

Product datasheet

Rabbit specific HRP/DAB Detection IHC Kit ab64261

★★★★★ [5 Abreviews](#) [162 References](#) [3 Images](#)

Overview

Product name	Rabbit specific HRP/DAB Detection IHC Kit
Product overview	<p>Rabbit specific HRP/DAB IHC Detection Kit ab64261 is a complete kit to detect a primary antibody raised in rabbit, using HRP-labeled-streptavidin and a biotinylated anti-rabbit 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">- 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.
Notes	<p>For alternative streptavidin-biotin IHC detection kits see:</p> <ul style="list-style-type: none">- Anti-Mouse / anti-Rabbit HRP-DAB IHC kit ab64264- Anti-Mouse HRP-DAB IHC Kit ab64259 <p>For high sensitivity HRP micro-polymer IHC detection kits see:</p> <ul style="list-style-type: none">- Anti-Mouse / anti-Rabbit HRP/DAB Micro-polymer IHC Kit ab236466- Anti-Rabbit HRP-DAB Micro-polymer IHC Kit ab236469 <p>Alternatively, see HRP-AEC IHC detection kits.</p> <p>Find more kits and reagents for antigen retrieval, blocking, signal amplification, visualization, counterstaining, and mounting in the IHC kits and reagents guide.</p>
Tested applications	Suitable for: IHC-P, IHC-Fr

Properties

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

Components	15 ml	15 ml

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

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

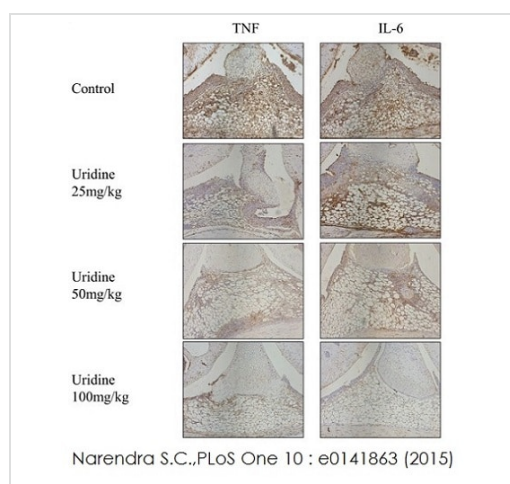
Applications

The Abpromise guarantee Our **Abpromise guarantee** covers the use of ab64261 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.
IHC-Fr		Use at an assay dependent concentration.

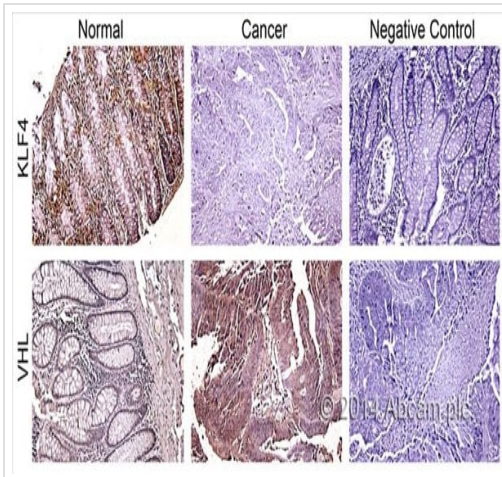
Images



Immunohistochemical analysis staining TNF or IL6 in mouse tissue sections. Samples are incubated in murine TNF ([ab9739](#)) or IL6 primary antibody. Bound primary antibodies were detected using IHC detection kit ([ab64261](#)). Immunohistochemistry was used to determine the effect of uridine on synovial expression of TNF and IL6.

Immunohistochemistry - ab64261

Image from Narendra S.C., PLoS One 10(10), Fig 5a.
doi: 10.1371/journal.pone.0141863 Reproduced under the Creative Commons license
<http://creativecommons.org/licenses/by/4.0/>



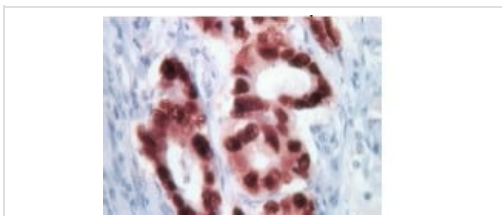
Immunohistochemistry (Formalin/PFA-fixed paraffin-embedded sections) - Anti Rabbit HRP/DAB

Detection Kit (ab64261)

This image is courtesy of an anonymous Abreview

Immunohistochemical analysis of KLF4 and pVHL expression in colon cancer tissues. Sections from colon cancer and adjacent tissues were analyzed by immunostaining against KLF4 and pVHL. Staining without primary antibody served as negative control.

Tissue sections were de-waxed with xylene and rehydrated through gradient ethanol into water. For antigen retrieval, sections were heated in citrate buffer (pH6.0) for 10 min at 95°C in a microwave oven. After cooling to room temperature, the sections were then digested with 0.05% trypsin for 10 min at 37°C. Endogenous peroxidase activity was quenched with 0.3% H₂O₂ in methanol for 30 min at room temperature. After PBS washes, nonspecific antibody binding was blocked by pre-incubating slides with 10% normal goat non-immune serum at 37° for 30 min. After blotting off the blocking serum, sections were incubated with primary antibody against LKF4 as well as primary antibody against VHL at 4° overnight. After PBS washes again, sections were incubated with biotinylated secondary antibody at 1:200 dilution for 30 min at room temperature. After incubating with Vectastain ABC reagent (Vector Laboratories, Inc., Burlingame, CA) for 30 min at room temperature, the sections were developed with Anti Rabbit HRP/DAB Detection Kit (ab64261). Sections were washed in running tap water and lightly counterstained with hematoxylin, followed by dehydration and coverslip mounting. Negative controls were obtained by omitting the primary antibody.



Immunohistochemistry (Formalin/PFA-fixed paraffin-embedded sections) - Anti Rabbit HRP/DAB

Detection Kit (ab64261)

Human colon carcinoma fixed in 10% NBF for 24 hrs and stained with anti p53 antibody using Anti Rabbit HRP/DAB Detection Kit (ab64261).

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