

Human Galectin-3BP ELISA Kit ab213784

1 References 1 Image

Overview

Product name Human Galectin-3BP ELISA Kit

Detection method Colorimetric

Precision Intra-assay

Sample	n	Mean	SD	CV%
1	16	312pg/ml	18.72	= 6%
2	16	1090pg/ml	63.22	= 5.8%
3	16	4158pg/ml	282.74	= 6.8%

Inter-assay

Sample	n	Mean	SD	CV%
1	24	326pg/ml	20.86	= 6.4%
2	24	1154pg/ml	84.24	= 7.3%
3	24	3768pg/ml	271.29	= 7.2%

Sample type Cell culture supernatant, Serum, Hep Plasma

Assay type Sandwich (quantitative)

Sensitivity < 10 pg/ml

Range 156 pg/ml - 10000 pg/ml

Assay time 3h 30m

Assay duration Multiple steps standard assay

Species reactivity Reacts with: Human

Product overview The Human Galectin-3BP Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213784) is designed for the quantitative measurement of Human Galectin-3BP in cell culture supernatants, serum and plasma (heparin).

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

A monoclonal antibody from mouse specific for Galectin-3BP has been pre-coated onto 96-well plates. Standards (NSO; V19-D585) and test samples are added to the wells, a biotinylated detection monoclonal antibody from mouse specific for Galectin-3BP 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-3BP amount of sample captured in plate.

Notes Galectin-3-binding protein is a protein that in humans is encoded by the LGALS3BP gene. The galectins are a family of beta-galactoside-binding proteins implicated in modulating cell–cell and cell–matrix interactions. Using fluorescence *in situ* hybridization, the full length 90K cDNA has been localized to chromosome 17q25. The native protein binds specifically to a human macrophage associated lectin known as Mac-2 and also binds to galectin 1.

Platform Pre-coated microplate (12 x 8 well strips)

Properties

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

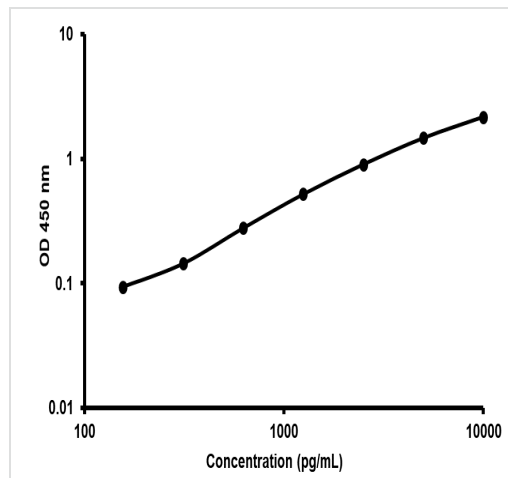
Components	Identifier	1 x 96 tests	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml	1 x 12ml
Adhesive Plate Seal		4 units	4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml	1 x 12ml
Anti-Human Galectin-3BP coated Microplate (12 x 8 wells)		1 unit	1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl	1 x 100µl
Biotinylated anti- Human Galectin-3BP antibody		1 x 100µl	1 x 100µl
Lyophilized recombinant Human Galectin-3BP standard		2 vials	2 vials
Sample Diluent Buffer	Green Cap	1 x 30ml	1 x 30ml
TMB Color Developing Agent	Black Cap	1 x 10ml	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml	1 x 10ml
Wash Buffer (25X)		1 x 20ml	1 x 20ml

Function Promotes intergrin-mediated cell adhesion. May stimulate host defense against viruses and tumor cells.

Tissue specificity Ubiquitous. Detected in body fluids such as semen, milk, serum, tears, saliva and urine. Expressed by keratinocytes and fibroblasts.

Sequence similarities Contains 1 BACK (BTB/Kelch associated) domain.
Contains 1 BTB (POZ) domain.
Contains 1 SRCR domain.

Images



Human Galectin-3BP ELISA Kit (ab213784) Standard Curve.

Human Galectin-3BP ELISA Kit (ab213784)

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