

Human Galectin-4 ELISA Kit ab213785

1 References 画像数 1

製品の概要

製品名	Human Galectin-4 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
	1	16	209pg/ml	11.7
	2	16	1291pg/ml	98.11
	3	16	4066pg/ml	223.63
				CV%
				= 7.6%
				= 7.6%
				= 5.5%
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	1	24	215pg/ml	12.04
	2	24	1324pg/ml	108.56
	3	24	4394pg/ml	320.76
				CV%
				= 5.6%
				= 8.2%
				= 7.3%
サンプルの種類	Cell culture supernatant, Serum, Cell Lysate, Hep Plasma, EDTA Plasma			
アッセイタイプ	Sandwich (quantitative)			
検出感度	< 10 pg/ml			
検出範囲	31.2 pg/ml - 2000 pg/ml			
全工程の試験時間	3h 30m			
ステップ	Multiple steps standard assay			
種交差性	交差種: Human			
製品の概要	The Human Galectin-4 Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213785) is designed for the quantitative measurement of Human Galectin-4 in cell culture supernatants, cell lysates, serum and plasma (heparin, EDTA).			

The ELISA kit is based on standard sandwich enzyme-linked immune-sorbent assay technology.

A monoclonal antibody from mouse specific for Galectin-4 has been pre-coated onto 96-well plates. Standards (Expression system for standard: *E.coli*; Immunogen sequence: A2-I323) and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for Galectin-4 is added subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates are washed away with PBS or TBS buffer. HRP substrate TMB is used to visualize HRP enzymatic reaction. TMB is catalyzed by HRP to produce a blue color product that changed into yellow after adding acidic stop solution. The density of yellow is proportional to the Human Galectin-4 amount of sample captured in plate.

特記事項

Galectin-4, also known as LGALS4 or GAL4, is a protein that in humans is encoded by the LGALS4 gene. It is mapped to 19q13.2. Galectins are a family of animal lectins defined by an affinity for beta galactoside-containing saccharides and by shared sequence elements. Galectin-4 is expressed as a 38-kD protein in the human colon adenocarcinoma T84 cell line. Galectin-4 is an S-type lectin that is strongly under expressed in colorectal cancer, and this gene plays a role in cell adhesion. It has been found that decreased Galectin-4 expression occurs early in colorectal carcinogenesis.

試験プラットフォーム

Pre-coated microplate (12 x 8 well strips)

製品の特性

保存方法

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

内容	ラベル	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml
Adhesive Plate Seal		4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml
Anti-Human Galectin-4 coated Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl
Biotinylated anti-Human Galectin-4 antibody		1 x 100µl
Lyophilized recombinant Human Galectin-4 standard		2 vials
Sample Diluent Buffer	Green Cap	1 x 30ml
TMB Color Developing Agent	Black Cap	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml
Wash Buffer (25X)		1 x 20ml

機能

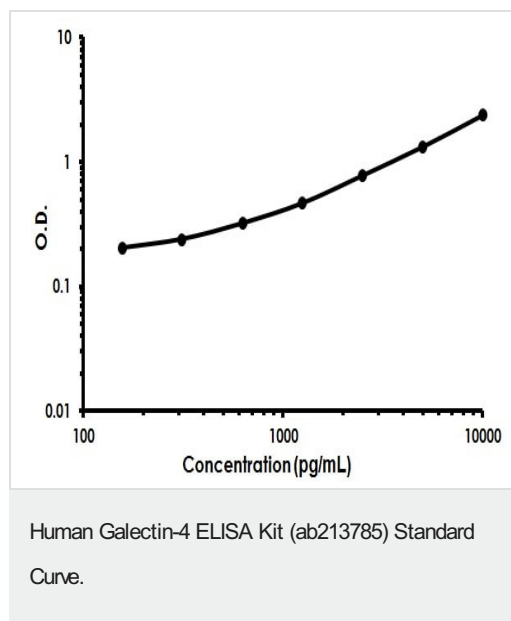
Galectin that binds lactose and a related range of sugars. May be involved in the assembly of adherens junctions.

配列類似性

Contains 2 galectin domains.

ドメイン

Contains two homologous but distinct carbohydrate-binding domains.



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