

DoubleStain IHC Kit: G&Rt on human/mouse tissue (DAB & AP/Red) ab183280

Overview

Product name	DoubleStain IHC Kit: G&Rt on human/mouse tissue (DAB & AP/Red)
Assay type	Polymer Multiplex
Product overview	<p>Abcam's DoubleStain IHC kit (ab183280) is designed to use with user supplied goat and rat primary antibodies to detect two distinct antigens on human or mouse tissue or cell samples. . This kit has been tested on paraffin embedded tissue, however it can also be used on frozen specimens or freshly prepared monolayer cell smears.</p> <p>Double staining is one of most common methods used in immunohistostaining allowing the detection of two distinct antigens in a single tissue. Abcam's DoubleStain IHC Kit (ab183280) supplies the user with two polymer enzyme conjugates: HRP polymer anti-Goat IgG and AP polymer anti-Rat IgG and two distinct substrates/chromogens, DAB (brown) and Permanent Red (red). Users will sequentially apply the two enzyme conjugates onto the specimen. When two proteins are present a brown/red color will develop depending on the presence and location of the antigens the two colors should be distinct. If only the anti-goat antigen is present only the DAB brown chromogen will be seen and if only the Rat antigen is present only the Permanent Red chromogen will be visible. The Kit is a non-biotin system avoiding endogenous biotin non-specific binding.</p>
Tested applications	Suitable for: IHC-Fr, IHC-P

Properties

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

Components	18 ml	6 ml	60 ml
Aqueous Mounting Medium	1 x 18ml	1 x 7ml	1 x 70ml
DAB Chromogen (20x)	1 x 2ml	1 x 1.5ml	1 x 3.5ml
DAB Substrate	1 x 30ml	1 x 12ml	1 x 70ml
DoubleStain Blocker	1 x 18ml	1 x 6ml	1 x 60ml
Goat HRP Polymer	1 x 18ml	1 x 6ml	1 x 60ml

Components	18 ml	6 ml	60 ml
Permanent Red Activator (5x)	1 x 7.2ml	1 x 3ml	1 x 14ml
Permanent Red Chromogen (100x)	1 x 360µl	1 x 150µl	1 x 0.7ml
Permanent Red Substrate	1 x 36ml	1 x 15ml	1 x 70ml
Rat AP Polymer	1 x 18ml	1 x 6ml	1 x 60ml
Rat Primer	1 x 18ml	1 x 6ml	1 x 60ml

Applications

The Abpromise guarantee Our **Abpromise guarantee** covers the use of ab183280 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-Fr		Use at an assay dependent concentration.
IHC-P		Use at an assay dependent concentration.

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