# 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**

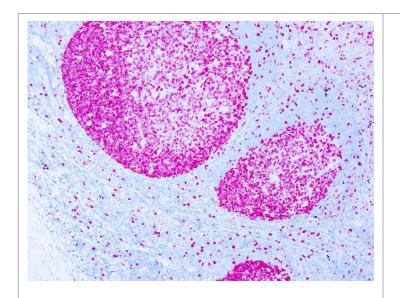
Order Online »

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 Dilutions/Conditions** Immunohistochemistry (paraffin sections) (1:25-1:50)

Suggested dilutions/conditions may not be available for all

applications.

Optimal conditions must be determined individually for

each application.

Species Reactivity Human

Last modified: June 6, 2025



info-

eu@enzolifesciences.com