

## Product Information

### eLuminol™ Protein Gel Stain, 1000X

Catalog Number	Packaging Size
P003A	0.5 mL
P003B	1 mL

#### Storage upon receipt:

- 2-25°C
- Protect from light

Ex/Em: 300, 460/600 nm

### 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 0.5 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.

### Staining Protocol

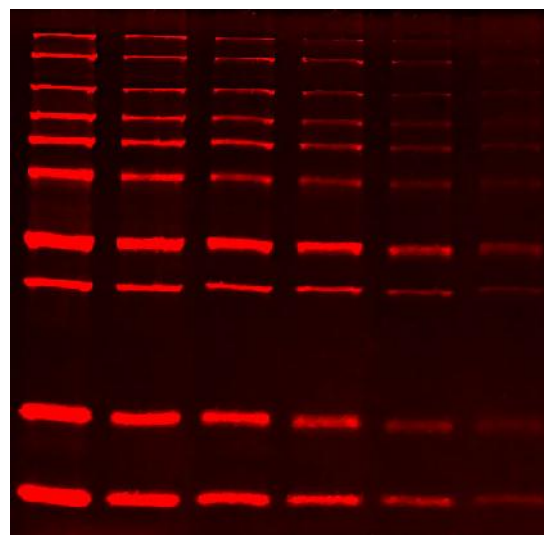
**Note:** The protocol is optimized for standard 1 mm thick, 8 cm × 8 cm SDS-PAGE minigels. Larger or thicker gels require additional volumes of reagents or longer incubation times.

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.

### Using eLuminol™ Stain as a Post-Stain (Optional)

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

### Related Products

Catalog No.	Product
P001	Coomassie® Blue Fast Stain, 1X, 1L