

Product datasheet

TripleStain IHC Kit: M&R&G on human tissue (DAB, AP/Red & HRP/Green) ab183290

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

Overview

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

Assay type Polymer Multiplex

Product overview Abcam's TripleStain IHC Kit (ab183290) is designed to be used with user-supplied one goat, one mouse and one rabbit primary antibodies to detect three distinct antigens on a single human tissue or cell samples. This kit has been tested on paraffin embedded tissue; however it may also be used on frozen or freshly prepared monolayer 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 (ab183290) supplies polymer enzyme conjugates: Goat HRP Polymer, Rabbit AP Polymer and Mouse HRP Polymer with three chromogens, DAB (brown); Permanent Red (red); and Emerald (green). Abcam's TripleStain IHC Kit (ab183290) is a non-biotin system, avoiding non-specific binding caused by endogenous biotin. This kit has been optimized to have no cross reaction when detecting three different primary antibodies using our unique blocking system. Simplified steps allow the user to complete triple staining within 3 hours (without antigen retrieval) or 4 hours (with antigen retrieval). The well tested protocol allows the user to permanently mount slides with a coverslip.

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

Properties

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

Components	60 ml	18 ml	6 ml	72 ml	240 ml
DAB Chromogen (20x)	1 x 6ml	1 x 2ml	1 x 1.5ml	1 x 2ml	1 x 6ml
DAB Substrate	1 x 120ml	1 x 30ml	1 x 12ml	2 x 18ml	1 x 120ml
Emerald Chromogen	1 x 60ml	1 x 18ml	1 x 6ml	2 x 18ml	1 x 120ml
Goat HRP Polymer	1 x 60ml	1 x 18ml	1 x 6ml	1 x 18ml	1 x 60ml

Components	60 ml	18 ml	6 ml	72 ml	240 ml
Mouse HRP Polymer	1 x 60ml	1 x 18ml	1 x 6ml	2 x 18ml	1 x 120ml
Non-aqueous Mounting Medium	0 x 0ml	1 x 18ml	1 x 6ml	2 x 18ml	2 x 30ml
Permanent Red Activator (5x)	1 x 24ml	1 x 7.2ml	1 x 3ml	1 x 7.2ml	2 x 12ml
Permanent Red Chromogen (100x)	1 x 1.2ml	1 x 360µl	1 x 150µl	1 x 360µl	1 x 1.2ml
Permanent Red Substrate	1 x 120ml	1 x 36ml	1 x 15ml	2 x 18ml	1 x 120ml
Rabbit AP Polymer	1 x 60ml	1 x 18ml	1 x 6ml	1 x 18ml	1 x 60ml
Triple Stain Blocker	1 x 60ml	1 x 18ml	1 x 6ml	2 x 18ml	1 x 120ml

Applications

The Abpromise guarantee Our **Abpromise guarantee** covers the use of ab183290 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.
IHC-Fr		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