

Mouse Galectin 1 ELISA Kit ab119595

3 References 画像数 1

製品の概要

製品名	Mouse Galectin 1 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%
	1	16	343pg/ml	22.29	= 6.5%
	2	16	1747pg/ml	132.77	= 7.6%
	3	16	4774pg/ml	300.76	= 8.5%
	Inter-Assay (日差再現性)				
	サンプル	N	平均値	SD	CV%
	1	24	318pg/ml	20.67	= 6.5%
	2	24	1697pg/ml	144.24	= 8.5%
	3	24	4684pg/ml	304.46	= 6.5%
サンプルの種類	Cell culture supernatant, Serum, Plasma, Cell Lysate				
アッセイタイプ	Sandwich (quantitative)				
検出感度	< 5 pg/ml				
検出範囲	156 pg/ml - 10000 pg/ml				
ステップ	Multiple steps standard assay				
種交差性	交差種: Mouse				
製品の概要	Abcam's mouse Galectin 1 <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed				

buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates are washed away with PBS or TBS buffer. TMB is then used to visualize the HRP enzymatic reaction. TMB is catalyzed by HRP 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 mouse Galectin 1 amount of sample captured in plate.

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

#### 機能

May regulate apoptosis, cell proliferation and cell differentiation. Binds beta-galactoside and a wide array of complex carbohydrates. Inhibits CD45 protein phosphatase activity and therefore the dephosphorylation of Lyn kinase.

#### 組織特異性

Expressed in placenta, maternal decidua and fetal membranes. Within placenta, expressed in trophoblasts, stromal cells, villous endothelium, syncytiotrophoblast apical membrane and villous stroma. Within fetal membranes, expressed in amnion, chorioamniotic mesenchyma and chorion (at protein level). Expressed in cardiac, smooth, and skeletal muscle, neurons, thymus, kidney and hematopoietic cells.

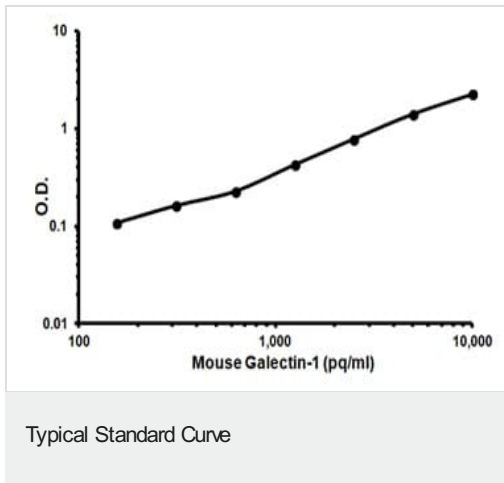
#### 配列類似性

Contains 1 galectin domain.

#### 細胞内局在

Secreted > extracellular space > extracellular matrix.

#### 画像



Representative Standard Curve using ab119595

**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