



3050 Spruce Street  
Saint Louis, Missouri 63103 USA  
Telephone 800-325-5832 • (314) 771-5765  
Fax (314) 286-7828  
email: techserv@sia.com  
sigma-aldrich.com

## Product Information

### Chemiluminescent Peroxidase Substrate

Product Codes **CPS-1, CPS-1-30, CPS-1-60, CPS-1-120, and CPS-1-300**

Storage Temperature 2-8 °C

## TECHNICAL BULLETIN

### Product Description

Sigma's Chemiluminescent Peroxidase Substrate can be used for the highly sensitive detection of peroxidase labeled material in a variety of Western blotting applications. This substrate is an enhanced luminol product with a stabilized peroxide buffer solution that provides picogram sensitivity with minimal background interference.

### Components

The Chemiluminescent Peroxidase Substrate is available in 4 package sizes each containing the Chemiluminescent Reagent (Product Code C 9107) and the Chemiluminescent Reaction Buffer (Product Code C 9232).

| Package Size | C 9107 | C 9232 |
|--------------|--------|--------|
| 30 ml        | 10 ml  | 20 ml  |
| 60 ml        | 20 ml  | 40 ml  |
| 120 ml       | 40 ml  | 80 ml  |
| 300 ml       | 100 ml | 200 ml |

### Precautions and Disclaimer

This product is for laboratory research use only. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

### Preparation Instructions

Prepare the Working Solution by mixing 1 part of the Chemiluminescent Reagent (Product Code C 9107) with 2 parts of the Chemiluminescent Reaction Buffer (Product Code C 9232). Mix well and protect from light. It is recommended to use 0.043 to 0.125 ml per cm<sup>2</sup> of membrane. For extended signal duration, a 1:1 ratio of Chemiluminescent Reagent to Chemiluminescent Reaction Buffer may be used.

### Storage/Stability

It is recommended to store the components at 2-8 °C. The components are stable for a minimum of 18 months when stored in the original container and protected from light. The Working Solution is stable for several hours at room temperature when protected from light.

### Procedure

Sigma's Chemiluminescent Peroxidase Substrate is very sensitive and great care must be taken to optimize the individual assay components (antibodies, conjugates, etc). In a Western blot, an optimized system is needed to minimize background reactivity associated with nonspecific immunochemical interactions. The following is a general guideline for the use of this product. The protocol starts with a transferred membrane.

### Notes:

- For optimal results, individual assay components must be optimized for minimal background and maximal signal.
- This product is designed for use only in Western blotting.
- All steps below should be performed with slight agitation on a rocker or an orbital shaker such that the membrane is freely floating.
- All incubations should be performed at room temperature.
- Gloves must be worn when working with the membrane to avoid contamination.
- Azide inhibits horseradish peroxidase (HRP) and should not be used as a buffer preservative for assay components.

1. Remove membrane from Western blotting apparatus and wash membrane for 1 minute in either Tris-buffered Saline with TWEEN<sup>®</sup> 20 (TBST, Product Code T 9039) or phosphate buffered saline with TWEEN 20 (PBST, Product Code P 3563). Note that either a TBS or PBS system can be used for Western blotting.
2. Block membrane in appropriate blocking agent for 30 minutes. Western Blocker Solution (Product Code W 0138) is recommended for high sensitivity detection.
3. Add primary antibody to the blocking agent. The final concentration of primary antibody in this solution can range from 0.2-20 µg/ml.
4. Incubate membrane with the primary antibody solution for at least 30 minutes.
5. Wash with TBST or PBST for 1 minute.
6. Remove TBST or PBST and add at least 10 ml of appropriate blocking agent to the membrane. Add secondary antibody; a 1:50,000 to 1:500,000 dilution in blocking agent may be used.
7. Incubate the membrane with the secondary antibody solution for 30 minutes.
8. Remove blocking solution and wash membrane 5 times for 5 minutes each with TBST or PBST.
9. Remove the membrane from the wash buffer and drain any excess liquid from the membrane. Keep the membrane damp; do not let the membrane dry out.
10. Place the membrane on a flat sheet of plastic wrap (or on any clean plastic surface).
11. Develop the blots with the Working Solution for 5 minutes.
12. Drain excess substrate and place in holder or plastic wrap.
13. Expose BioMax light film to the blot. Exposure times range from 30 seconds to 10 minutes. It is best to do a quick exposure of 10 to 30 seconds to determine what exposure time is needed. If the signal is too intense even at the short exposure times, let the signal decay from 1 to 8 hours and then re-expose the film.

#### Related Products

| Product Name             | Package Size | Product Code |
|--------------------------|--------------|--------------|
| TBS                      | 10 packets   | T 6664       |
| PBS                      | 10 packets   | P 3813       |
| Western Blocker Solution | 400 ml       | W 0138       |
| TBS + 3% milk            | 10 packets   | T 8793       |
| PBS + 3% milk            | 10 packets   | P 2194       |
| PBS + 5% milk            | 10 packets   | P 4739       |
| TBS + TWEEN 20           | 10 packets   | T 9039       |
| PBS + TWEEN 20           | 10 packets   | P 3563       |
| Anti-Mouse HRP Antibody  | 2 ml         | A 9044       |

TWEEN is a registered trademark of the ICI Group.

RBG/MKS/MAM 1/04

## Troubleshooting Guide

| Problem Type                                                  | Cause                                                            | Solution                                                                                                                              |
|---------------------------------------------------------------|------------------------------------------------------------------|---------------------------------------------------------------------------------------------------------------------------------------|
| Too much background signal observed.                          | Not enough wash steps were performed at the end of the blotting. | Double the number of washing steps.                                                                                                   |
|                                                               | Too much primary antibody used.                                  | Lower the amount of primary antibody used and wash with TBST for 5 minutes instead of 1 minute after the primary antibody incubation. |
|                                                               | Too much secondary antibody used.                                | Lower the amount of secondary antibody used.                                                                                          |
| Image is reversed on film (dark background and light bands).  | Too much secondary antibody used.                                | Lower the amount of secondary antibody used.                                                                                          |
| Bands on membrane have brown or yellow tone.                  | Too much secondary antibody used.                                | Lower the amount of secondary antibody used.                                                                                          |
| Nonspecific bands show up on membrane.                        | Too much primary antibody used.                                  | Lower the amount of primary antibody used and wash with TBST for 5 minutes instead of 1 minute after the primary antibody incubation. |
|                                                               | Too much secondary antibody used.                                | Lower the amount of secondary antibody used.                                                                                          |
| Membrane is stippled.                                         | Secondary antibody has some aggregate formation.                 | Filter secondary antibody.                                                                                                            |
| No signal is seen with chemiluminescent reaction on membrane. | Protein levels are too low for detection.                        | Increase exposure time of film and increase level of protein loads.                                                                   |
|                                                               | Not enough primary antibody used.                                | Use more primary antibody.                                                                                                            |
|                                                               | Not enough secondary antibody used.                              | Use more secondary antibody.                                                                                                          |

Sigma brand products are sold through Sigma-Aldrich, Inc.

Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.