

Rabbit specific IHC polymer detection kit HRP/DAB ab209101

[17 References](#) [8 图像](#)

概述

产品名称	兔specific IHC polymer detection试剂盒HRP/DAB
产品概述	Rabbit specific IHC polymer detection kit HRP/DAB (ab209101) is a sensitive immunoenzymatic system for detecting antibodies bound to antigens within tissue sections by light microscopy. Signal amplification is achieved using an indirect polymer detection system. As Rabbit specific IHC polymer detection kit HRP/DAB does not use avidin or biotin, nonspecific staining from endogenous avidin-biotin activity is eliminated and there is no need for additional steps to block endogenous biotin. The primary antibody specific to an antigen on the tissue section is detected by primary antibody Amplifier. The HRP polymer is added to enable detection with signal amplification. Finally, antibody-bound antigen sites are visualized by adding an appropriate substrate/chromogen (not provided).
经测试应用	适用于: IHC

性能

存放说明 Store at +4°C. Please refer to protocols.

组件	15 ml
Amplifier	1 x 15ml
Detector	1 x 15ml

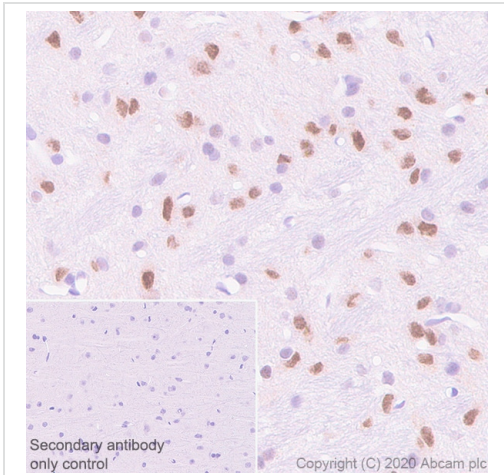
应用

The Abpromise guarantee **Abpromise™**承诺保证使用ab209101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		Use at an assay dependent concentration.

图片

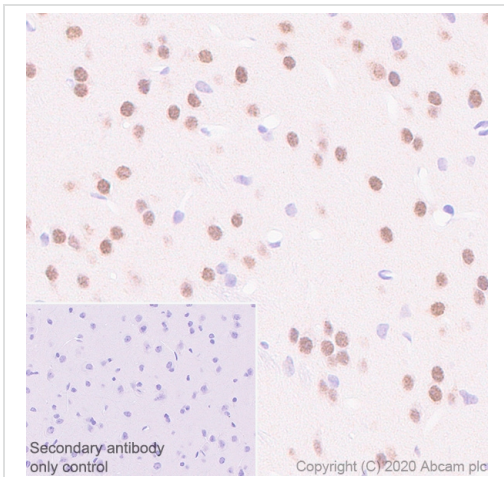


Immunohistochemistry paraffin embedded sections -
Rabbit specific IHC polymer detection kit HRP/DAB
(ab209101)

Immunohistochemical analysis of paraffin-embedded Rat cerebrum tissue labeling USP22 with **ab195289** at 1/1000 dilution followed by a ready to use Rabbit specific IHC polymer detection kit HRP/DAB (ab209101). Nuclear staining on rat cerebrum. The section was incubated with **ab195289** for 30 mins at room temperature. The immunostaining was performed on a Leica Biosystems BOND[®] RX instrument. Counterstained with Hematoxylin.

Secondary antibody only control: Secondary antibody is a ready to use Rabbit specific IHC polymer detection kit HRP/DAB (ab209101).

Heat mediated antigen retrieval with Citrate buffer (pH 6.0, epitope retrieval solution 1) for 20 mins

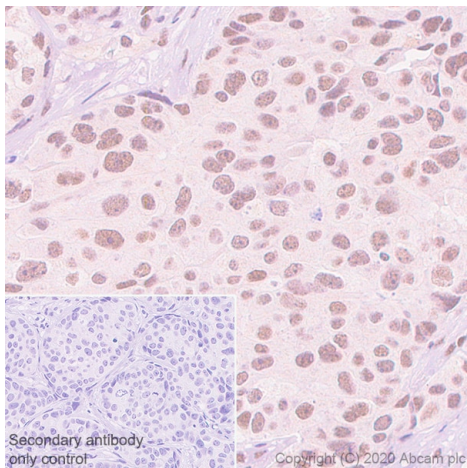


Immunohistochemistry paraffin embedded sections -
Rabbit specific IHC polymer detection kit HRP/DAB
(ab209101)

Immunohistochemical analysis of paraffin-embedded Mouse cerebrum tissue labeling USP22 with **ab195289** at 1/1000 dilution followed by a ready to use Rabbit specific IHC polymer detection kit HRP/DAB (ab209101). Nuclear staining on mouse cerebrum. The section was incubated with **ab195289** for 30 mins at room temperature. The immunostaining was performed on a Leica Biosystems BOND[®] RX instrument. Counterstained with Hematoxylin.

Secondary antibody only control: Secondary antibody is a ready to use Rabbit specific IHC polymer detection kit HRP/DAB (ab209101).

Heat mediated antigen retrieval with Citrate buffer (pH 6.0, epitope retrieval solution 1) for 20 mins

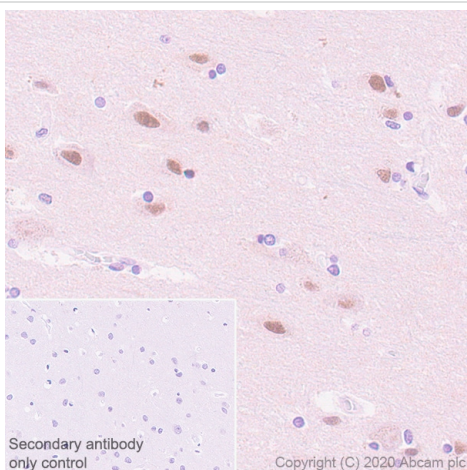


Immunohistochemistry paraffin embedded sections -
Rabbit specific IHC polymer detection kit HRP/DAB
(ab209101)

Immunohistochemical analysis of paraffin-embedded Human breast carcinoma tissue labeling USP22 with **ab195289** at 1/100 dilution followed by a ready to use Rabbit specific IHC polymer detection kit HRP/DAB (ab209101). Nuclear staining on human breast carcinoma. The section was incubated with **ab195289** for 30 mins at room temperature. The immunostaining was performed on a Leica Biosystems BOND[®] RX instrument. Counterstained with Hematoxylin.

Secondary antibody only control: Secondary antibody is a ready to use Rabbit specific IHC polymer detection kit HRP/DAB (ab209101).

Heat mediated antigen retrieval with Tris-EDTA buffer (pH 9.0, epitope retrieval solution2) for 20 mins

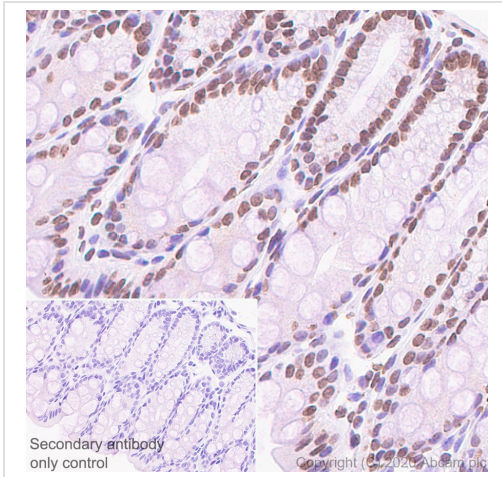


Immunohistochemistry paraffin embedded sections -
Rabbit specific IHC polymer detection kit HRP/DAB
(ab209101)

Immunohistochemical analysis of paraffin-embedded Human cerebrum tissue labeling USP22 with **ab195289** at 1/100 dilution followed by a ready to use Rabbit specific IHC polymer detection kit HRP/DAB (ab209101). Nuclear staining on human cerebrum. The section was incubated with **ab195289** for 30 mins at room temperature. The immunostaining was performed on a Leica Biosystems BOND[®] RX instrument. Counterstained with Hematoxylin.

Secondary antibody only control: Secondary antibody is a ready to use Rabbit specific IHC polymer detection kit HRP/DAB (ab209101).

Heat mediated antigen retrieval with Tris-EDTA buffer (pH 9.0, epitope retrieval solution2) for 20 mins

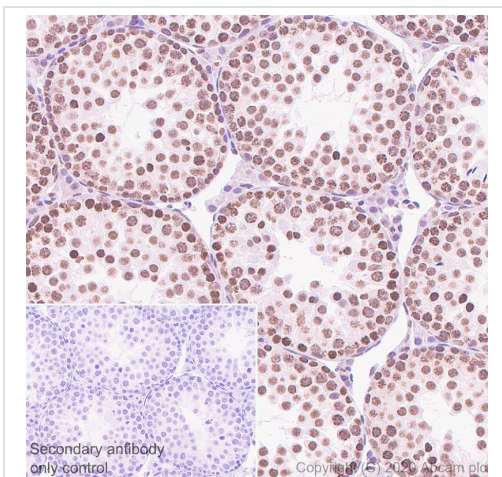


Immunohistochemistry paraffin embedded sections -
Rabbit specific IHC polymer detection kit HRP/DAB
(ab209101)

Immunohistochemical analysis of paraffin-embedded Mouse colon tissue labeling HSF1 with **ab52757** at 1/1000 dilution followed by a ready to use Rabbit specific IHC polymer detection kit HRP/DAB (ab209101). Nuclear staining on mouse colon. The section was incubated with **ab52757** for 30 mins at room temperature. The immunostaining was performed on a Leica Biosystems BOND® RX instrument Counterstained with Hematoxylin.

Secondary antibody only control: Secondary antibody is a ready to use Rabbit specific IHC polymer detection kit HRP/DAB (ab209101).

Heat mediated antigen retrieval with Citrate buffer (pH 6.0, epitope retrieval solution 1) for 20 mins

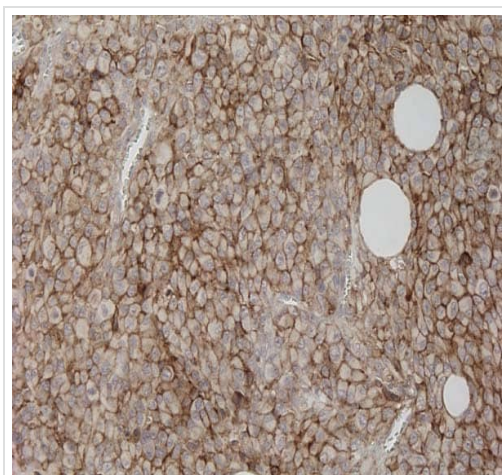


Immunohistochemistry paraffin embedded sections -
Rabbit specific IHC polymer detection kit HRP/DAB
(ab209101)

Immunohistochemical analysis of paraffin-embedded Mouse testis tissue labeling HSF1 with **ab52757** at 1/1000 dilution followed by a ready to use Rabbit specific IHC polymer detection kit HRP/DAB (ab209101). Nuclear staining on mouse testis. The section was incubated with **ab52757** for 30 mins at room temperature. The immunostaining was performed on a Leica Biosystems BOND® RX instrument Counterstained with Hematoxylin.

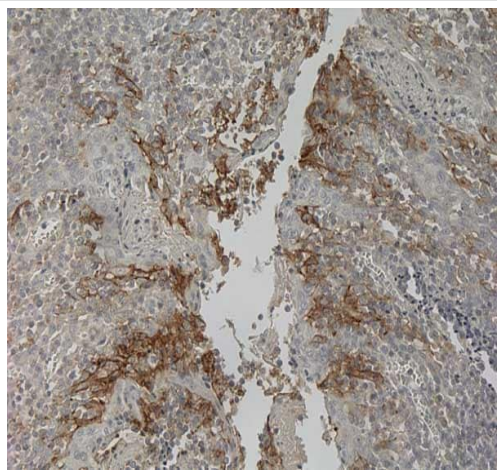
Secondary antibody only control: Secondary antibody is a ready to use Rabbit specific IHC polymer detection kit HRP/DAB (ab209101).

Heat mediated antigen retrieval with Citrate buffer (pH 6.0, epitope retrieval solution 1) for 20 mins



ab209101 Melanoma

Immunohistochemistry of human melanoma tissue (formalin/PFA-fixed paraffin embedded sections) stained with Anti-PD-L1 antibody [28-8] (**ab205921**) and detected with Rabbit specific IHC polymer detection kit HRP/DAB (ab209101).



ab209101 Tonsil

Immunohistochemistry of human tonsil tissue (formalin/PFA-fixed paraffin embedded sections) stained with Anti-PD-L1 antibody [28-8] ([ab205921](#)) and detected with Rabbit specific IHC polymer detection kit HRP/DAB ([ab209101](#)).

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