

StayBlue/AP Plus ab176915

1 Image

Overview

| | |
|------------------|--|
| Product name | StayBlue/AP Plus |
| Assay time | 0h 20m |
| Product overview | StayBlue/AP Plus is a substrate chromogen system designed to be used for either IHC or ISH when using alkaline phosphate detection. StayBlue/AP Plus produces a blue color. It is insoluble in alcohol and xylene substitutes (both aliphatic hydrocarbon and citrus based) therefore sections can be dehydrated in alcohol, cleared in xylene substitute and permanently mounted. This chromogen substrate system may be used for both automation and manual use. |

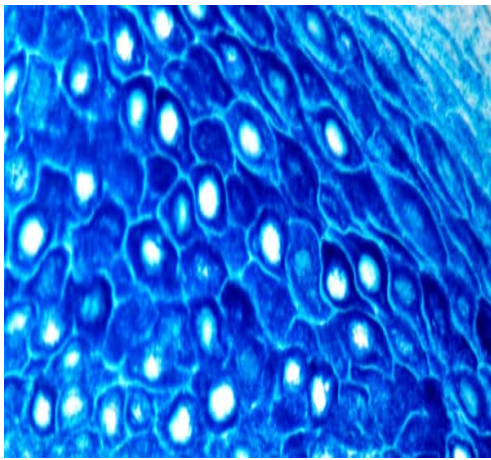
| | |
|-------|---|
| Notes | Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances. It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses. |
|-------|---|

Properties

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

| Components | 30 ml | 110 ml |
|-----------------------------------|----------|-----------|
| Empty Mixing Dropping Bottle | 1 x 30ml | 1 x 110ml |
| StayBlue/AP Plus Chromogen | 1 x 1ml | 1 x 3ml |
| StayBlue/AP Plus Substrate Buffer | 1 x 30ml | 1 x 110ml |

Images



Formalin-fixed; paraffin-embedded Human tonsil tissue stained with HMW CK antibody labeled with StayBlue/AP Plus (ab176915) produces a strong blue color.

Immunohistochemistry (Formalin/PFA-fixed paraffin-embedded sections) - StayBlue/AP Plus (Alcohol and Xylene Substitute Compatible) (ab176915)

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