

Mouse and Rabbit Specific HRP Detection IHC Kit ab93677

10 References 画像数 1

製品の概要

製品名	Mouse and Rabbit Specific HRP Detection IHC Kit
製品の概要	Mouse and Rabbit Specific HRP Detection IHC Kit ab93677 contains complet reagents for a labeled streptavidin-biotin immunoenzymatic antigen detection system. This technique involves the sequential incubation of the specimen with an unconjugated primary antibody specific to the target antigen, a biotinylated secondary antibody which reacts with the primary antibody, enzyme-labeled streptavidin, and substrate-chromogen.
	Reacts with Mouse and Rabbit IgG (H+L)
特記事項	ab93677 does not contain chromogen or substrate.
アプリケーション	適用あり: IHC-P

製品の特性

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

内容	60 ml
Biotinylated Goat Anti-Polyvalent	1 x 60ml
Protein block	1 x 60ml
Streptavidin Peroxidase	1 x 60ml

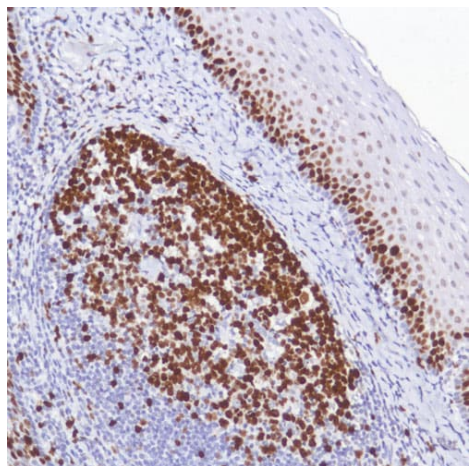
関連性 Polyvalent HRP is a labeled streptavidin-biotin immunoenzymatic antigen detection system. Horseradish peroxidase (HRP) readily combines with hydrogen peroxide (H₂O₂) and the resultant complex can oxidize a wide variety of chromogenic substrates.

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ab93677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-P		Use at an assay dependent dilution.

画像



Immunohistochemical analysis of Human Tonsil tissue labeling Ki-67 using the ab93677 anti-Mouse and Rabbit specific HRP Detection IHC kit.

Anti-Mouse and Rabbit Specific HRP Detection IHC Kit

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