

Human Platelet Glycoprotein IIb + IIIa ELISA Kit

ab108851

[6 References](#) [1 Image](#)

Overview

Product name	Human Platelet Glycoprotein IIb + IIIa ELISA Kit				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Overall				4.4%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Overall				8.2%
Sample type	Plasma, Platelets, Cell Lysate				
Assay type	Sandwich (quantitative)				
Sensitivity	= 0.5 ng/ml				
Range	1.25 ng/ml - 40 ng/ml				
Recovery	101 %				
Assay time	4h 00m				
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Human				
Product overview	Abcam's Platelet Glycoprotein IIb + IIIa Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Platelet Glycoprotein IIb+IIIa concentrations in Human platelets, platelet-rich plasma and cell culture lysate.				

A Platelet Glycoprotein IIb + IIIa specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a Platelet Glycoprotein IIb+IIIa 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 Platelet Glycoprotein IIb + IIIa 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

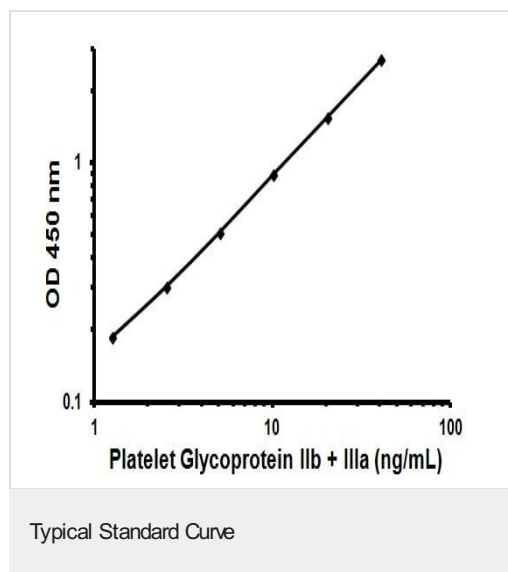
Properties

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 Platelet Glycoprotein IIb + IIIa Antibody	1 x 120µl
Chromogen Substrate	1 x 7ml
Platelet Glycoprotein IIb + IIIa Microplate (12 x 8 well strips)	1 unit
Platelet Glycoprotein IIb + IIIa Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11ml

Images



Representative Standard Curve using ab108851.

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