

Version 2 Last updated 8 April 2020

How to use Matched Antibody Pair Kit

This product is for research use only and is not intended for diagnostic use.

Table of Contents

1. Overview	1
2. Precautions	2
3. Storage and Stability	2
4. Limitations	2
5. Materials Required, Not Supplied	3
6. Reagent Preparation	5
7. Standard Preparation	7
8. Assay Procedure	8
9. Calculations	10
10. Troubleshooting	11
11. Notes	12

1. Overview

Matched antibody pair kits include a capture and a biotinylated detector antibody pair and a calibrated protein standard.

Matched antibody pair kits and reagents deliver consistent, specific, and sensitive results.

- Batch-to-batch consistency: only recombinant monoclonal antibodies are used in our matched antibody pairs.
- Specificity: antibody pairs are screened in plasma and serum to ensure specificity in complex samples.
- Sensitivity: benchmarked against commercially available antibody pairs to ensure equivalent or superior performance compared with the competition.

Additional buffers and plates are required for the assay but not provided with the kit. An accessory pack can be purchased separately which includes buffer reagents and plates required to perform 10 x 96-well plate sandwich ELISAs (ab210905).

2. Precautions

Please read these instructions carefully prior to beginning the assay.

- Some components in this kit contain ProClin® which may cause an allergic skin reaction or respiratory irritation.
- The Stop Solution suggested for use with this kit is a concentrated acid solution and should be used with caution and adequate personal protective equipment (PPE).
- The TMB substrate suggested for use with this kit may cause skin, eye, and respiratory irritation. Wear PPE. Wash hands thoroughly after handling.
- Please review the SDS on the Abcam website prior to use.
- Capture antibody contains sodium azide, which may react with lead and copper plumbing to form explosive metallic azides. Do not dispose the leftover component directly into sink.

3. Storage and Stability

Store kit at -20°C immediately upon receipt.

4. Limitations

- The supplied kit is intended for research use only. Not for use in diagnostic procedures.
- This kit contains sufficient materials to perform multiple sandwich ELISA, provided the following conditions are met:
 - The reagents are prepared as described in the booklet.
 - The assay is run as described in the Assay Procedure.
 - The recommended reagents and solutions are used.

5. Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully perform this assay:

Recommended reagent, solution and consumables:

- Matched Antibody Pair Kit ELISA Accessory Pack (10 x 96-well plates) (ab210905): containing Nunc™ MaxiSorp™ 96-well microplates, plate seals, Coating Buffer, 10X Wash Buffer, 10X Blocking Buffer, TMB and Stop Solution.

The reagent and solutions listed above may also be purchased separately:

- Nunc™ MaxiSorp™ 96-well plates (ab210903)
- Coating buffer 1X (ab210899)
35 mM NaHCO₃, 15 mM Na₂CO₃, pH 9.6
- Blocking buffer 10X (ab210904)
Dilute in PBS to 1X: 1% BSA*, 0.05% Tween® 20, in 1X PBS, pH 7.2 - 7.4
*We recommend ≥96% purity BSA as less pure BSA can increase background
- Wash Buffer 10X (ab206977)
Dilute in water to 1X: 0.05% Tween® 20 in 1X PBS
- TMB substrate (ab210902)
- Stop solution (ab210900): 4.9% orthophosphoric acid
- 10X Phosphate Buffered Saline (PBS) (ab128983)
Dilute in water to 1X: 0.14 M NaCl, 0.003 M KCl, 0.002 M KH₂PO₄, 0.01 M Na₂HPO₄
- Streptavidin-HRP Solution (ab210901), 50 µg/mL
Recommend diluting to 0.01-0.05 µg/mL in 1X Blocking Buffer
- Microcentrifuge tubes for dilution of standards
- Double distilled water (ddH₂O)
- Optional: Protease Inhibitor Cocktail (ab65621)
- Optional: BCA Protein Quantification Kit (ab102536)

Required Equipment:

- Microplate reader capable of measuring absorbance at 450 nm.
- Multi-channel and single-channel pipettes.

Recommended Equipment

- Plate shaker or rocker.

6. Reagent Preparation

- Equilibrate all reagents to room temperature (18-25°C) prior to use.
- Allow all reagents to sit for a minimum of 15 minutes with gentle shaking after the initial reconstitution. Mix well before use.
- Dilute reagents to 1X working concentrations, if required before use.
- Reagent dilutions should be prepared and used immediately.
- Prepare only as much reagents as is needed on the day of the experiment.
- Antibody buffer information: The Capture antibody is provided in a glycerol free formulation and the Detector antibody is provided in a buffer that contains glycerol.

6.1 Capture Antibody

Capture Antibody is provided at 1 mg/mL. Dilute the Capture Antibody to the suggested working concentration of 2 µg/mL in Coating Buffer (ab210899). Add 50 µL per well.

- Δ Note:** For best results in your application, optimization of the concentration of capture antibody in Coating Buffer may be required.

6.2 Detector Antibody

Detector Antibody is provided at 0.25 mg/mL. Dilute the biotin-labeled Detector Antibody to the suggested working concentration of 0.5 µg/mL in Blocking Buffer or other appropriate diluent. Add 50 µL per well.

- Δ Note:** For best results in your application, optimization of the concentration of detector antibody in 1X Blocking Buffer may be required.

6.3 Plate Preparation

- 6.3.1 Add 50 μL of 2 $\mu\text{g}/\text{mL}$ Capture antibody to each well of a 96-well of a high bind microplate (we recommend Nunc™ Maxisorp™ 96-well plate (ab210903)).
- 6.3.2 Seal the plate with a plate seal. Incubate the plate either overnight at 4°C or for 2 hours at room temperature on a plate rocker or shaker.
- 6.3.3 Wash plate three times with 350 μL of the recommended 1X Wash Buffer (ab206977). Remove liquid completely from last wash by tapping the plate vigorously against a pad of absorbent towels.
- 6.3.4 Reduce non-specific binding by adding 300 μL of 1X Blocking Buffer (ab210904) to each well, seal the plate and incubate either overnight at 4°C or for 2 hours at room temperature.
- 6.3.5 Repeat the wash procedure in step 6.3.3.
- 6.3.6 Plate should be used immediately after blocking.

7. Standard Preparation

Prepare serially diluted standards immediately prior to use. Always prepare a fresh set of positive controls for every use.

The following section describes the preparation of a two-fold diluted standard curve for duplicate measurements (recommended).

- 7.1 Reconstitute the protein standard sample by adding 100 μL ddH₂O water. Gently mix at room temperature for 10 minutes to ensure that the protein is completely in solution. This is the **Stock Standard** Solution. Unused reconstituted protein standard should be aliquoted and stored at -80°C .

Δ Note: Refer to the vial label to see the quantity of protein standard provided.

- 7.2 Label eight tubes, Standards #1– 8.

- 7.3 To prepare Standard #1, dilute an aliquot of the Stock Standard in 1X Blocking Buffer to the highest concentration specified in the product datasheet. A seven-point standard curve using 2-fold serial dilutions in 1X Blocking Buffer is recommended.

Δ Note: Each well will require 50 μL of standard. Prepare enough standard dilutions to allow for duplicate readings.

Example of Stock standard and standard #1 preparation

Label states that 10 ng protein standard is provided.

Reconstitute in 100 μL H₂O. Stock Standard concentration = 100 ng/mL

Product states required Standard #1 concentration should be 1 ng/mL (i.e. 1000 pg/mL)

Dilute Stock standard 100X by adding 10 μL of Stock Standard to 1 mL of 1X Blocking Buffer for Standard #1

Create rest of the two-fold dilution series in 1X Blocking Buffer.

- 7.4 Standard #8 is the Blank control (buffer only) and contains no standard protein.

8. Assay Procedure

- Equilibrate all materials and prepared reagents to room temperature prior to use.
 - We recommend that you assay all standards, controls and samples in duplicate.
 - Prepare all reagents and standards as directed in the previous sections.
- 8.1 Dilute the experimental sample with 1X Blocking Buffer. Dilute the sample so that the resulting concentration is within the dynamic range of the assay. Multiple sample dilutions using 1:2 dilution series is advised if the concentration of the target protein is unknown.
 - 8.2 Add 50 μ L diluted standard and samples to each well. Seal the plate and incubate for 2 hours at room temperature on a plate shaker set to 400 rpm.
 - 8.3 Wash plate three times with 350 μ L of 1X Wash Buffer (ab206977). Remove liquid completely from last wash by tapping the plate vigorously against a pad of absorbent towels.
 - 8.4 Dilute Detector Antibody from stock concentration of 0.25 mg/mL to the suggested working concentration of 0.5 μ g/mL in 1X Blocking Buffer (ab210904) or other appropriate diluent.
 - 8.5 Add 50 μ L of diluted Detector antibody to each well. Seal the plate and incubate for 1 hour at room temperature on a plate shaker set to 400 rpm.
- Δ Note:** For best results, the concentration of detector antibody in the working solution may require optimization.
(See <http://www.abcam.com/protocols/elisa-troubleshooting-tips> for more information.)
- 8.6 Repeat wash step as described as in step 8.3.
 - 8.7 Dilute HRP-Streptavidin solution (ab210901) to 0.05 μ g/mL in 1X Blocking buffer. Add 50 μ L of diluted HRP-Streptavidin solution to each well. Seal the plate and incubate for 1 hour at room temperature on a plate shaker set to 400 rpm.

Δ **Note:** For best results, the concentration of diluted HRP streptavidin solution may require optimization.

8.8 Repeat wash step as described in step 8.3.

8.9 Add 100 μL of TMB Substrate to each well and incubate for **up to 20 minutes** in the dark on a plate shaker set to 400 rpm.

Δ **Note:** For best results, the incubation time requires optimization.

8.10 Before add Stop Solution, plate can be read kinetically at 600nm to monitor proper incubation time.

8.11 Add 100 μL of Stop Solution to each well. Shake plate on a plate shaker at 400 rpm for 1 minute to mix.

(See <http://www.abcam.com/protocols/elisa-troubleshooting-tips> for more information)

8.12 Measure the endpoint of the plate at 450 nm.

8.13 Analyze the data as described below.

9. Calculations

Calculate the average absorbance value for the blank control (zero) standards. Subtract the average blank control standard absorbance value from all other absorbance values.

Create a standard curve by plotting the average blank control subtracted absorbance value for each standard concentration (y-axis) against the target protein concentration (x-axis) of the standard. Use graphing software to draw the best smooth curve through these points to construct the standard curve. Note: Most microplate reader software or graphing software will plot these values and fit a curve to the data. A four-parameter curve fit (4PL) is often the best choice; however, other algorithms (e.g. linear, semi-log, log/log, 5-parameter logistic) can also be tested to determine if it provides a better curve fit to the standard values.

See our webinar www.abcam.com/kits/calculate-accurate-protein-concentrations for more information on calculating ELISA raw data.

See our ELISA optimization guide: www.abcam.com/kits/technical-tips-for-elisa-and-multiplex-immunoassay-development to design your own sandwich ELISA.

10. Troubleshooting

Problem	Reason	Solution
Poor standard curve	Inaccurate Pipetting	Check pipettes
	Improper standard dilution	Prior to opening, briefly spin the stock standard tube and dissolve the powder thoroughly by gentle mixing
Low Signal	Incubation times too brief	Ensure sufficient incubation times; increase to 2 or 3-hour standard/sample incubation
	Inadequate reagent volumes or improper dilution	Check pipettes and ensure correct preparation
	Incubation times with TMB too brief	Ensure sufficient incubation time until blue color develops prior addition of Stop solution
Large CV	Plate is insufficiently washed	Review manual for proper wash technique. If using a plate washer, check all ports for obstructions.
	Contaminated wash buffer	Prepare fresh wash buffer
Low sensitivity	Improper storage of the ELISA kit	Aliquot and store your reconstituted standards at -80°C, avoid freeze/thaw. Store antibodies at -20°C. Keep TMB substrate solution protected from light.

11. Notes

Nunc™ MaxiSorp™ are unregistered trademarks owned by Thermo Fisher Scientific.
TWEEN® is a registered trademark of Croda International PLC
PROCLIN® is a registered trademark of The Dow Chemical Company or an affiliated company of
Dow

Technical Support

Copyright © 2020 Abcam, All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

www.abcam.com/contactus

www.abcam.cn/contactus (China)

www.abcam.co.jp/contactus (Japan)