

ProLite™ FAST Blue Protein Gel Stain *200 Gels*

Catalog number: 18002
Unit size: Set

Component	Storage	Amount (Cat No. 18002)
Component A: ProLite™ FAST Blue Protein Gel Stain	Freeze (< -15 °C), Minimize light exposure	1 Set (200 Gels)
Component B: Staining Buffer	Freeze (< -15 °C), Minimize light exposure	1 Bottle (10 mL)

OVERVIEW

ProLite™ FAST Blue Protein Gel Stain is a ready-to-use solution. It is used to rapidly stain proteins on SDS-PAGE gels. The proteins stained with the ProLite™ FAST Blue Protein Gel Stain can be directly observed visually during or after electrophoresis. No post-staining, washing and tedious destaining are required. The staining can be easily recorded with a phone camera, eliminating the need for a sophisticated instrument. To stain a protein, one simply needs to add the dye to the protein solution, followed by brief incubation at room temperature. The staining does not require any additional steps, e.g., post-gel staining steps. It might be the quickest protein stain available.

KEY PARAMETERS

Gel Imager

Emission	Coomassie Blue Filter Set
Excitation	Coomassie Blue Filter Set

PREPARATION OF STOCK SOLUTIONS

1. Prepare a stock solution by adding 500 µL of DMSO into the vial of ProLite™ FAST Blue Protein Gel Stain (Component A).

Note: Any unused stock solution should be divided into single-use aliquots and stored at ≤-20 °C for two weeks. Avoid repeated freeze-thaw cycles and minimize light exposure.

Note: The 500 µL stock solution can effectively stain up to 200 gels, with each gel containing 10 wells.

PREPARATION OF WORKING SOLUTION

1. Prepare a working solution by adding 50 µL of the ProLite™ FAST Blue Protein Gel Stain stock solution to 1 mL of Staining buffer (Component B).

Note: Any unused working solution should be divided into single-use aliquots and stored at ≤-20 °C for two weeks. Avoid repeated freeze-thaw cycles and minimize light exposure.

SAMPLE EXPERIMENTAL PROTOCOL

1. Add 0.25 µL of the ProLite™ FAST Blue Protein Gel Stain stock solution to 10 µL of the protein lysate.

Note: The dye volume can be adjusted based on the total volume and protein concentration used in the lysate.

2. Incubate the mixture at room temperature for 15 to 60 minutes, protected from light.
3. Add the loading dye to the protein lysate as per the manufacturer's guidelines.
4. Load the samples onto the gel and run.

5. Following the completion of the gel run, the samples should be visibly present within the gel.

EXAMPLE DATA ANALYSIS AND FIGURES

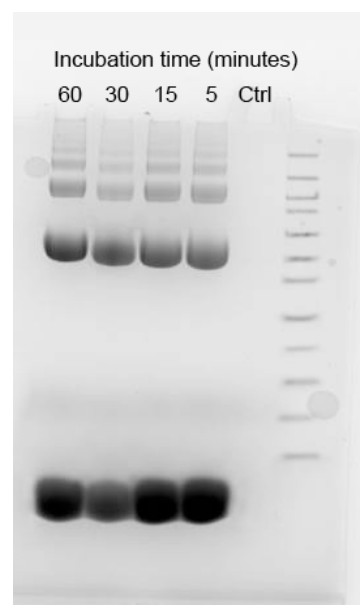


Figure 1. Bovine serum albumin (BSA) was stained with ProLite™ FAST Blue Protein Gel Stain (Cat No. 18002) for 60, 30, 15, and 5 minutes, and then loaded on a 4-12% Bis-Tris gel. A control sample was also included in the experiment, where no dye was added. The image was captured using a Coomassie Blue filter set.

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.