

Human Galectin-4 ELISA Kit ab213785

[1 References](#) [1 Image](#)

Overview

Product name Human Galectin-4 ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

| Sample | n | Mean | SD | CV% |
|--------|----|-----------|--------|--------|
| 1 | 16 | 209pg/ml | 11.7 | = 7.6% |
| 2 | 16 | 1291pg/ml | 98.11 | = 7.6% |
| 3 | 16 | 4066pg/ml | 223.63 | = 5.5% |

Inter-assay

| Sample | n | Mean | SD | CV% |
|--------|----|-----------|--------|--------|
| 1 | 24 | 215pg/ml | 12.04 | = 5.6% |
| 2 | 24 | 1324pg/ml | 108.56 | = 8.2% |
| 3 | 24 | 4394pg/ml | 320.76 | = 7.3% |

Sample type

Cell culture supernatant, Serum, Cell Lysate, Hep Plasma, EDTA Plasma

Assay type

Sandwich (quantitative)

Sensitivity

< 10 pg/ml

Range

31.2 pg/ml - 2000 pg/ml

Assay time

3h 30m

Assay duration

Multiple steps standard assay

Species reactivity

Reacts with: Human

Product overview

The Human Galectin-4 Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213785) is designed for the quantitative measurement of Human Galectin-4 in cell culture supernatants, cell lysates, serum and plasma (heparin, EDTA).

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

A monoclonal antibody from mouse specific for Galectin-4 has been pre-coated onto 96-well plates. Standards (Expression system for standard: *E.coli*; Immunogen sequence: A2-I323) and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for Galectin-4 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-4 amount of sample captured in plate.

Notes Galectin-4, also known as LGALS4 or GAL4, is a protein that in humans is encoded by the LGALS4 gene. It is mapped to 19q13.2. Galectins are a family of animal lectins defined by an affinity for beta galactoside-containing saccharides and by shared sequence elements. Galectin-4 is expressed as a 38-kD protein in the human colon adenocarcinoma T84 cell line. Galectin-4 is an S-type lectin that is strongly under expressed in colorectal cancer, and this gene plays a role in cell adhesion. It has been found that decreased Galectin-4 expression occurs early in colorectal carcinogenesis.

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

Properties

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

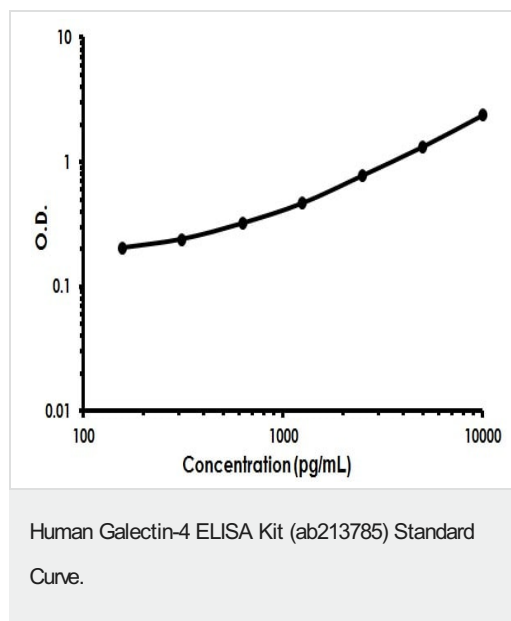
| Components | Identifier | 1 x 96 tests |
|--|------------|--------------|
| ABC Diluent Buffer | Blue Cap | 1 x 12ml |
| Adhesive Plate Seal | | 4 units |
| Antibody Diluent Buffer | Green Cap | 1 x 12ml |
| Anti-Human Galectin-4 coated Microplate (12 x 8 wells) | | 1 unit |
| Avidin-Biotin-Peroxidase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

Function Galectin that binds lactose and a related range of sugars. May be involved in the assembly of adherens junctions.

Sequence similarities Contains 2 galectin domains.

Domain Contains two homologous but distinct carbohydrate-binding domains.

Images



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