



Product datasheet

Chemiluminescent HRP Substrate (membrane) ab5801

2 References

概述

产品名称	Chemiluminescent HRP底物(membrane)
产品概述	A 2 component chemiluminescent substrate for the detection of horseradish peroxidase (HRP) in western blotting.
说明	<p>The Abcam Chemiluminescent HRP Substrate (membrane) is very sensitive and great care must be taken to optimize the individual assay components (antibodies, conjugates, solid phase, etc.) to minimize background reactivity associated with nonspecific immunochemical reactivities</p> <p>Prepare 1x substrate by mixing 1 ml of the chemiluminescent reagent with 2 ml of the chemiluminescent reaction buffer. Mix well and protect from light. More or less of the 1x substrate can be prepared using the same ratio.</p> <p>Stability: Stable for a minimum of 18 months when stored at 2-8°C in the original container and protected from light. The 1x diluted reagent is stable for several hours at room temperature when protected from light.</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	1 kit
luminol reagent.	1 x 20ml
reaction buffer.	2 x 20ml

应用

The Abpromise guarantee **Abpromise™**承诺保证使用ab5801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. Incubate membrane in chemiluminescent substrate for 1 minute then rinse off.

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