

ECL Western Blotting Substrate Kit ab65623

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

概述

产品名称	ECL Western Blotting底物试剂盒
产品概述	<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">- 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 <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>
说明	<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 cm² of membrane.</p>
经测试应用	适用于: WB

性能

存放说明 Store at +4°C. Please refer to protocols.

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

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

应用

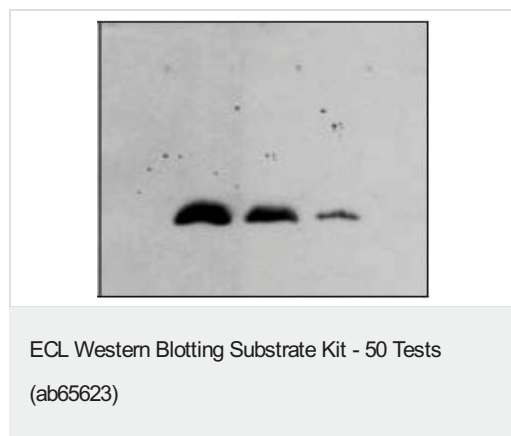
The Abpromise guarantee

Abpromise™承诺保证使用ab65623于以下的经测试应用

“应用说明”部分 下显示的仅为推荐的起始稀释度;实际最佳的稀释度/浓度应由使用者检定。

应用	Ab评论	说明
WB		Use at an assay dependent dilution.

图片



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.cn/abpromise> or contact our technical team.

Terms and conditions

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