

Ki-67 monoclonal antibody (SP6)

Ki-67 is a nuclear protein associated with cell proliferation. It is expressed during all active phases of the cell cycle but not in resting cells, making it a key marker for assessing growth fractions in tissues and tumors.

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

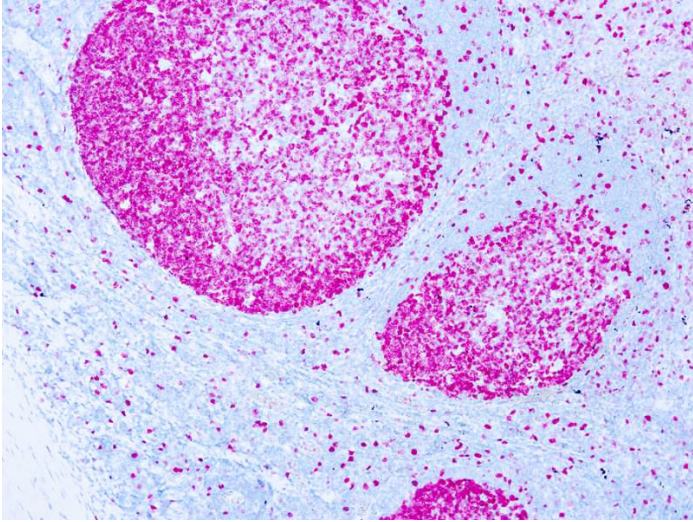
Ordering Information

[Order Online »](#)

ENZ-ABS1082-0500	500µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Formalin fixed paraffin embedded human tonsil stained with Ki-67 (SP6) antibody

Handling & Storage

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 (PS)

Application Notes Pretreatment – EDTA Buffer pH 8.0. Incubation & Temperature 30-60 min @ RT

Clone SP6

Formulation Contains sodium azide as a preservative.

Host Rabbit

Immunogen Synthetic peptide corresponding to the C-terminus of human Ki-67.

Isotype IgG

Positive Control Tonsil

Purity Detail Purified immunoglobulin.

Recommendation Immunohistochemistry (paraffin sections) (1:25-1:50)

Dilutions/Conditions Suggested dilutions/conditions may not be available for all applications.
Optimal conditions must be determined individually for each application.

Species Reactivity 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)