

Human beta 2 Microglobulin ELISA Kit ab108885

[5 References](#) [4 图像](#)

概述

产品名称 人beta 2 Microglobulin ELISA试剂盒

检测方法 Colorimetric

精确度 批次内

样品	n	Mean	SD	CV%
Overall				5.3%

批次间

样品	n	Mean	SD	CV%
Overall				10.3%

样品类型 Cell culture supernatant, Saliva, Milk, Urine, Serum, Plasma

检测类型 Sandwich (quantitative)

灵敏度 = 33 pg/ml

范围 0.19 ng/ml - 12.5 ng/ml

回收率 97 %

检测时间 4h 00m

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

产品概述 Abcam's beta 2 Microglobulin *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of beta 2 Microglobulin levels in plasma, serum, milk, saliva, urine, CSF and cell culture supernatants.

A beta 2 Microglobulin specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a beta 2 Microglobulin 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 beta 2 Microglobulin captured in plate.

Get results in 90 minutes with Human Beta-2-Microglobulin ELISA Kit ([ab181423](#)) from our SimpleStep ELISA® range.

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 Biotinylated beta 2 Microglobulin Antibody	1 x 60µl
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent N Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
beta 2 Microglobulin Microplate (12 x 8 well strips)	1 unit
beta 2 Microglobulin Standard	2 vials
Chromogen Substrate	1 x 7ml
Sealing Tapes	3 units
Stop Solution	1 x 11ml

功能 Component of the class I major histocompatibility complex (MHC). Involved in the presentation of peptide antigens to the immune system.

疾病相关 Defects in B2M are the cause of hypercatabolic hypoproteinemia (HYCATHYP) [MIM:241600]. Affected individuals show marked reduction in serum concentrations of immunoglobulin and albumin, probably due to rapid degradation.
Note=Beta-2-microglobulin may adopt the fibrillar configuration of amyloid in certain pathologic states. The capacity to assemble into amyloid fibrils is concentration dependent. Persistently high beta(2)-microglobulin serum levels lead to amyloidosis in patients on long-term hemodialysis.

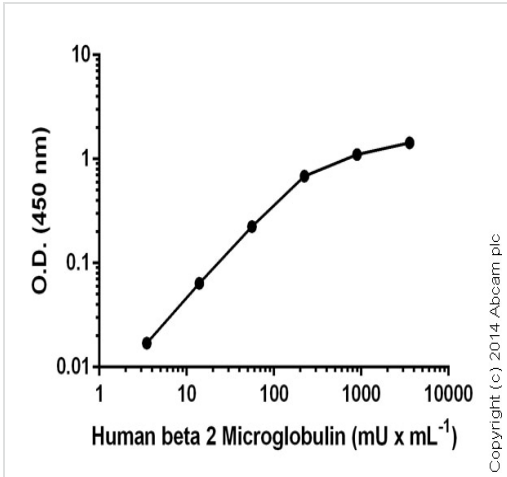
序列相似性 Belongs to the beta-2-microglobulin family.

Contains 1 Ig-like C1-type (immunoglobulin-like) domain.

翻译后修饰 Glycation of Ile-21 is observed in long-term hemodialysis patients.

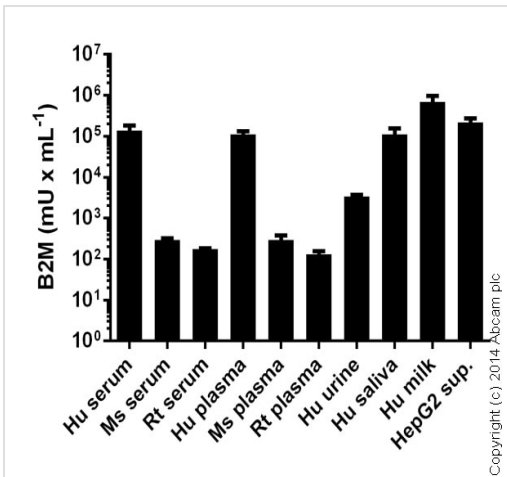
细胞定位 Secreted. Detected in serum and urine.

图片



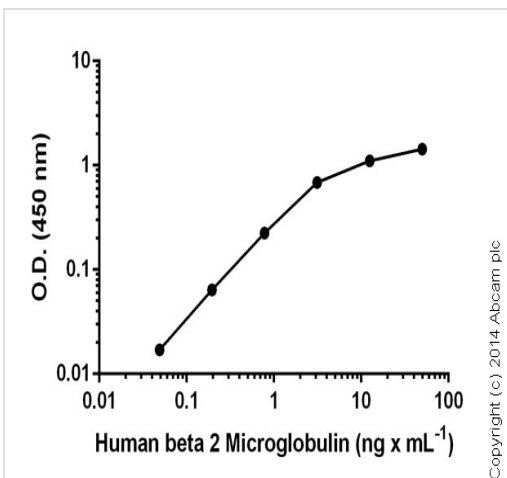
Standard curve: mean of duplicates (+/- SD) with background reads subtracted

Sandwich ELISA - beta 2 Microglobulin Human
ELISA Kit (ab108885)



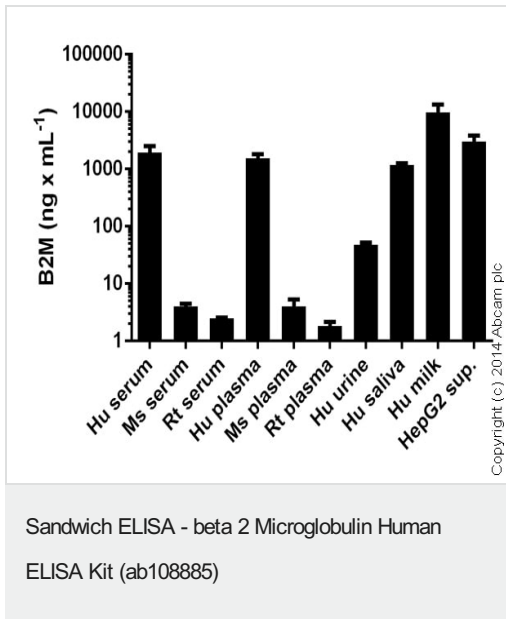
Beta 2 Microglobulin measured in biological fluids and cell culture supernatants showing quantity (mU) per mL of tested sample. Human serum and plasma samples diluted 1000-8000 fold. Human milk diluted 4000-32000 fold. Human urine diluted 100-200 fold. Human saliva diluted 400-1600 fold. Mouse and rat samples diluted 10 fold. HepG2 supernatant diluted 1000 fold.

Sandwich ELISA - beta 2 Microglobulin Human
ELISA Kit (ab108885)



Standard curve: mean of duplicates (+/- SD) with background reads subtracted

Sandwich ELISA - beta 2 Microglobulin Human
ELISA Kit (ab108885)



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