



Product datasheet

Human on human IHC kit (HRP/DAB) ab214749

1 References

製品の概要

製品名	Human on human IHC kit (HRP/DAB)
種交差性	交差種: Human
製品の概要	<p>Abcam's Human on human IHC kit (HRP/DAB) (ab214749) is designed for staining human primary antibodies on human tissues without background staining.</p> <p>Antigen detection using a primary antibody of the same species as the test tissue yields high background when an indirect detection method is used. This severely limits the ability to screen human antibodies on human tissues. Abcam's Human on human IHC kit (HRP/DAB) (ab214749) is a ready to use system that results in excellent sensitivity and high specificity using a simple protocol. Furthermore, as the protocol uses a polymer-based, biotin-free system, it does not require blocking of endogenous biotin.</p>
アプリケーション	適用あり: IHC

製品の特性

保存方法	Store at +4°C. Please refer to protocols.
------	---

内容	6 ml	18 ml	110 ml
20X DAB Chromogen	1 x 1.5ml	1 x 2ml	1 x 6ml
5X Quenching Buffer	1 x 1.5ml	2 x 2.3ml	2 x 13ml
Blocking Buffer A	1 x 6ml	1 x 18ml	1 x 110ml
Blocking Buffer B	1 x 6ml	1 x 18ml	1 x 110ml
DAB Substrate Buffer	1 x 12ml	2 x 15ml	1 x 120ml
Human HRP Polymer	1 x 6ml	1 x 18ml	1 x 110ml
Human Primer	1 x 6ml	1 x 18ml	1 x 110ml

アプリケーション

The Abpromise guarantee Abpromise保証は、次のテスト済みアプリケーションにおけるab214749の使用に適用されます
アプリケーションノートには、推奨の開始希釈率がありますが、適切な希釈率につきましてはご検討ください。

アプリケーション	Abreviews	特記事項
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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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