

NBP1-48335 Protocol

SDS-Page protocol for Galectin-8 Protein (NBP1-48335)

SDS-PAGE (NBP1-48335): https://www.novusbio.com/products/galectin-8-protein_nbp1-48335

1. Mix equal volumes of human blood and Alsever's solution (pH 7.0).
(Alsever's solution: NaCl 0.42 g, Sodium citric acid 0.8g, Citric acid 0.055 g, D-glucose 2.05g in DW100 ml) <p/>
2. Centrifuge at 15000rpm for 10 minutes and wash four times with PBS. <p/>
3. Dilute packed cells in a 0.5 mg/ml trypsin-EDTA solution to give 4% red cell suspension. <p/>
4. Incubate for 1h at 37C and wash four times with PBS. <p/>
5. Dilute packed cells in PBS to give 4% red cell suspension. <p/>
6. Load 50ul of 0.5%BSA-in-0.15M-NaCl solution and 25ul of 4%-Red-Cell-in-PBS in u shaped wells. <p/>
7. Add 25ul of serial diluted galectin protein in PBS to each well plate. (Round bottom 96 well plate) <p/>
8. Incubate for 30min at room temperature to observe visible agglutination.