

Product datasheet

Human AACT Matched Antibody Pair Kit ab221445

1 Image

Overview

<b>Product name</b>	Human AACT Matched Antibody Pair Kit
<b>Detection method</b>	Colorimetric
<b>Assay type</b>	ELISA set
<b>Range</b>	500 pg/ml - 32000 pg/ml
<b>Species reactivity</b>	<b>Reacts with:</b> Human
<b>Product overview</b>	Matched Antibody Pair Kits include a capture and a biotinylated detector antibody pair, along with a calibrated protein standard, suitable for sandwich ELISA. The Matched Antibody Pair Kit can be used to quantify native and recombinant human AACT.

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. An accessory pack can be purchased which includes buffer reagents required to perform 10 x 96-well plate sandwich ELISAs ([ab210905](#)).

For additional information on the performance of the antibody pair used in this kit, please see our equivalent SimpleStep ELISA® ([ab217779](#)) which uses the same antibody pair.

**To receive an electronic copy of the Certificate of Analysis, please send an [email](#) with "CoA for matched antibody pair kit" in the subject line and the desired product number and lot number in the body of the email.**

<b>Tested applications</b>	<b>Suitable for:</b> ELISA
<b>Platform</b>	Reagents

## Properties

**Storage instructions** Store at -20°C. Please refer to protocols.

Components	10 x 96 tests	2 x 96 tests
Human AACT Capture Antibody	1 x 100µg	1 x 20µg
Human AACT Detector Antibody	1 x 25µg	1 x 5µg
Human AACT Lyophilized Protein	1 vial	1 vial

**Function** Although its physiological function is unclear, it can inhibit neutrophil cathepsin G and mast cell chymase, both of which can convert angiotensin-1 to the active angiotensin-2.

**Tissue specificity** Plasma. Synthesized in the liver. Like the related alpha-1-antitrypsin, its concentration increases in the acute phase of inflammation or infection. Found in the amyloid plaques from the hippocampus of Alzheimer disease brains.

**Involvement in disease** Defects in SERPINA3 may be a cause of chronic obstructive pulmonary disease (COPD) [MIM:107280].

**Sequence similarities** Belongs to the serpin family.

**Domain** The reactive center loop (RCL) extends out from the body of the protein and directs binding to the target protease. The protease cleaves the serpin at the reactive site within the RCL, establishing a covalent linkage between the carboxyl group of the serpin reactive site and the serine hydroxyl of the protease. The resulting inactive serpin-protease complex is highly stable.

**Cellular localization** Secreted.

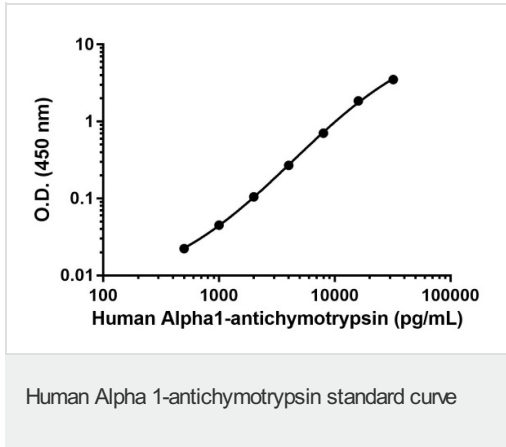
## Applications

Our [Abpromise guarantee](#) covers the use of **ab221445** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
ELISA		Use at an assay dependent concentration.

## Images



Standard calibration curve. Background subtracted values graphed.

**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.com/abpromise> or contact our technical team.

### Terms and conditions

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors