

TripleStain IHC Kit: R&R&M on human tissue (DAB, AP/Red & Green/HRP) ab183288

Overview

Product name	TripleStain IHC Kit: R&R&M on human tissue (DAB, AP/Red & Green/HRP)
Assay type	Polymer Multiplex
Product overview	Abcam's TripleStain IHC Kit (ab183288) is designed to be used with one mouse and two rabbit primary antibodies, supplied by the user, to detect three distinct antigens on human tissue or cell samples. This kit has been tested on paraffin embedded tissue specimens, however, it can also be used for frozen tissue or cell smears.

Triple staining uses traditional and non-traditional methods in immunohistostaining to reveal three distinct antigens and their co-expression on a single tissue. Abcam's TripleStain IHC Kit (ab183288) supplies polymer enzyme conjugates: Mouse HRP Polymer, Rabbit AP Polymer and Rabbit HRP Polymer with three chromogens, DAB (brown); Permanent Red (red); and Emerald (green). Abcam's TripleStain IHC Kit (ab183288) is a non-biotin system, avoiding non-specific binding caused by endogenous biotin. This kit has been optimized to have no cross detection when detecting two primary antibodies from the same host species using our unique blocking system. Simplified steps allow users to complete triple staining within 5 hours (without antigen retrieval) or 6 hours (with antigen retrieval). The well tested protocol provides the user with the ability to permanently mount slides with a coverslip.

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

Properties

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

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

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

Applications

The Abpromise guarantee Our **Abpromise guarantee** covers the use of ab183288 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