

DoubleStain CL IHC Kit: G&Rt on human/mouse tissue (HRP/Green & AP/Red) ab183282

Overview

Product name	DoubleStain CL IHC Kit: G&Rt on human/mouse tissue (HRP/Green & AP/Red)
Assay type	Polymer Multiplex
Product overview	<p>Abcam's Doublestain IHC Kit (ab183282) is designed to be used with user supplied goat and rat primary antibodies to detect two distinct antigens on human or mouse tissue or cell samples. The kit has been tested on paraffin-embedded human and mouse tissues.</p> <p>Double staining is one of most common methods used in immunohistostaining allowing for the detection of two distinct antigens in a single tissue. Abcam's Doublestain IHC Kit (ab183282) supplies the user with two polymer enzyme conjugates: Goat AP Polymer and Rat HRP Polymer with two distinct substrates/chromogens, Emerald and Permanent Red. With this protocol the goat and rat primary antibodies will be applied to the tissue at the same time but the user will sequentially apply the polymer enzyme conjugate to the specimen. When two proteins are present in the same location a blue or purple color will develop depending on the quantity of each antigen screened. If only the anti-goat antigen is present only the Permanent Red chromogen will be visible and if only the Rat antigen is present then only the Emerald chromogen will be seen. Abcam's Doublestain IHC Kit (ab183282) is a non-biotin system avoiding endogenous biotin non-specific binding.</p>
Tested applications	Suitable for: IHC-P, IHC-Fr

Properties

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

Components	18 ml	6 ml	60 ml
DoubleStain Blocker	1 x 18ml	1 x 6ml	1 x 60ml
Emerald Chromogen	1 x 18ml	1 x 6ml	1 x 70ml
Goat AP Polymer	1 x 18ml	1 x 6ml	1 x 60ml
Non-aqueous Mounting Medium	1 x 18ml	1 x 6ml	0 x 0ml

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 HRP 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 ab183282 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 concentration.
IHC-Fr		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