

SimpleStep ELISA® Custom ELISA Kit ab270552

SimpleStep ELISA

1 References [画像数 3](#)

製品の概要

| | |
|-------|-----------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| 製品名 | SimpleStep ELISA® Custom ELISA Kit |
| 製品の概要 | <p>SimpleStep ELISA® Custom ELISA Kit (ab270552) is the latest cutting-edge ELISA kit with a built-in system for antibody labeling, enabling the rapid and easy development of in-house ELISAs using any antibody pair, while eliminating the need to perform plate coating procedures. It is powered by the combination of a proprietary technologies, enabling you to reduce the amount of capture antibody required, measure different targets on the same plate with different assay antibody pairs, and streamline your ELISA to a single-wash format if desired.</p> <p>The capture antibody is quickly and easily conjugated with the peptide supplied within the kit. Once added to the plate, the interaction between the antibody and the peptide immobilizes the immune-complex to the plate.</p> <p>The detection antibody is similarly conjugated with HRP using the HRP conjugation reagent, also supplied in the kit. Thus, with minimal hands-on time, you can set up both of your antibodies to ensure the best performance of your ELISA.</p> <p>This kit has ready-to-go assay plates, pre-coated with antibody, which are specific for the peptide.</p> <p>This product is manufactured by Expedeon, an Abcam company. It was previously called CaptSure DIY ELISA Kit (kit sufficient for 5 x 96 well plates). Expedeon product code 6300020 is the same as the 5 x 96 tests size of this product.</p> |

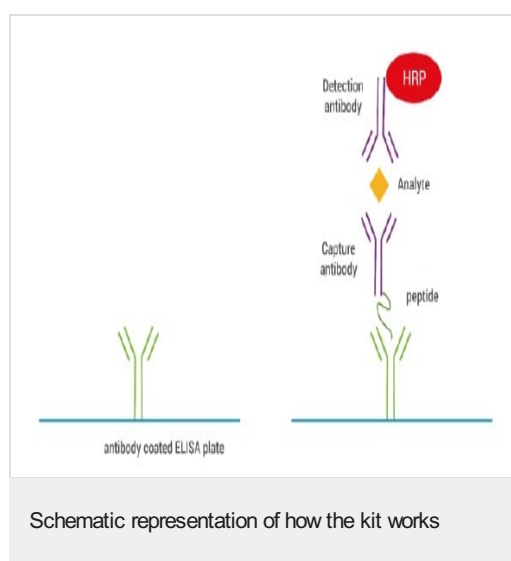
| | |
|------|--------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| 特記事項 | <p>Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.</p> <p>It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.</p> |
|------|--------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|

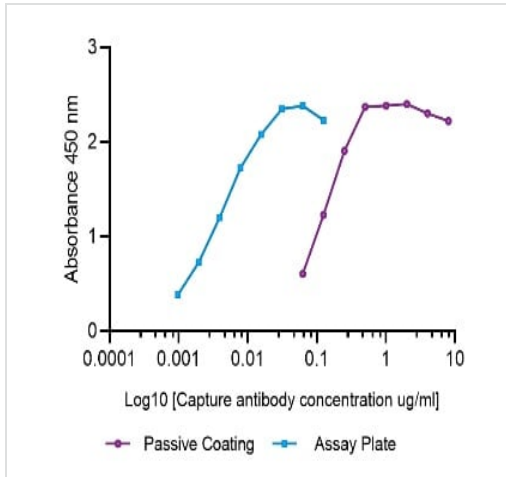
製品の特性

| | |
|------|----------------------------|
| 保存方法 | Please refer to protocols. |
|------|----------------------------|

| 内容 | 5 x 96 tests | 5 x 96 tests |
|----------------------------------------------------|------------------|------------------|
| ab275264 - Assay Plates (ab275264) | 1 x 5 x 96 tests | 1 x 5 x 96 tests |
| Capture Modifier reagent (ab270690) | 1 vial | 1 vial |
| Capture Peptide (ab270690) | 3 vials | 3 vials |
| Capture Quencher reagent (ab270690) | 1 vial | 1 vial |
| Detection Reagent (ab270697) | 1 unit | 1 unit |
| HRP mix | 3 x 100µg | 3 x 100µg |
| HRP-Labeled Antibody Dilution Buffer (ab270697) | 1 unit | 1 unit |
| Lysis Extraction Buffer (ab270697) | 1 unit | 1 unit |
| Modifier reagent | 1 x 200µl | 1 x 200µl |
| Peptide tagged Antibody Dilution Buffer (ab270697) | 1 unit | 1 unit |
| Quencher reagent | 1 x 200µl | 1 x 200µl |
| Sample Dilution Buffer (ab270697) | 2 units | 2 units |
| Stop Solution (ab270697) | 1 unit | 1 unit |
| Wash Buffer (ab270697) | 1 unit | 1 unit |

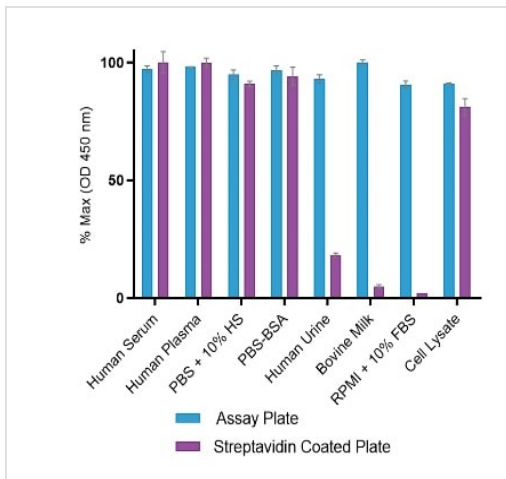
画像





SimpleStep ELISA® Custom ELISA Kit and conventional passive coating approach were compared. With ab270552, 10x less capture antibody can be used without affecting the assay performance.

Example data



Other solution-phase binding ELISA formats typically use biotin to immobilize immunocomplexes. These can constrain the sample types that can be used with the assay. In contrast, ab270552 has been specifically designed to be used with most sample matrices without interference.

Example data

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