

Human beta 2 Microglobulin ELISA Kit ab108885

5 References 画像数 4

製品の概要

製品名	Human beta 2 Microglobulin ELISA Kit			
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%
				5.3%
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	Overall			CV%
				10.3%
サンプルの種類	Cell culture supernatant, Saliva, Milk, Urine, Serum, Plasma			
アッセイタイプ	Sandwich (quantitative)			
検出感度	= 33 pg/ml			
検出範囲	0.19 ng/ml - 12.5 ng/ml			
添加回収試験	97 %			
全工程の試験時間	4h 00m			
ステップ	Multiple steps standard assay			
種交差性	交差種: Human			
製品の概要	Abcam's beta 2 Microglobulin <i>in vitro</i> 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.

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 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

機能

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

関連疾患

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.

配列類似性

Belongs to the beta-2-microglobulin family.

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

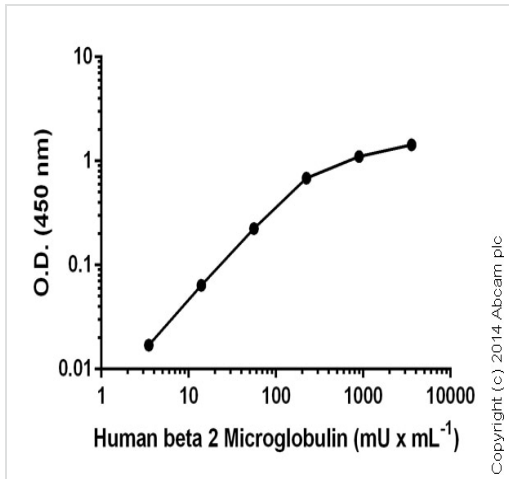
翻訳後修飾

Glycation of Ile-21 is observed in long-term hemodialysis patients.

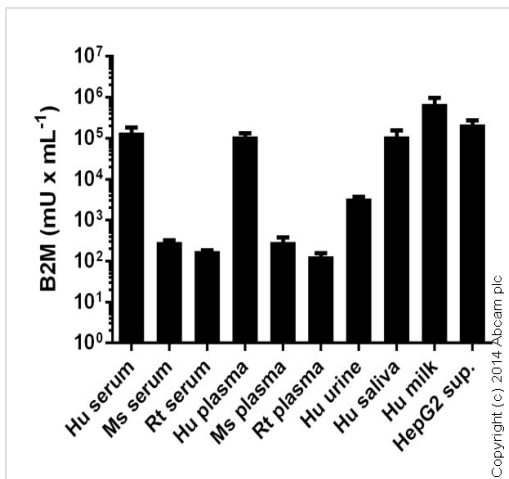
細胞内局在

Secreted. Detected in serum and urine.

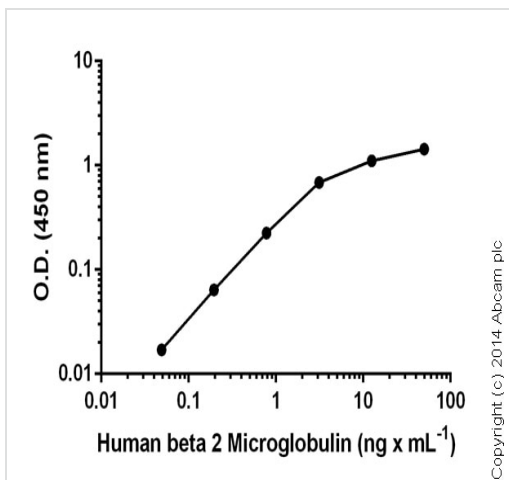
画像



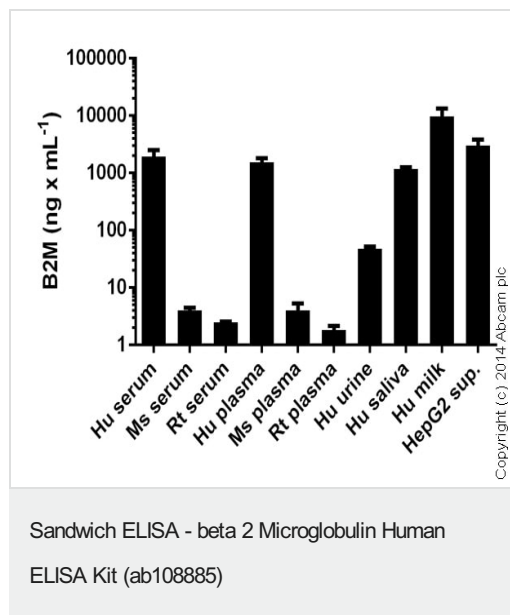
Sandwich ELISA - beta 2 Microglobulin Human
ELISA Kit (ab108885)



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



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.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