

# 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 »](#)

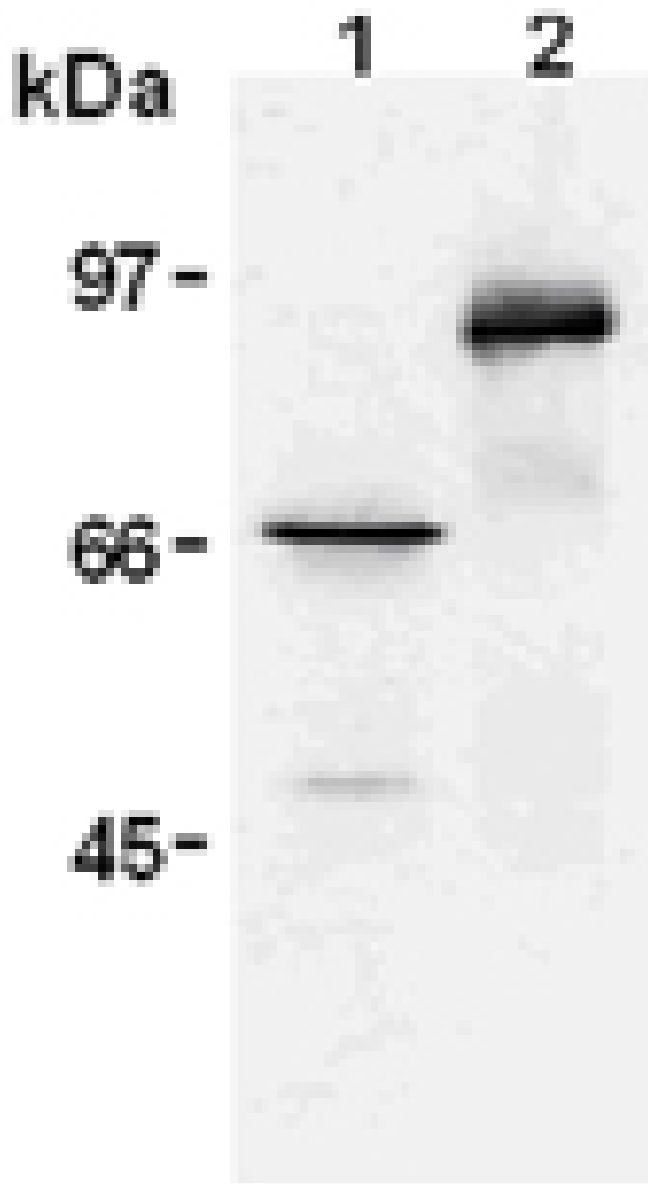
## Ordering Information

[Order Online »](#)

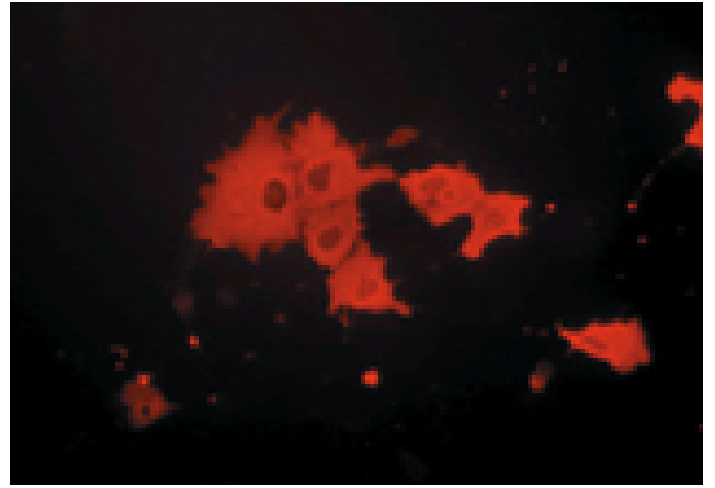
ADI-MSA-110-F	200µg
---------------	-------

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 <a href="#">Worry-Free Guarantee</a>

Last modified: May 29, 2024



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-fr@enzolifesciences.com](mailto:info-fr@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)