

## Product datasheet

# TripleStain IHC Kit: M&R&Rt on rodent tissue (DAB, HRP/Green & AP/Red) ab183297

★★★★★ 1 Abreviews

### Overview

<b>Product name</b>	TripleStain IHC Kit: M&R&Rt on rodent tissue (DAB, HRP/Green & AP/Red)
<b>Assay type</b>	Polymer Multiplex
<b>Product overview</b>	The TripleStain IHC kit (ab183297) is designed to be used with user supplied mouse, rat and rabbit primary antibodies to detect three distinct antigens on mouse tissue and cell samples or human tissue. This kit has been tested on paraffin embedded tissue, however it can also be used on frozen tissue or cell smears.

Triple staining is a common method used in immunohistochemistry for the detection of three distinct antigens and their co-expression in a single tissue. Abcam's TripleStain IHC Kit supplies polymer enzyme conjugates: polymer-HRP anti-mouse IgG, polymer-AP anti-rat IgG, and polymer-HRP anti-rabbit IgG with three substrates/chromogens; DAB (brown), Emerald (green), and Permanent Red (Red). The TripleStain Kit is a non-biotin system avoiding non-specific background due to endogenous biotin binding. This kit has been optimized to have no cross detection when detecting two primary antibodies from the mouse and rat host species using a unique blocking system. Simplified steps allow users to complete triple staining within 5 hours (without antigen retrieval) or 6-7 hours (with antigen retrieval). This protocol also includes a method to dehydrate, clear and permanently mount slides with coverslip.

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

### Properties

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

Components	Identifier	24 ml	72 ml	240 ml
—				
Antibody Blocker (40x)		2 x 15ml	1 x 50ml	1 x 100ml
DAB Chromogen (20x)		1 x 1.5ml	1 x 2ml	1 x 6ml
DAB Substrate		1 x 15ml	2 x 18ml	1 x 120ml

Components	Identifier	24 ml	72 ml	240 ml
Emerald Chromogen		1 x 15ml	2 x 18ml	1 x 120ml
Mouse HRP Polymer		1 x 12ml	2 x 18ml	1 x 120ml
Non-aqueous Mounting Medium		1 x 12ml	2 x 18ml	
Permanent Red Activator (5x)		1 x 3ml	1 x 7.2ml	2 x 12ml
Permanent Red Chromogen (100x)		1 x 150µl	1 x 360µl	1 x 1.2ml
Permanent Red Substrate		1 x 15ml	2 x 18ml	1 x 120ml
Plastic pipette		1 unit	1 unit	1 unit
Rabbit HRP Polymer		1 x 6ml	1 x 18ml	1 x 60ml
Rat AP Polymer		1 x 6ml	1 x 18ml	1 x 60ml
Rat Primer		1 x 12ml	2 x 18ml	1 x 120ml

## Applications

Our [Abpromise guarantee](#) covers the use of **ab183297** 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