

Mouse specific HRP/DAB Detection IHC Kit ab64259

★★★★★ [2 Abreviews](#) [42 References](#) [画像数 2](#)

製品の概要

製品名

Mouse specific HRP/DAB Detection IHC Kit

製品の概要

Mouse specific HRP/DAB IHC Detection Kit ab64259 is a complete kit to detect a primary antibody raised in mouse, using HRP-labeled-streptavidin and a biotinylated anti-mouse secondary antibody.

It is supplied with a full protocol and with peroxidase and protein blocking reagents, biotinylated secondary antibody, streptavidin-HRP, and DAB substrate.

In the method:

- a sample is incubated with an unconjugated primary antibody specific to the target antigen
- a biotinylated secondary antibody then binds to the primary antibody
- HRP-labeled streptavidin then binds to the secondary antibody
- the HRP produces a brown colored substance at the site of primary antibody binding by reacting with DAB.

特記事項

For alternative streptavidin-biotin IHC detection kits see:

- Anti-Mouse / anti-Rabbit HRP-DAB IHC kit [ab64264](#)
- Anti-Rabbit HRP-DAB IHC Kit [ab64261](#)

For high sensitivity HRP micro-polymer IHC detection kits see:

- Anti-Mouse / anti-Rabbit HRP/DAB Micro-polymer IHC Kit [ab236466](#)
- Anti-Rabbit HRP-DAB Micro-polymer IHC Kit [ab236469](#)

Alternatively, see [HRP-AEC IHC detection 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](#).

アプリケーション

適用あり: IHC-P

製品の特性

保存方法

Store at +4°C. Please refer to protocols.

内容	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

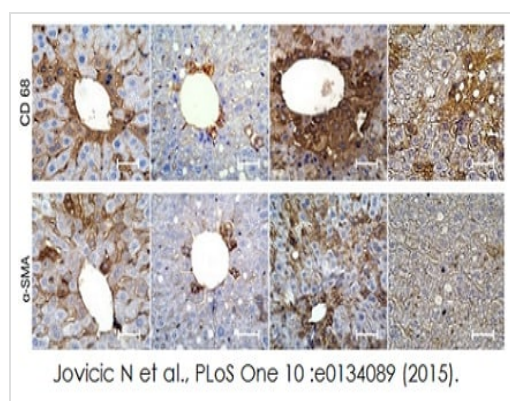
関連性 HRP/DAB detection kit is a immunoenzymatic antigen detection system for immunohistochemistry techniques.

アプリケーション

The Abpromise guarantee Abpromise保証は、次のテスト済みアプリケーションにおけるab64259の使用に適用されます
 アプリケーションノートには、推奨の開始希釈率がありますが、適切な希釈率につきましてはご検討ください。

アプリケーション	Abreviews	特記事項
IHC-P		Use at an assay dependent dilution.

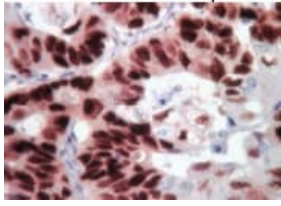
画像



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.co.jp/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors