

Human GASP-1 ELISA Kit ab267602

1 Image

Overview

Product name	Human GASP-1 ELISA Kit				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 10%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 12%
Sample type	Serum, Plasma, Cell culture media				
Assay type	Sandwich (quantitative)				
Sensitivity	20 pg/ml				
Range	20.48 pg/ml - 5000 pg/ml				
Recovery	Sample specific recovery				
	Sample type		Average %		Range
	Serum		108		92% - 127%
	Plasma		101		89% - 123%
	Cell culture media		94		80% - 108%
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Human				
Product overview	Human GASP-1 ELISA Kit is designed for the quantitative determination of GASP-1 in cell culture supernatants, plasma and serum samples.				

This assay employs an antibody specific for human GASP-1 coated on a 96-well plate. Standards and samples are pipetted into the wells and GASP-1 present in a sample is bound to the wells by

the immobilized antibody. The wells are washed and biotinylated anti-human GASP-1 antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of GASP-1 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Cross Reactivity: This ELISA kit shows no cross-reactivity with the following cytokines tested: human ACE-2, aFGF, Albumin, AMICA, Angiopoietin-4, BAFF, CA19-9, CD163, Clusterin, CRTAM, CXCL14, Cystatin C, Decorin, Dkk-3, DLL1, Fetuin A, FOLR1, Furin, GASP-2, GCSF R, HAI-2, IL-17 RB, IL-27, LAG-3, LDL R, Pepsinogen 1, RANK, RBP4, SOST, Syndecan-1, TACI, TFPI, Thrombospondin 1, TRAIL R2, TRANCE, Troponin I, uPA, VECadherin, WISP-1.

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
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent	1 x 15ml
Anti-Human GASP-1 coated Microplate (12 x 8 wells)	1 unit
Biotinylated Anti-Human GASP-1 Detection Antibody	2 vials
Human GASP-1 Standard (Lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB Substrate Solution	1 x 12ml

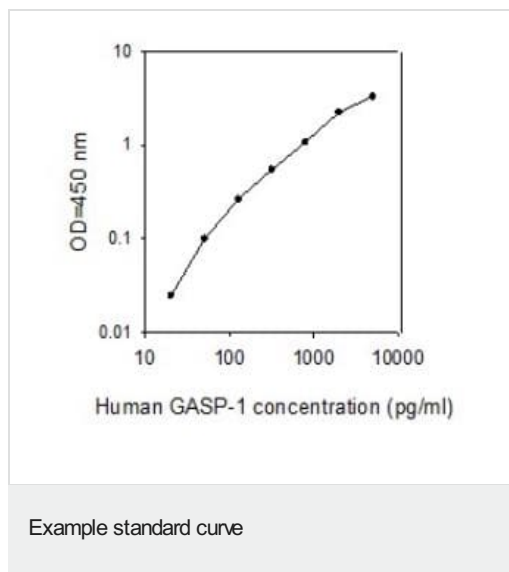
Function May modulate lysosomal sorting and functional down-regulation of a variety of G-protein coupled receptors. Targets receptors for degradation in lysosomes.

Tissue specificity Expressed in the brain, with lower expression in medulla, spinal cord and substantia nigra.

Sequence similarities Belongs to the GPRASP family.

Cellular localization Cytoplasm.

Images



Human GASP-1 Standard curve.

These standard curves are for demonstration only. A standard curve must be run with each assay

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