

## Product datasheet

# Mouse specific HRP/DAB Detection IHC Kit **ab64259**

★★★★★ [2 Abreviews](#) [42 References](#) [2 Images](#)

### Overview

<b>Product name</b>	Mouse specific HRP/DAB Detection IHC Kit
<b>Product overview</b>	<p>Mouse specific HRP/DAB IHC Detection Kit <b>ab64259</b> is a complete kit to detect a primary antibody raised in mouse, using HRP-labeled-streptavidin and a biotinylated anti-mouse secondary antibody.</p> <p>It is supplied with a full protocol and with peroxidase and protein blocking reagents, biotinylated secondary antibody, streptavidin-HRP, and DAB substrate.</p> <p>In the method:</p> <ul style="list-style-type: none"> <li>- a sample is incubated with an unconjugated primary antibody specific to the target antigen</li> <li>- a biotinylated secondary antibody then binds to the primary antibody</li> <li>- HRP-labeled streptavidin then binds to the secondary antibody</li> <li>- the HRP produces a brown colored substance at the site of primary antibody binding by reacting with DAB.</li> </ul>

<b>Notes</b>	<p>For alternative streptavidin-biotin IHC detection kits see:</p> <ul style="list-style-type: none"> <li>- Anti-Mouse / anti-Rabbit HRP-DAB IHC kit <a href="#">ab64264</a></li> <li>- Anti-Rabbit HRP-DAB IHC Kit <a href="#">ab64261</a></li> </ul> <p>For high sensitivity HRP micro-polymer IHC detection kits see:</p> <ul style="list-style-type: none"> <li>- Anti-Mouse / anti-Rabbit HRP/DAB Micro-polymer IHC Kit <a href="#">ab236466</a></li> <li>- Anti-Rabbit HRP-DAB Micro-polymer IHC Kit <a href="#">ab236469</a></li> </ul> <p>Alternatively, see <a href="#">HRP-AEC IHC detection kits</a>.</p> <p>For staining mouse tissues with mouse antibodies, use mouse-on-mouse polymer kit <a href="#">ab127055</a>.</p> <p>Find more kits and reagents for antigen retrieval, blocking, signal amplification, visualization, counterstaining, and mounting in the <a href="#">IHC kits and reagents guide</a>.</p>
--------------	---

<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
50x DAB Chromogen	1 x 0.5ml
Biotinylated goat anti-mouse IgG(H+L) (ab64255)	1 x 15ml
DAB substrate	1 x 15ml
Hydrogen Peroxide Block	1 x 15ml
Protein block	1 x 15ml
Streptavidin Peroxidase	1 x 15ml

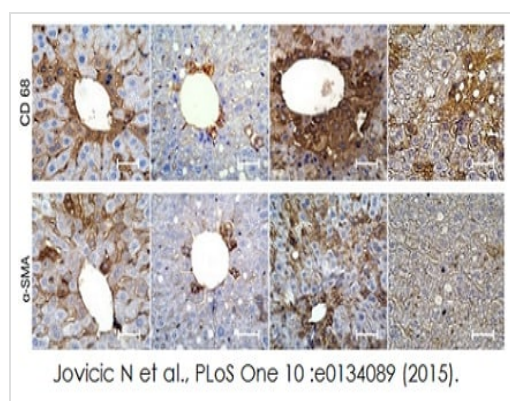
**Relevance** HRP/DAB detection kit is a immunoenzymatic antigen detection system for immunohistochemistry techniques.

## Applications

**The Abpromise guarantee** Our **Abpromise guarantee** covers the use of ab64259 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 dilution.

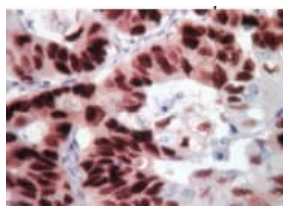
## Images



Immunohistochemical analysis staining SMA or CD68 in paraffin embedded mouse liver tissue sections Samples were incubated with anti-α- SMA (**ab7817**) or anti-CD68 (**ab49777**). A Mouse specific HRP/DAB Detection IHC Kit (AB64259 was used as a stain and Mayer's hematoxylin was used as a counterstain. Top panel of images staining CD68 and bottom panel staining α-SMA.

### Immunohistochemistry - ab64259

Jovicic N et al., PLoS One 10(7). Fig 6D & E. DOI: 10.1371/journal.pone.0134089. Reproduced under the Creative Commons license <https://creativecommons.org/licenses/by/4.0/>



Human colon carcinoma stained with anti p53 antibody using Anti Mouse HRP/DAB Detection Kit (ab64259).

Immunohistochemistry (Formalin/PFA-fixed paraffin-embedded sections) - Anti Mouse HRP/DAB Detection Kit (ab64259)

**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