

# Mouse and Rabbit Specific HRP Detection IHC Kit ab93677

[10 References](#) [1 图像](#)

### 概述

#### 产品名称

小鼠and兔Specific HRP Detection IHC试剂盒

#### 产品概述

Mouse and Rabbit Specific HRP Detection IHC Kit ab93677 contains complet reagents for a labeled streptavidin-biotin immunoenzymatic antigen detection system. This technique involves the sequential incubation of the specimen with an unconjugated primary antibody specific to the target antigen, a biotinylated secondary antibody which reacts with the primary antibody, enzyme-labeled streptavidin, and substrate-chromogen.

#### 说明

Reacts with Mouse and Rabbit IgG (H+L)

ab93677 **does not** contain chromogen or substrate.

#### 经测试应用

适用于: IHC-P

### 性能

#### 存放说明

Store at +4°C. Please refer to protocols.

组件	60 ml
Biotinylated Goat Anti-Polyvalent	1 x 60ml
Protein block	1 x 60ml
Streptavidin Peroxidase	1 x 60ml

#### 相关性

Polyvalent HRP is a labeled streptavidin-biotin immunoenzymatic antigen detection system. Horseradish peroxidase (HRP) readily combines with hydrogen peroxide (H<sub>2</sub>O<sub>2</sub>) and the resultant complex can oxidize a wide variety of chromogenic substrates.

### 应用

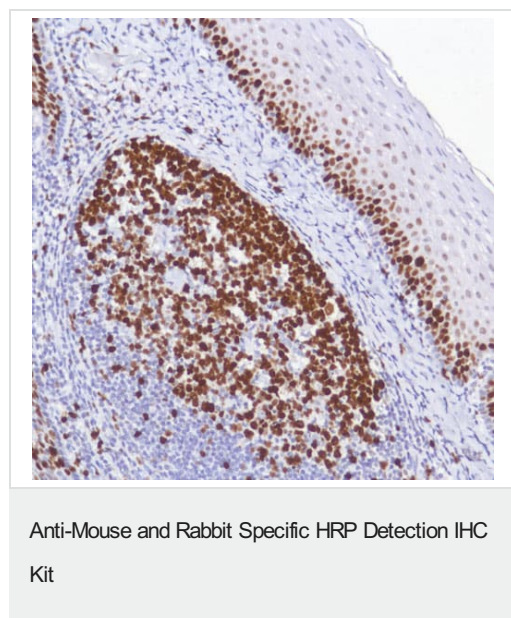
#### The Abpromise guarantee

**Abpromise™**承诺保证使用ab93677于以下的经测试应用

“应用说明”部分 下显示的仅为推荐的起始稀释度;实际最佳的稀释度/浓度应由使用者检定。

应用	Ab评论	说明
IHC-P		Use at an assay dependent dilution.

## 图片



Immunohistochemical analysis of Human Tonsil tissue labeling Ki-67 using the ab93677 anti-Mouse and Rabbit specific HRP Detection IHC kit.

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

## Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.cn/abpromise> or contact our technical team.

## Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors