

DoubleStain IHC Kit: R&Rt on human/mouse tissue (Green/HRP & AP/Red) ab183285

★★★★★ [1 Abreviews](#) [2 References](#)

概述

产品名称 DoubleStain IHC试剂盒: R&Rt on人/小鼠组织(Green/HRP & AP/Red)

检测类型 Polymer Multiplex

产品概述 Abcam's Doublestain IHC Kit (ab183285) is designed for use with user supplied rabbit and rat primary antibodies to detect two distinct antigens on human and mouse tissue or cell samples. This kit has been tested on paraffin embedded tissue, however, it can also be used to stain frozen specimens or freshly prepared monolayer cell smears.

Double staining is a common method used in immunohistostaining, allowing for the detection of two distinct antigens in a single tissue. Abcam's Doublestain IHC Kit (ab183285) supplies the user with two polymer enzyme conjugates: Rat HRP Polymer (minimal cross reaction to mouse) and Rabbit AP Polymer with two distinct substrates/chromogens, Emerald and Permanent Red. Emerald chromogen reacts with the Rat HRP polymer conjugate to produce a green color. Permanent Red reacts with the Rabbit AP polymer to produce a red color. When two proteins are co-localized a blue/ purple color will develop depending on which antigen is stronger. If only the rat antigen is present only the Emerald chromogen will be present and if the rabbit antigen is present only the Permanent Red chromogen will be present. Abcam's Doublestain IHC Kit (ab183285) is a non-biotin system avoiding the extra steps involved in blocking non-specific binding due to endogenous biotin.

经测试应用 **适用于:** IHC-Fr, IHC-P

性能

存放说明 Store at +4°C. Please refer to protocols.

组件	36 ml	60 ml	6 ml	120 ml
Emerald Chromogen	2 x 18ml	1 x 60ml	1 x 6ml	1 x 120ml
Non-aqueous Mounting Medium	4 x 9ml	0 x 0ml	1 x 6ml	0 x 0ml
Permanent Red Activator (5x)	1 x 7.2ml	1 x 24ml	1 x 3ml	2 x 12ml

组件	36 ml	60 ml	6 ml	120 ml
Permanent Red Chromogen (100x)	1 x 360µl	1 x 1.2ml	1 x 150µl	1 x 1.2ml
Permanent Red Substrate	2 x 18ml	1 x 120ml	1 x 15ml	1 x 120ml
Rabbit AP Polymer	1 x 18ml	1 x 60ml	1 x 6ml	1 x 60ml
Rat HRP Polymer	1 x 18ml	1 x 60ml	1 x 6ml	1 x 60ml

应用

The Abpromise guarantee **Abpromise™**承诺保证使用ab183285于以下的经测试应用

“应用说明”部分 下显示的仅为推荐的起始稀释度;实际最佳的稀释度/浓度应由使用者检定。

应用	Ab评论	说明
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.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors