

ECL Western Blotting Substrate Kit ab65623

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

Overview

Product name ECL Western Blotting Substrate Kit

Product overview 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.

Our ECL kits include this kit ab65623 and our high sensitivity ECL substrate kits:

- High Sensitivity [ECL Substrate Kit ab133406](#) to detect 23pg-187ng of protein per band
- Very High Sensitivity [ECL Substrate Kit ab133408](#) to detect 4.6pg-4.7ng of protein per band
- Ultra High Sensitivity [ECL Substrate Kit ab133409](#) to detect 1.2pg-2ng of protein per band

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.

Notes 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.

Reagents provided are sufficient for 400 cm² of membrane.

Tested applications **Suitable for:** WB

Properties

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

Components	50 tests	500 tests
Reagent A	1 x 25ml	1 x 250ml
Reagent B	1 x 25ml	1 x 250ml

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

Applications

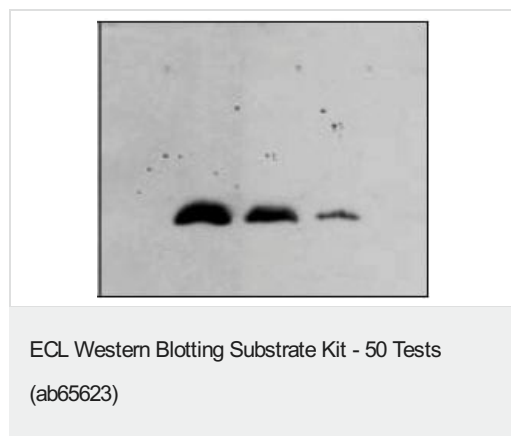
The Abpromise guarantee

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

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