



Product datasheet

Human SAP ELISA Kit (APCS) ab137970

★★★★★ [1 Abreviews](#) [2 References](#) [2 Images](#)

Overview

Product name Human SAP ELISA Kit (APCS)

Detection method Colorimetric

Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Overall				= 4.4%

	Inter-assay				
	Sample	n	Mean	SD	CV%
	Overall				= 9.2%

Sample type Cell culture supernatant, Urine, Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity > 24 pg/ml

Range 0.078 ng/ml - 5 ng/ml

Recovery 98 %

	Sample specific recovery		
	Sample type	Average %	Range
	Serum		97% - 104%
	Plasma		88% - 109%

Assay time 4h 00m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Human SAP (APCS) *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of human SAP in plasma, serum, urine, and cell culture supernatants.

A SAP specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a SAP specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Conjugate is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize Streptavidin-Peroxidase enzymatic reaction. TMB is catalyzed by Streptavidin-Peroxidase to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the amount of SAP captured in plate.

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

Platform Microplate

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
20X Wash Buffer Concentrate	2 x 30ml
50X Biotinylated Human SAP Antibody	1 x 120µl
Chromogen Substrate	1 x 7ml
SAP Microplate (12 x 8 well strips)	1 unit
SAP Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11ml

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.

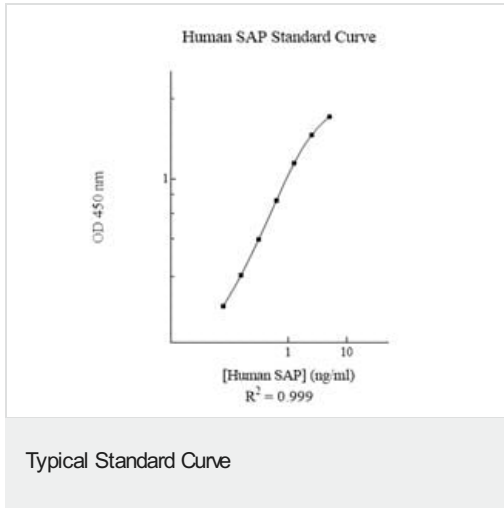
Involvement in disease Note=SAP is a precursor of amyloid component P which is found in basement membrane and associated with amyloid deposits.

Sequence similarities Belongs to the pentaxin family.
Contains 1 pentaxin domain.

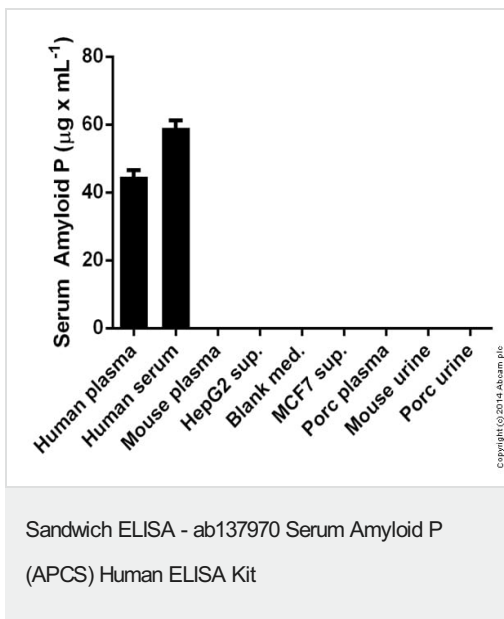
Post-translational modifications N-glycosylated with a complex biantennary oligosaccharide chain with a sialic acid at the end (disialo-SAP). Monosialo-SAP as well as asialo-SAP are also detected (PubMed:15174148).

Cellular localization Secreted.

Images



Representative Standard Curve using ab137970.



SAP measured in various samples showing quantity (microgram) per mL of tested sample

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