

Mouse and Rabbit Specific HRP/DAB Detection IHC kit ab64264

★★★★★ [3 Abreviews](#) [174 References](#) [2 图像](#)

概述

产品名称

小鼠和兔Specific HRP/DAB Detection IHC试剂盒

产品概述

Mouse and Rabbit specific HRP/DAB Detection IHC Kit ab64264 is a complete kit to detect a primary antibody raised in either mouse or rabbit, using HRP-labeled-streptavidin and a biotinylated anti-rabbit/anti-mouse secondary antibody.

It is supplied with a full protocol and with peroxidase and protein blocking reagents, biotinylated secondary antibody, streptavidin-HRP, and DAB substrate.

In the method:

- a sample is incubated with an unconjugated primary antibody specific to the target antigen
- a biotinylated secondary antibody then binds to the primary antibody
- HRP-labeled streptavidin then binds to the secondary antibody
- the HRP produces a brown colored substance at the site of primary antibody binding by reacting with DAB.

说明

For alternative streptavidin-biotin IHC detection kits see:

- Anti-Rabbit HRP-DAB IHC kit [ab64261](#)
- Anti-Mouse HRP-DAB IHC Kit [ab64259](#)

For high sensitivity HRP micro-polymer IHC detection kits see:

- Anti-Mouse / anti-Rabbit HRP/DAB Micro-polymer IHC Kit [ab236466](#)
- Anti-Rabbit HRP-DAB Micro-polymer IHC Kit [ab236469](#)

Alternatively, see [HRP-AEC IHC detection kits](#).

For staining mouse tissues with mouse antibodies, use mouse-on-mouse polymer kit [ab127055](#).

Find more kits and reagents for antigen retrieval, blocking, signal amplification, visualization, counterstaining, and mounting in the [IHC kits and reagents guide](#).

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	15 ml
50x DAB Chromogen	1 x 0.5ml
Biotinylated Goat Anti-Polyvalent	1 x 15ml
DAB substrate	1 x 15ml
Hydrogen Peroxide Block	1 x 15ml
Protein block	1 x 15ml
Streptavidin Peroxidase	1 x 15ml

相关性

HRP/DAB detection kit is a immunoenzymatic antigen detection system for immunohistochemistry techniques.

应用

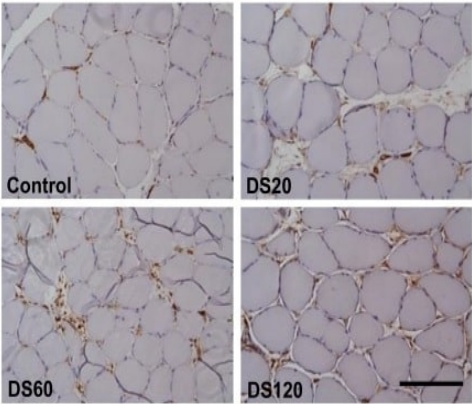
The Abpromise guarantee

Abpromise™承诺保证使用ab64264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



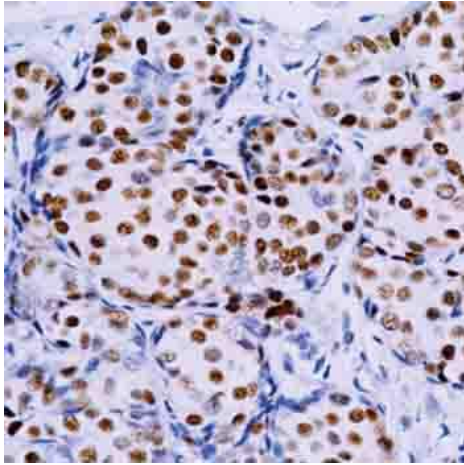
Control DS20 DS60 DS120

Yu S.H., PLoS One (9):e114649 (2014)

Immunohistochemistry - ab64264

Yu S.H., PLoS One 9 (12), Fig 7c. doi: 10.1371/journal.pone.0114649 Reproduced under the Creative Commons license <http://creativecommons.org/licenses/by/4.0/>

Immunohistochemical analysis staining CD163 in mouse soleus muscle section. Samples were incubated in rabbit CD163 antibody at 1:450 dilution and visualized using mouse and rabbit specific HRP/DAB detection IHC Kit (ab64264).



Immunohistochemistry (Formalin/PFA-fixed paraffin-embedded sections) - Anti Mouse & Rabbit HRP/DAB Detection Kit (ab64264)

Staining of human estrogen receptor in breast carcinoma tissue using **ab16660** - Rabbit monoclonal [SP1] to Estrogen Receptor, and **ab64246** - Mouse & Rabbit HRP/DAB Detection 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