

Product datasheet

StayBlack/HRP ab179917

1 Image

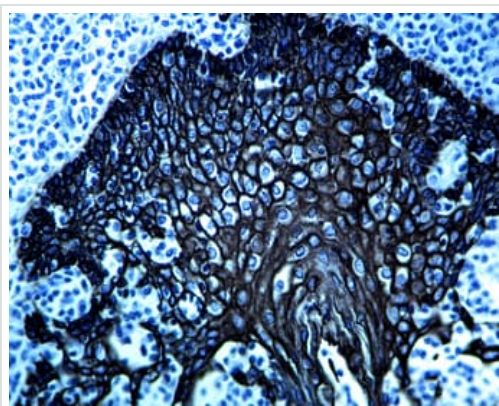
Overview

Product name	StayBlack/HRP
Assay time	0h 10m
Product overview	StayBlack/HRP is a substrate chromogen system designed to be used for either IHC or ISH when using horseradish peroxidase detection. StayBlack/HRP produces a black color. This chromogen substrate system may be used for both automation and manual use.

Properties

Components	30 ml	110 ml
Empty Mixing Dropping Bottle	1 unit	1 unit
StayBlack/HRP Chromogen	1 x 1ml	1 x 3ml
StayBlack/HRP Substrate Buffer	1 x 30ml	1 x 110ml

Images



Immunohistochemistry (Formalin/PFA-fixed paraffin-embedded sections) - StayBlack/HRP (ab179917)

Formalin-fixed; paraffin-embedded Human tonsil tissue stained with HMW Cytokeratin antibody ([ab850](#)) labeled with StayBlack/HRP (ab179917) produces a strong black color.

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