

Mouse Galectin-3 ELISA Kit (Colorimetric)

Catalog Number EA-9009

(For Research Use Only)

Introduction

Galectin-3 is a member of the lectin family and contains a carbohydrate-recognition-binding domain that binds to βgalactoside. Due to its ability to bind to β-galactoside, galectin-3 plays an important role in cell-cell adhesion, cell-matrix interactions, macrophage activation, angiogenesis, metastasis, apoptosis. Given galectin-3's broad biological functionality, it has been demonstrated to be involved in a number of diseases including cancer, fibrosis, heart disease, and stroke. Overexpression of galectin-3 is commonly seen in cancerous conditions and can affect cancer cell growth and differentiation, chemoattraction, apoptosis, and metastasis. Because of this, galectin-3 is increasingly being used as a diagnostic marker for different cancers.

Principle of the assay

The wells of the plate are coated with capture antibody specific to galectin-3. In this assay, the test sample initially reacts with the solid phase capture antibody, resulting in galectin-3 being bound to the well. The wells are then washed to remove unbound proteins, and biotin-linked antibodies are added to bind to the immobilized galectin-3. After washing away the unbound antibodies, Streptavidin-HRP conjugate is added to form a complex with the antibody-bound galectin-3. After incubation, the wells are washed to remove unbound Streptavidin-HRP conjugate. TMB substrate is then added and forms a blue color when the HRP-linked antibodies are detected. The reaction is then terminated with Stop Solution, which changes the color from blue to yellow. The galectin-3 concentration in each well is directly proportional to its color intensity and can be quantified by measuring its optical density at 450 nm (OD450) in a microplate reader.

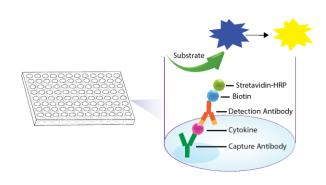


Diagram of Mouse Galectin-3 ELISA

Materials provided with the kit

Component	Qty	Store at
One clear plate coated with antibody against galectin-3	1	4°C
Biotin-labeled anti-galectin-3 antibody	200 µL	-20°C
Streptavidin-HRP conjugate	50 µL	4°C
1x Diluent buffer	40 mL	4°C
5x Assay wash buffer	40 mL	4°C
Substrate	10 mL	4°C
Stop solution	5 mL	4°C

Material required but not provided

- Microplate reader
- Distilled H2O

Reagent preparation before starting experiment

- Dilute the 5x Assay wash buffer to 1x buffer
 40 ml 5x Assay wash buffer
 160 ml ddH2O
- Dilute 50 times of biotin labeled antibody with 1X Diluent buffer.

(AVOID FREEZE/THAW OF ANTIBODY MIX)

• Dilute 200 times of streptavidin-HRP with 1x Diluent buffer.

Sample preparation before starting experiment

- For **cell culture medium samples**, add 100 ul directly to the well or dilute 2-fold with 1X Diluent buffer.
- For **cell lysate samples**, use cell lysis buffer (Catalog# EA-0001). Follow protocol on Cell Lysate Buffer User Manual on our website.
- For serum or plasma samples, we recommend a 1:10 to 1:20 dilution with 1X diluent buffer. When serum-containing conditional media is required, be sure to use serum as control.

Assay procedure

- 1. Take the plate from the aluminized bag. Seal the unused wells with a film.
- Prepare 2.5 ml sample and add 100 μl of sample per well to one section and incubate for 2 hours at room temperature with gentle shaking.
 Optional: If you want to have a blank reading, you

can design one well as a blank well by adding diluent buffer instead of your sample.

- 3. Invert the plate over an appropriate container and expel the contents forcibly. Wash the plate by adding 200 μ l of 1x Assay wash buffer. Repeat the washing process two times for a total of three washes. Complete removal of liquid at each wash by firmly tapping the plate against a pile of clean paper towels.
- 4. Add $100 \ \mu l$ of diluted biotin-labeled antibody to each well and incubate for 1 hour at room temperature with gentle shaking.
- 5. Repeat the aspiration/wash as in step 3.
- Add 100 µl of diluted streptavidin-HRP conjugate to each well and incubate for 45 min at room temperature with gentle shaking.
- 7. Repeat the aspiration/wash as in step 3.
- 8. Add 100 μ l substrate to each well and incubate for 30-40 minutes at least.

Note: Substrate incubation time may vary due to different antibodies reactivity. Stronger signals (Strong blue color) could be stopped early after 5 minutes. Weaker signals should be incubated for 10-30 minutes.

- Add 50 µl of Stop solution to each well. The color in the wells should change from blue to yellow.
- 10. Determine the optical density of each well with a microplate reader at 450 nm within 30 minutes.