

Product datasheet

DoubleStain IHC Kit: M&R on human tissue (BCIP & AEC) ab210060

Overview

Product name	DoubleStain IHC Kit: M&R on human tissue (BCIP & AEC)
Sample type	Tissue
Product overview	<p>Abcam's DoubleStain IHC Kit: M&R on human tissue (BCIP & AEC) (ab210060) is designed to detect two distinct antigens on human tissue or cell samples using a simplified protocol with user-supplied mouse and rabbit antibodies. This kit has been tested in paraffin embedded tissues. However, this kit can be used to stain frozen specimen and/or freshly prepared monolayer cell smears.</p>

Double staining is a common method used in immunohistochemistry to evaluate two distinct antigens in a single tissue. This kit contains a HRP-Polymer anti-Mouse IgG, an AP-Polymer anti-Rabbit IgG, and two distinct chromogens. The AEC chromogen (red color) is used with HRP Polymer anti-Mouse IgG and the BCIP/NBT chromogen (purple/blue color) is used with AP Polymer anti-Rabbit IgG. DoubleStain IHC Kit: M&R on human tissue (BCIP & AEC) is a non-biotin system that avoids the extra steps involved in blocking non-specific binding due to endogenous biotin.

Tested applications	Suitable for: IHC
----------------------------	--------------------------

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 Buffer (20X)	1 x 1ml	1 x 2ml	1 x 6ml
Rabbit AP polymer	1 x 6ml	1 x 18ml	1 x 60ml
BCIP/NBT Solution	1 x 12ml	2 x 18ml	1 x 120ml

Components	12 ml	36 ml	120 ml
Mouse HRP Polymer	1 x 6ml	1 x 18ml	1 x 60ml
Hydrogen Peroxide (20x)	1 x 1ml	1 x 2ml	1 x 6ml
Aqueous Mounting Medium	1 x 12ml	2 x 18ml	1 x 120ml

Applications

The Abpromise guarantee Our [Abpromise guarantee](#) covers the use of ab210060 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		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