

Product datasheet

DoubleStain IHC Kit: M&Rt on mouse tissue (BCIP & AEC) ab183278

Overview

Product name	DoubleStain IHC Kit: M&Rt on mouse tissue (BCIP & AEC)
Assay type	Polymer Multiplex
Product overview	<p>Abcam's DoubleStain IHC Kit (ab183278) is designed to use with user supplied mouse and rat primary antibody to detect two distinct antigens on mouse tissue or cell samples. This kit has been tested on paraffin embedded tissue, however it can also be used on frozen specimens or freshly prepared monolayer cell smears. This kit is designed not to give background with most mouse strains however there may be some, especially when using frozen tissue, that require additional blocking.</p> <p>Double staining is one of most common methods used in immunohistostaining allowing the detection of two distinct antigens in a single tissue. Abcam's DoubleStain IHC Kit (ab183278) supplies the user with two polymer enzyme conjugates: Mouse HRP Polymer and Rat AP Polymer and two distinct substrates/chromogens, AEC (Red color, used with the Mouse HRP Polymer) and BCIP/NBT (Purple color, used with the Rat AP Polymer). A Primer step is used to increase specificity of antibody staining. Both enzyme conjugates are applied to the specimen at the same time and mixed on the slide. This kit offers simplified steps making for a quicker and easier protocol than those used in a sequential procedure. Abcam's DoubleStain IHC Kit (ab183278) is a non-biotin system that avoids endogenous biotin non-specific binding.</p>

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

Properties

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

Components	12 ml	36 ml	120 ml
AEC Chromogen (20x)	1 x 2ml	1 x 4ml	1 x 12ml
AEC Substrate (20x)	1 x 1ml	1 x 2ml	1 x 6ml
Aqueous Mounting Medium	1 x 15ml	2 x 18ml	1 x 120ml
Hydrogen Peroxide (20x)	1 x 1ml	1 x 2ml	1 x 6ml

Components	12 ml	36 ml	120 ml
Mouse HRP (AEC) Polymer	1 x 6ml	1 x 18ml	1 x 60ml
Mouse Primer	1 x 12ml	2 x 18ml	1 x 120ml
MRt Block A	1 x 6ml	1 x 18ml	1 x 60ml
MRt Block B	1 x 6ml	1 x 18ml	1 x 60ml
Rat AP Polymer	1 x 6ml	1 x 18ml	1 x 60ml
Ready to Use BCIP/NBT	1 x 15ml	2 x 18ml	1 x 120ml

Applications

Our [Abpromise guarantee](#) covers the use of **ab183278** 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