

Human CD44 ELISA Kit ab45912

7 References 画像数 2

製品の概要

製品名	Human CD44 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	CV%
	Sample A	8	9.85ng/ml	1.55	1.52%
	Sample B	8	10.5ng/ml	0.055	4.76%
	Inter-Assay (日差再現性)				
	サンプル	N	平均値	SD	CV%
	Sample A	50	9.95ng/ml	0.307	3.08%
	Sample B	50	1.07ng/ml	0.105	9.34%
サンプルの種類	Cell culture supernatant, Serum, Plasma				
アッセイタイプ	Sandwich (quantitative)				
検出感度	< 0.113 ng/ml				
検出範囲	0.31 ng/ml - 10 ng/ml				
全工程の試験時間	2h 15m				
ステップ	Multiple steps standard assay				
種交差性	交差種: Human				
製品の概要	Abcam's CD44 Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for				

A monoclonal antibody specific for CD44 has been coated onto the wells of the microtiter strips provided. Samples, including standards of known CD44 concentrations, control specimens or unknowns are pipetted into these wells. During the first incubation, the standards or samples and a biotinylated monoclonal antibody specific for CD44 are simultaneously incubated. After washing, the enzyme Streptavidin-HRP, that binds the biotinylated antibody is added, incubated and washed. A TMB substrate solution is added which acts on the bound enzyme to induce a

colored reaction product. The intensity of this colored product is directly proportional to the concentration of CD44 present in the samples.

This kit will recognize both endogenous and recombinant Human CD44.

試験プラットフォーム

Microplate

製品の特性

保存方法

Store at +4°C. Please refer to protocols.

内容	ラベル	2 x 96 tests	1 x 96 tests
10X Standard Diluent Buffer	Black	1 x 25ml	1 x 15ml
200X Wash Buffer	White	2 x 10ml	1 x 10ml
Biotinylated Antibody Diluent	Red	1 x 13ml	1 x 7ml
Biotinylated anti-CD44 antibody	Red	2 x 0.4ml	1 x 0.4ml
CD44 Control	Silver	4 vials	2 vials
CD44 Microplate (12 x 8 well strips)		2 units	1 unit
CD44 Standard	Yellow	4 vials	2 vials
Chromogen TMB Substrate Solution		1 x 24ml	1 x 11ml
HRP Diluent	Red	1 x 23ml	1 x 12ml
Plastic Plate Covers		4 units	2 units
Stop Reagent	Black	2 x 11ml	1 x 11ml
Streptavidin-HRP		4 x 5µl	2 x 5µl

機能

Receptor for hyaluronic acid (HA). Mediates cell-cell and cell-matrix interactions through its affinity for HA, and possibly also through its affinity for other ligands such as osteopontin, collagens, and matrix metalloproteinases (MMPs). Adhesion with HA plays an important role in cell migration, tumor growth and progression. Also involved in lymphocyte activation, recirculation and homing, and in hematopoiesis. Altered expression or dysfunction causes numerous pathogenic phenotypes. Great protein heterogeneity due to numerous alternative splicing and post-translational modification events.

組織特異性

Isoform 10 (epithelial isoform) is expressed by cells of epithelium and highly expressed by carcinomas. Expression is repressed in neuroblastoma cells.

配列類似性

Contains 1 Link domain.

ドメイン

The lectin-like LINK domain is responsible for hyaluronan binding.

翻訳後修飾

Proteolytically cleaved in the extracellular matrix by specific proteinases (possibly MMPs) in several cell lines and tumors.

N-glycosylated.

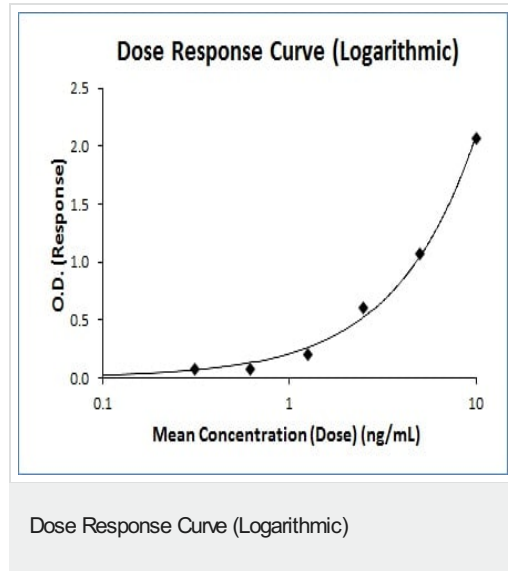
O-glycosylated; contains more-or-less-sulfated chondroitin sulfate glycans, whose number may

affect the accessibility of specific proteinases to their cleavage site(s).
 Phosphorylated; activation of PKC results in the dephosphorylation of Ser-706 (constitutive phosphorylation site), and the phosphorylation of Ser-672.

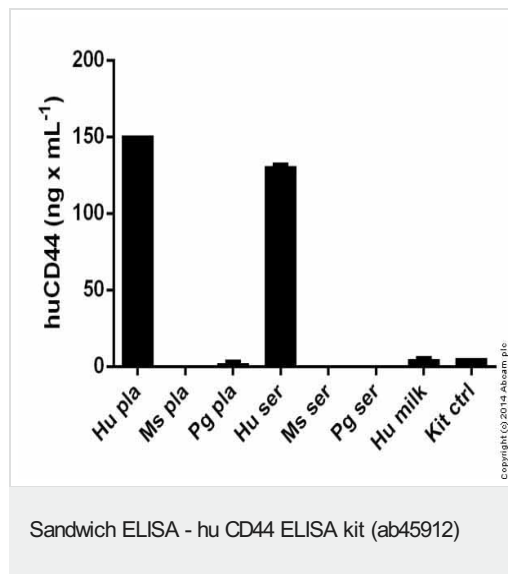
細胞内局在

Membrane.

画像



Representative Dose Response Curve (Logarithmic) using ab45912.



huCD44 measured in biofluids showing quantity (ng) per mL of tested sample

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