

Human Alpha 1 acid Glycoprotein / AGP ELISA Kit (AGP) ab108854

8 References 2 Images

Overview

Product name	Human Alpha 1 acid Glycoprotein / AGP ELISA Kit (AGP)			
Detection method	Colorimetric			
Precision	Intra-assay			
	Sample	n	Mean	SD
	Overall			5.3%
	Inter-assay			
	Sample	n	Mean	SD
	Overall			11.1%
Sample type	Serum, Plasma			
Assay type	Competitive			
Sensitivity	= 94 ng/ml			
Range	0.25 µg/ml - 2 µg/ml			
Recovery	= 98 %			
Assay time	3h 00m			
Assay duration	Multiple steps standard assay			
Species reactivity	Reacts with: Human			
Product overview	Abcam's Alpha 1 acid Glycoprotein / AGP Human <i>in vitro</i> competitive ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Alpha 1 acid Glycoprotein / AGP in plasma and serum.			

An Alpha 1 acid Glycoprotein / AGP specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently biotinylated Alpha 1 acid Glycoprotein / AGP 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 inversely proportional to the amount of Alpha 1 acid Glycoprotein / AGP 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
1X Biotinylated Human Alpha 1 acid Glycoprotein / AGP (Lyophilized)	1 vial
20X Wash Buffer Concentrate	1 x 30ml
Alpha 1 acid Glycoprotein / AGP Microplate (12 x 8 well strips)	1 unit
Alpha 1 acid Glycoprotein / AGP Standard	1 vial
Chromogen Substrate	1 x 7ml
Sealing Tapes	3 units
Stop Solution	1 x 11ml

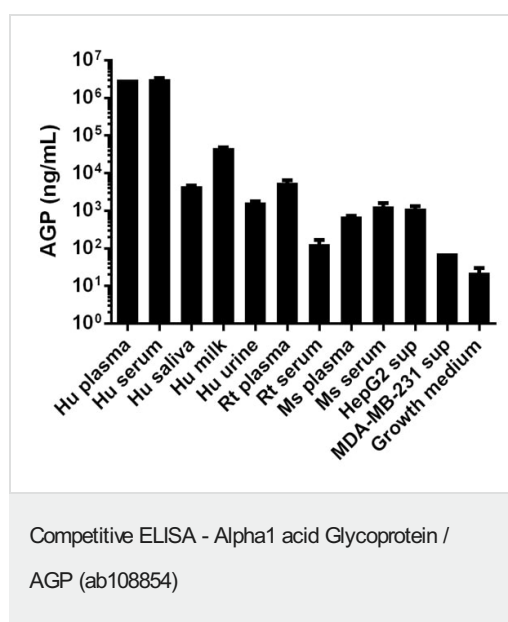
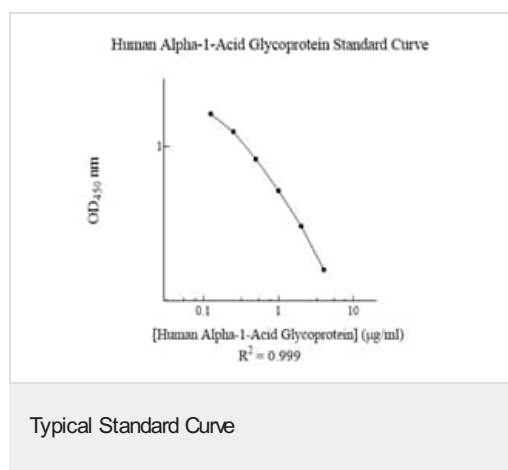
Function Appears to function in modulating the activity of the immune system during the acute-phase reaction.

Tissue specificity Expressed by the liver and secreted in plasma.

Sequence similarities Belongs to the calycin superfamily. Lipocalin family.

Cellular localization Secreted.

Images



Alpha1 acid Glycoprotein / AGP measured in various biological fluids showing quantity (ng) per mL of sample tested

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