

ECL Western Blotting Substrate Kit ab65623

★★★★★ [1 Abreviews](#) [110 References](#) [画像数 1](#)

製品の概要

製品名 ECL Western Blotting Substrate Kit

製品の概要 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.

特記事項 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.

アプリケーション **適用あり:** WB

製品の特性

保存方法 Store at +4°C. Please refer to protocols.

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

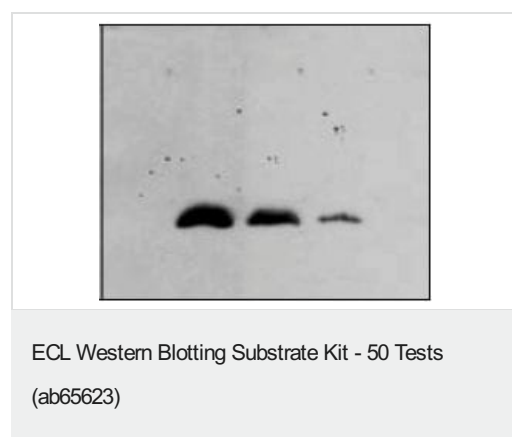
関連性 ECL (enhanced luminol-based chemiluminescent) detection is a sensitive, nonradioactive method for easy detection of horseradish peroxidase (HRP) on western blots.

アプリケーション

The Abpromise guarantee Abpromise保証は、 次のテスト済みアプリケーションにおけるab65623の使用に適用されます
アプリケーションノートには、推奨の開始希釈率がありますが、適切な希釈率につきましてはご検討ください。

アプリケーション	Abreviews	特記事項
WB		Use at an assay dependent dilution.

画像



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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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