

## Product datasheet

# Rabbit specific HRP/DAB Detection IHC Detection Kit - Micro-polymer ab236469

★★★★★ [2 Abreviews](#) [15 References](#) [5 Images](#)

### Overview

<b>Product name</b>	Rabbit specific HRP/DAB Detection 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 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.

**Notes** The previous version of this kit, the EXPOSE Rabbit Specific HRP/DAB Detection IHC Kit, was used by researchers to stain over 75,000 slides since 2012 (based on volumes of reagent sold).

#### Other IHC detection products

For alternative kits, see anti-Mouse / anti-Rabbit HRP-DAB Micro-polymer IHC Kit [ab236466](#), and [HRP-AEC micro-polymer IHC kits](#).

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

**Tested applications** **Suitable for:** IHC-P

### Properties

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

Components	60 ml	15 ml	15 ml	60 ml
50x DAB Chromogen	1 x 2ml	1 x 0.5ml	1 x 0.5ml	1 x 2ml

Components	60 ml	15 ml	15 ml	60 ml
DAB substrate	1 x 60ml	1 x 15ml	1 x 15ml	1 x 60ml
Goat anti-rabbit HRP Conjugate	1 x 60ml	1 x 15ml	1 x 15ml	1 x 60ml
Hydrogen Peroxide Block	1 x 60ml	1 x 15ml	1 x 15ml	1 x 60ml
Protein block	1 x 60ml	1 x 15ml	1 x 15ml	1 x 60ml

**Relevance**

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

**Applications**

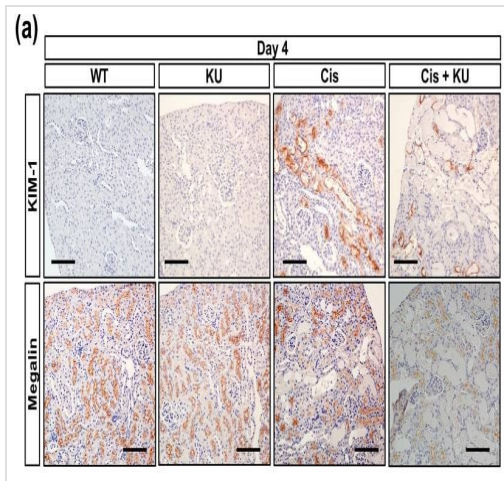
**The Abpromise guarantee**

Our **Abpromise guarantee** covers the use of ab236469 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.

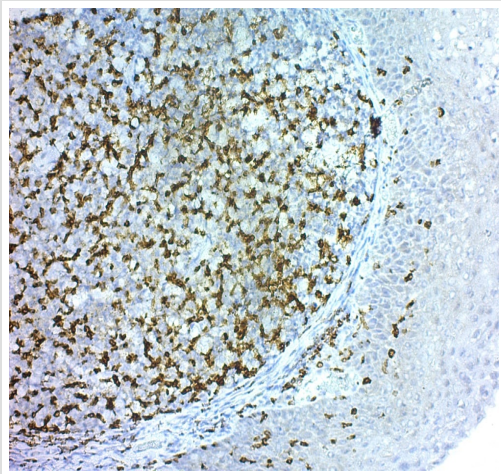
**Images**



ATM inhibition upregulated proinflammatory and profibrotic signaling in cisplatin nephropathy. a) Immunostaining of kidney sections for Kim-1 and megalin (4 days after treatment).

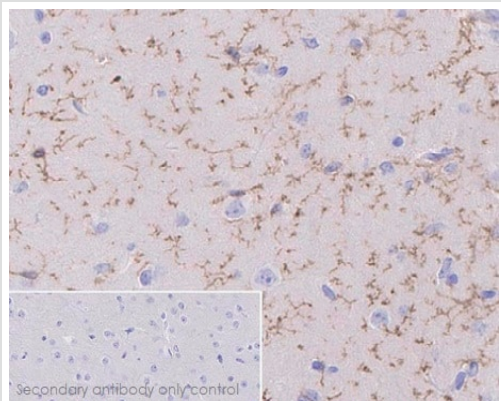
Immunohistochemistry (Formalin/PFA-fixed paraffin-embedded sections) - Rabbit specific HRP/DAB Detection IHC Detection Kit - Micro-polymer (ab236469)

Image from Uehara et al., Sci Rep., 10(1):4441; doi: 10.1038/s41598-020-61456-7. Reproduced under the Creative Commons license <https://creativecommons.org/licenses/by/4.0/>



ab236469 Rabbit Specific HRP-DAB IHC Detection Kit - Micro-Polymer on Human Tonsil staining CD3

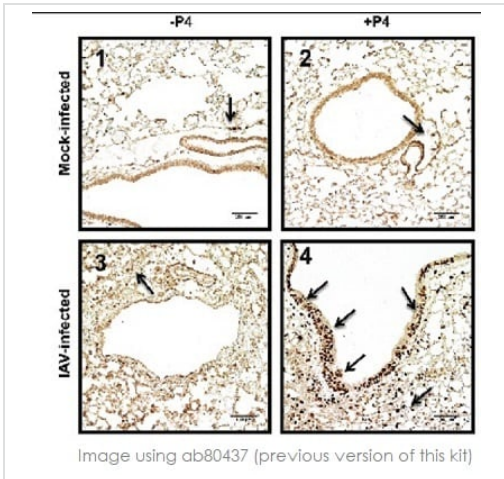
Immunohistochemistry (Formalin/PFA-fixed paraffin-embedded sections) - Rabbit specific HRP/DAB Detection IHC Detection Kit - Micro-polymer (ab236469)



**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.

Image produced during antibody validation of TMEM antibody ab209064 in IHC using the same HRP micro-polymer secondary antibody as this kit

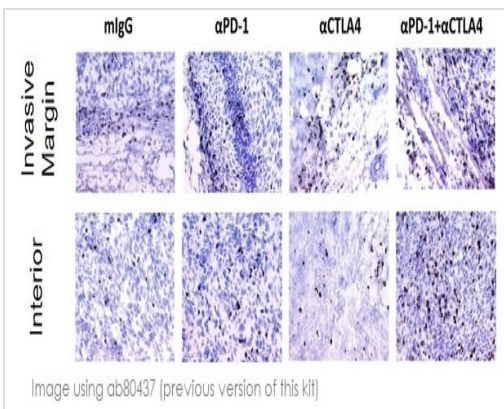
Immunohistochemistry (Formalin/PFA-fixed paraffin-embedded sections)



Immunohistochemical analysis of mouse lung tissue labeling Ki67 detected with **ab80437** (this kit is identical to **ab80437**, except that it uses the HRP micro-polymer secondary antibody used by Abcam's R&D team for antibody validation in IHC). Tissue sections were embedded in paraffin and were counterstained with hematoxylin.

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

Hall OJ et al. PLoS Pathog. 12:e1005840 (2016)  
 Reproduced under the Creative Commons license  
<http://creativecommons.org/licenses/by/4.0/>



Immunohistochemical (frozen) analysis of MC38 tumor sections labeling CD3 with **ab16669** at 1/400 dilution. Sections were fixed with acetone, treated with peroxidase block to quench endogenous peroxidase, and then further blocked with a 10% goat serum and 5% BSA solution. CD3 positive T cells were detected using **ab80437** (this kit is identical to **ab80437**, except that it uses the HRP micro-polymer antibody used by Abcam's R&D team for antibody validation in IHC). Invasive margin (top) shows sections derived from the periphery of the tumor and interior (bottom) shows sections from within the tumor.

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

Selby MJ et al. PLoS One (2016)

**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