

### Mouse SAP ELISA Kit (PTX2) ab213888

[1 Image](#)

#### Overview

**Product name** Mouse SAP ELISA Kit (PTX2)

**Detection method** Colorimetric

#### Precision

Intra-assay

Sample	n	Mean	SD	CV%
1	16	3.608ng/ml	0.144	= 4%
2	16	12.778ng/ml	0.715	= 5.6%
3	16	47.613ng/ml	2.809	= 5.9%

Inter-assay

Sample	n	Mean	SD	CV%
1	24	3.726ng/ml	0.176	= 4.7%
2	24	13.449ng/ml	0.981	= 7.3%
3	24	43.062ng/ml	2.626	= 6.1%

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

**Assay type** Sandwich (quantitative)

**Sensitivity** < 50 pg/ml

**Range** 1.56 ng/ml - 100 ng/ml

**Assay time** 3h 30m

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Mouse

**Product overview** The Mouse SAP Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213888) is designed for the quantitative measurement of Mouse SAP in cell culture supernatants, cell lysates, serum and plasma (heparin, EDTA).

The ELISA kit is based on standard sandwich enzyme-linked immunosorbent assay technology. A

monoclonal antibody from rat specific for SAP has been pre-coated onto 96-well plates. Standards (Expression system for standard: NSO; Immunogen sequence: Q21- D224) and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for SAP is added subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex was added and unbound conjugates were washed away with PBS or TBS buffer. HRP substrate TMB was used to visualize HRP enzymatic reaction. TMB was catalyzed by HRP to produce a blue color product that changed into yellow after adding acidic stop solution. The density of yellow is proportional to the Mouse Serum Amyloid P / SAP amount of sample captured in plate.

**Notes** SAP also known as PTX2 or APCS, is the identical serum form of amyloid P component (AP), a 25kDa pentameric protein first identified as the pentagonal constituent of *in vivo* pathological deposits called "amyloid". It belongs to the pentraxins family, characterized by calcium dependent ligand binding and distinctive flattened beta-jellyroll structure similar to that of the legume lectins. This gene is mapped to 1q23.2. The binding of the encoded protein to proteins in the pathological amyloid cross-beta fold suggests its possible role as a chaperone. This protein is also thought to control the degradation of chromatin. It has been demonstrated that this protein binds to apoptotic cells at an early stage, which raises the possibility that it is involved in dealing with apoptotic cells *in vivo*.

**Platform** Pre-coated microplate (12 x 8 well strips)

**Properties**

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

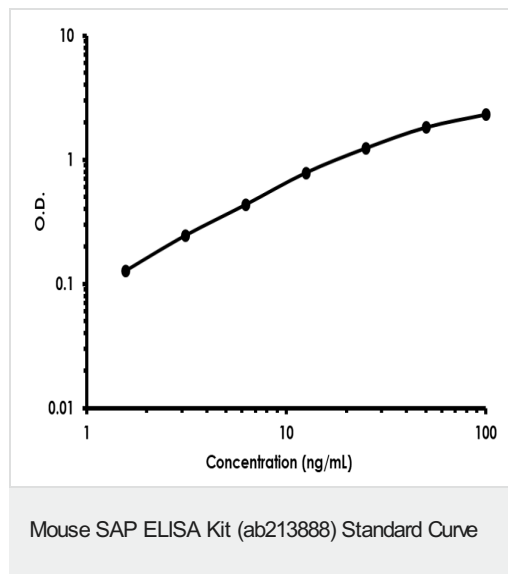
Components	Identifier	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml
Adhesive Plate Seal		4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml
Anti-Mouse SAP coated Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl
Biotinylated anti- Mouse SAP antibody		1 x 100µl
Lyophilized recombinant Mouse SAP standard		2 vials
Sample Diluent Buffer	Green Cap	1 x 30ml
TMB Color Developing Agent	Black Cap	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml
Wash Buffer (25X)		1 x 20ml

**Function** Can interact with DNA and histones and may scavenge nuclear material released from damaged circulating cells. May also function as a calcium-dependent lectin.

**Tissue specificity** Found in serum and urine.

<b>Involvement in disease</b>	Note=SAP is a precursor of amyloid component P which is found in basement membrane and associated with amyloid deposits.
<b>Sequence similarities</b>	Belongs to the pentaxin family. Contains 1 pentaxin domain.
<b>Post-translational modifications</b>	N-glycosylated with a complex biantennary oligosaccharide chain with a sialic acid at the end (disialo-SAP). Monosialo-SAP as well as asialo-SAP are also detected (PubMed:15174148).
<b>Cellular localization</b>	Secreted.

## Images



Mouse Serum Amyloid P / SAP ELISA Kit (ab213888) 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