

Product datasheet

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

★★★★★ 2 Abreviews 7 References 5 Images

Overview

Product name	Rabbit specific HRP/DAB Detection IHC Detection Kit - Micro-polymer
Detection method	Colorimetric
Assay type	Micro-Polymer
Product overview	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.

This kit is an exact copy of EXPOSE Rabbit Specific HRP/DAB Detection IHC Kit [ab80437](#), except that it uses the HRP micro-polymer antibody used during the validation of Abcam's antibodies in IHC. This HRP micro-polymer antibody was identified by our R&D team as giving the most sensitive and specific staining.

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	15 ml	60 ml
50x DAB Chromogen	1 x 0.5ml	1 x 2ml
DAB substrate	1 x 15ml	1 x 60ml
Goat anti-rabbit HRP Conjugate	1 x 15ml	1 x 60ml
Hydrogen Peroxide Block	1 x 15ml	1 x 60ml
Protein block	1 x 15ml	1 x 60ml

Relevance

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

Applications

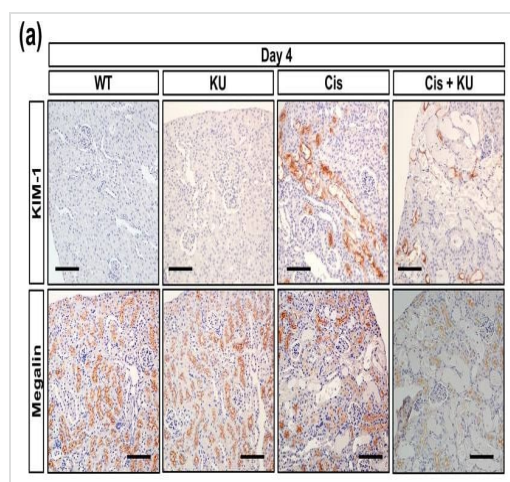
The Abpromise guarantee

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.

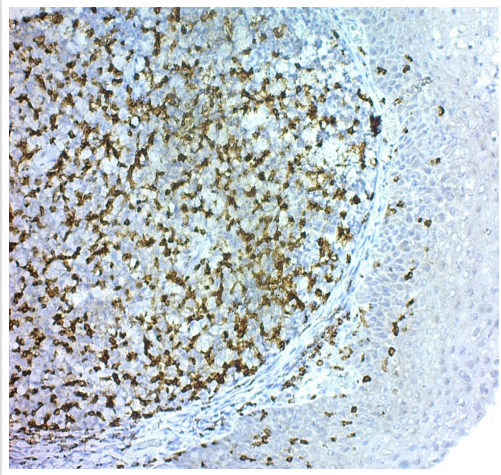
Images



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/>



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

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

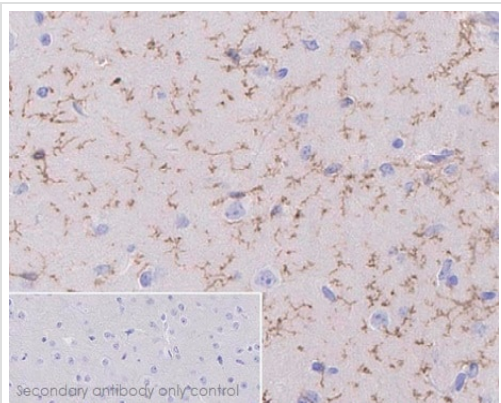
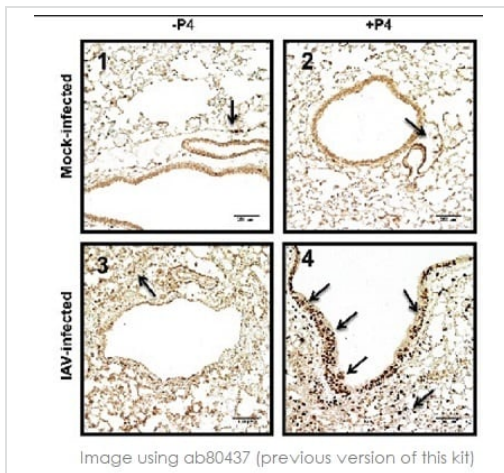


Image produced during antibody validation of TMEM antibody ab209064 in IHC using the same HRP micro-polymer secondary antibody as this kit

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

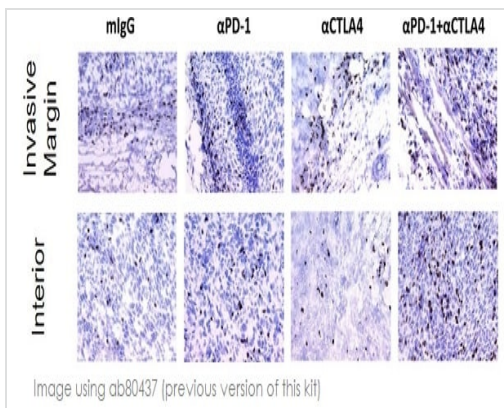
ab209064 at 1:2000 staining TMEM119 antibody in mouse cerebrum tissue by immunohistochemistry (FFPE). Immunohistochemical analysis of paraffin-embedded mouse cerebrum tissue labeling TMEM119 with ab209064 at 1/2000 dilution followed by the same HRP micro-polymer goat anti-rabbit secondary antibody used 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.



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