

Polyubiquitinated conjugates recombinant monoclonal antibody (UBQM)

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

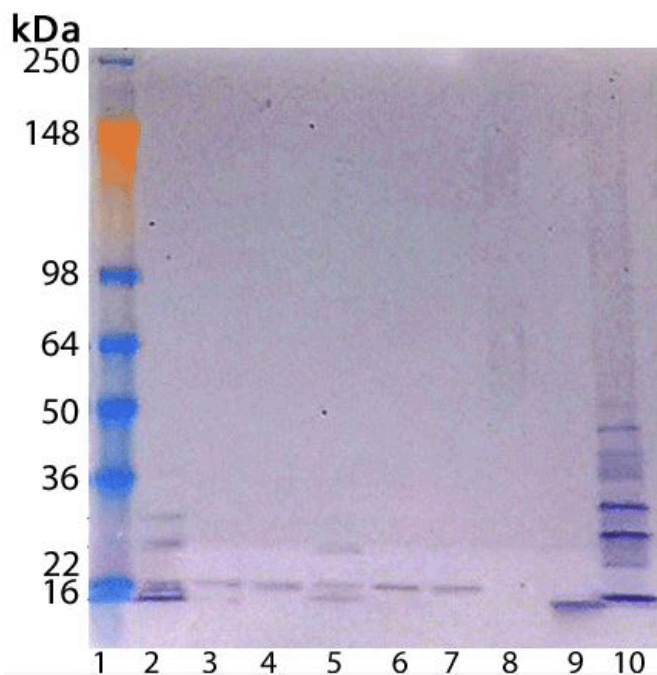
Ordering Information

[Order Online »](#)

ENZ-ABS842-0100	100µg
ENZ-ABS842-0500	500µg

Manuals, SDS & CofA

[View Online »](#)



Western Blot analysis of ENZ-ABS842 detecting ubiquitin chains: Lane 1: MW Marker, Lane 2-7: E2 ubiquitinated conjugates (Prod. No. BML-UW9050, BML-UW9060), Lane 9: [K6] mutant Ubiquitin, Lane 10: Ubiquitin chains (Ub2-7).

Handling & Storage

Use/Stability	Stable for at least 6 months after receipt when stored as recommended.
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, WB
Clone	UBQM
Formulation	Liquid. In PBS containing 0.09% sodium azide.
Host	Mouse
Immunogen	Poly-ubiquitinated lysozyme.
Isotype	IgG1k
Purity Detail	Protein G 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 polyubiquitinated conjugates
UniProt ID	P0CG47 (UBB), P0CG48 (UBC), P62979 (RPS27A), P62987 (UBA52)

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)

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