

Mono- and polyubiquitinated conjugates recombinant monoclonal antibody

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

(UBCJ)
Citations: 38

[View Online »](#)

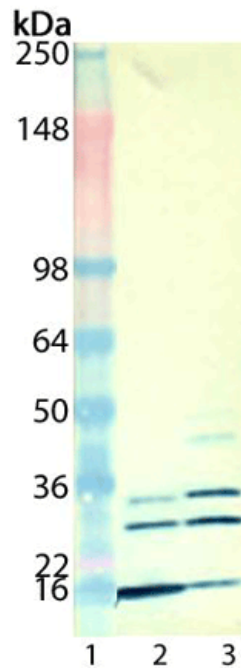
Ordering Information

[Order Online »](#)

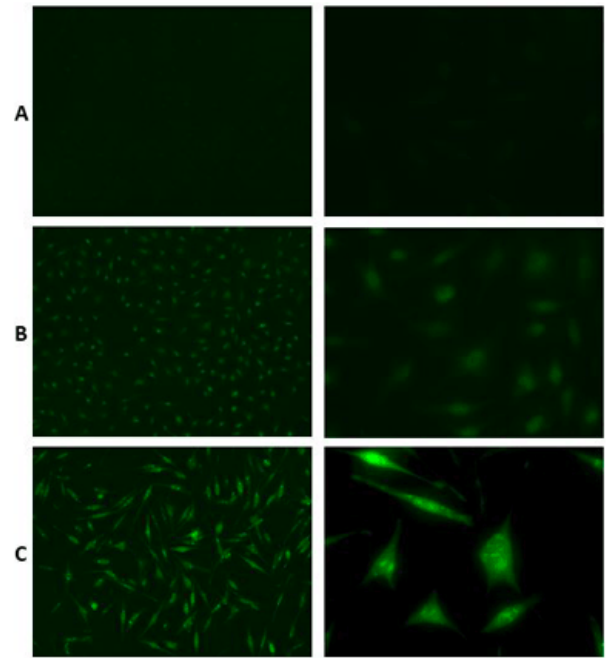
ENZ-ABS840-0100	100µg
ENZ-ABS840-0500	500µg

Manuals, SDS & CofA

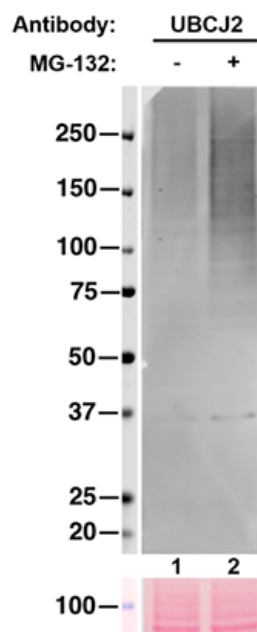
[View Online »](#)



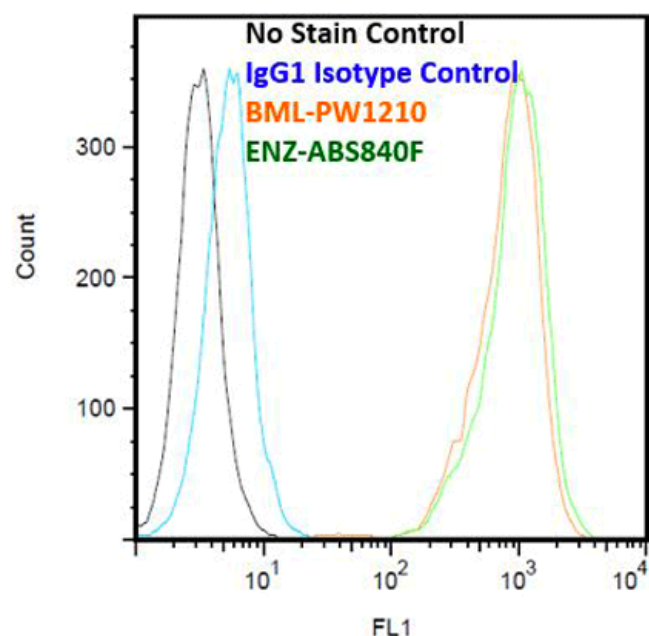
Western Blot analysis of ENZ-ABS840HRP detecting ubiquitin chains: Lane 1: MW Marker, Lane 2: K48-linked, Lane 3: K63-linked.



Immunocytochemistry (ICC) analysis of ubiquitinated protein conjugates using FITC labeled mono- and polyubiquitinated conjugates mAb (UBCJ2), (Prod. No. ENZ-ABS840F): HeLa cells, pre-treated with proteasome inhibitor MG-132 (Prod. No. BML-PI102, or ENZ-51035) at 5 μ M for 12 hours prior to analysis with antibody at 1:250 dilution. Panel A: isotype control (Prod. No. ADI-SAB-600FI); Panel B: untreated; Panel C: proteasome inhibitor-treated HeLa cells. Left: 20X and Right: 63X amplified.



Western Blot analysis of ubiquitinated conjugates in whole cell lysates using the Mono- and polyubiquitinated conjugates mAb (UBCJ2) (Prod. No. ENZ-ABS840). HeLa cells were treated for 12 hours with DMSO (vehicle, lane 1) or proteasome inhibitor MG-132 at 25 μ M (lane 2). Membrane was incubated O.N. (4°C) with 1 μ g/mL of mAb clone UBCJ2. Detection was done with a goat anti-mouse secondary antibody. Ponceau-S staining was performed as loading control (bottom panel). Image is a courtesy of Dr Jacob Seeler, Nuclear Organization and Oncogenesis Unit, Pasteur Institute, Paris (FR).



Flow cytometry analysis of 106 Jurkat cells stained using Mono- and polyubiquitinated conjugates mAb (UBCJ2) (FITC conjugate), (Prod. No. ENZ-ABS840F), at a concentration of 10 μ g/ml, side-by-side comparison to Mono- and polyubiquitinated conjugates mAb (FK2) (fluorescein labeled), (Prod. No. BML-PW1210).

Handling & Storage

Handling	Store unopened vial at -20°C. Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	ELISA, Flow Cytometry, ICC, WB
Clone	UBCJ2
Formulation	Liquid. In PBS containing 0.09% sodium azide.
Host	Mouse
Immunogen	Poly-ubiquitinylated lysozyme.
Isotype	IgG1κ
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	ELISA (1:1000)Immunocytochemistry (1:250)Flow cytometry (1:100)Western Blot (1:1000)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.
Species Reactivity	Species independent
Specificity	Recognizes mono- and polyubiquitinylated protein conjugates in a wide range of species.
UniProt ID	P0CG47 (UBB), P0CG48 (UBC)
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)