

Mouse Galectin 1 ELISA Kit ab119595

3 References 1 Image

Overview

Product name	Mouse Galectin 1 ELISA Kit				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	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				
	Sample	n	Mean	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%
Sample type	Cell culture supernatant, Serum, Plasma, Cell Lysate				
Assay type	Sandwich (quantitative)				
Sensitivity	< 5 pg/ml				
Range	156 pg/ml - 10000 pg/ml				
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Mouse				
Product overview	<p>Abcam’s mouse Galectin 1 <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of mouse Galectin 1 in cell culture supernatants, serum and plasma (heparin, EDTA).</p> <p>A Galectin 1 specific rat monoclonal antibody has been precoated onto 96-well plates. Standards and test samples are added to the wells and incubated. A biotinylated detection polyclonal antibody from goat specific for Galectin 1 is then added followed by washing with PBS or TBS</p>				

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.

Platform Microplate

Properties

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

Components	Identifier	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

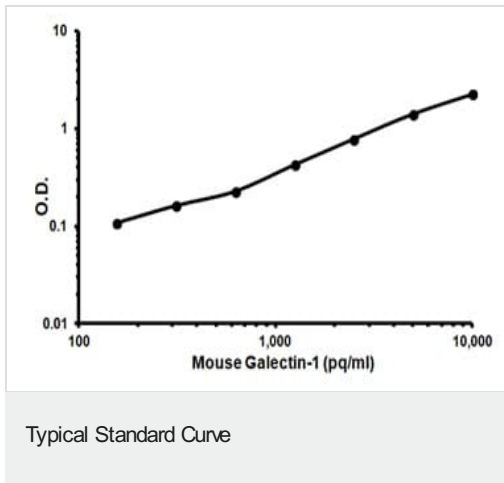
Function May regulate apoptosis, cell proliferation and cell differentiation. Binds beta-galactoside and a wide array of complex carbohydrates. Inhibits CD45 protein phosphatase activity and therefore the dephosphorylation of Lyn kinase.

Tissue specificity 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.

Sequence similarities Contains 1 galectin domain.

Cellular localization Secreted > extracellular space > extracellular matrix.

Images



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.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors