

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.

Isotype IgG1

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 ubiquitinated 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](#).



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)