



Orders: [orders@novusbio.com](mailto:orders@novusbio.com)

Support: [technical@novusbio.com](mailto:technical@novusbio.com)

Web: [www.novusbio.com](http://www.novusbio.com)

Protocols, Publications, Related Products, Reviews and more:

[www.novusbio.com/NBP1-30270](http://www.novusbio.com/NBP1-30270)

## NBP1-30270 Protocol

### SDS-Page protocol for Galectin-4 Biologically Active Protein (NBP1-30270)

SDS-PAGE (NBP1-30270): [https://www.novusbio.com/products/galectin-4-biologically-active-protein\\_nbp1-30270](https://www.novusbio.com/products/galectin-4-biologically-active-protein_nbp1-30270)

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 DW 100 ml) <br/>
2. Centrifuge at 15,000rpm for 10 minutes and wash four times with PBS.<br/>
3. Dilute packed cells in a 0.5 mg/ml trypsin-EDTA solution to give 4% red cell suspension. <br/>
4. Incubate for 1hr at 37C and wash four times with PBS.<br/>
5. Dilute packed cells in PBS to give 4% red cell suspension.<br/>
6. Load 50ul of 0.5%BSA-in-0.15M-NaCl solution and 25ul of 4%-Red-Cell-in-PBS in u shaped wells. <br/>
7. Add 25ul of serial diluted galectin protein in PBS to each well plate. (Round bottom 96 well plate)<br/>
8. Incubate for 30min at room temperature to observe visible agglutination.