

Human GC Globulin ELISA Kit (Vitamin D Binding Protein) ab108853

5 References 画像数 1

製品の概要

製品名	Human GC Globulin ELISA Kit (Vitamin D Binding Protein)			
検出方法	Colorimetric			
再現性	Intra-Assay (同時再現性)			
	サンプル	N	平均値	SD
	Overall			CV%
				6.2%
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	Overall			CV%
				9.9%
サンプルの種類	Serum, Plasma			
アッセイタイプ	Competitive			
検出感度	> 0.06 µg/ml			
検出範囲	0.098 µg/ml - 100 µg/ml			
添加回収試験	97 %			
全工程の試験時間	3h 00m			
ステップ	Multiple steps standard assay			
種交差性	交差種: Human			
製品の概要	Abcam's Human GC Globulin (Vitamin D Binding Protein) <i>in vitro</i> competitive ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of GC Globulin in plasma, serum, saliva, milk and cell-culture supernatants.			

A GC Globulin specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently biotinylated GC Globulin 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 GC Globulin captured in plate.

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

試験プラットフォーム

Microplate

製品の特性

保存方法

Store at -20°C. Please refer to protocols.

内容	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent M Concentrate	1 x 30ml
20X Wash Buffer Concentrate	1 x 30ml
2X Biotinylated Human GC Globulin (Lyophilized)	1 vial
Chromogen Substrate	1 x 7ml
GC Globulin Microplate (12 x 8 well strips)	1 unit
GC Globulin Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11ml

機能

Multifunctional protein found in plasma, ascitic fluid, cerebrospinal fluid, and urine and on the surface of many cell types. In plasma, it carries the vitamin D sterols and prevents polymerization of actin by binding its monomers. DBP associates with membrane-bound immunoglobulin on the surface of B-lymphocytes and with IgG Fc receptor on the membranes of T-lymphocytes.

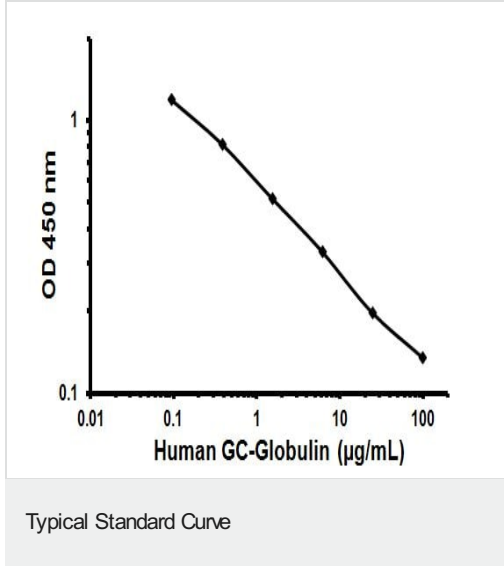
配列類似性

Belongs to the ALB/AFP/VDB family.
Contains 3 albumin domains.

細胞内局在

Secreted.

画像



Representative Standard Curve using ab108853.

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors