

Myc-tag monoclonal antibody (PL14)

Epitope tagging is a widely accepted technique that fuses an epitope peptide to a certain protein as a marker for gene expression. With this technique, gene expression can be easily monitored on Western blotting, immunoprecipitation and immunofluorescence utilizing an antibody that recognizes such an epitope. The Myc-Tag (EQKLISEEDL) corresponds to the partial peptide sequence of the human c-myc gene product.

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

Citations: 3

[View Online »](#)

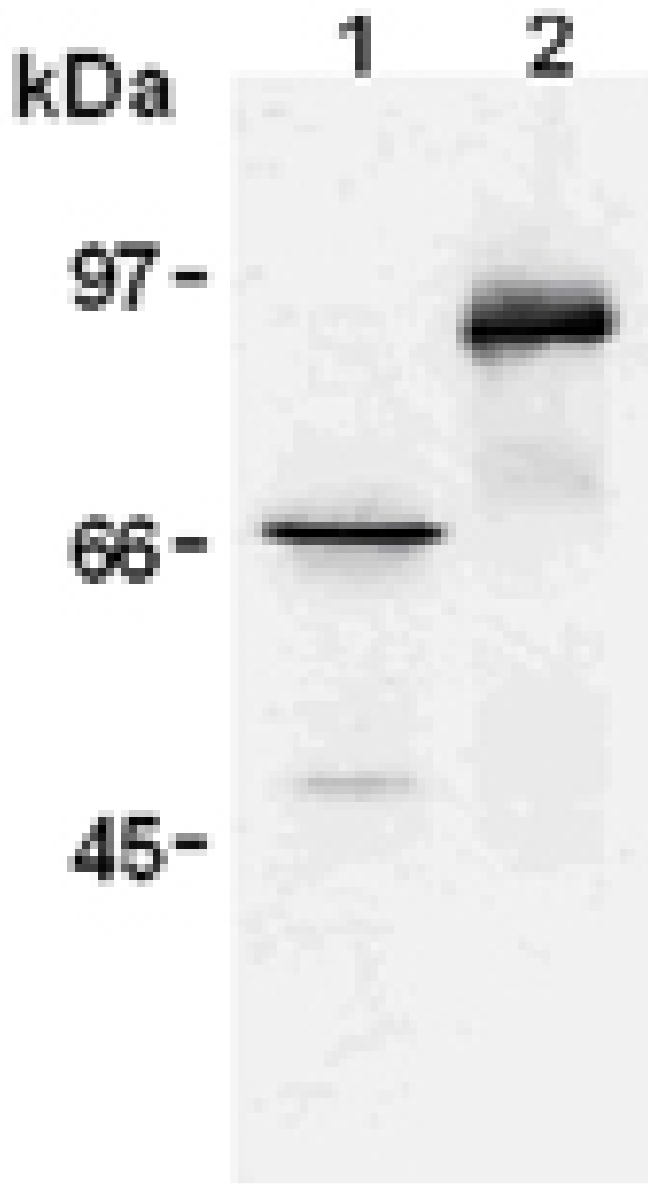
Ordering Information

[Order Online »](#)

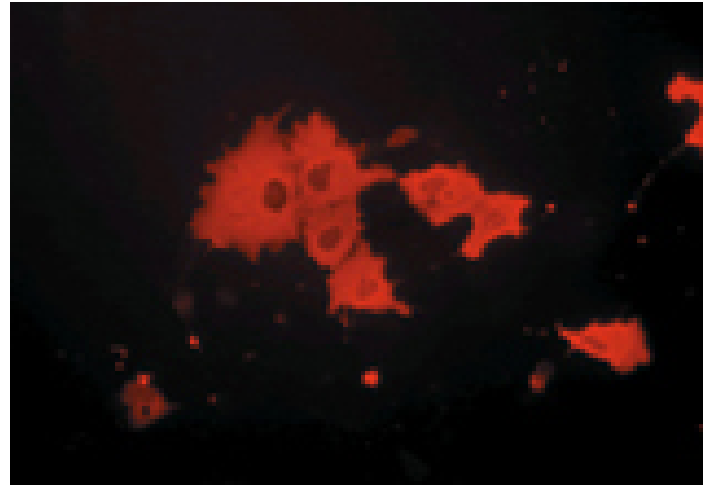
ADI-MSA-110-F	200µg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of 293T cells transfected with 6Myc/Heparanase (lane 1) and 6Myc/cdc25 (lane 2), probed with Myc-tag mAb (PL14).



Immunocytochemistry analysis of monkey Cos-7 cells transfected with 6myc-hCAP1, stained with Myc-tag mAb (PL14).

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ICC, IP, WB

Clone PL14

Formulation Liquid. In PBS, pH 7.2, containing 50% glycerol. No preservative has been added.

Host Mouse

Immunogen 6myc-tag GST fusion protein.

Isotype IgG1

Purity Detail Protein A affinity purified.

Recommendation
Dilutions/Conditions Immunocytochemistry (2µg/ml)Immunoprecipitation (5µg/ml)Western Blot (1µg/ml, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Specificity Recognizes the Myc-tag of Myc-tagged recombinant proteins.

UniProt ID P01106 (human)

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)