

Product datasheet

Human Glypican 1 ELISA Kit ab267650

[1 Image](#)

Overview

Product name	Human Glypican 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	Cell culture supernatant, Serum, Plasma				
Assay type	Sandwich (qualitative)				
Sensitivity	0.15 ng/ml				
Range	0.164 ng/ml - 40 ng/ml				
Recovery	Sample specific recovery				
	Sample type	Average %	Range		
	Cell culture supernatant	96.14	86% - 102%		
	Serum	101.4	91% - 112%		
	Plasma	95.11	85% - 102%		
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Human				
Product overview	Human Glypican 1 ELISA Kit is designed for the quantitative determination of Glypican 1 in cell culture supernatants, plasma and serum samples.				

This assay employs an antibody specific for human Glypican 1 coated on a 96-well plate.

Standards and samples are pipetted into the wells and Glypican 1 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-human Glypican 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 Glypican 1 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

This ELISA kit shows no cross-reactivity with the following cytokines tested: human ALK-1, B7-H2, BLAME, BMP-8, CD28, Common beta Chain, Contactin-1, Desmoglein-1, Desmoglein-3, EDAR, EphA1, EphB6, Ephrin-B3, Epregrulin, FGF-12, FGF17, FOLR2, Galectin-8, Glypican 5, GHR, IFN-gamma R1, IL-22 R alpha 1, IL-22BP, IL-23 R, IL-31 RA, IL-7 R alpha, Integrin alpha 5, MDM2, Nectin-1, NKp30, Nogo, Receptor, Notch-3, OSM R beta, Prolactin R, RELT, Ryk, Semaphorin 6D, Semaphorin 7A, Siglec-11.

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 Glypican 1 coated Microplate (12 x 8 wells)	1 unit
Biotinylated Anti-Human Glypican 1 Detection Antibody	2 vials
Human Glypican 1 Standard (Lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB Substrate Solution	1 x 12ml

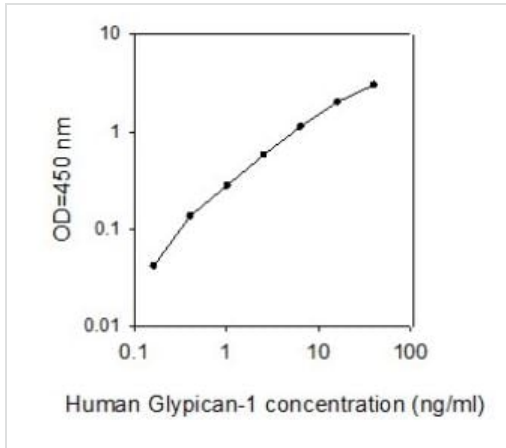
Function Cell surface proteoglycan that bears heparan sulfate.

Sequence similarities Belongs to the glypican family.

Post-translational modifications This cell-associated glypican is further processed to give rise to a medium-released species.

Cellular localization Cell membrane and Secreted > extracellular space.

Images



This standard curve is for demonstration only. A standard curve must be run with each assay.

Example 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