

## Product datasheet

# DoubleStain IHC Kit: R&Rt on human/mouse tissue (Green/HRP & AP/Red) ab183285

★★★★★ [1 Abreviews](#) [2 References](#)

### Overview

<b>Product name</b>	DoubleStain IHC Kit: R&Rt on human/mouse tissue (Green/HRP & AP/Red)
<b>Assay type</b>	Polymer Multiplex
<b>Product overview</b>	Abcam's Doublestain IHC Kit (ab183285) is designed for use with user supplied rabbit and rat primary antibodies to detect two distinct antigens on human and mouse tissue or cell samples. This kit has been tested on paraffin embedded tissue, however, it can also be used to stain frozen specimens or freshly prepared monolayer cell smears.

Double staining is a common method used in immunohistostaining, allowing for the detection of two distinct antigens in a single tissue. Abcam's Doublestain IHC Kit (ab183285) supplies the user with two polymer enzyme conjugates: Rat HRP Polymer (minimal cross reaction to mouse) and Rabbit AP Polymer with two distinct substrates/chromogens, Emerald and Permanent Red. Emerald chromogen reacts with the Rat HRP polymer conjugate to produce a green color. Permanent Red reacts with the Rabbit AP polymer to produce a red color. When two proteins are co-localized a blue/ purple color will develop depending on which antigen is stronger. If only the rat antigen is present only the Emerald chromogen will be present and if the rabbit antigen is present only the Permanent Red chromogen will be present. Abcam's Doublestain IHC Kit (ab183285) is a non-biotin system avoiding the extra steps involved in blocking non-specific binding due to endogenous biotin.

**Tested applications**                      **Suitable for:** IHC-Fr, IHC-P

### Properties

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

Components	36 ml	60 ml	6 ml	120 ml
Emerald Chromogen	2 x 18ml	1 x 60ml	1 x 6ml	1 x 120ml
Non-aqueous Mounting Medium	4 x 9ml	0 x 0ml	1 x 6ml	0 x 0ml
Permanent Red Activator (5x)	1 x 7.2ml	1 x 24ml	1 x 3ml	2 x 12ml

Components	36 ml	60 ml	6 ml	120 ml
Permanent Red Chromogen (100x)	1 x 360µl	1 x 1.2ml	1 x 150µl	1 x 1.2ml
Permanent Red Substrate	2 x 18ml	1 x 120ml	1 x 15ml	1 x 120ml
Rabbit AP Polymer	1 x 18ml	1 x 60ml	1 x 6ml	1 x 60ml
Rat HRP Polymer	1 x 18ml	1 x 60ml	1 x 6ml	1 x 60ml

## Applications

**The Abpromise guarantee** Our **Abpromise guarantee** covers the use of ab183285 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