

Human Galectin 7 ELISA Kit (GAL7) ab100523

2 Images

Overview

Product name	Human Galectin 7 ELISA Kit (GAL7)
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 20 pg/ml
Range	13.72 pg/ml - 10000 pg/ml
Recovery	> 90 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	96.55	87% - 105%
Serum	107.1	95% - 115%
Plasma	91.82	70% - 125%

Assay duration Multiple steps standard assay

Species reactivity Reacts with: Human

Product overview Abcam's Galectin 7 (GAL7) Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human Galectin 7 in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for Human Galectin 7 coated on a 96-well plate. Standards and samples are pipetted into the wells and Galectin 7 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human Galectin 7 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 7 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Notes Optimization may be required with urine samples

Platform Microplate

Properties

Storage instructions

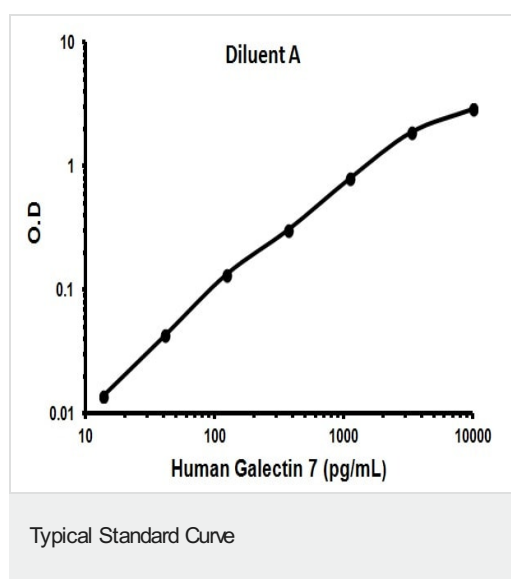
Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Human Galectin 7	2 vials
Galectin 7 Microplate (12 x 8 well strips)	1 unit
Galectin 7 Standard Human (recombinant) (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

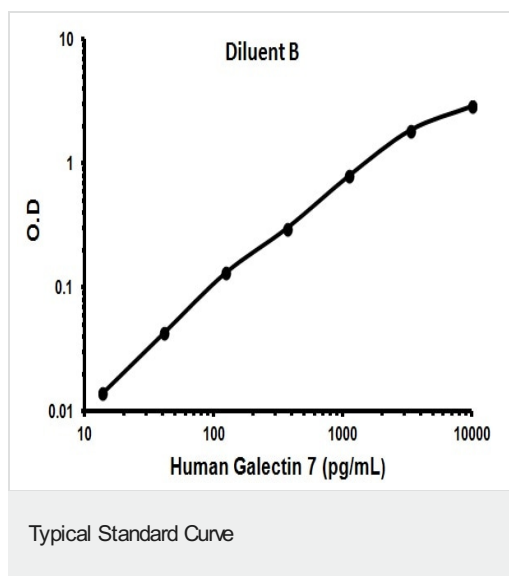
Relevance

Galectins belong to a family of related beta-galactoside-binding lectins, also referred to as S-type or S-Lac lectins. Members of this family have been implicated in a variety of functions, including growth regulation, cell adhesion, migration, neoplastic transformation, and immune responses. Galectin 7 is thought to be involved in cell-cell and/or cell-matrix interactions necessary for normal growth control. Galectin 7 has been implicated in apoptosis and is believed to upregulate MMP9 and as such be involved in the progression of tumorigenesis.

Images



Representative Standard Curve using ab100523.



Representative Standard Curve using ab100523.

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