

# EXPOSE Rabbit Specific AP (red) Detection IHC Kit ab94737

**概述**

<b>产品名称</b>	EXPOSE兔Specific AP (red) Detection IHC试剂盒
<b>检测方法</b>	Colorimetric
<b>检测类型</b>	Micro-Polymer
<b>产品概述</b>	This detection kit offers a red alternative to the brown of the DAB product for antigen detection. This technique involves the sequential incubation of the specimen with an unconjugated rabbit primary antibody specific to the target antigen, a secondary antibody- AP conjugate which reacts with the primary antibody, and AP-compatible chromogen (Fast Red), so that a detectable signal is generated from converting Naphthol Phosphate substrate to a red colored product at the site where the antibody portion of the conjugate is bound to its target.
<b>说明</b>	Handle Enhancer, Naphthol Phosphate and Fast Red Chromogen with care.
<b>经测试应用</b>	<b>适用于:</b> IHC-P

**性能**

**存放说明** Store at +4°C. Please refer to protocols.

Components	
Protein block	1 x 125ml
Co-factor (Enhancer)	1 x 125ml
Naphthol Phosphate	1 x 125ml
Fast Red Chromogen	1 x 125ml
AP Conjugate GRI	1 x 125ml
Protein block	1 x 60ml
Co-factor (Enhancer)	1 x 60ml
Naphthol Phosphate	1 x 60ml
Fast Red Chromogen	1 x 60ml

Components	
AP Conjugate GRI	1 x 60ml
Protein block	1 x 15ml
Co-factor (Enhancer)	1 x 15ml
Naphthol Phosphate	1 x 15ml
Fast Red Chromogen	1 x 15ml
AP Conjugate GRI	1 x 15ml

**相关性** Alkaline phosphatase (AP) is a hydrolase enzyme responsible for removing phosphate groups from many types of molecules, including nucleotides, proteins, and alkaloids. Alkaline phosphatase conjugate is used in detection systems for immunohistochemistry techniques.

## 应用

Our [Abpromise guarantee](#) covers the use of **ab94737** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

应用	Ab评论	说明
IHC-P		Use at an assay dependent dilution.

**Please note:** All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

## 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 <http://www.abcam.cn/abpromise> or contact our technical team.

## Terms and conditions

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