

Product datasheet

Fast Red Substrate Kit ab64254

[13 References](#) [2 Images](#)

Overview

Product name Fast Red Substrate Kit

Product overview 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.

Notes This product was previously called Liquid Fast Red Substrate Kit.

Streptavidin - Alkaline phosphatase ab246833 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 [IHC kits and reagents guide](#).

Tested applications **Suitable for:** IHC-P, WB

Properties

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

Storage buffer pH: 8.50
Constituents: 20% Fast Red TR salt, 20% Naphthol AS-MX phosphate, Tris buffer

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

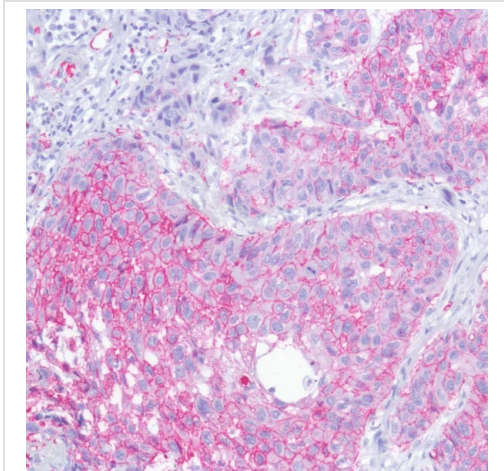
Relevance Fast-Red is a widely used chromogen which produces a red reaction product in the presence of alkaline phosphatase enzyme.

Applications

The Abpromise guarantee Our [Abpromise guarantee](#) covers the use of ab64254 in the following tested applications. The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

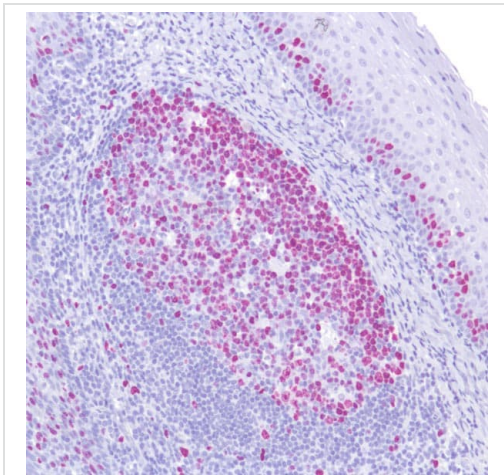
Application	Abreviews	Notes
IHC-P		Use at an assay dependent dilution.
WB		Use at an assay dependent dilution.

Images



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.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