

ab286894 – UltraBrite Yellow IHC chromogen (HRP)

For staining tissue for immunohistochemistry or ISH.
For research use only – not intended for diagnostic use.

For overview, typical data and additional information please visit:

<http://www.abcam.com/ab286894>

Storage and Stability:

Store at +4°C.

Materials Supplied

Item	Quantity	Storage Condition
UltraBrite Yellow IHC Chromogen (HRP) Substrate buffer	30 mL	+4°C
UltraBrite Yellow IHC Chromogen (HRP) Chromogen	1 mL	+4°C
Empty Mixing Bottle	1 Unit	+4°C

Reagent Preparation

- Aliquot 1 mL of UltraBrite™ yellow IHC chromogen (HRP) substrate Buffer in a mixing bottle.
- Add one drop (~20 µL) of concentrated UltraBrite yellow IHC chromogen (HRP) chromogen solution.
- Replace tip, mix, and allow solution to reach room temperature before using.

Δ Note: The UltraBrite yellow IHC chromogen (HRP) chromogen-substrate working solution is light sensitive and should be kept away from light as much as possible. Working solution is stable for up to 1 day, for optimal results prepare fresh reagent.

Assay Protocol

Staining Procedures:

– Following peroxidase incubation, wash tissue sections with wash buffer, then follow protocol of choice (working solution incubation time: 10 – 20 mins):

Pre-Mix Working Solution (Automation): UltraBrite yellow IHC chromogen (HRP) has a working solution stability of at least 1 day and can be loaded directly onto instrument as a single solution. Reduce exposure to light to achieve optimal staining. Working solution is applied directly to slide.

On Board Mixing (Automation): Instruments that have on-board mixing capability can load the chromogen and substrate-buffer components independently. Working solution is made mixing reagents 1:50 using on-board mixing station before application to slide.

Manual Use: Mix substrate-chromogen and buffer in a 1:50 ratio and apply directly to slide.

Counterstain:

Counterstain with Hematoxylin or Nuclear Fast Red for good contrast. Wash with distilled or deionized H₂O followed by immuno wash buffer.

Mounting:

Slides should be air dried (do not dehydrate in alcohol or xylene). After rinsing off counterstain in distilled or de-ionized H₂O, leave slides on benchtop for at least 20 minutes to air dry, then permanently mount.

Technical Support

Copyright © 2025 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

www.abcam.com/contactus

www.abcam.cn/contactus (China)

www.abcam.co.jp/contactus (Japan)