

Phosphate Buffered Saline (PBS) Casein ELISA Reagent ab171532

1 References

概述

产品名称	Phosphate Buffered Saline (PBS) Casein ELISA试剂盒
经测试应用	适用于: ELISA 不适用于: IHC-Fr or IHC-P
常规说明	<p>Abcam's Phosphate Buffered Saline (PBS) Casein ELISA Reagent consists of many phosphoserine groups that function to bind Ca^{2+}. The solution is ready for dilution to specific application requirements. ab171532 contributes to the minimization of non-specific interactions of the secondary reagents with each other and with the primary solid phase binding sites. The high blocking efficiency of casein is due to its content of small molecular weight proteins. Phosphate Buffered Saline (PBS), pH 7.4, is a poor buffer above pH 7.5. Phosphate buffer should not be used in assays where competition for phosphate groups or complex formation with a metal ion is essential for the enzyme activation. Phosphate ions will inhibit carboxypeptidase, carboxylase, urease, muscle diaminease, formase and phosphoglucomutase.</p> <p>Storage Instructions:</p> <p>This product should be clear and colorless in appearance. To minimize the risk of contamination, a fresh working solution should be prepared each day or store unused working solutions at the recommended storage temperature. If proper technique is not performed the possibility of introducing contaminants is always possible.</p> <p>Product Use:</p> <p>ab171532 should be diluted with reagent quality water. To produce a 1X working solution, dilute 1 mL of Phosphate Buffered Saline (PBS) Casein ELISA Reagent in 4 mL reagent quality water. To use as a blocking reagent, treat target surface by covering completely with 1X solution. To use as a diluent, dilute antigen/antibody to desired concentration in 1X solution and treat surface by covering completely. When using 1X solution for coating plates, best results are obtained following an overnight incubation at +4°C.</p>

性能

形式	Liquid
存放说明	Store at room temperature.
存储溶液	pH: 7 Constituent: Casein

应用

The Abpromise guarantee

Abpromise™承诺保证使用ab171532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
ELISA		Use at an assay dependent concentration.

应用说明

Is unsuitable for IHC-Fr or IHC-P.

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