

## Product datasheet

# DoubleStain IHC Kit: M&R on human tissue (DAB & AP/Red) ab210059

★★★★★ [1 Abreviews](#) [10 References](#)

### Overview

<b>Product name</b>	DoubleStain IHC Kit: M&R on human tissue (DAB & AP/Red)
<b>Sample type</b>	Tissue
<b>Species reactivity</b>	<b>Reacts with:</b> Human
<b>Product overview</b>	<p>Abcam's DoubleStain IHC Kit: M&amp;R on human tissue (DAB &amp; AP/Red) (ab210059) 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>

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, an AP-Polymer anti-Rabbit IgG, and two distinct chromogens. The DAB chromogen (brown color) is used with the HRP-Polymer anti-Mouse IgG and Permanent Red (red color) is used with the AP-Polymer anti-Rabbit IgG. DoubleStain IHC Kit: M&R on human tissue (DAB & AP/Red) is a non-biotin system that avoids the extra steps involved in blocking non-specific binding due to endogenous biotin.

**Tested applications** **Suitable for:** IHC

### Properties

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

Components	36 ml	60 ml	6 ml	12 ml
Aqueous Mounting Medium	2 x 18ml	1 x 60ml	1 x 6ml	1 x 12ml
DAB Chromogen (20X)	1 x 2ml	1 x 6ml	1 x 1.5ml	1 x 1.5ml
DAB Substrate	2 x 18ml	1 x 120ml	1 x 12ml	1 x 12ml
Mouse HRP polymer	1 x 18ml	1 x 60ml	1 x 6ml	1 x 6ml
Permanent Red Activator (5X)	1 x 7.2ml	1 x 24ml	1 x 3ml	1 x 3ml

Components	36 ml	60 ml	6 ml	12 ml
Permanent Red Chromogen (100X)	1 x 360µl	1 x 1.2ml	1 x 150µl	1 x 150µl
Permanent Red Substrate (RTU)	2 x 18ml	1 x 120ml	1 x 15ml	1 x 15ml
Rabbit AP polymer	1 x 18ml	1 x 60ml	1 x 6ml	1 x 6ml

## Applications

**The Abpromise guarantee** Our **Abpromise guarantee** covers the use of ab210059 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