

Product datasheet

TripleStain IHC kit: M&M&R on rodent tissue (DAB, HRP/Green & AP/Red) ab183292

Overview

Product name	TripleStain IHC kit: M&M&R on rodent tissue (DAB, HRP/Green & AP/Red)
Assay type	Polymer Multiplex
Product overview	The TripleStain IHC kit (ab183292) is designed to use with two mouse antibodies and one rabbit antibody (user supplied) to detect three distinct antigens on a single mouse/rat tissue or cell sample. This kit has been tested in paraffin-embedded formalin-fixed tissues. However, it can be used on frozen specimens and freshly prepared monolayer cell smears.

Triple staining is a common method used in immunohistochemistry for the detection of three distinct antigens in a single tissue. Abcam's TripleStain IHC Kit supplies polymer enzyme conjugates: polymer-HRP anti-mouse IgG, polymer-AP anti-mouse IgG, and polymer-HRP anti-rabbit IgG with three substrates/chromogens; DAB (brown), Emerald (green), and Permanent Red (Red). DAB (Brown) and Permanent Red (red). The TripleStain Kit is a non-biotin system avoiding non-specific background due to endogenous biotin binding. A Primer step is used to increase specificity of antibody staining. This kit has been optimized to have no cross detection when detecting two primary antibodies from the same host species using a unique blocking system.

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

Properties

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

Components	72 ml	60 ml	6 ml
Antibody Blocker (40x)	1 x 50ml	1 x 100ml	1 x 30ml
Blocker A	2 x 18ml	1 x 60ml	1 x 6ml
Blocker B	2 x 18ml	1 x 60ml	1 x 6ml
DAB Chromogen (20x)	1 x 2ml	1 x 6ml	1 x 1.5ml
DAB Substrate	2 x 18ml	1 x 120ml	1 x 12ml

Components	72 ml	60 ml	6 ml
Emerald Chromogen	2 x 18ml	1 x 60ml	1 x 6ml
Mouse AP Polymer	1 x 18ml	1 x 60ml	1 x 6ml
Mouse HRP Polymer	2 x 18ml	1 x 60ml	1 x 6ml
Mouse Primer	2 x 18ml	1 x 120ml	1 x 6ml
Non-aqueous Mounting Medium	2 x 18ml	0 x 0ml	1 x 6ml
Permanent Red Activator (5x)	1 x 7.2ml	1 x 24ml	1 x 3ml
Permanent Red Chromogen (100x)	1 x 360µl	1 x 1.2ml	1 x 150µl
Permanent Red Substrate	2 x 18ml	1 x 120ml	1 x 15ml
Rabbit HRP Polymer	1 x 18ml	1 x 60ml	1 x 6ml

Applications

The Abpromise guarantee Our **Abpromise guarantee** covers the use of ab183292 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