

ECL Substrate Kit (High Sensitivity) ab133406

★★★★★ [5 Abreviews](#) [46 References](#) [2 图像](#)

概述

产品名称

ECL底物试剂盒(High Sensitivity)

产品概述

High Sensitivity ECL Substrate Kit ab133406 is designed for the detection of proteins with 23pg-187ng of protein per band. It uses an enhanced chemiluminescent substrate for western blotting that was developed for film imaging and is also compatible with CCD imaging. The high sensitivity ECL substrate produces a strong signal with very low background. Additionally, the ECL signal is long lasting, allowing repeated exposures without fear of losing data.

Our ECL kits include our popular [ECL Substrate Kit ab65623](#) and our high sensitivity ECL substrate kits:

- this kit (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.

This product was previously called Optiblot ECL Detect Kit (23pg-187ng).

说明

Use:

- 200ml kit for 2000cm² membrane
- 500ml kit for 5000cm² membrane

The primary antibody can often be diluted 5 to 10 fold more than usual when using this ECL substrate. A typical primary antibody dilution range using the substrate is 1/5000 - 1/20,000, with a typical secondary antibody dilution range of 1/20,000 - 1/100,000. Some optimisation may be required.

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	200 ml	20 ml	500 ml
Luminol/Enhancer Solution	1 x 100ml	1 x 10ml	1 x 250ml
Peroxide Chemiluminescent Detection Reagent	1 x 100ml	1 x 10ml	1 x 250ml

应用

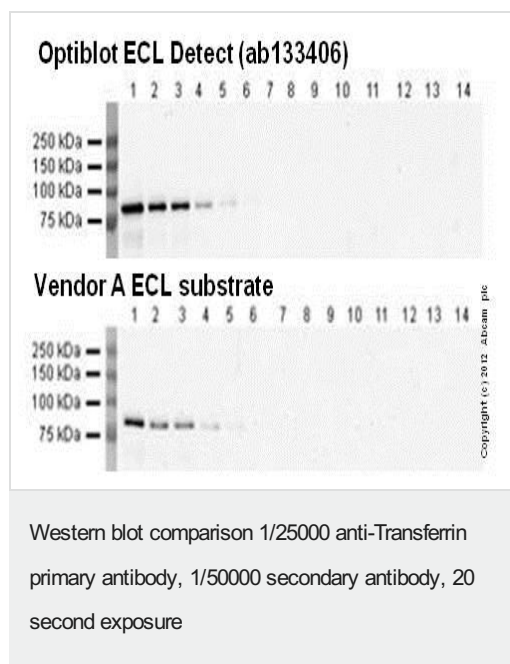
The Abpromise guarantee

Abpromise™承诺保证使用ab133406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 concentration.

图片

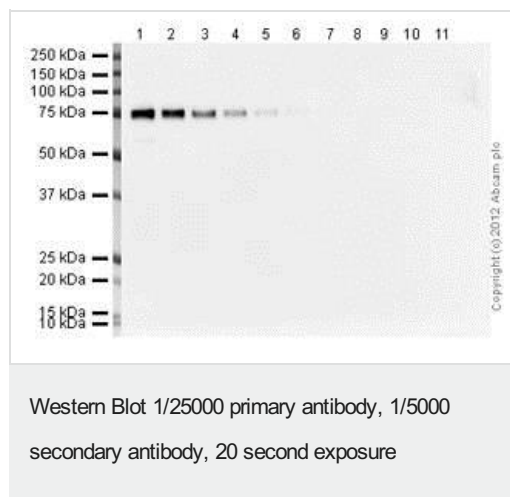


Anti-Transferrin antibody (**ab1223**) at **1/25000** dilution. Lanes 1-14 Transferrin protein (**ab91435**), Loading dilution (Lanes 1-8): 12.5, 6.2, 3.1, 1.6, 0.8, 0.3, 0.19µg protein.

Secondary

HRP conjugated polyclonal to Rabbit IgG (**ab97080**) at **1/50000** developed using Optiblot ECL Detect and Vendor A ECL substrate.

Exposure time: 20 seconds



Each lane contains the following amount of COX2 recombinant protein (**ab58868**): 1) 100ng 2) 50ng 3) 25ng 4) 12.5ng 5) 6.25ng 6) 3.13ng 7) 1.56ng 8) 780pg 9) 390pg 10) 195pg 11) 98pg. The blot was probed with rabbit anti-COX2 antibody (**ab15191**) at 1/25000 dilution and with a rabbit secondary (**ab97080**) at 1/5000 dilution. The blot was developed with 20ml of Optiblot ECL Detect.

Exposure time: 20 seconds

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