

## Product datasheet

# AEC Substrate Kit ab64252

4 References 1 Image

### Overview

---

<b>Product name</b>	AEC Substrate Kit
<b>Description</b>	AEC Substrate Kit
<b>Tested applications</b>	<b>Suitable for:</b> IHC-P
<b>General notes</b>	AEC substrate kit for use in IHC using HRP / peroxidase detection. Ready to use formulation. Previously called AEC Substrate System (Ready to Use).

#### IHC protocol suitable for use with AEC Substrate Kit ab64252:

1. Deparaffinize and rehydrate formalin-fixed paraffin-embedded tissue section.
2. Use appropriate [antigen retrieval buffer or enzyme](#) (primary antibody dependent) to treat sections. Wash 3 times in buffer.
3. Add enough [hydrogen peroxide blocking solution](#) to cover the sections. Incubate for 10 minutes. Wash 2 times in buffer. If necessary, use [avidin biotin blocking](#).
4. Apply [protein block](#) (or [normal serum](#) from same species as secondary antibody) and incubate for 5 minutes at room temperature to block nonspecific background staining. Wash once in buffer.
5. Apply primary antibody in [antibody diluent](#) and incubate.
6. Wash 4 times in buffer. Incubate slide with [biotinylated secondary antibody](#) (or [HRP polymer secondary antibody](#) and skip step 7). Wash 4 times in buffer.
7. Apply [streptavidin-HRP](#) and incubate for 10 minutes at room temperature.
8. Rinse 4 times in buffer. **Place slide in DAB substrate or AEC Substrate ab64252 and incubate until desired color is achieved (1-10 mins, recommended 10 mins for AEC).** Rinse 4 times in buffer.
9. Add enough drops of [hematoxylin](#) to cover the section. Incubate for 1 minute.
10. Rinse 7-8 times in tap water. Add [mounting medium](#) to cover the section.

Find complete IHC kits, and reagents for antigen retrieval, blocking, signal amplification, visualization, counterstaining, and mounting in the [IHC kits and reagents guide](#).

### Properties

---

<b>Form</b>	Liquid
-------------	--------

**Storage instructions** Store at +4°C.

**Storage buffer** Constituents: 0.03% Hydrogen peroxide, 3% DMF stabilizers

**Relevance** AEC (3-Amino-9-Ethylcarbazole) is a widely used chromogen for immunohistochemical staining. AEC produces a red end product that is soluble in alcohol and must be used with an aqueous counterstain and mounting media. AEC produces a clean, very intense stain that is free of unwanted background.

## Applications

Our [Abpromise guarantee](#) covers the use of **ab64252** 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.

## Images



Human colon carcinoma stained with anti-Keratin 20 antibody using AEC Substrate System (Ready to Use) (ab64252).

Immunohistochemistry (Formalin/PFA-fixed paraffin-embedded sections) - AEC Substrate System (Ready to Use) (ab64252)

**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