

Optiblot SDS-PAGE Sample Preparation Kit ab133414

画像数 1

製品の概要

製品名	Optiblot SDS-PAGE Sample Preparation Kit
製品の概要	<p>The Optiblot SDS-PAGE Sample Preparation Kit (ab133414) provides a means to quickly concentrate protein samples, and separate them from buffers that interfere with electrophoresis.</p> <p>The fast, efficient protocol generates samples ready to load on a gel in less than ten minutes, much more quickly than can be achieved with alternate methods such as dialysis, or acetone or TCA precipitation. Additionally, the Optiblot SDS-PAGE Sample Preparation protocol is easily scaled up, allowing multiple samples to be prepared in parallel.</p>
アプリケーション	適用あり: Purification

製品の特性

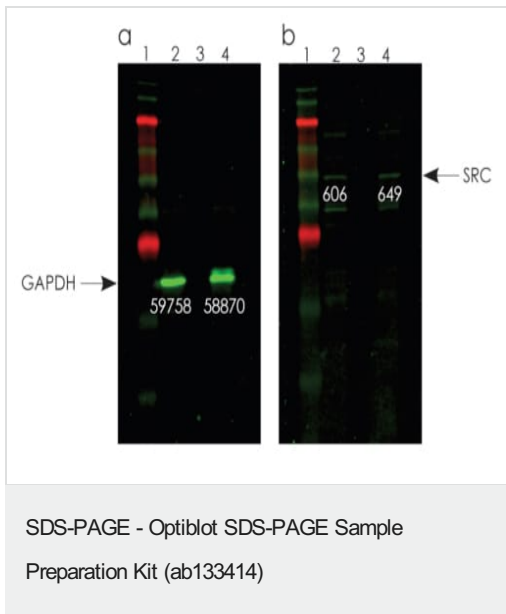
内容	10 tests	25 tests
Extraction loading buffer	1 x 1ml	1 x 1ml
Protein filters	10 units	25 units
Resin	1 x 500ml	2 x 500ml

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ab133414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
Purification		Use at an assay dependent concentration.

画像



Resin concentration is compatible with Western blot

detection. Western blots were created with samples of A431 cell extract (panels a and b, lane 2), diluted cell extract (lane 3) and the diluted extract concentrated using the Optiblot SDS-PAGE Sample Preparation Kit protocol (lane 4). The blots were detected using the Optiblot Fluorescent Western Blot Kit ([ab133410](#)). No bands were detectable in the diluted cell extract when stained for either GAPDH (panel a, lane 3) or SRC (panel b, lane 3). However, after resin concentration, bands are easily visualized for both proteins (lanes 4). Band intensities are indicated.

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