

Product Information

Phos-Tag™ Phosphoprotein Gel Stain

Catalog Number	Packaging Size
P005A	500 mL

Storage upon receipt:

- 2-25°C
- Protect from light

Ex/Em: 550/580 nm

Product Description

Phos-Tag™ Phosphoprotein Gel Stain is a high sensitive fluorescent stain designed for selectively detecting phosphoproteins in polyacrylamide gels. This stain contains a phos-tag™ group, which allows direct, in-gel detection of phosphate groups attached to tyrosine, serine, or threonine residues, without the need for antibodies or radioisotopes. The stain can be used with standard SDS-polyacrylamide gels or with 2-D gels. **Phos-Tag™ Phosphoprotein Gel Stain** has the following advantages:

- **High sensitivity.** Detect as little as 1 ng phosphoprotein.
- **Simple and fast staining.**
- **Compatibility with standard laboratory equipment.**
- **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.

Sample Preparation

A delipidated and desalted sample is essential for adequate separation of proteins by electrophoresis and subsequent staining by Phos-Tag™ Phosphoprotein Gel Stain.

1. For a 150 µL sample (~150-300 µg of protein), add 600 µL of methanol and mix well by vortexing.
2. Add 150 µL of chloroform and mix well by vortexing.
3. Add 450 µL of ultrapure water and mix well by vortexing.
4. Centrifuge at ~12,000 rpm for 5 min.
5. Discard the upper phase, keeping the white precipitation disc that forms between the upper and lower phases.
6. Add 450 µL of methanol and mix well by vortexing.
7. Centrifuge at ~12,000 rpm for 5 minutes.
8. Discard the supernatant and dry the pellet in a vacuum centrifuge for 10 minutes.
9. Resuspend the pellet in standard 1X sample buffer for electrophoresis.

Staining Protocol

Note: The protocol is optimized for standard 1 mm thick, 8 cm x 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. **Fix** gel with 100 mL of fix solution (50% methanol, 10% acetic acid), and agitate on an orbital shaker for 30 min. Repeat one more time with 100 mL fresh fix solution.
3. **Wash** the gel in 100 mL of ultrapure water with gentle agitation for 10 minutes. Repeat this step twice, for a total of three washes.
4. **Stain** the gel with enough **Phos-Tag™ Phosphoprotein Gel Stain** (40-60 mL) to cover the gel, and agitate on an orbital shaker for 60-90 min.
5. **Destain** the gel with **Phos-Tag™ Phosphoprotein Destain Solution** (P005B) with gentle agitation for 30 minutes. Repeat this procedure two more times.
6. **Wash** the gel twice with ultrapure water for 5 minutes per wash. If the background is high or irregular, the gel may be left in the second wash for 20-30 minutes and re-imaged.
7. **Image** gel with a 300 nm UV transilluminator, blue-light transilluminator or a laser scanner.

Protocol Quick Reference

	Reagent	Protocol
Fix	50% methanol, 10% acetic acid	100 mL, 30 min
		100 mL, 30 min
Wash	Ultrapure water	100 mL, 10 min
		100 mL, 10 min
		100 mL, 10 min
Stain	Phos-Tag™ Phosphoprotein Gel Stain	40-60 mL 60-90 minutes.
Destain	Phos-Tag™ Phosphoprotein Destain Solution	60 mL, 30 min
		60 mL, 30 min
		60 mL, 30 min
Wash	Ultrapure water	100 mL, 5 min
		100 mL, 5 min

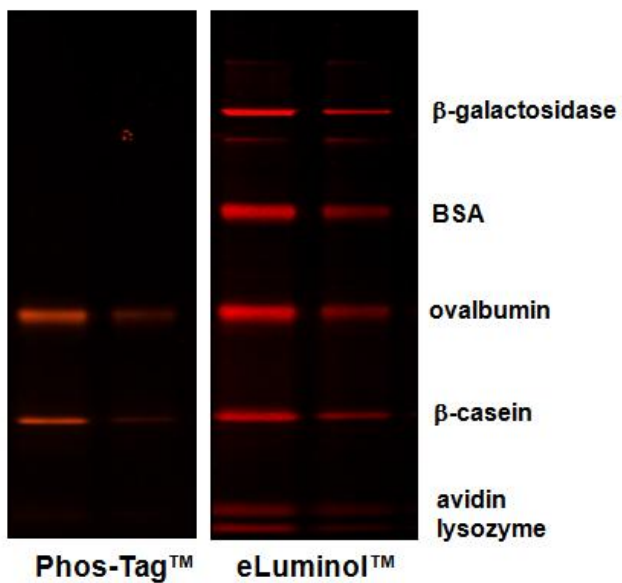
Staining the Gel for Total Protein

After staining with **Phos-Tag™ Phosphoprotein Gel Stain**, the gel can be stained with a total-protein 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 eLuminol™ Protein Gel Stain solution (40-60 mL). Microwave 45 seconds, and agitate on an orbital shaker for 15 min. Repeat microwave 45 seconds, and agitate on an orbital shaker for another 15 min.
4. **Wash** gel with 100 mL wash solution (10% methanol, 7% acetic acid) for 30 min.
5. **Image** gel with a 300 nm UV transilluminator, blue-light transilluminator or a laser scanner.

Related Products

Catalog No.	Product
P003A	eLuminol™ Protein Gel Stain, 0.5 mL
P003B	eLuminol™ Protein Gel Stain, 1 mL
P005B	Phos-Tag™ Phosphoprotein Destain Solution, 500mL



Protein gel stain results with Phos-Tag™ Phosphoprotein Gel Stain, followed by eLuminol™ Protein Gel Stain.