

Product datasheet

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.

Tested applications **Suitable for:** Sandwich ELISA

Platform Microplate

Properties

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

Components	1 x 96 tests
20X Wash Buffer Concentrate	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

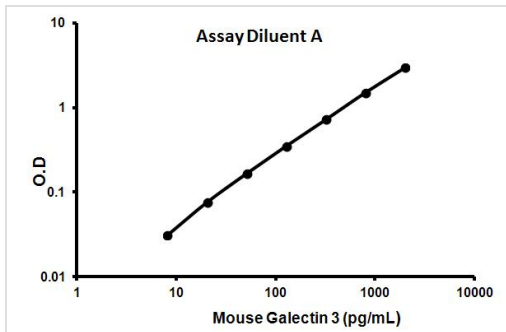
Applications

Our [Abpromise guarantee](#) covers the use of **ab113343** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

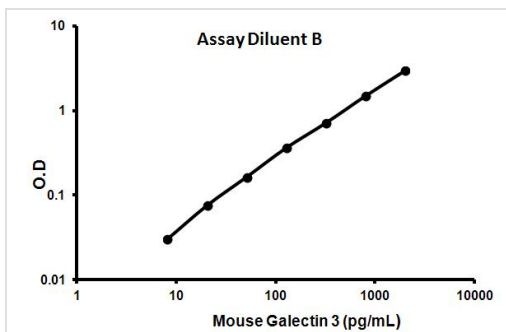
Application	Abreviews	Notes
Sandwich ELISA		Use at an assay dependent concentration.

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 AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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