

# Antibody Diluent (HAMA Blocker) for ELISA ab193969

SimpleStep ELISA

7 References 画像数 1

製品の概要

**製品名** Antibody Diluent (HAMA Blocker) for ELISA

**製品の概要** Antibody Diluent (HAMA Blocker) for ELISA ab193969 is used for sandwich antibody pair based assays that are susceptible to heterophilic antibody interference. This diluent is specially formulated to quantitate your proteins of interest in complex sample types, such as Human serum or Human plasma. Our formulation is shown to alleviate antibody interference due to heterophilic antibodies such as Human Anti-Mouse Antibody (HAMA). Heterophilic antibodies are a common problem in immunoassays such as ELISA, often causing false positive or false negative results. Abcam’s Antibody Diluent (HAMA Blocker) for ELISA is shown to eliminate heterophilic/HAMA interferences without compromising true signal quantitation.

The Antibody Diluent (HAMA Blocker) for ELISA solution is also known as Antibody Diluent CPI when included in our SimpleStep ELISA™ kits.

**特記事項** **DO NOT FREEZE.** Buffer should be stored at +4°C.

Instructions for use

Antibody Diluent (HAMA Blocker) for ELISA comes as a ready to use solution.

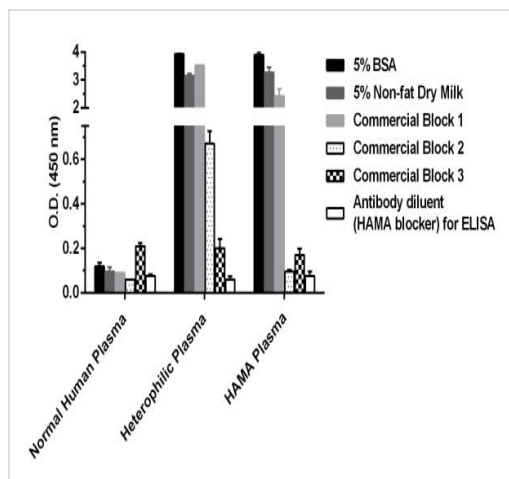
Samples should be diluted to within the desired working range of your assay in Antibody Diluent (HAMA Blocker) for ELISA.

製品の特性

**保存方法** Store at +4°C. Please refer to protocols.

内容	12 ml	50 ml
Antibody Diluent CPI - HAMA Blocker (ab193969)	1 x 12ml	1 x 50ml

画像



Example of Heterophilic Antibody Interference in Human Plasma

Example of Heterophilic Antibody Interference in Human Plasma. False positive signals are detected in a sandwich ELISA assay in Heterophilic positive and HAMA positive plasma. This interference was not observed in a control normal human plasma sample. When using both standard blocking buffers and commercially formulated heterophilic blocking buffers (1, 2, 3). Abcam's Serum and Plasma Assay Diluent outperformed its competitors by eliminating the heterophilic interference or false positive signal. Mean raw absorbance values  $\pm$  SD are graphed.

Note on above data:

Commercial Block 1 – strong false positive signal detected in both heterophilic and HAMA positive plasmas.

Commercial Block 2 – Moderate false positive signal detected in heterophilic positive plasma.

Commercial Block 3 – Elevated background (OD  $\sim$  0.2) in all human plasma samples. This high of a background signal negatively affects sensitivity of the assay.

Our Serum and Plasma Assay Diluent – Completely eliminated the heterophilic interference in all human plasma samples tested. (OD  $\sim$  0.08)

**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