# Ubiquitin monoclonal antibody (P4D1) (HRP conjugate)

Ubiquitin modification of cellular proteins has been shown to be involved with a wide range of biochemical processes including proteasomal degradation, signal transduction, DNA repair, endocytosis and autophagy.

This antibody is covered by our Worry-Free Guarantee.

Citations: 9

View Online »

# **Ordering Information**

**Order Online** »

BML-PW0935-0025	25µl
BML-PW0935-0100	100µl

Manuals, SDS & CofA

View Online »

## **Handling & Storage**

**Handling** Avoid freeze/thaw cycles.

Short Term Storage +4°C

Long Term Storage +4°C

Shipping Blue Ice

### Regulatory Status RUO - Research Use Only

### **Product Details**

**Application** WB

Clone P4D1

**Formulation** Liquid. In PBS containing 0.01% sodium azide.

**Host** Mouse

**Immunogen** Denatured bovine ubiquitin.

lsotype lgG1

Purity Detail Ion exchange purified.

**Recommendation Dilutions/Conditions**Western Blot (1:1,000)Suggested dilutions/conditions may

not be available for all applications. Optimal conditions must be determined individually for each application.

Source Purified from tissue culture supernatant. HRP conjugated

chemically in vitro.

Species Reactivity Species independent

**Specificity** Recognizes mono- and poly-ubiquitin protein conjugates,

free polyubiquitin chains and free ubiquitin.

### Technical Info / Product Notes

P4D1 antibody was developed using a denatured glutaraldehyde cross-linked ubiquitin IgG complex as the immunogen. Monoclonal antibody was purified from tissue culture supernatant by ion exchange chromatography. The general ubiquitin reactivity of the monoclonal antibody has been demonstrated by Western blotting using synthetic ubiquitin chains, single isopeptide-linked polyubiquitin chains, and with endogenously ubiquitinylated species in cell lysates.

**UniProt ID** 

P0CG53 (UBB), P0CH28 (UBC), P62992 (RPS27A), P63048 (UBA52)

**Worry-free Guarantee** 

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-

uk@enzolifesciences.com