

Product datasheet

Monkey Galectin-3 ELISA Kit (LGALS3) ab269556

Recombinant SimpleStep ELISA

[5 Images](#)

Overview

Product name Monkey Galectin-3 ELISA Kit (LGALS3)

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Serum	8			3.2%

Inter-assay

Sample	n	Mean	SD	CV%
Serum	3			2.6%

Sample type

Cell culture supernatant, Serum, Cell culture media, Hep Plasma, EDTA Plasma, Cit plasma

Assay type

Sandwich (quantitative)

Sensitivity

115 pg/ml

Range

208.9 pg/ml - 6000 pg/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	87	83% - 90%
Serum	105	100% - 110%
Cell culture media	83	81% - 85%
Hep Plasma	109	108% - 110%
EDTA Plasma	106	102% - 110%
Cit plasma	117	116% - 118%

Assay time	1h 30m
Assay duration	One step assay
Species reactivity	Reacts with: Monkey
Product overview	Monkey Galectin-3 ELISA Kit (LGALS3) (ab269556) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of Galectin-3 (LGALS3) protein in edta plasma, hep plasma, serum, cell culture media, cit plasma, and cell culture supernatant. It uses our proprietary SimpleStep ELISA® technology. Quantitate Monkey Galectin-3 (LGALS3) with 115 pg/ml sensitivity.

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate ([ab203359](#)) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

Notes	Galectin-3 (Gal-3, Galactoside-binding protein, Mac-2 antigen) is a member of the galectin family of carbohydrate-binding proteins. Galectin-3 is active intracellularly and extracellularly. It is widely expressed, including by fibroblasts and macrophages. Intracellularly, Galectin-3 can function in the nucleus as a pre-mRNA splicing factor and in the cytoplasm it has roles in endocytosis, exocytosis and apoptosis. Extracellularly, Galectin-3 is often pro-inflammatory and is involved in innate immunity. Galectin-3 can be monomeric or form higher order homo and hetero oligomers.
--------------	---

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

Properties

Storage instructions	Store at +4°C. Please refer to protocols.
-----------------------------	---

Components	1 x 96 tests
10X Monkey Galectin-3 (LGALS3) Capture Antibody	1 x 600µl
10X Monkey Galectin-3 (LGALS3) Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent 4BI	1 x 6ml
Monkey Galectin-3 (LGALS3) Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit

Components	1 x 96 tests
Sample Diluent NS (ab193972)	2 x 50ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

Function

Galactose-specific lectin which binds IgE. May mediate with the alpha-3, beta-1 integrin the stimulation by CSPG4 of endothelial cells migration. Together with DMBT1, required for terminal differentiation of columnar epithelial cells during early embryogenesis.

Tissue specificity

A major expression is found in the colonic epithelium. It is also abundant in the activated macrophages.

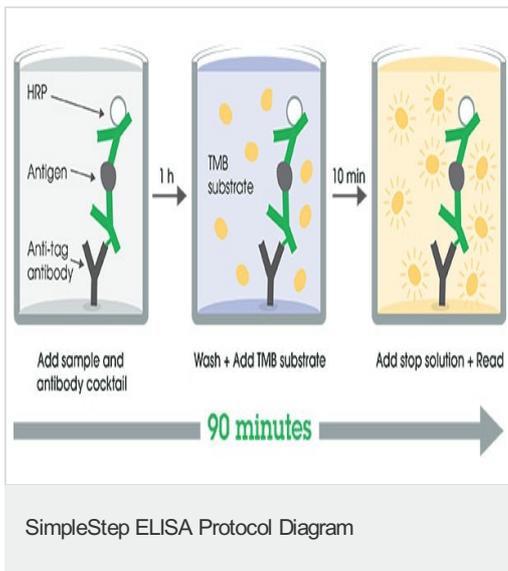
Sequence similarities

Contains 1 galectin domain.

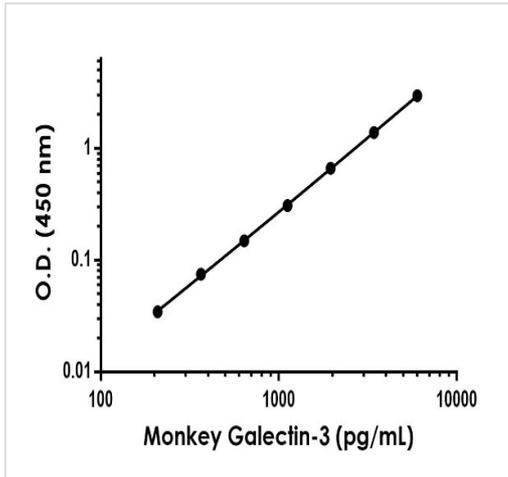
Cellular localization

Nucleus. Cytoplasmic in adenomas and carcinomas. May be secreted by a non-classical secretory pathway and associate with the cell surface.

Images

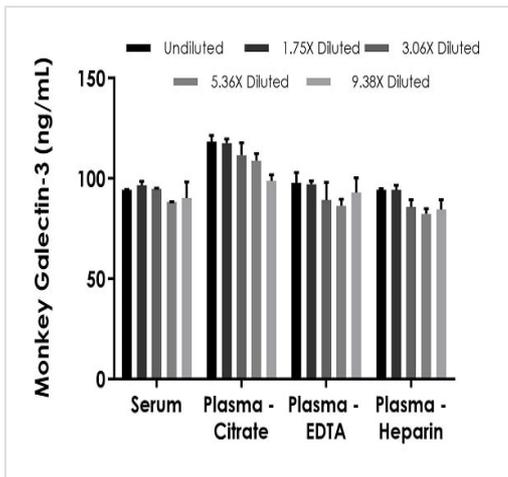


SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



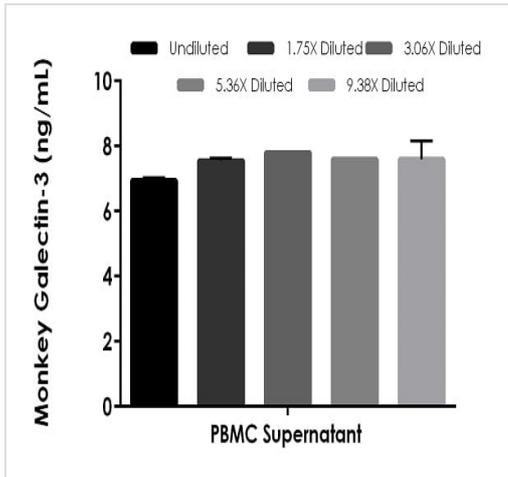
The Galectin-3 standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.

Example of monkey Galectin-3 standard curve in Sample Diluent NS.



The concentrations of Galectin-3 were measured in duplicates, interpolated from the Galectin-3 standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 4%, plasma (citrate) 4%, plasma (EDTA) 4% and plasma (heparin) 4%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean Galectin-3 concentration was determined to be 93 ng/mL in neat serum, 111 ng/mL in neat plasma (citrate), 93 ng/mL in neat plasma (EDTA) and 88 ng/mL in neat plasma (heparin).

Interpolated concentrations of native Galectin-3 in Rhesus Macaque monkey serum and plasma samples.



Interpolated concentrations of native Galectin-3 in Rhesus Macaque monkey PBMC cell culture supernatant cultured for 24 hours.

The concentrations of Galectin-3 were measured in duplicates, interpolated from the Galectin-3 standard curves and corrected for sample dilution. Undiluted samples are as follows: PBMC supernatant 57%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean Galectin-3 concentration was determined to be 7.5 ng/mL in PBMC supernatant.

Powered by
recombinant antibodies

Research with confidence
Consistent and reproducible results

Long-term and scalable supply
Recombinant technology

Success from the first experiment
Confirmed specificity

Ethical standards compliant
Animal-free production

Sandwich ELISA - Monkey Galectin-3 ELISA Kit
(LGALS3) (ab269556)

To learn more about the advantages of recombinant antibodies see [here](#).

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