

## Product datasheet

# Human Annexin A1 ELISA Kit ab222868

[1 Image](#)

### Overview

**Product name** Human Annexin A1 ELISA Kit

**Detection method** Colorimetric

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
Plasma 1	20			5%
Plasma 2	20			5.3%
Plasma 3	20			5.9%

Inter-assay

Sample	n	Mean	SD	CV%
Plasma 1	20			9.5%
Plasma 2	20			9.7%
Plasma 3	20			10.4%

**Sample type** Cell culture supernatant, Serum, Plasma, Cell Lysate

**Assay type** Sandwich (quantitative)

**Sensitivity** 0.07 ng/ml

**Range** 0.313 ng/ml - 20 ng/ml

**Recovery** 97 %

**Assay time** 4h 00m

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Human

**Does not react with:** Rabbit, Cow

**Product overview** ab222868 Human Annexin A1 ELISA Kit is designed for the quantitative measurement of Annexin A1 in plasma, serum, cell lysate and cell culture samples.

The kit employs a quantitative sandwich enzyme immunoassay technique that measures human Annexin A1 in approximately 4 hours. A polyclonal antibody specific for human Annexin A1 has been pre-coated onto a 96-well microplate with removable strips. Annexin A1 in standards and samples is sandwiched by the immobilized antibody and biotinylated polyclonal antibody specific for Annexin A1, which is recognized by a streptavidin-peroxidase conjugate. All unbound material is washed away and a peroxidase enzyme substrate is added. The color development is stopped and the intensity of the color is measured.

**The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.**

**Notes**

Annexin A1, also known as lipocortin I, annexin I, chromobindin-9, and calpactin II, belongs to the annexin family of calcium-dependent phospholipid-binding proteins. The annexin proteins are structurally characterized by a core domain containing a conserved 70 amino acid sequence repeated four or eight times. This core is preceded by a unique N-terminal domain that varies in sequence and length and is thought to be responsible for specifying different annexin functions. Annexin A1 has an apparent relative molecular mass of 40 kDa with 346 amino acids and is preferentially located on the cytosolic face of the plasma membrane. It mimics the effect of steroids, mediates anti-inflammatory activity and is a potent inhibitor of phospholipase A2 activity. Annexin A1 regulates neutrophil extravasation by interacting with the formyl peptide receptor and triggers different signaling pathways.

**Platform**

Pre-coated microplate (12 x 8 well strips)

**Properties**

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

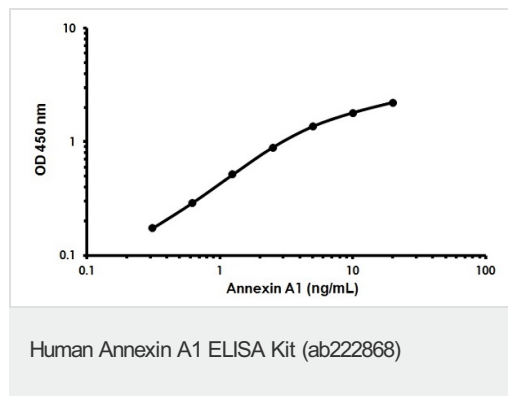
Components	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent N Concentrate	1 x 30ml
1X Standard Diluent	1 x 2ml
20X Wash Buffer Concentrate	2 x 30ml
30x Biotinylated Human Annexin A1	1 x 180µl
Anti-Human Annexin A1 coated Microplate (12 x 8 wells)	1 unit
Chromogen Substrate	1 x 8ml
Human Annexin A1 Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 12ml

**Function**

Calcium/phospholipid-binding protein which promotes membrane fusion and is involved in exocytosis. This protein regulates phospholipase A2 activity. It seems to bind from two to four calcium ions with high affinity.

<b>Sequence similarities</b>	Belongs to the annexin family. Contains 4 annexin repeats.
<b>Domain</b>	A pair of annexin repeats may form one binding site for calcium and phospholipid.
<b>Post-translational modifications</b>	Phosphorylated by protein kinase C, epidermal growth factor receptor/kinase and TRPM7. Phosphorylation results in loss of the inhibitory activity.
<b>Cellular localization</b>	Nucleus. Cytoplasm. Cell projection > cilium. Basolateral cell membrane. Found in the cilium, nucleus and basolateral cell membrane of ciliated cells in the tracheal endothelium (By similarity). Found in the cytoplasm of type II pneumocytes and alveolar macrophages.

## Images



Example of human Annexin A1 standard curve.

**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