

Product datasheet

DoubleStain IHC kit: G&M on human tissue (BCIP & AEC) ab183271

Overview

Product name	DoubleStain IHC kit: G&M on human tissue (BCIP & AEC)
Assay type	Polymer Multiplex
Product overview	The DoubleStain IHC kit (ab183271) is designed to use with user supplied goat and mouse antibodies to detect two distinct antigens on human tissue or cell samples. This kit has been tested in paraffin tissues, however it can also be used on frozen specimens and freshly prepared monolayer cell smears.

Double staining is a common method used in immunohistochemistry for the detection of two distinct antigens in a single tissue. Abcam's DoubleStain IHC Kit supplies two polymer enzyme conjugates: AP polymer anti-Goat IgG and HRP(AEC) polymer anti-Mouse IgG with two distinct substrates/chromogens, BCIP/NBT (Purple) and AEC (Red). Users will apply the two enzyme conjugates onto the specimen sequentially. If only the anti-goat antigen is present only the BCIP/NBT(Purple) will be present and if the mouse antigen is present only the AEC(Red) will be present. Abcam's DoubleStain IHC Kit is a non-biotin system avoiding endogenous biotin non-specific binding.

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

Properties

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

Components	120 ml	18 ml	6 ml
AEC Chromogen (20x)	1 x 8ml	1 x 4ml	1 x 2ml
AEC Substrate (20x)	1 x 4ml	1 x 2ml	1 x 1ml
Aqueous Mounting Medium	1 x 70ml	1 x 18ml	1 x 6ml
DoubleStain Blocker	1 x 60ml	1 x 18ml	1 x 18ml
Goat AP Polymer	1 x 60ml	1 x 18ml	1 x 6ml

Components	120 ml	18 ml	6 ml
Hydrogen Peroxide (20x)	1 x 4ml	1 x 2ml	1 x 2ml
Mouse HRP (AEC) Polymer	1 x 60ml	1 x 18ml	1 x 6ml
Ready to Use BCIP/NBT	1 x 120ml	1 x 18ml	1 x 6ml

Applications

The Abpromise guarantee Our **Abpromise guarantee** covers the use of ab183271 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-Fr		Use at an assay dependent concentration.
IHC-P		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