

# ECL Western Blotting Substrate Kit ab65623

★★★★★ [1 Abreviews](#) [110 References](#) [1 Image](#)

### Overview

<b>Product name</b>	ECL Western Blotting Substrate Kit
<b>Product overview</b>	<p>The ECL Western Blotting Substrate is a highly sensitive, nonradioactive, enhanced luminol-based chemiluminescent substrate for easy detection of horseradish peroxidase (HRP) on immunoblots. The ECL Western Blotting Substrate offers superior signal and clear background. Blots can be repeatedly exposed to X-ray film to obtain optimal results or stripped of the immunodetection reagents and re-probed.</p> <p>Our ECL kits include this kit ab65623 and our high sensitivity ECL substrate kits:</p> <ul style="list-style-type: none"> <li>- High Sensitivity <a href="#">ECL Substrate Kit ab133406</a> to detect 23pg-187ng of protein per band</li> <li>- Very High Sensitivity <a href="#">ECL Substrate Kit ab133408</a> to detect 4.6pg-4.7ng of protein per band</li> <li>- Ultra High Sensitivity <a href="#">ECL Substrate Kit ab133409</a> to detect 1.2pg-2ng of protein per band</li> </ul> <p>Detection ranges in pg and ng stated above should be used for guidance only as detection range is dependent on the molecular weight of a protein.</p>
<b>Notes</b>	<p>This product is manufactured by BioVision, an Abcam company and was previously called K820 ECL Western Blotting Substrate Kit. K820-500 is the same size as the 500 test size of ab65623.</p> <p>Reagents provided are sufficient for 400 cm<sup>2</sup> of membrane.</p>
<b>Tested applications</b>	<b>Suitable for:</b> WB

### Properties

**Storage instructions** Store at +4°C. Please refer to protocols.

Components	50 tests
ECL Reagent A	1 x 25ml
ECL Reagent B	1 x 25ml

**Relevance** ECL (enhanced luminol-based chemiluminescent) detection is a sensitive, nonradioactive method for easy detection of horseradish peroxidase (HRP) on western blots.

## Applications

### The Abpromise guarantee

Our **Abpromise guarantee** covers the use of ab65623 in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
WB		Use at an assay dependent dilution.

## Images



ECL Western Blotting Substrate

ECL Western Blotting Substrate Kit - 50 Tests  
(ab65623)

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

### Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

### Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors