

46117 Landing Parkway, Fremont, CA 94538 U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352

www.biotium.com

## **PRODUCT AND SAFETY DATA SHEET**

**PRODUCT NAME: Ponceau S Solution** CATALOG # 22001 **PROPERTIES:** Color & Form Red solution N/A **Solubility** Absorption/Emission N/A Extinction N/A Coefficient STORAGE AND Store at room temperature or 4 °C. HANDLING: **APPLICATION:** Ponceau S Solution is used for the detection of proteins on PVDF and nitrocellulose membranes. This stain is not recommended with nylon membranes. Ponceau S staining is reversible to allow further immunological detection. The limit of detection for this stain is about 250 ng of transferred protein. Ponceau S is a negative stain which binds to the positively charged amino groups of the protein. PVDF membranes must be rinsed in methanol then rinsed in water just before staining. Nitrocellulose membranes can be stained directly. Typically, membranes are incubated for 5-10 min in the stain, then rinsed in water until the background is clean. The stain can be completely removed from protein bands by continued washing in water or buffer. The Ponceau S Solution contains 0.1% Ponceau S (w/v) and 5% acetic acid. **TOXICITY:** Not established. Not listed by NTP, IARC or OSHA. FIRST AID: Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

**Disclaimer:** *Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.*