



Product datasheet

DoubleStain Co-Localization IHC Kit: M&M on rodent tissue (Green/HRP & AP/Red) ab183275

Overview

Product name	DoubleStain Co-Localization IHC Kit: M&M on rodent tissue (Green/HRP & AP/Red)
Assay type	Polymer Multiplex
Product overview	<p>Abcam's Doublestain IHC Kit (ab183275) is designed to use with two user supplied mouse antibodies to detect two distinct antigens on mouse and rat tissue or cell samples. The advantage of this kit series is that it will allow you to visualize when two proteins are co localized by producing a third color blue purple. Specimens can be frozen or paraffin embedded, or freshly prepared monolayer cell smears.</p> <p>Double staining is a common method used in immunohistochemistry that allows for detection of two distinct antigens in a single tissue. This kit uses an HRP or AP polymer based technology combined with a proprietary blocking buffer system that achieves ultra sensitivity with no background or cross reactivity. Abcam's Doublestain IHC Kit (ab183275) supplies the user with primer system to enhance the two polymer enzyme conjugates; Mouse HRP Polymer and Mouse AP Polymer with two distinct substrates/chromogens, Permanent Red and Emerald. Permanent Red reacts with Mouse AP Polymer conjugate to produce a red color. Emerald chromogen reacts with Mouse HRP Polymer conjugate to produce a green color. However when the chromogens are produced in the same place the color appears blue to purple in color. Abcam's Doublestain IHC Kit (ab183275) is a non-biotin system that avoids the extra steps involved in blocking non-specific binding due to endogenous biotin.</p>
Tested applications	Suitable for: IHC-P, IHC-Fr

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 125ml
Blocker A	1 x 6ml	1 x 18ml	1 x 60ml
Blocker B	1 x 6ml	1 x 18ml	1 x 60ml

Components	6 ml	18 ml	60 ml
Emerald Chromogen	1 x 6ml	1 x 18ml	1 x 60ml
Mouse AP Polymer	1 x 6ml	1 x 18ml	1 x 60ml
Mouse HRP Polymer	1 x 6ml	1 x 18ml	1 x 60ml
Mouse Primer	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 1.4ml	1 x 3.6ml	1 x 12ml
Permanent Red Chromogen (100x)	1 x 70µl	1 x 180µl	1 x 600µl
Permanent Red Substrate	1 x 15ml	1 x 36ml	1 x 120ml

Applications

The Abpromise guarantee Our **Abpromise guarantee** covers the use of ab183275 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