

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

[5 References](#) [1 图像](#)

概述

产品名称 人GC Globulin ELISA试剂盒(Vitamin D Binding蛋白)

检测方法 Colorimetric

精确度 批次内

样品	n	Mean	SD	CV%
Overall				6.2%

批次间

样品	n	Mean	SD	CV%
Overall				9.9%

样品类型 Serum, Plasma

检测类型 Competitive

灵敏度 > 0.06 µg/ml

范围 0.098 µg/ml - 100 µg/ml

回收率 97 %

检测时间 3h 00m

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

产品概述 Abcam's Human GC Globulin (Vitamin D Binding Protein) *in vitro* 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.

平台 Microplate

性能

存放说明 Store at -20°C. Please refer to protocols.

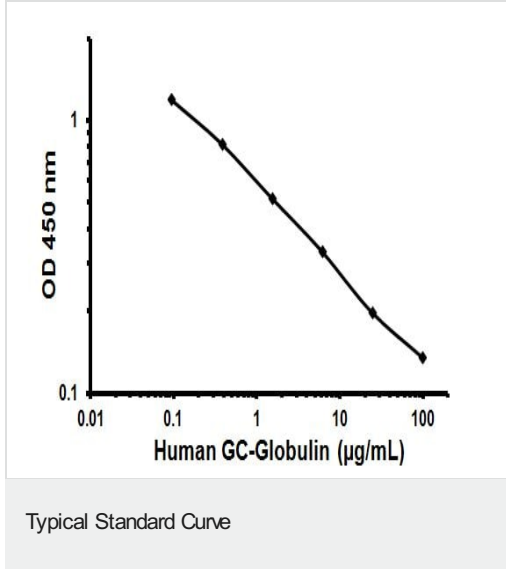
组件	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent M Concentrate	1 x 30ml
20X Wash Buffer Concentrate	1 x 30ml
2X Biotinylated Human GC Globulin (Lyophilized)	1 vial
Chromogen Substrate	1 x 7ml
GC Globulin Microplate (12 x 8 well strips)	1 unit
GC Globulin Standard	1 vial
Sealing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

功能 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.

序列相似性 Belongs to the ALB/AFP/VDB family.
Contains 3 albumin domains.

细胞定位 Secreted.

图片



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.cn/abpromise> or contact our technical team.

Terms and conditions

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