

Fast Red Substrate Kit ab64254

13 References 画像数 2

製品の概要

製品名	Fast Red Substrate Kit
製品の概要	Fast Red Substrate kit ab64254 for immunohistochemical staining using alkaline phosphatase-based detection. Fast Red can be used as chromogen for immunohistochemical staining. In the presence of alkaline phosphatase enzyme, Fast Red produces a red reaction product that can be seen using either brightfield or fluorescent microscopy.
特記事項	This product was previously called Liquid Fast Red Substrate Kit. <u>Streptavidin - Alkaline phosphatase ab246833</u> is suitable for use with Fast red substrate. Find more kits and reagents for antigen retrieval, blocking, signal amplification, visualization, counterstaining, and mounting in the <u>IHC kits and reagents guide</u> .
アプリケーション	適用あり: IHC-P, WB

製品の特性

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

内容	125 ml	60 ml
75x Fast-Red Chromogen	1 x 2ml	1 x 1ml
Liquid Fast Red Substrate	1 x 125ml	1 x 60ml

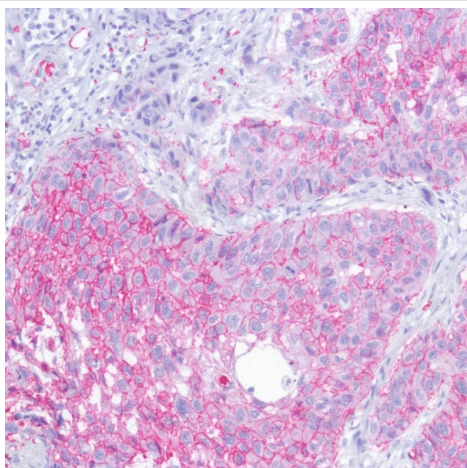
関連性 Fast-Red is a widely used chromogen which produces a red reaction product in the presence of alkaline phosphatase enzyme.

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ab64254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。

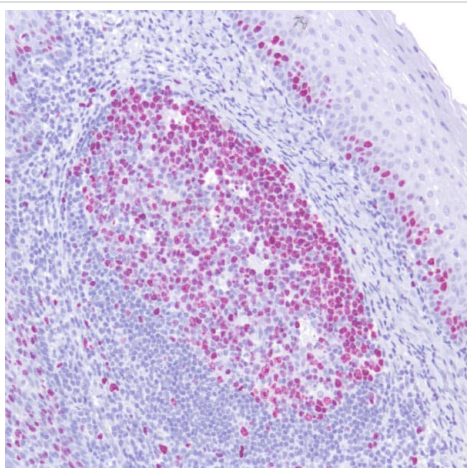
アプリケーション	Abreviews	特記事項
IHC-P		Use at an assay dependent dilution.
WB		Use at an assay dependent dilution.

画像



IHC of human esophageal carcinoma

Immunohistochemical staining of human esophageal carcinoma with **ab40084** at a dilution of 1/200 and detected using ab64254.



IHC of human tonsil

Immunohistochemical staining of human tonsil with **ab16667** at a dilution of 1/200 and detected using ab64254.

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