

## ab64212 Endogenous Avidin/Biotin Blocking Kit

For the blocking of endogenous biotin during immunohistochemistry.

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

### Overview

In the demonstration of tissue antigens by Immunohistochemistry, the evaluation of specific staining is sometimes impeded by the presence of endogenous biotin, which may lead to an erroneous interpretation or false positive results. This blocking system is intended for use in immunohistochemical staining to block nonspecific staining due to endogenous biotin during the IHC detection of antigens.

### Precautions

Please observe safe laboratory practice and consult the safety datasheet.

For general guidelines, precautions, limitations on the use of our assay kits and general assay troubleshooting tips, particularly for first time users, please consult our guide:

[www.abcam.com/assaykitguidelines](http://www.abcam.com/assaykitguidelines)

For typical data produced using the assay, please see the assay kit datasheet on our website.

### Storage and Stability:

Store kit at 4°C immediately on receipt and check below for storage for individual components. Kit can be stored for 1 year from receipt, if components have not been reconstituted.

Aliquot components in working volumes before storing at the recommended temperature.

Avoid repeated freeze-thaws of reagents.

### Materials Supplied:

Item	Quantity	Storage temperature (before prep)	Storage temperature (after prep)
Avidin Block	15 mL	4°C	N/A
Biotin Block	15 mL	4°C	N/A

### Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully perform this assay:

- Buffer of choice for washing tissue sections

### Reagent Preparation

**Avidin and Biotin Block:** Ready to use as supplied.

### Assay Procedure

#### **Blocking/staining protocol:**

1. Incubate tissue sections with avidin block for 10-15 minutes.
2. Wash 2 times with buffer.
3. Incubate tissue sections with biotin block for 10-15 minutes.
4. Wash 2 times with buffer.
5. Apply primary antibody to the tissue sections.

### Troubleshooting

General troubleshooting points are found at [www.abcam.com/assaykitguidelines](http://www.abcam.com/assaykitguidelines)

### Technical Support

Copyright © 2020 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/contactus](http://www.abcam.com/contactus)

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

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