

Product datasheet

Mouse Specific HRP/AEC (ABC) Detection IHC kit ab64258

[1 Image](#)

Overview

Product name Mouse Specific HRP/AEC (ABC) Detection IHC kit

Product overview Mouse specific HRP/AEC (ABC) IHC Detection Kit ab64258 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 AEC 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 colored substance at the site of primary antibody binding by reacting with AEC.

Notes

For alternative AEC streptavidin-biotin IHC detection kits see:

- anti-Mouse / anti-Rabbit HRP-AEC IHC kit [ab93705](#)
- Anti-Rabbit HRP-AEC IHC Kit [ab64260](#)

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

- Anti-Mouse / anti-Rabbit HRP-AEC Micro-polymer IHC Kit [ab236467](#)
- Anti-Rabbit HRP-AEC Micro-polymer IHC Kit [ab236468](#)

Alternatively, see [HRP-DAB IHC detection 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
AEC Single Solution	1 x 15ml
Biotinylated goat anti-mouse IgG(H+L) (ab64255)	1 x 15ml
Hydrogen Peroxide Block	1 x 15ml
Protein block	1 x 15ml
Streptavidin Peroxidase	1 x 15ml

Relevance

An HRP/AEC Detection Kit is a labeled immunoenzymatic antigen detection system. Horseradish peroxidase (HRP) readily combines with hydrogen peroxide (H₂O₂) and the resultant complex can oxidize a wide variety of chromogenic substrates. AEC is 3-amino-9-ethyl-carbazole, a chromogenic substrate.

Applications

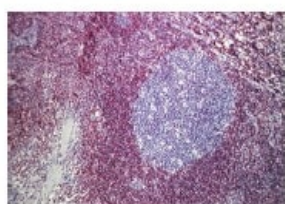
The Abpromise guarantee

Our **Abpromise guarantee** covers the use of ab64258 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 of Bcl-2 expression in Human tonsil, using ab64258 Detection Kit.

Immunohistochemistry (Formalin/PFA-fixed paraffin-embedded sections) - Anti-Mouse/HRP/AEC Detection Kit (ab64258)

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