

# Ki-67 monoclonal antibody (Ready-to-Use)

## Ready-to-use monoclonal antibody to Ki-67 for use with IHC applications

Ki-67 is a nuclear, non-histone protein that is expressed only during active phases of the cell cycle (G1, S, G2 and M), but not in the resting phases (G0 and G1 early phase). Although the antigen has also been associated with ribosomal RNA transcription, it is strongly linked to cell proliferation and has thus been indicated as an effective marker in grading the proliferation rate of tumors, including those of the brain, breast, cervix, and prostate.

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

### Ordering Information

[Order Online »](#)

ENZ-ABS682-0015	15ml
-----------------	------

### Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	When stored correctly, the antibody is stable until the date indicated on the label. To ensure proper stability of the antibody after each run, replace the cap and immediately place the bottle in a refrigerator in an upright position.
<b>Handling</b>	Avoid freeze/thaw cycles.
<b>Short Term Storage</b>	+4°C
<b>Long Term Storage</b>	+4°C
<b>Shipping</b>	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	MKI67
<b>Application</b>	IHC
<b>Formulation</b>	Liquid. Ready-to-use solution in PBS containing stabilizers and preservatives.
<b>Host</b>	Mouse
<b>Immunogen</b>	Recombinant Ki-67
<b>Isotype</b>	IgG
<b>Purity Detail</b>	Affinity purified.
<b>Recommendation Dilutions/Conditions</b>	Immunohistochemistry (RTU)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
<b>Species Reactivity</b>	Human
<b>UniProt ID</b>	P46013

### 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)