


Ponceau S Staining Solution

Catalog Numbers A40000279 and A40000278

Pub. No. MAN0025939 Rev. A.0

 **WARNING!** Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from [thermofisher.com/support](https://www.thermofisher.com/support).

Product description

Thermo Scientific™ Ponceau S Staining Solution (Cat. No. [A40000279](#), [A40000278](#)) is a ready-to-use membrane stain for evaluating protein transfer efficiency or for use in total protein normalization for western blot analysis. The stain has minimal nonspecific interactions with nitrocellulose or PVDF membranes, providing a reliable method for visualization of membrane-bound proteins. Staining with Ponceau S Staining Solution results in pink-reddish protein bands that are easily photographed and reversed in less than 5 minutes. Subsequent immunodetection is unaffected, as the stain does not alter membrane-bound proteins.

Contents and storage

Cat. No.	Amount	Storage
A40000279	500 mL	Room temperature
A40000278	1,000 mL	

Procedural guidelines

- Do not dilute before use. Ponceau S Staining Solution is supplied in a ready-to-use format.
- Ensure that membranes are completely immersed in staining solution throughout the incubation.
- Ensure that all staining solution is removed before blocking membranes for fluorescent western blot detection.

Stain nitrocellulose or PVDF membranes with Ponceau S Staining Solution

1. Wash the membrane in ultrapure water 3 times for 1 minute each time with gentle agitation to remove all transfer buffer.
2. Immerse the membrane in Ponceau S Staining Solution, then incubate with gentle agitation for 5–15 minutes.
3. Carefully wash the membrane in ultrapure water for 30–90 seconds or until the desired staining intensity is achieved.

IMPORTANT! To avoid uneven staining, do not pour water directly onto the membrane surface.

4. (Optional) Destain the membrane—Wash with a 0.1% NaOH solution for 1–2 minutes, then rinse in ultrapure water.

Related products

Product	Source
Cell lysis buffers	thermofisher.com/celllysis
Precast protein gels	thermofisher.com/proteingels
Western blotting reagents and accessories	thermofisher.com/westernblot
Western blot imaging and analysis	thermofisher.com/westernimaging

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Revision history: Pub. No. MAN0025939

Revision	Date	Description
A.0	11 November 2021	New manual for Ponceau S Staining Solution.

Important Licensing Information: This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2021 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.