

Product datasheet

TripleStain IHC Kit: M&M&R on Human tissue (DAB, AP/Red & HRP/Green) ab183286

★★★★★ [1 Abreviews](#) [5 References](#) [1 Image](#)

Overview

Product name	TripleStain IHC Kit: M&M&R on Human tissue (DAB, AP/Red & HRP/Green)
Assay type	Polymer Multiplex
Product overview	<p>Abcam's TripleStain IHC kit (ab183286) is designed to use with user supplied two mouse primary antibodies and one rabbit primary antibody to detect three distinct antigens on a single human tissue or cell samples. This kit has been tested on paraffin embedded tissue only; however it may also be used on frozen or freshly prepared monolayer cell smears.</p> <p>Triple staining uses traditional and non-traditional methods in immunohistostaining to reveal three distinct antigens and their co-expression on a single tissue. This kit supplies polymer enzyme conjugates: Rabbit HRP Polymer, Mouse HRP Polymer and Mouse AP Polymer with three chromogens, DAB Chromogen (brown); Permanent Red Chromogen (red); and Emerald Chromogen (green). Abcam's TripleStain IHC kit (ab183286) 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.</p>
Tested applications	Suitable for: IHC-P

Properties

Storage instructionsStore at +4°C. Please refer to protocols.

Components	60 ml	6 ml	18 ml	24 ml
Antibody Blocker (40x)	1 x 100ml	1 x 30ml	1 x 50ml	1 x 30ml
Blocker A	1 x 60ml	1 x 6ml	1 x 18ml	1 x 12ml
Blocker B	1 x 60ml	1 x 6ml	1 x 18ml	1 x 12ml
DAB Chromogen (20x)	1 x 6ml	1 x 1.5ml	1 x 2ml	1 x 1.5ml

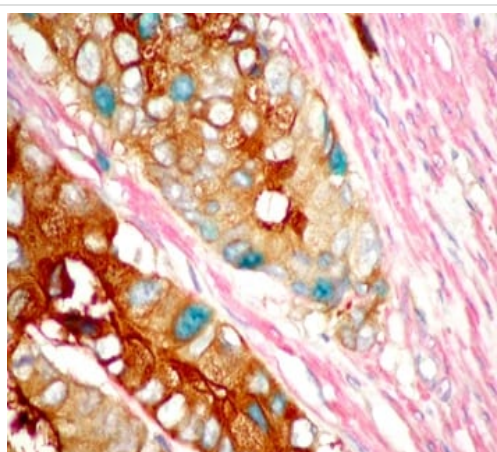
Components	60 ml	6 ml	18 ml	24 ml
DAB Substrate	1 x 120ml	1 x 15ml	1 x 36ml	1 x 15ml
Emerald Chromogen	1 x 60ml	1 x 6ml	1 x 18ml	1 x 15ml
Mouse AP Polymer	1 x 60ml	1 x 6ml	1 x 18ml	1 x 6ml
Mouse HRP Polymer	1 x 60ml	1 x 6ml	1 x 18ml	1 x 12ml
Non-aqueous Mounting Medium	0 x 0ml	1 x 6ml	1 x 18ml	1 x 3ml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

Applications

The Abpromise guarantee Our **Abpromise guarantee** covers the use of ab183286 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-P		Use at an assay dependent concentration.

Images



Immunohistochemistry (Paraffin-embedded sections)

Immunohistochemical analysis of paraffin embedded human colon cancer tissue. Mouse anti-smooth muscle actin (red), rabbit anti-CEA (brown), mouse anti-PCNA (emerald).

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