

## Product datasheet

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

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

### Overview

---

|                         |   |
|-------------------------|---|
| <b>Product name</b>     | Mouse and Rabbit Specific HRP/DAB IHC Detection Kit - Micro-polymer   |
| <b>Detection method</b> | Colorimetric  |
| <b>Assay type</b>       | Micro-Polymer   |
| <b>Product overview</b> | 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.

|              |   |
|--------------|---|
| <b>Notes</b> | 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](#).

|                            |                            |
|----------------------------|----------------------------|
| <b>Tested applications</b> | <b>Suitable for:</b> IHC-P |
|----------------------------|----------------------------|

### Properties

---

|                             |   |
|-----------------------------|---|
| <b>Storage instructions</b> | Store at +4°C. Please refer to protocols. |
|-----------------------------|---|

| Components                            | 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 |

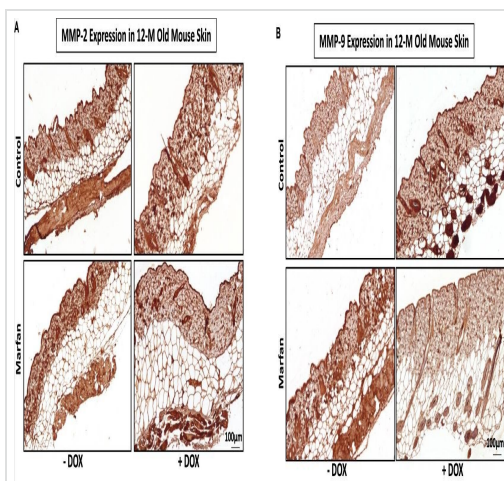
**Relevance** HRP/DAB detection kit is a immunoenzymatic antigen detection system for immunohistochemistry techniques.

## Applications

**The Abpromise guarantee** Our **Abpromise guarantee** covers the use of ab236466 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 concentration. |

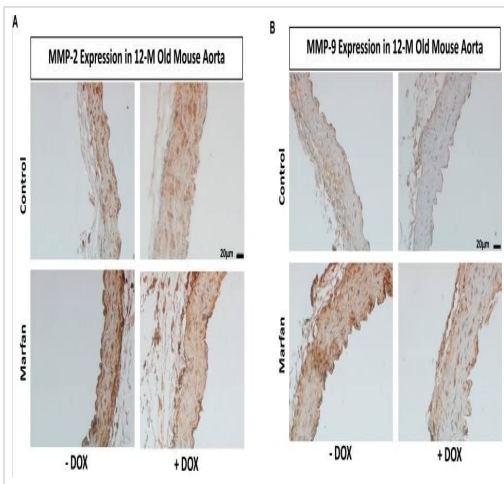
## Images



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

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

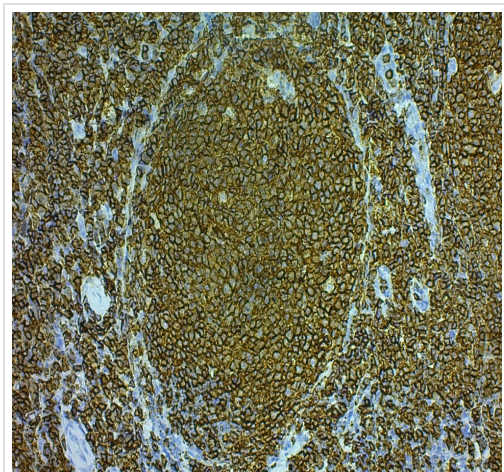
Effect of doxycycline (DOX) treatment on the expression of MMP-2 and MMP-9 in the skin. (A) The upper panel shows representative MMP-2 staining of skin from CTRL and MFS mice at 12 months of age with or without DOX, and the lower panel shows bar graphs representing the expression of MMP-2 in the skin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

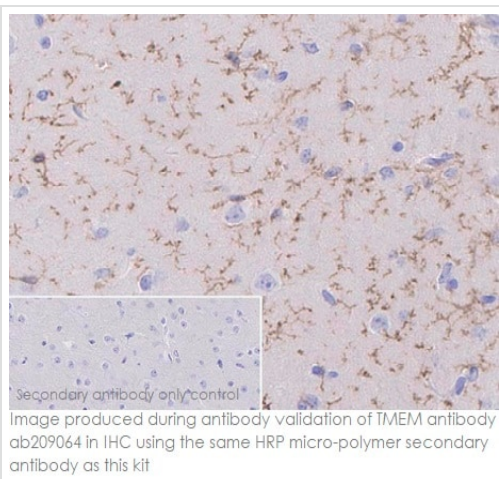
Image from Arash et al., Sci Rep; 10(1):7154. doi: 10.1038/s41598-020-64071-8.

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.



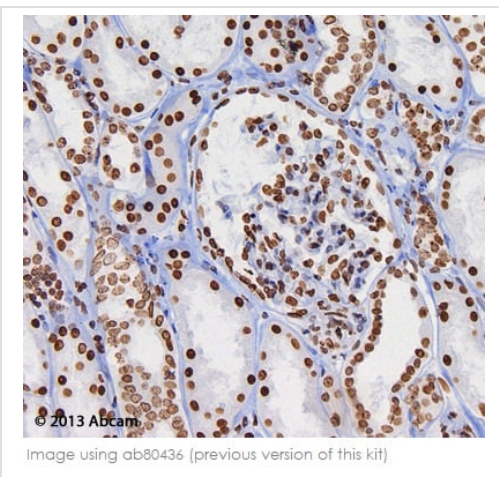
Immunohistochemistry (Formalin/PFA-fixed paraffin-embedded sections) - Mouse and Rabbit Specific HRP/DAB IHC Detection Kit - Micro-polymer (ab236466)

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



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.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