

Mouse Galectin 3 ELISA Kit ab113343

[2 Images](#)

Overview

Product name	Mouse Galectin 3 ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 6 pg/ml
Range	8.19 pg/ml - 2000 pg/ml
Recovery	> 100 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	124.3	116% - 131%
Serum	111.8	101% - 120%
Plasma	99.08	90% - 107%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Mouse

Product overview Abcam's Galectin 3 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Mouse Galectin 3 in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for Mouse Galectin 3 coated on a 96- well plate. Standards and samples are pipetted into the wells and Galectin 3 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Mouse Galectin 3 antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of Galectin 3 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Platform Microplate

Properties

Storage instructions

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

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
400X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Mouse Galectin 3	2 vials
Galectin 3 Microplate (12 x 8 wells)	1 unit
Recombinant Mouse Galectin 3 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

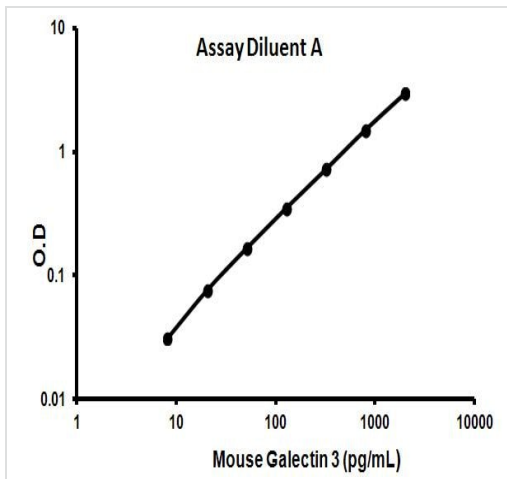
Relevance

Galactose-specific lectin which binds IgE. May mediate with the alpha-3, beta-1 integrin the stimulation by CSPG4 of endothelial cells migration. Together with DMBT1, required for terminal differentiation of columnar epithelial cells during early embryogenesis. By similarity. In the nucleus: acts as a pre-mRNA splicing factor. Involved in acute inflammatory responses including neutrophil activation and adhesion, chemoattraction of monocytes macrophages, opsonization of apoptotic neutrophils, and activation of mast cells. Probably forms homo- or heterodimers. Interacts with DMBT1. Forms a complex with the ITGA3, ITGB1 and CSPG4. Interacts with LGALS3BP, LYPD3, CYHR1 and UACA. A major expression is found in the colonic epithelium. It is also abundant in the activated macrophages. Contains 1 galectin domain.

Cellular localization

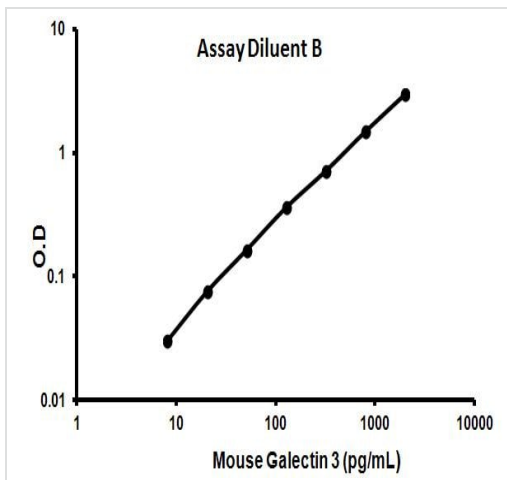
Cytoplasmic, Nuclear and Secreted

Images



Representative Standard Curve using ab113343

Representative Standard Curve



Representative Standard Curve using ab113343

Representative Standard Curve

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