

ProLite™ Blue 660 Protein Assay Reagent

Catalog number: 18003
Unit size: 100 mL

Component	Storage	Amount (Cat No. 18003)
ProLite™ Blue 660 Protein Assay Reagent	Refrigerated (2-8 °C), Minimize light exposure	100 mL

OVERVIEW

ProLite™ Blue 660 Protein Assay Reagent provides a simple, mix-and-read solution for quantifying total protein concentrations in solution using absorbance at 660 nm. The assay is compatible with BSA and other common protein standards, and performs reliably across a broad dynamic range. Optimized for microplate-based workflows, the assay requires only 10 µL of sample and a 5-minute incubation period, offering a convenient format for routine protein quantification in biochemical and life science research.

AT A GLANCE

1. Prepare a BSA standards.
2. Add 10 µL of standards and samples into the wells of a microplate.
3. Add 150 µL of ProLite™ Blue 660 Protein Assay Reagent to each well, mix gently, and incubate for 5 minutes at room temperature.
4. Measure absorbance at 660 nm and calculate protein concentrations.

KEY PARAMETERS

Absorbance microplate reader

Absorbance	660 nm
Recommended plate	Clear bottom

SAMPLE EXPERIMENTAL PROTOCOL

Microplate procedure (15.6 to 1000 µg/mL)

1. Prepare BSA standard solutions: Add 50 µL of 2 mg/mL BSA Standard (not provided) to 50 µL of PBS (not provided) to generate 1000 µg/mL BSA standard solution. Then use 1:2 serial dilutions in PBS to get serially diluted BSA standards to achieve 1000 to 15.6 µg/mL).
Note: It is necessary to create a standard curve during each assay.
2. Add 10 µL of each replicate of standard, unknown samples in appropriate microplate well.
3. Add 150 µL of the ProLite™ Blue 660 Protein Assay Reagent to each well and mix gently.
4. Cover plate and incubate at room temperature for 5 minutes.
5. Measure the absorbance of the standards and unknown samples at 660 nm.
6. Prepare a standard curve by plotting the average bBlank-corrected 660nm measurement for each BSA standard vs. its concentration in µg/mL. Use the standard curve to determine the protein concentration of each unknown sample.

IMPORTANT NOTES

1. Certain detergents (such as tween 20 at more than 10% or Triton X-100 at more than 1%) at very high concentrations may interfere with the performance of ProLite™ Blue 660 Protein Assay Reagent.
2. Precipitate may form in the ProLite™ Blue 660 Protein Assay Reagent over time. This may be more noticeable if the product is exposed to air for extended periods of time (3 to 4 months). Minimal precipitation will not impact performance and can be

removed by centrifugation.

EXAMPLE DATA ANALYSIS AND FIGURES

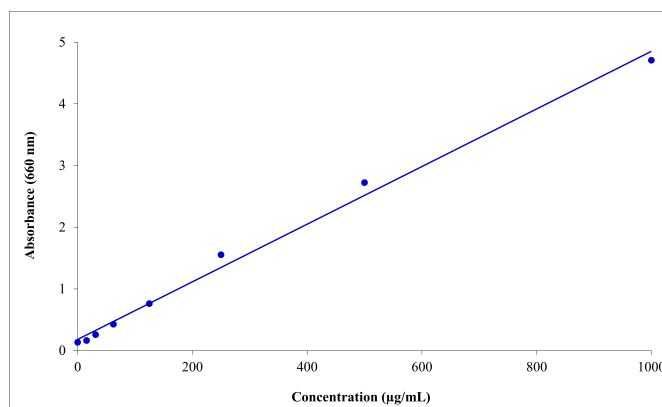


Figure 1. BSA dose responses was measured with ProLite™ Blue 660 Protein Assay Reagent using a clear bottom 96-well plate.

DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.