

Product datasheet

DoubleStain IHC Kit: M&R on human tissue (HRP/Green & AP/Red) ab210061

★★★★★ [1 Abreviews](#) [3 References](#)

Overview

Product name	DoubleStain IHC Kit: M&R on human tissue (HRP/Green & AP/Red)
Sample type	Tissue
Species reactivity	Reacts with: Human
Product overview	<p>Abcam's DoubleStain IHC Kit: M&R on human tissue (HRP/Green & AP/Red) (ab210061) is designed to detect two distinct antigens on human tissue or cell samples using a simplified protocol with user-supplied mouse and rabbit antibodies. This kit has been tested in paraffin embedded tissues. However, this kit can be used to stain frozen specimen and/or freshly prepared monolayer cell smears.</p> <p>Double staining is a common method used in immunohistochemistry to evaluate two distinct antigens in a single tissue. This kit contains a HRP-Polymer anti-Mouse IgG, AP-Polymer anti-Rabbit IgG, and two distinct chromogens, Emerald (green) and Permanent Red (red). User may apply the two enzyme conjugates onto the specimen at the same time and mix them on the slide. Simplified steps offer user much faster and quicker protocol than a sequential procedure. A second advantage is it allows the researcher to visualize when two proteins are co-localized because of the color change when the chromogens overlap that can be semi-quantitative. For example, if the area of co-localization stains blue, the antigen indicated by Emerald is expressed at higher concentration in the cell and if the color is purple, the antigen indicated by Permanent-Red is expressed at higher concentrations DoubleStain IHC Kit: M&R on human tissue (HRP/Green & AP/Red) is a non-biotin system that avoids endogenous biotin non-specific binding.</p> <p>The Organic Mounting Medium is not supplied with the 120 mL kit.</p> <p>The Permanent Red working solution is not stable. It must be used (freshly) immediately after it is made.</p>
Tested applications	Suitable for: IHC

Properties

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

Components	18 ml	6 ml	60 ml	12 ml
Emerald Chromogen	1 x 18ml	1 x 6ml	1 x 60ml	1 x 15ml
Mouse HRP Polymer	1 x 18ml	1 x 6ml	1 x 60ml	1 x 6ml
Organic Mounting Medium	1 x 18ml	1 x 6ml	0 x 0ml	1 x 12ml
Permanent Red Activator (5X)	1 x 7.2ml	1 x 3ml	1 x 24ml	1 x 3ml
Permanent Red Chromogen (100X)	1 x 360µl	1 x 150µl	1 x 1.2ml	1 x 150µl
Permanent Red Substrate (RTU)	1 x 36ml	1 x 15ml	1 x 120ml	1 x 15ml
Rabbit AP polymer	1 x 18ml	1 x 6ml	1 x 60ml	1 x 6ml

Applications

The Abpromise guarantee Our **Abpromise guarantee** covers the use of ab210061 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		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