

Polyubiquitin (K63-linkage-specific) monoclonal antibody (HWA4C4)

Modification of proteins by addition of K⁶³-linked polyubiquitin chains is implicated in a variety of cellular events, including DNA repair, signal transduction and receptor endocytosis.

A monoclonal antibody (mAb) has been generated that specifically recognizes K63-linked polyubiquitin chains, but not any other isopeptide-linked (K⁶, K¹¹, K²⁷, K²⁹, K³³, or K⁴⁸) polyubiquitin chain or monoubiquitin. The specificity of this K⁶³-linked ubiquitin-specific monoclonal antibody was demonstrated using synthetic ubiquitin chains, single isopeptide-linked polyubiquitin chains, and with endogenously ubiquitinated species in both cell lysates and intact cells.

The hybridoma secreting the antibody [clone HWA4C4] was generated by fusion of splenocytes from Balb/c mice which had received repeated immunisation with a synthetic peptide encompassing the ubiquitin K⁶³-isopeptide motif. Immunoglobulin was purified from tissue culture supernatant by ion exchange chromatography.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 44

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-PW0600-0025	25µl
BML-PW0600-0100	100µl

Manuals, SDS & CofA

[View Online »](#)

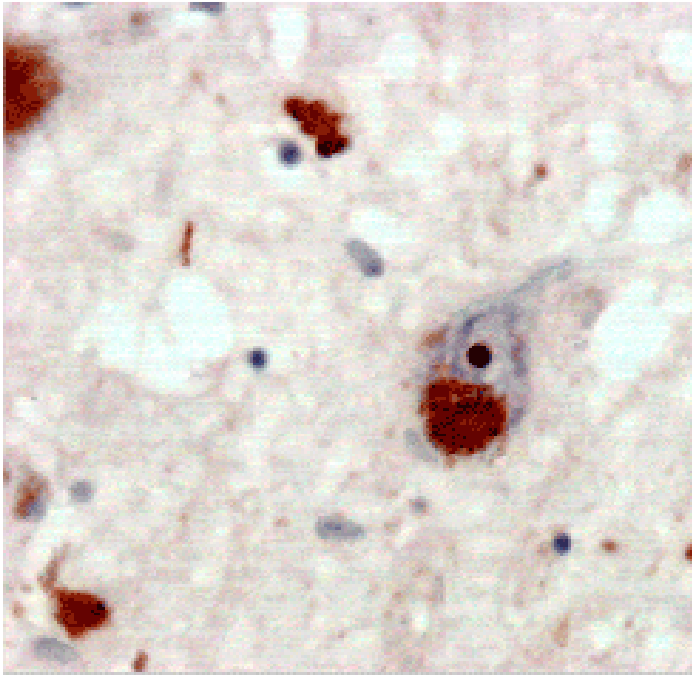


Figure 3: Formalin-fixed paraffin-embedded sections of hippocampus from human Alzheimer's disease brain stained with MAb to Polyubiquitin (K⁶³-linkage-specific) (HWA4C4) at 1:100 dilution with biotinylated secondary antibody and DAB detection.

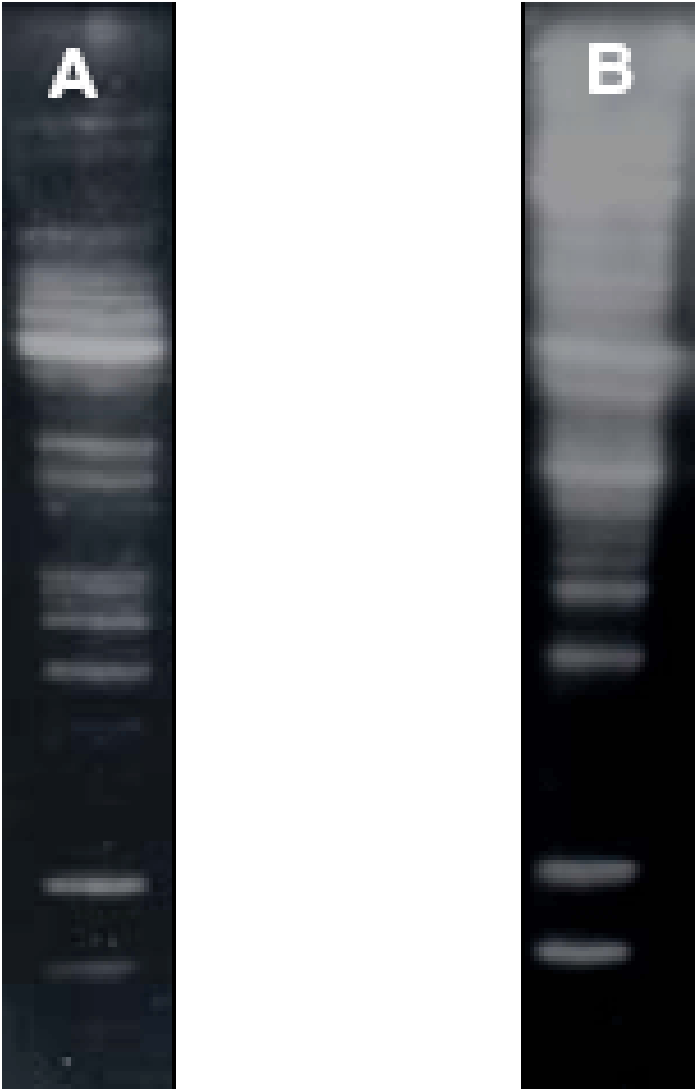


Figure 2: Western blot of HeLa S100 Fraction (BML-SW8750, 5µg) probed with MAb to Polyubiquitin (K⁶³-linkage-specific) (HWA4C4) (BML-PW0600) (Lane A) and MAb to Polyubiquitinated Conjugates (FK1) (BML-PW8805) (Lane B) at 1:1'000 dilution.

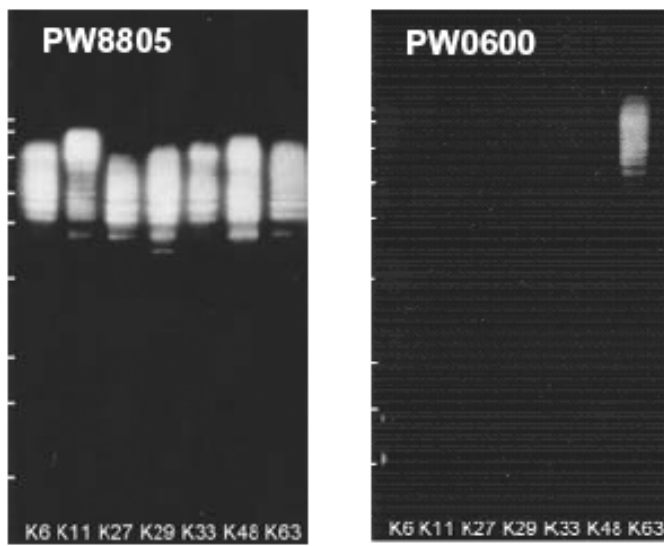


Figure 1: Western blot of single isopeptide-linkage polyubiquitinated conjugates K⁶ (BML-UW0615), K¹¹ (BML-UW0620), K²⁷ (BML-UW0625), K²⁹ (BML-UW0630), K³³ (BML-UW0635), K⁴⁸ (BML-UW0640), K⁶³ (BML-UW0645), 1µg each, using MAb to Polyubiquitinated Conjugates (FK1) (BML-PW8805) and MAb to Polyubiquitin (K⁶³-linkage-specific) (HWA4C4) (BML-PW0600) at 1:1'000 dilution.

Handling & Storage

Handling Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ELISA, ICC, IHC, WB

Clone HWA4C4

Crossreactivity Does not cross-react with any other isopeptide-linked (K6, K11, K27, K29, K33, or K48) polyubiquitinated species.

Formulation Liquid. In PBS containing 0.09% sodium azide.

Host Mouse

Immunogen Synthetic peptide containing the ubiquitin K63-isopeptide motif.

Isotype IgG2a

Recommendation Dilutions/Conditions Immunocytochemistry (1:100)Immunohistochemistry (1:100)Western Blot (1:500 – 1:1,000)Specificity for K63 of mAb to Polyubiquitin (K63-linkage-specific) (HWA4C4) is temperature sensitive. Western blot membranes must not be autoclaved. Excessive heat should be avoided in all applications. K63-linked polyubiquitin modified proteins comprise only a proportion of total cellular polyubiquitinated species. Increased sample loading may therefore be required for analysis by Western blot. Do not use milk as blocking solution. Use 1% BSA in an appropriate buffer instead.Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Source Purified from hybridoma tissue culture supernatant.

Species Reactivity Species independent

Specificity	Recognizes K ⁶³ -linked polyubiquitin in wide range of species.
UniProt ID	P0CG47 (UBB), P0CG48 (UBC)
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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