

Product datasheet

TripleStain IHC Kit: M&R&Rt on rodent tissue (DAB, DAB/Ni & AP/Red) ab183298

1 References

Overview

Product name	TripleStain IHC Kit: M&R&Rt on rodent tissue (DAB, DAB/Ni & AP/Red)
Assay type	Polymer Multiplex
Product overview	The TripleStain IHC Kit (ab183298) 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 (**ab183198**) supplies polymer enzyme conjugates: mouse HRP polymer, Rabbit HRP polymer, and Rat AP-polymer with three substrates/chromogens; DAB (brown), DAB-Ni (black), 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 the user to complete triple staining within 5 hours (without antigen retrieval) or 6 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	60 ml	6 ml	18 ml
Antibody Blocker (40x)	1 x 100ml	1 x 30ml	1 x 50ml
Aqueous Mounting Medium	1 x 60ml	1 x 6ml	1 x 18ml
DAB Chromogen (20x)	1 x 12ml	1 x 2ml	1 x 4ml
DAB Substrate	1 x 120ml	1 x 12ml	1 x 36ml

Components	60 ml	6 ml	18 ml
DAB-Ni Substrate (20x)	1 x 6ml	1 x 1ml	1 x 2ml
Hydrogen Peroxide (20x)	1 x 6ml	1 x 1ml	1 x 2ml
Mouse HRP Polymer	1 x 60ml	1 x 6ml	1 x 18ml
Nickel Solution (7x)	1 x 18ml	1 x 3ml	1 x 6ml
Permanent Red Activator (5x)	1 x 24ml	1 x 3ml	1 x 7.2ml
Permanent Red Chromogen (100x)	1 x 1.2ml	1 x 150µl	1 x 360µl
Permanent Red Substrate	1 x 120ml	1 x 15ml	1 x 36ml
Rabbit HRP Polymer	1 x 60ml	1 x 6ml	1 x 18ml
Rat AP Polymer	1 x 60ml	1 x 6ml	1 x 18ml
Rat Primer	1 x 60ml	1 x 6ml	1 x 18ml

Applications

The Abpromise guarantee Our **Abpromise guarantee** covers the use of ab183298 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