

Product datasheet

Mouse and Rabbit Specific HRP Plus (ABC) Detection IHC Kit (Ready To Use) ab93697

6 References 1 Image

Overview

Product name	Mouse and Rabbit Specific HRP Plus (ABC) Detection IHC Kit (Ready To Use)
Product overview	The reagents in this kit constitute a labeled streptavidin-biotin immunoenzymatic antigen detection system with increased sensitivity. This technique involves the sequential incubation of the specimen with an unconjugated primary antibody specific to the target antigen, a biotinylated secondary antibody which reacts with the primary antibody, enzyme-labeled streptavidin, and substrate-chromogen.
	Reacts with Mouse and Rabbit IgG (H+L)
Tested applications	Suitable for: IHC-P

Properties

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

Components	60 ml	125 ml
Biotinylated Goat Anti-Polyvalent Plus	1 x 60ml	1 x 125ml
Protein block	1 x 60ml	1 x 125ml
Streptavidin Peroxidase Plus	1 x 60ml	1 x 125ml

Relevance Polyvalent HRP is a labeled streptavidin-biotin 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.

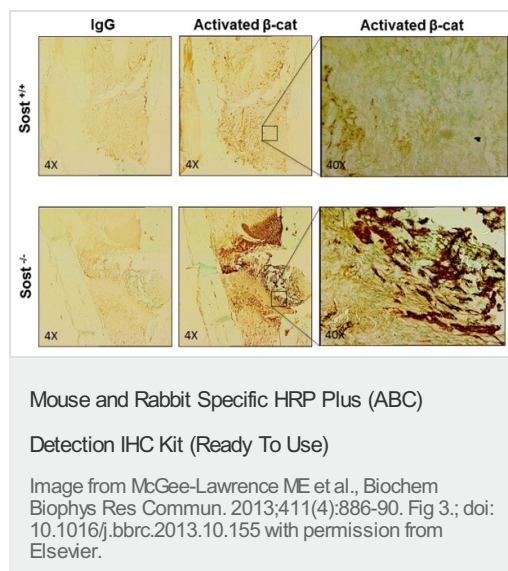
Applications

Our [Abpromise guarantee](#) covers the use of **ab93697** 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



Immunohistochemistry analysis of mouse femur tissue sections. Femurs were fixed in 10% neutral buffered formalin and decalcified for 7 days in 15% EDTA; complete decalcification was confirmed by X-ray. Decalcified bones were embedded in paraffin and sectioned longitudinally through the defect at a thickness of 8 microns. Immunohistochemical staining was performed with monoclonal anti-activated beta-catenin (un-phosphorylated on Ser33/37 and Thr41), or an IgG isotype control. Chromogens were developed using ab93697, Mouse and Rabbit Specific HRP Plus (ABC) Detection IHC Kit (Ready to Use), followed by 3'-3'-diaminobenzidine (DAB), and then counterstained with fast green. Beta-catenin levels are elevated in defects of *Sost*^{-/-} mice.

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