

Ki-67 monoclonal antibody

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](#).

Citations: 1

[View Online »](#)

Ordering Information

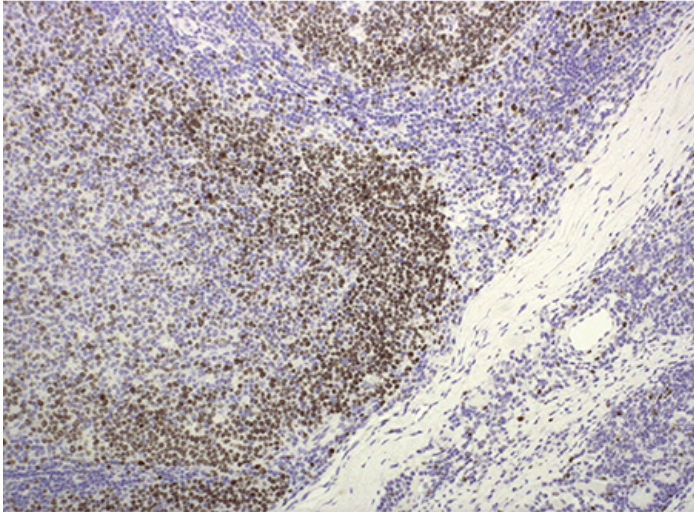
[Order Online »](#)

ENZ-ABS678-1000	1ml
ENZ-ABS678-0100	100µl

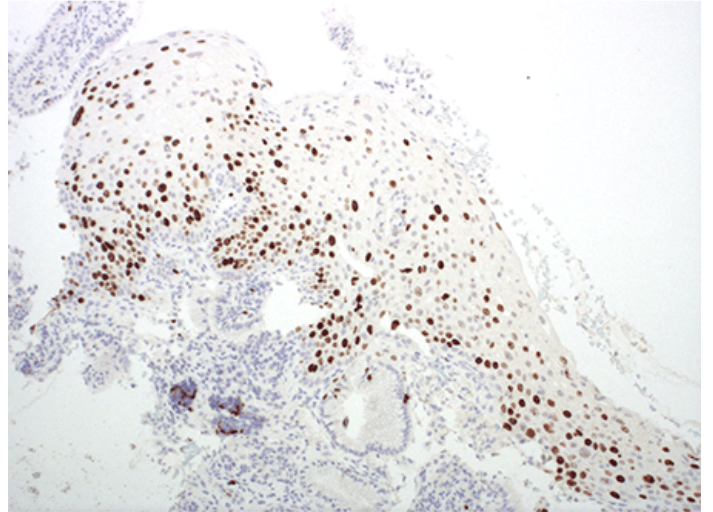
Manuals, SDS & CofA

[View Online »](#)

- Strong, clear staining with minimal background
- Cost-effective option compared to competitor's antibody



Cervical tissue incubated with Ki-67 monoclonal antibody and then stained with POLYVIEW® PLUS (anti-Mouse) HRP-DAB Kit (ENZ-KIT160).



Cervical tissue incubated with Ki-67 monoclonal antibody and then stained with POLYVIEW® PLUS (anti-Mouse) HRP-DAB Kit (ENZ-KIT160).

Handling & Storage

Use/Stability	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.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	MKI67
Application	IHC
Application Notes	For IHC applications: Dilute the concentrate as appropriate in Antibody Blocking Buffer (ENZ-ACC108), and then apply 150 µL to the slide, as indicated in the user's IHC protocol.
Formulation	Liquid. In Tris buffer, pH 7.5, containing 1% BSA and <0.1% sodium azide.
Host	Mouse
Immunogen	Recombinant Ki-67.
Isotype	IgG
Purity Detail	Affinity-purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (1:100-1:400)Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	P46013

Worry-free Guarantee

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

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-
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)