

## 5 References 画像数 2

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

| 内容                                    | 15 ml    | 125 ml    |
|---------------------------------------|----------|-----------|
| Goat anti-rabbit HRP Conjugate        | 1 x 15ml | 1 x 125ml |
| Hydrogen Peroxide Block               | 1 x 15ml | 1 x 125ml |
| Mouse Specifying Reagent (Complement) | 1 x 15ml | 1 x 125ml |
| Protein block                         | 1 x 15ml | 1 x 125ml |

**関連性** Biotin-Free Polyvalent AEC detection is a biotin-free immunoenzymatic antigen detection system. AEC is 3-amino-9-ethyl-carbazole, a chromogenic substrate.

## 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ab236467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 concentration. |

## 画像

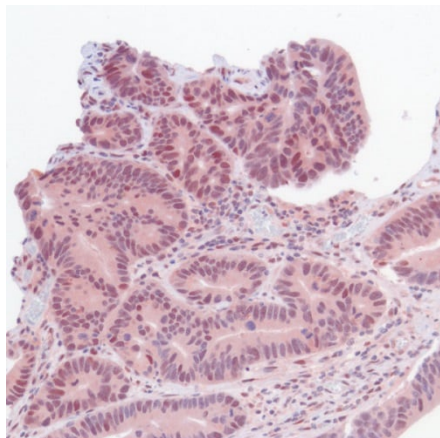


Image using ab93686 (previous version of this kit)

Immunohistochemical staining of p53 in human colon adenocarcinoma with anti-p53 antibody at a dilution of 1/400 and detected using [ab93686](#) (this kit is identical to [ab93686](#), except that it uses the same HRP micro-polymer antibody used by Abcam's R&D team for antibody validation in IHC).

Immunohistochemistry (Formalin/PFA-fixed paraffin-embedded sections)

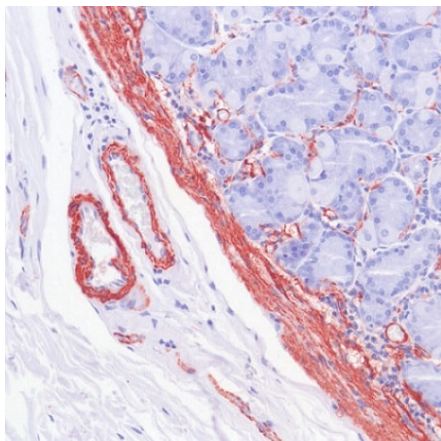


Image using ab93686 (previous version of this kit)

Immunohistochemistry (Formalin/PFA-fixed paraffin-embedded sections)

Immunohistochemical staining of smooth muscle actin in human small intestine with **ab7817** at a dilution of 1/200 and detected using **ab93686** (this kit is identical to **ab93686**, except that it uses the HRP micro-polymer secondary antibody used by Abcam's R&D team for antibody validation in IHC).

**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