



## **eLuminol™ Protein Gel Stain, 1000X**

Catalog Number	Packaging Size
<b>P003A</b>	<b>0.5 mL</b>
<b>P003B</b>	<b>1 mL</b>

### **Storage upon receipt:**

2-25°C

Protect from light

**Ex/Em:** 300, 460/600 nm

## **User Manual**

GeneCopoeia, Inc.  
9620 Medical Center Drive, #101  
Rockville, MD 20850  
USA

301-762-0888

866-360-9531

[inquiry@genecopoeia.com](mailto:inquiry@genecopoeia.com)

[www.genecopoeia.com](http://www.genecopoeia.com)

## I. Product Description

**eLuminol Protein Gel Stain** is a high sensitive fluorescent stain designed for detecting proteins in polyacrylamide gels. Compared to traditional Coomassie® stains, **eLuminol Protein Gel Stain** has the following advantages:

- **High sensitivity.** eLuminol can detect as little as 0.2 ng protein.
- **Simple and fast staining.** After electrophoresis, the gel is simply stained, and washed. Take about 90 min.
- **Compatibility with standard laboratory equipment.** Stained protein can be viewed using a 300 nm UV transilluminator, blue light transilluminator or a laser scanner.
- **Wide linear detection range.** At least three orders of magnitude.
- **Compatible with downstream analysis:** Compatible with MS and sequencing.
- **Stable:** Stable at room temperature for 1 year.

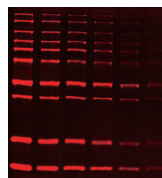
## II. Staining Protocol

**Note:** The protocol is optimized for standard 1 mm thick, 8 cm × 8 cm SDS-PAGE minigels. Larger or thicker gels require additional volumes of reagents or longer incubation times.

1. **Run** gel as usual according to your standard protocol.
2. **Prepare** 1X eLuminol stain solution by diluting 1,000X eLuminol Protein Gel Stain with a combination of water, methanol and acetic acid. For 1 mL of 1,000X eLuminol Stain, add 600 mL of water, 300 mL of methanol, and 100 mL of acetic acid. Store the 1X eLuminol Stain solution in a plastic bottle at room temperature or at 4 °C protected from light.
3. **Stain** gel with 1X eLuminol Stain solution (50~80 mL) at room temperature for 90 min with shaking.
4. **Wash** gel with 100 mL wash solution (10% methanol, 7% acetic acid) for 20 min with shaking.
5. **Image** gel with a 300 nm UV transilluminator, blue light transilluminator or a laser scanner.

## III. Using eLuminol Stain as a Post-Stain

1. Image the gel following staining with the first gel stain.
2. Rinse the gel with ultrapure water for 5 minutes. Repeat one more time.
3. Incubate gel with 1X eLuminol stain solution (50~80 mL) at room temperature for 90 min with shaking..
4. Wash gel with 100 mL wash solution (10% methanol, 7% acetic acid) for 20 min with shaking.
5. Image gel with a 300 nm UV transilluminator, blue light transilluminator or a laser scanner.



eLuminol Protein Gel Stain

## IV. Related Products

Catalog Number	Product
P001A	Coomassie Blue Fast Stain, 1X, 250 mL
P001B	Coomassie Blue Fast Stain, 1X, 500 mL
P001C	Coomassie Blue Fast Stain, 1X, 1000 mL