

# ab282569 Western Blot Stripping Buffer

Western Blot Stripping Buffer (ab282569) uniformly removes antibodies from Western blots so that the blots may be reprobed with different antibodies. Western Blot Stripping Buffer is compatible with nitrocellulose and PVDF membranes.

View [ab282569](#)

Western Blot Stripping Buffer datasheet:

<https://www.abcam.com/ab282569>

(use <https://www.abcam.cn/ab282569> for China,  
or <https://www.abcam.co.jp/ab282569> for Japan)

This product is for research use only and is not intended for diagnostic use.

## Materials Supplied and Storage

- Upon receipt store at room temperature.
- Supplied as 500 ml.

## Important Notes

- Blots may be stored in PBS or TBS at 4°C until the stripping procedure is performed. DO NOT STORE BLOTS IN DRY FORM.
- Probe the Western blot for the lowest expressed antigen first, as stripping the blot may reduce signal intensity.

## Stripping Protocol

Allow all reagents to warm to room temperature before use.

- Rinse blot with 20 mL of water for 5 minutes at room temperature with agitation.
- Strip blot with 20 mL of Western Blot Stripping Buffer for 5 minutes at room temperature with agitation.
- Wash blot 3 times with 20 mL 1X PBS-T; 5 minutes per wash at room temperature with agitation.
- Block membrane and carry out antibody incubations.

***Δ Note:*** Probe the Western blot for the lowest expressed antigen first, as stripping the blot may reduce signal intensity.

## Technical Support

Copyright © 2021 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print | Version 1

**For all technical or commercial enquiries please go to:**

[www.abcam.com/contactus](http://www.abcam.com/contactus)

[www.abcam.cn/contactus](http://www.abcam.cn/contactus) (China)

[www.abcam.co.jp/contactus](http://www.abcam.co.jp/contactus) (Japan)