# 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 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: 13

View Online »

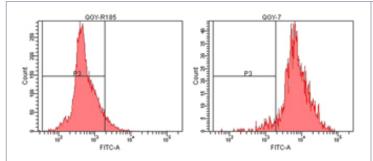
# **Ordering Information**

**Order Online** »

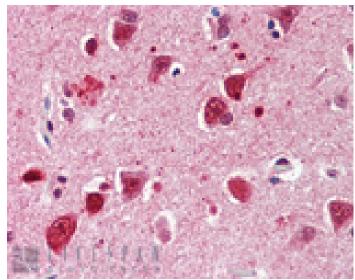
ADI-SPA-203-D	50µg
ADI-SPA-203-F	200μg

Manuals, SDS & CofA

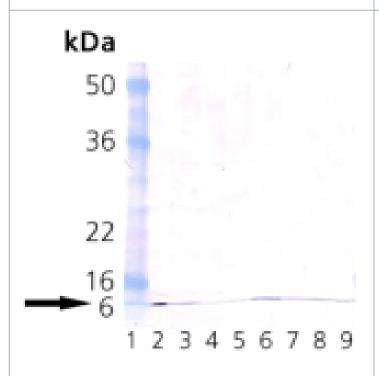
View Online »



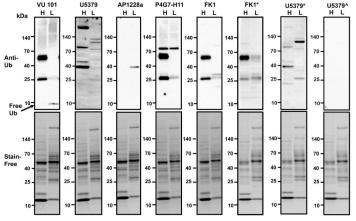
Flow cytometry analysis of human hepatoma QGY cells analyzed by flow cytometry using isotype control antibody (left) or Ubiquitin Monoclonal Antibody (P4G7-H11) (right) at 10µg/mL



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 µg 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. Stainfree 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.

lsotype lgG1

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



info-

eu@enzolifesciences.com