

Version 2b: Last updated 6 August 2025

# **ab16667**

## Protocol for Ki-67

Clone ID: SP6 Host Species: Rabbit monoclonal

This product is for research use only and is not intended for diagnostic use.

# Protocol

## 1. Deparaffinization/Rehydration

- 1.1. Heat slides in an oven at 63 °C - 65 °C for about 1 hour.
- 1.2. Deparaffinize/hydrate using the following series of washes: twice Xylene rinses (5-10 min each), followed by twice anhydrous ethanol rinses (5 min each), twice 95% ethanol rinses (5 min each), 85% ethanol rinse (3 min), 75% ethanol rinse (3 min), and finally thrice distilled water washes.

## 2. Antigen Retrieval

This is recommended Heat Induced Epitope Retrieval (HIER) using Decloaking Chamber/Pressure Cooker.

- 2.1. Add 500 ml of distilled water to Decloaker/Pressure Cooker.
- 2.2. Immerse slides into staining dish containing Antigen Retrieval Solution: **Heat mediated antigen retrieval using ab93678 (Citrate buffer, pH 6.0)**. Place staining dish into Decloaking chamber.
- 2.3. Program to run for 30 s at 125 °C or 15-30 min at 110 °C.
- 2.4. Let it cool down to room temperature (10-30 min).
- 2.5. Removes slides and rinse in distilled water.
- 2.6. Proceed to Staining step.

## 3. Staining

- 3.1. Wash slides with distilled water thrice (3 min each if on a shaker).
- 3.2. Inactivate endogenous peroxidase by covering tissue with 3% hydrogen peroxide for 5-10 min.

- 3.3. Wash slides three times with TBST (3 min each on a shaker).
- 3.4. Block slides with the blocking solution for 30-60 min.
- 3.5. Dilute primary antibody in primary antibody diluent per recommendation on datasheet.
- 3.6. Apply primary antibody to each section and **incubate overnight in the humidified chamber (2~8 °C)**.
- 3.7. Wash slides three times with TBST (3 min each on a shaker).
- 3.8. Apply to each section secondary HRP-conjugated anti-rabbit antibody diluted in the blocking solution per manufacturer's recommendation; incubate for 30 min at room temperature.
- 3.9. Wash slides three times with TBST (5 min each on a shaker).
- 3.10. Add freshly prepared DAB substrate to the sections and incubate until stain develops (10-60 s, according to the instruction).
- 3.11. Rinse sections with distilled water.
- 3.12. Counterstain with Hematoxylin (1-2 min, according to instruction).
- 3.13. Rinse sections with water.
- 3.14. Dehydrate samples using 75% ethanol rinse (3 min), 95% ethanol rinse (3 min), twice anhydrous ethanol rinses (3 min each), followed by twice rinses with Xylene or HistoClear (5 min each).
- 3.15. Dry slides and mount the coverslips onto slides using permanent mounting medium.

## Technical Support

Copyright © 2025 Abcam, All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

**For all technical or commercial enquiries please go to:**

[www.abcam.com/en-us/contact-us](http://www.abcam.com/en-us/contact-us)

[www.abcam.cn/contactus](http://www.abcam.cn/contactus) (China)

[www.abcam.co.jp/contactus](http://www.abcam.co.jp/contactus) (Japan)