

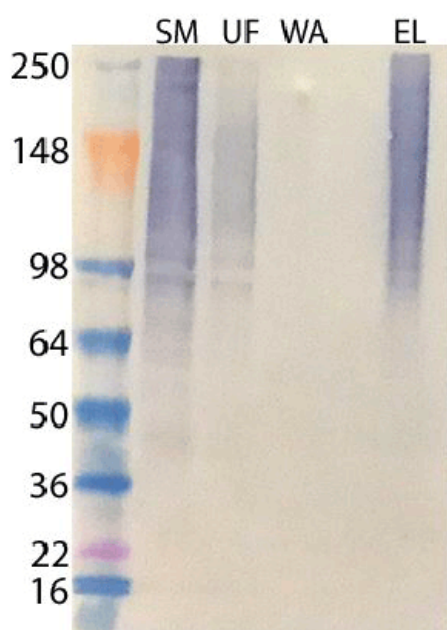
Mono- and polyubiquitinated conjugates recombinant monoclonal antibody (UBC12) (HRP conjugate)

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

Ordering Information [Order Online »](#)

ENZ-ABS840HRP-0025	25µg
ENZ-ABS840HRP-0100	100µg

Manuals, SDS & CofA [View Online »](#)



Western Blot analysis of ENZ-ABS840HRP detecting ubiquitinated protein conjugates captured in UBIQAPTURE-Q® kit (Prod. No. BML-UW8995): SM = Starting Material; UF = Unbound Fraction; WA = Washing Fraction; EL = Elution Fraction.

Handling & Storage

Handling	Do not freeze. Protect from light.
Short Term Storage	+4°C
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	ELISA, WB
Clone	UBCJ2
Formulation	Liquid. In PBS.
Host	Mouse
Immunogen	Poly-ubiquitinated lysozyme.
Isotype	IgG1κ
Purity Detail	Protein A affinity 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.
Species Reactivity	Species independent
Specificity	Recognizes mono- and polyubiquitinated protein conjugates in a wide range of species.
UniProt ID	P0CG47 (UBB), P0CG48 (UBC)



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

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

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

France
Phone: +33 1 42 44 55 55
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

Worry Free Guarantee This antibody is covered by our **Worry Free Guarantee**