



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

Ponceau S Solution

A red visible dye that rapidly stains proteins on western blotting membranes for easy visual inspection of protein transfer.



Product attributes

Product Description

Ponceau S is a classic visible red protein dye. It is commonly used during western blotting for easy visual inspection of protein transfer to PVDF and nitrocellulose membranes.

- Stains proteins on membranes pink or light red
- Limit of detection of about 250 ng of transferred protein
- Reversible and compatible with subsequent western blotting
- Solution contains 0.1% Ponceau S (w/v) and 5% acetic acid

Ponceau S staining is fast, easy and does not interfere with downstream western blotting. However, it is not very sensitive and staining can be uneven. To measure transferred protein quantitatively in fluorescent western blotting, we recommend the [VersaBlot™ Total Protein Normalization Kits](#). For safe and sensitive staining of proteins in gels, we recommend our [One-Step Protein Gel Stains](#).

See our [Western Blotting](#) Technology page to learn about more products for optimal western blots.

References

Analytical Chemistry (2016), doi:10.1021/acs.analchem.6b00699

This datasheet was generated on June 11, 2026 at 08:10:07 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/ponceau-s-solution/>