

Human Galectin-9 / Gal-9 ELISA Kit ab213786

1 References 1 Image

Overview

Product name Human Galectin-9 / Gal-9 ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
1	16	168pg/ml	7.9	= 4.7%
2	16	731pg/ml	51.9	= 7.1%
3	16	3580pg/ml	218.38	= 6.1%

Inter-assay

Sample	n	Mean	SD	CV%
1	24	177pg/ml	9.55	= 5.4%
2	24	792pg/ml	66.52	= 8.4%
3	24	3920pg/ml	274.4	= 7%

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

Assay type Sandwich (quantitative)

Sensitivity < 10 pg/ml

Range 93.8 pg/ml - 6000 pg/ml

Assay time 3h 30m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview The Human Galectin-9 / Gal-9 Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213786) is designed for the quantitative detection of Human Galectin-9 in cell culture supernatants, cell lysates, serum and plasma (heparin, EDTA).

The ELISA kit is based on standard sandwich enzyme-linked immunosorbent assay technology. A

polyclonal antibody from goat specific for Galectin-9 has been precoated onto 96-well plates. Standards (Expression system for standard: E.coli; Immunogen sequence: M1-T355) and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for Galectin-9 is added subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex was added and unbound conjugates were washed away with PBS or TBS buffer. HRP substrate TMB was used to visualize HRP enzymatic reaction. TMB was 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-9 amount of sample captured in plate.

Notes Galectin-9, also called HUAT or LGALS9A is a protein that in humans is encoded by the LGALS9 gene. This gene is mapped to 17q11.2. The galectins are a family of beta-galactoside-binding proteins implicated in modulating cell-cell and cell-matrix interactions. The protein encoded by this gene is an S-type lectin. It is overexpressed in Hodgkin's disease tissue and might participate in the interaction between the H and RS cells with their surrounding cells and might thus, play a role in the pathogenesis of this disease and/or its associated immunodeficiency. The protein has N- and C- terminal carbohydrate-binding domains connected by a link peptide.

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

Properties

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

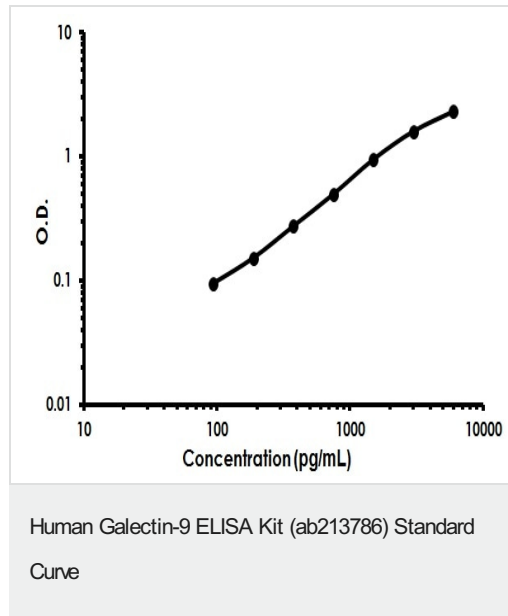
Components	Identifier	1 x 96 tests	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml	1 x 12ml
Adhesive Plate Seal		4 units	4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml	1 x 12ml
Anti-Human Galectin-9 coated Microplate (12 x 8 wells)		1 unit	1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl	1 x 100µl
Biotinylated anti-Human Galectin-9 antibody		1 x 100µl	1 x 100µl
Lyophilized recombinant Human Galectin-9 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 Binds galactosides. Has high affinity for the Forssman pentasaccharide. May play a role in thymocyte-epithelial interactions relevant to the biology of the thymus. Inhibits cell proliferation. It is a ligand for HAVCR2/TIM3. Induces T-helper type 1 lymphocyte (Th1) death. Isoform Short acts as an eosinophil chemoattractant.

Tissue specificity Peripheral blood leukocytes and lymphatic tissues. Overexpressed in Hodgkin disease tissue.

Sequence similarities	Contains 2 galectin domains.
Domain	Contains two homologous but distinct carbohydrate-binding domains.
Cellular localization	Cytoplasm. Secreted. May also be secreted by a non-classical secretory pathway.

Images



Human Galectin-9 / Gal-9 ELISA Kit (ab213786) 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