

## Product datasheet

# Mouse and Rabbit Specific HRP Detection IHC Kit ab93677

[10 References](#) [1 Image](#)

### Overview

<b>Product name</b>	Mouse and Rabbit Specific HRP Detection IHC Kit
<b>Product overview</b>	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.  Reacts with Mouse and Rabbit IgG (H+L)
<b>Notes</b>	ab93677 <b>does not</b> contain chromogen or substrate.
<b>Tested applications</b>	<b>Suitable for:</b> IHC-P

### Properties

**Storage instructions** Store at +4°C. Please refer to protocols.

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

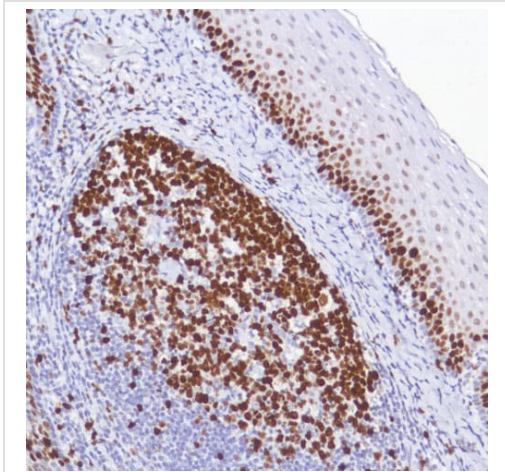
**Relevance** Polyvalent HRP is a labeled streptavidin-biotin immunoenzymatic antigen detection system. Horseradish peroxidase (HRP) readily combines with hydrogen peroxide (H<sub>2</sub>O<sub>2</sub>) and the resultant complex can oxidize a wide variety of chromogenic substrates.

### Applications

**The Abpromise guarantee** Our **Abpromise guarantee** covers the use of ab93677 in the following tested applications. The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
IHC-P		Use at an assay dependent dilution.

## Images



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

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.com/abpromise> or contact our technical team.

## Terms and conditions

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