

---

## Product Information

### ECL Western Blotting Detection Kit

**Catalog Number: P007**

#### Kit Contents

**P007A: 50 mL of Solution A;**

**P007B: 50 mL of Solution B.**

<b>Storage upon receipt:</b>
------------------------------

- |  |
|--|
| <ul style="list-style-type: none"><li>• 4°C</li><li>• Protect from light</li></ul> |
|--|

### Product Description

The **ECL Western Blotting Detection Kit** is long-lasting and contains a highly sensitive, nonradioactive, enhanced luminol-based chemiluminescent substrate for the detection of horseradish peroxidase (HRP) conjugated antibodies on immunoblots. This kit contains 50ml of Solution A and 50ml of Solution B and is easy to use and economical. Mix solutions A and B in a 1:1 ratio right before the application. The mixture is then ready for detection of any immobilized specific antigens in chemiluminescent Western blots through horseradish peroxidase (HRP) labeled antibodies. The kit is designed for a high quality and long lasting glow.

### Protocol

The following procedure is recommended:

1. Wash membrane thoroughly before proceeding with ECL.
2. Immediately before use, mix solutions A and B in a 1:1 ratio.
3. Add the reagents to the blot, and incubate for 1 minute.
4. Pour out the solution, and wrap the blot in clear plastic.
5. Place the blot into a film cassette and expose to X-ray film in a dark-room setting immediately.