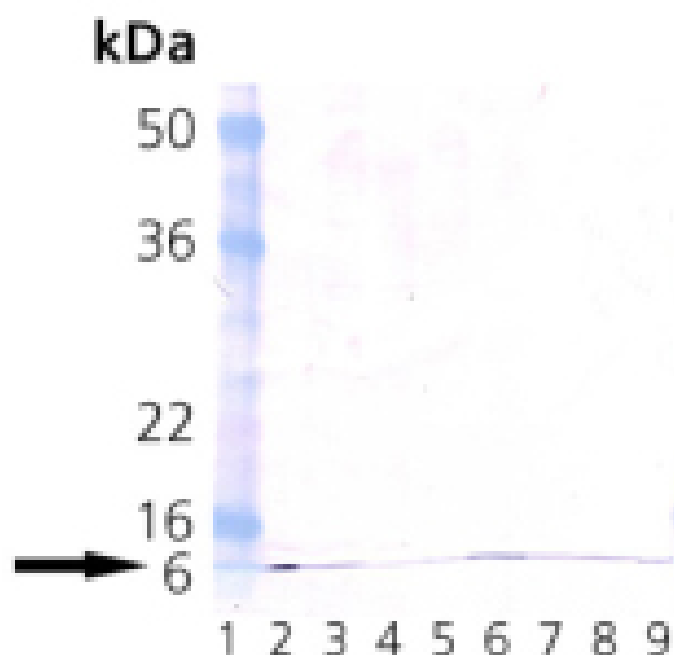
[View Online »](#)



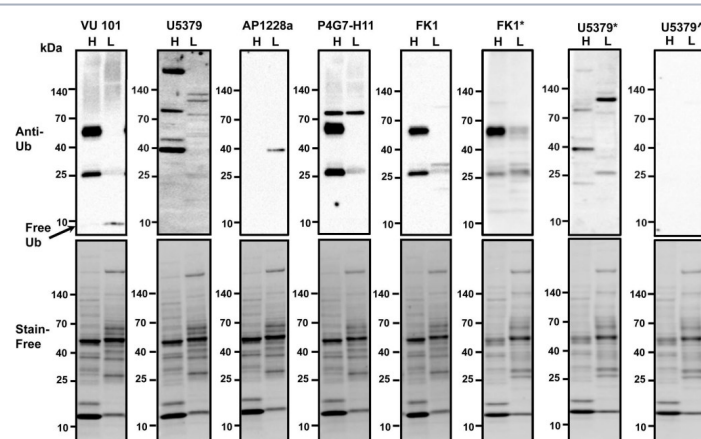
Flow cytometry analysis of human hepatoma QGY cells analyzed by flow cytometry using isotype control antibody (left) or Ubiquitin Monoclonal Antibody (P4G7-H11) (right) at 10 $\mu$ g/mL



Immunohistochemistry analysis of human brain, cortex tissue stained with Ubiquitin, mAb (P4G7-H11) at 5-10 $\mu$ 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  $\mu$ 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<sup>A</sup> 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



**Worry Free Guarantee** This antibody is covered by our **Worry Free Guarantee**

ENZO LIFE SCIENCES, European Sales Office Belgium, The Netherlands France Germany  
Phone: 800.942.0430 (ELS) AG Phone: +32 3 466 0420 info- info-  
usa@enzolifesciences.com info- fr@enzolifesciences.com de@enzolifesciences.com  
eu@enzolifesciences.com be@enzolifesciences.com

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
info-  
uk@enzolifesciences.com