

# Mouse and Rabbit Specific HRP/DAB IHC Detection Kit - Micro-polymer ab236466

★★★★★ [1 Abreviews](#) [47 References](#) [5 图像](#)

### 概述

产品名称	小鼠和兔Specific HRP/DAB IHC Detection试剂盒- Micro-polymer
检测方法	Colorimetric
检测类型	Micro-Polymer
产品概述	This IHC kit is used to detect mouse and rabbit primary antibodies in immunohistochemistry (IHC). It uses an HRP micro-polymer secondary antibody for highly sensitive detection.

HRP micro-polymer methods are more sensitive than the traditional streptavidin-biotin method, and they do not require a blocking step for endogenous biotin.

The kit includes ready-to-use hydrogen peroxide blocking reagent, protein blocking reagent, HRP micro-polymer conjugated goat anti-rabbit secondary antibody, and DAB chromogen/substrate. It includes an unconjugated rabbit anti-mouse antibody, which is used together with the HRP polymer anti-rabbit antibody to detect mouse primary antibodies.

说明	The previous version of this kit, the EXPOSE Mouse and Rabbit Specific HRP/DAB Detection IHC Kit, was used by researchers to stain over 150,000 slides since 2012 (based on volumes of reagent sold).
----	---

### Other IHC detection products

For alternative kits, see anti-Rabbit HRP-DAB Micro-polymer IHC Kit [ab236469](#), and [HRP-AEC micro-polymer IHC 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	60 ml	125 ml
50x DAB Chromogen	1 x 0.5ml	1 x 2ml	1 x 4ml
DAB substrate	1 x 15ml	1 x 60ml	1 x 125ml
Goat anti-rabbit HRP Conjugate	1 x 15ml	1 x 60ml	1 x 125ml
Hydrogen Peroxide Block	1 x 15ml	1 x 60ml	1 x 125ml
Mouse Specifying Reagent (Complement)	1 x 15ml	1 x 60ml	1 x 125ml
Protein block	1 x 15ml	1 x 60ml	1 x 125ml

相关性

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

应用

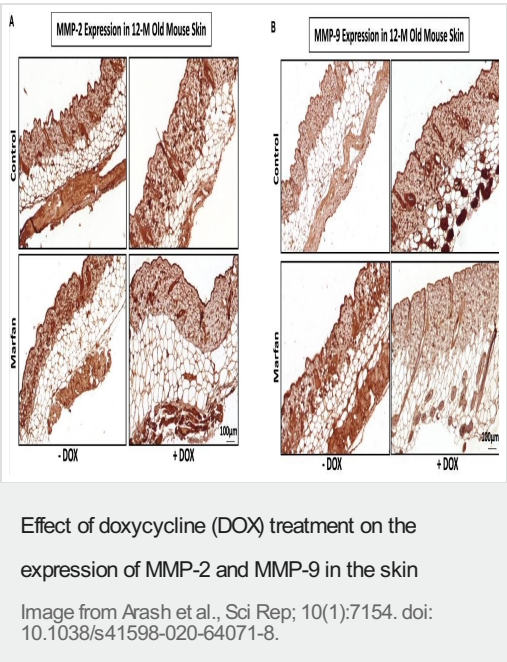
The Abpromise guarantee

Abpromise™承诺保证使用ab236466于以下的经测试应用

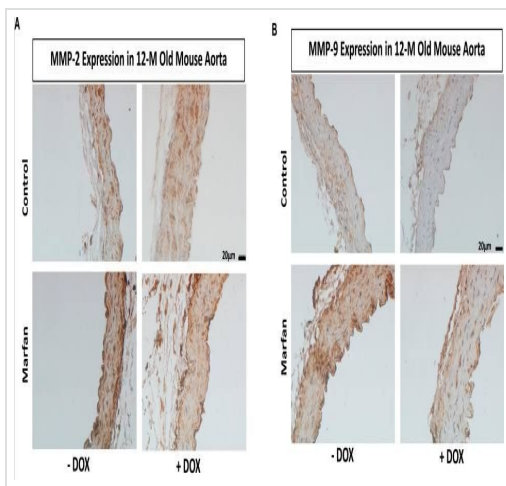
“应用说明”部分 下显示的仅为推荐的起始稀释度;实际最佳的稀释度/浓度应由使用者检定。

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

图片



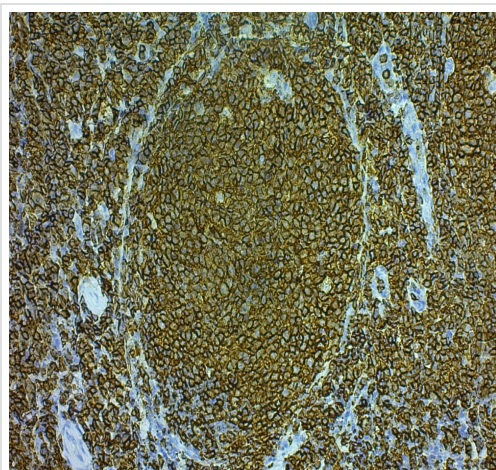
Effect of doxycycline (DOX) treatment on the expression of MMP-2 and MMP-9 in the skin. (A) The upper panel shows representative MMP-2 staining of skin from CTRL and MFS mice at 12 months of age with or without DOX, and the lower panel shows bar graphs representing the expression of MMP-2 in the skin of CTRL and MFS mice at 12 weeks of age with or without DOX. (B) The upper panel shows representative MMP-9 staining of skin from CTRL and MFS mice at 12 months of age with or without DOX treatment. The lower panel shows bar graphs representing the expression of MMP-9 in the skin of CTRL and MFS mice at 12 weeks of age treated with or without DOX.



Effect of doxycycline (DOX) treatment on the expression of MMP-2 and MMP-9

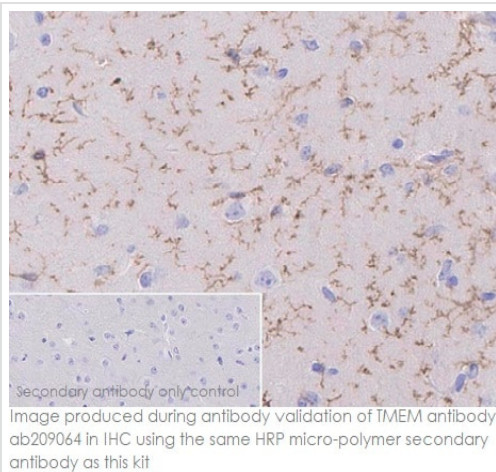
Image from Arash et al., Sci Rep; 10(1):7154. doi: 10.1038/s41598-020-64071-8.

Effect of doxycycline (DOX) treatment on the expression of MMP-2 and MMP-9 in the aorta. Immunohistochemistry staining images and bar graphs representing the expression of (A) MMP-2 and (B) MMP-9 in the aortic roots of CTRL and MFS mice at 12 weeks of age treated with or without DOX.



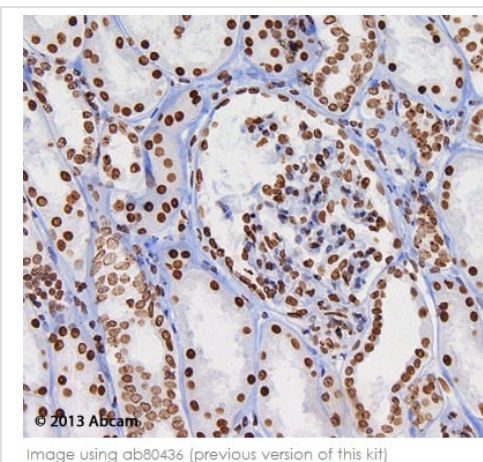
ab236466 Mouse and Rabbit Specific HRP-DAB Detection IHC Kit on Human Tonsil staining CD20

Immunohistochemistry (Formalin/PFA-fixed paraffin-embedded sections) - Mouse and Rabbit Specific HRP/DAB IHC Detection Kit - Micro-polymer (ab236466)



Immunohistochemistry (Formalin/PFA-fixed paraffin-embedded sections)

**ab209064** at 1:2000 staining TMEM119 antibody in mouse cerebrum tissue by immunohistochemistry (FFPE). Immunohistochemical analysis of paraffin-embedded mouse cerebrum tissue labeling TMEM119 with **ab209064** at 1/2000 dilution followed by the same HRP micro-polymer goat anti-rabbit secondary antibody used in this kit. Positive staining on glial cells in mouse cerebrum is observed. Perform heat mediated antigen retrieval using **ab93684** (Tris/EDTA buffer, pH 9.0) before commencing with IHC staining protocol. Counter stained with hematoxylin.



Immunohistochemistry (Formalin/PFA-fixed paraffin-embedded sections)

**ab1220** staining human kidney sections by IHC-P using EXPOSE IHC detection kit **ab80436** (this kit is identical to **ab80436**, except that it uses the same HRP polymer secondary antibody used by Abcam's R&D team for antibody validation). Formalin fixed paraffin embedded tissue sections were pre-treated using heat mediated antigen retrieval (using a pressure cooker) with sodium citrate buffer (pH 6) for 30 minutes. The section was incubated with **ab1220**, 5 µg/ml, for 1 hour at room temperature. DAB was used as the chromogen and the section was counterstained with haematoxylin and mounted with DPX.

**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