

Tris Buffered Saline (TBS) Casein ELISA Reagent ab171533

概述

产品名称	Tris Buffered Saline (TBS) 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 Tris Buffered Saline (TBS) Casein ELISA Reagent consists of many phosphoserine groups that function to bind Ca^{2+}. The 10X solution is ready for dilution to specific application requirements. ab171533 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. Tris Buffered Saline (TBS), pH 7.6, is a poor buffer below pH 7.5. Tris has the disadvantages of its reactivity as a primary amine and its appreciable solubility in organic solvents, which can lead to an accumulation in the biological phase of a reaction system. To avoid complex formation with ionic species such as calcium and magnesium in blood, tris buffers are preferable over phosphate buffers.</p> <p>Storage Instructions: Store at room temperature. 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: ab171533 should be diluted 10X with reagent quality water. To produce a 1X working solution, dilute 1 mL of Tris Buffered Saline (TBS) Casein ELISA Reagent in 9 mL reagent quality water. Allow working solution to equilibrate to room temperature (25°C) prior to use. To use as a blocking reagent, treat target surface by covering completely with solution. To use working solution as a diluent, dilute antigen/antibody to desired concentration and treat surface by covering completely with solution. When solution is used for coating plates are obtained at +4°C.</p>

性能

形式	Liquid
存放说明	Please see notes section.
存储溶液	pH: 7 Constituent: Casein

应用

The Abpromise guarantee **Abpromise™**承诺保证使用ab171533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