

## Bradford Assay Protocol

A Bradford Assay is used to determine the concentration of a protein before or after modification. The Protein Assay Kit I (from BioRad) is recommended. A reference protocol is provided below.

### A. Sample Preparation

1. Prepare 2 mL of Bradford working solution by adding 400  $\mu$ L Bradford dye reagent to 1600  $\mu$ L ultrapure water (1:4 ratios).
2. Prepare bovine IgG standards (or other appropriate standard) and a blank in 1.5 mL tubes as follows:
  - a. Add 100  $\mu$ L 2 mg/mL bovine IgG standard to 300  $\mu$ L PBS (0.5 mg/mL standard)
  - b. Add 200  $\mu$ L 0.5 mg/ml standard to 200  $\mu$ L PBS (0.25 mg/mL standard)
  - c. Add 200  $\mu$ L 0.25 mg/mL standard to 200  $\mu$ L PBS (0.125 mg/mL standard)
  - d. Add 200  $\mu$ L 0.125 mg/mL standard to 200  $\mu$ L PBS (0.0625 mg/mL standard)
  - e. 100  $\mu$ L PBS (buffer blank)
3. Dilute the protein sample to approximately 0.25 mg/mL with PBS to fall within the standard curve. Note the dilution factor used.

### B. Well Loading

1. In a flat-bottom 96-well plate, prepare standards by pipetting 11  $\mu$ L of each standard (and the blank) into separate wells.
2. Add 10  $\mu$ L of protein sample to 3 separate wells.
3. Add 100  $\mu$ L of Bradford working solution to each well using a multi-channel pipet.
4. Seal the plate with adhesive tape and shake for 15 seconds using a plate reader to mix.

### C. Plate Reading

1. Incubate the plate at room temperature for 5-60 min.
2. Measure absorbance at 595 nm using pre-programmed BCA assay plate reader software.

## Materials Required

Reagents	Equipment
Bradford Reagent	96-Well Plate
10x PBS	Plate Reader
Protein Standards	Microcentrifuge Tubes