



Laminin (Mouse) ELISA Kit

Catalog Number KA0380

96 assays

Version: 05

Intended for research use only

www.abnova.com

Table of Contents

Introduction	3
Intended Use	3
Background	3
Principle of the Assay	3
General Information	4
Materials Supplied	4
Storage Instruction	4
Materials Required but Not Supplied	4
Precautions for Use	5
Assay Protocol	6
Reagent Preparation	6
Sample Preparation	6
Assay Procedure	7
Data Analysis	9
Calculation of Results	9
Performance Characteristics	9
Resources	10
References	10
Plate Layout	11

Introduction

Intended Use

For quantitative detection of mouse Laminin in cell culture supernates, serum and plasma (heparin, EDTA).

Background

Laminin is a large basement membrane glycoprotein composed of three subunits designated the A, B1, and B2.¹ Laminin has diverse biological functions, which include stimulating epithelial cell growth and differentiation.² The nucleotide sequence of human laminin A chain has an open reading frame encoding 3075-amino acids.¹ The human laminin A chain is at locus 18p11.3.³ The nucleotide sequence of the human laminin B1 reveals a 5358-base pair open reading frame that potentially codes for 1786 amino acids, including 20 amino acids of a presumptive signal peptide.² The gene for the human laminin-B1 chain has been localized to chromosome 7, band q31.⁴ The B2 chain consists of six distinct domains, including two domains with alpha-helical, coiled-coil structures, two domains with cysteine-rich homologous repeats, and two globular domains. The amino acid sequences of the B2 and B1 chains demonstrate considerable homology.⁵ The human laminin B2 chain gene maps to the long arm of chromosome 1 in the band q31.6 The standard product used in this kit is isolated from plasma.

Principle of the Assay

The Laminin (Mouse) ELISA Kit was based on standard sandwich enzyme-linked immune-sorbent assay technology. A monoclonal antibody from rat specific for Laminin has been precoated onto 96-well plates. Standards and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for Laminin 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 mouse Laminin amount of sample captured in plate.

General Information

Materials Supplied

List of component

Component	Amount
96-well plate precoated with anti- mouse Laminin antibody	96 (8x12) wells
Lyophilized mouse Laminin standard	10ng/tube x 2
Biotinylated anti-mouse Laminin antibody, dilution 1:100	130 μ l
Avidin-Biotin-Peroxidase Complex (ABC), dilution 1:100	130 μ l
Sample diluent buffer	30 ml
Antibody diluent buffer	12 ml
ABC diluent buffer	12 ml
TMB color developing agent	10 ml
TMB stop solution	10 ml

Storage Instruction

Store at 4°C for 6 months, at -20°C for 12 months. Avoid multiple freeze-thaw cycles.

Materials Required but Not Supplied

- ✓ Microplate reader in standard size.
- ✓ Automated plate washer.
- ✓ Adjustable pipettes and pipette tips. Multichannel pipettes are recommended in the condition of large amount of samples in the detection.
- ✓ Clean tubes and Eppendorf tubes.
- ✓ Washing buffer (neutral PBS or TBS).

Preparation of 0.01M TBS:

Add 1.2g Tris, 8.5g NaCl; 450 μ l of purified acetic acid or 700 μ l of concentrated hydrochloric acid to 1000ml H₂O and adjust pH to 7.2-7.6. Finally, adjust the total volume to 1L.

Preparation of 0.01 M PBS:

Add 8.5g sodium chloride, 1.4g Na₂HPO₄ and 0.2g NaH₂PO₄ to 1000ml distilled water and adjust pH to 7.2-7.6. Finally, adjust the total volume to 1L.

Precautions for Use

- ✓ To inspect the validity of experiment operation and the appropriateness of sample dilution proportion, pilot experiment using standards and a small number of samples is recommended.
- ✓ The TMB Color Developing agent is colorless and transparent before using, contact us freely if it is not the case.
- ✓ Before using Kit, spin tubes and bring down all components to bottom of tubes.
- ✓ Duplicate well assay is recommended for both standard and sample testing.
- ✓ Don't let 96-well plate dry, for dry plate will inactivate active components on plate.
- ✓ Don't reuse tips and tubes to avoid cross contamination.
- ✓ To avoid to use the reagents from different batches together.
- ✓ In order to avoid marginal effect of plate incubation due to temperature difference (reaction may be stronger in the marginal wells), it is suggested that the diluted ABC and TMB solution will be pre-warmed in 37°C for 30 min before using

Assay Protocol

Reagent Preparation

- ✓ Reconstitution of the mouse Laminin standard: Laminin standard solution should be prepared no more than 2 hours prior to the experiment. Two tubes of Laminin standard (10ng per tube) are included in each kit. Use one tube for each experiment.
 - 10,000pg/ml of mouse Laminin standard solution: Add 1 ml sample diluent buffer into one tube, keep the tube at room temperature for 10 min and mix thoroughly.
 - 5000pg/ml→156pg/ml of mouse Laminin standard solutions: Label 6 Eppendorf tubes with 5000pg/ml, 2500pg/ml, 1250pg/ml, 625pg/ml, 312pg/ml, 156pg/ml, respectively. Aliquot 0.3 ml of the sample diluent buffer into each tube. Add 0.3 ml of the above 10,000pg/ml Laminin standard solution into 1st tube and mix. Transfer 0.3 ml from 1st tube to 2nd tube and mix. Transfer 0.3 ml from 2nd tube to 3rd tube and mix, and so on.
- Note: The standard solutions are best used within 2 hours. The 10ng/ml standard solution should be stored at 4 °C for up to 12 hours, or at -20 °C for up to 48 hours. Avoid repeated freeze-thaw cycles.*
- ✓ Preparation of biotinylated anti-mouse Laminin antibody working solution: The solution should be prepared no more than 2 hours prior to the experiment.
 - The total volume should be: 0.1ml/well x (the number of wells). (Allowing 0.1-0.2 ml more than total volume).
 - Biotinylated anti-mouse Laminin antibody should be diluted in 1:100 with the antibody diluent buffer and mixed thoroughly. (i.e. Add 1µl Biotinylated anti-mouse Laminin antibody to 99µl antibody diluent buffer.)
- ✓ Preparation of Avidin-Biotin-Peroxidase Complex (ABC) working solution: The solution should be prepared no more than 1 hour prior to the experiment.
 - The total volume should be: 0.1ml/well x (the number of wells). (Allowing 0.1-0.2 ml more than total volume).
 - Avidin- Biotin-Peroxidase Complex (ABC) should be diluted in 1:100 with the ABC dilution buffer and mixed thoroughly. (i.e. Add 1µl ABC to 99µl ABC diluent buffer.)

Sample Preparation

- ✓ Sample Preparation and Storage

Store samples to be assayed within 24 hours at 2-8 °C. For long-term storage, aliquot and freeze samples at -20 °C. Avoid repeated freeze-thaw cycles.

- ✓ Cell culture supernate: Remove particulates by centrifugation, analyze immediately or aliquot and store at -20 °C.
- ✓ Serum: Allow the serum to clot in a serum separator tube (about 4 hours) at room temperature. Centrifuge at approximately 2000 X g for 15 min. Analyze the serum immediately or aliquot and store

frozen at -20 °C.

- ✓ Plasma: Collect plasma using EDTA or Heparin as an anticoagulant. Centrifuge for 15 min at 2000 x g within 30 min of collection. Analyze immediately or aliquot and store frozen at -20 °C. Citrate is not recommended as the anticoagulant.

- ✓ Sample Dilution Guideline
The user needs to estimate the concentration of the target protein in the sample and select a proper dilution factor so that the diluted target protein concentration falls near the middle of the linear regime in the standard curve. Dilute the sample using the provided diluent buffer. The following is a guideline for sample dilution. Several trials may be necessary in practice. The sample must be well mixed with the diluents buffer.
- ✓ High target protein concentration (100-1000ng/ml). The working dilution is 1:100. i.e. Add 1 µl sample into 99 µl sample diluent buffer.
- ✓ Medium target protein concentration (10-100ng/ml). The working dilution is 1:10. i.e. Add 10 µl sample into 90 µl sample diluent buffer.
- ✓ Low target protein concentration (156-10,000pg/ml). The working dilution is 1:2. i.e. Add 50 µl sample to 50 µl sample diluent buffer.
- ✓ Very Low target protein concentration (≤ 156 pg/ml). No dilution necessary, or the working dilution is 1:2.

Assay Procedure

The ABC working solution and TMB color developing agent must be kept warm at 37 °C for 30 min before use. When diluting samples and reagents, they must be mixed completely and evenly. Standard Laminin detection curve should be prepared for each experiment. The user will decide sample dilution fold by crude estimation of Laminin amount in samples.

1. Aliquot 0.1ml per well of the 10,000pg/ml, 5000pg/ml, 2500pg/ml, 1250pg/ml, 625pg/ml, 312pg/ml, 156pg/ml mouse Laminin standard solutions into the precoated 96-well plate. Add 0.1ml of the sample diluent buffer into the control well (Zero well). Add 0.1ml of each properly diluted sample of mouse cell culture supernates, serum and plasma (heparin, EDTA) to each empty well. See "Sample Dilution Guideline" above for details. We recommended that each mouse Laminin standard solution and each sample is measured in duplicate.
2. Seal the plate with the cover and incubate at 37 °C for 90 min.
3. Remove the cover, discard plate content, and blot the plate onto paper towels or other absorbent material. Do NOT let the wells completely dry at any time.
4. Add 0.1ml of biotinylated anti-mouse Laminin antibody working solution into each well and incubate the plate at 37 °C for 60 min.
5. Wash the plate three times with 0.01M TBS or 0.01M PBS, and each time let washing buffer stay in the wells for 1 min. Discard the washing buffer and blot the plate onto paper towels or other absorbent material. (Plate Washing Method: Discard the solution in the plate without touching the side walls. Blot

the plate onto paper towels or other absorbent material. Soak each well with at least 0.3 ml PBS or TBS buffer for 1~2 minutes. Repeat this process two additional times for a total of THREE washes. *Note: For automated washing, aspirate all wells and wash THREE times with PBS or TBS buffer, overfilling wells with PBS or TBS buffer. Blot the plate onto paper towels or other absorbent material.*)

6. Add 0.1ml of prepared ABC working solution into each well and incubate the plate at 37°C for 30 min.
7. Wash plate 5 times with 0.01M TBS or 0.01M PBS, and each time let washing buffer stay in the wells for 1-2 min. Discard the washing buffer and blot the plate onto paper towels or other absorbent material.
8. Add 90 µl of prepared TMB color developing agent into each well and incubate plate at 37°C in dark for 15-20 min (For reference only, the optimal incubation time should be determined by end user. And the shades of blue can be seen in the wells with the four most concentrated mouse Laminin standard solutions; the other wells show no obvious color).
9. Add 0.1ml of prepared TMB stop solution into each well. The color changes into yellow immediately.
10. Read the O.D. absorbance at 450nm in a microplate reader within 30 min after adding the stop solution.

Summary

1. Add samples and standards and incubate the plate at 37°C for 90 min. Do not wash.
2. Add biotinylated antibodies and incubate the plate at 37°C for 60 min. Wash plate 3 times with 0.01M TBS.
3. Add ABC working solution and incubate the plate at 37°C for 30 min. Wash plate 5 times with 0.01M TBS.
4. Add TMB color developing agent and incubate the plate at 37°C in dark for 15-20 min.
5. Add TMB stop solution and read.

Data Analysis

Calculation of Results

For calculation, (the relative O.D.450) = (the O.D.450 of each well) – (the O.D.450 of Zero well). The standard curve can be plotted as the relative O.D.450 of each standard solution (Y) vs. the respective concentration of the standard solution (X). The mouse Laminin concentration of the samples can be interpolated from the standard curve.

Note: if the samples measured were diluted, multiply the dilution factor to the concentrations from interpolation to obtain the concentration before dilution.

✓ Typical result

Typical Data Obtained from Mouse Laminin

Concentration	0.0pg/ml	156pg/ml	312pg/ml	625pg/ml	1250pg/ml	2500pg/ml	5000pg/ml	10,000pg/ml
O.D.	0.082	0.115	0.156	0.223	0.412	0.739	1.210	2.152

(TMB reaction incubate at 37°C for 17 min)

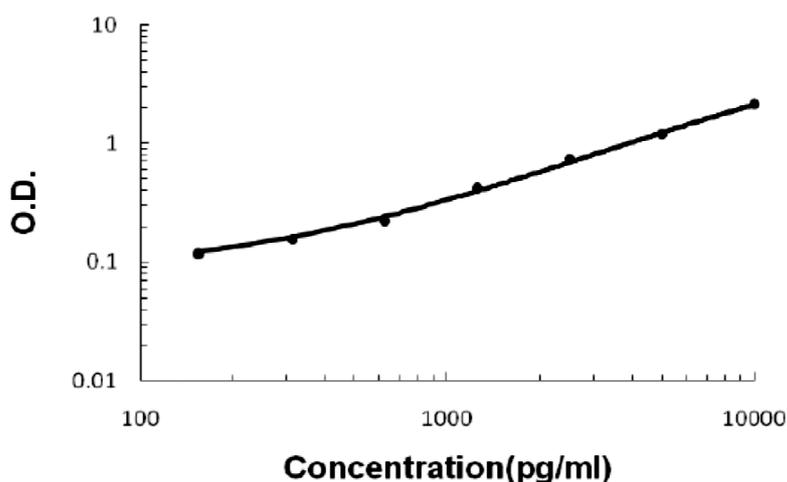


Figure 1: Typical Standard Curve for Laminin (Mouse) ELISA Kit.

Performance Characteristics

- ✓ Range
156pg/ml-10,000pg/ml
- ✓ Sensitivity
< 10pg/ml
- ✓ Specificity
Nature and recombinant mouse Laminin
- ✓ Cross-reactivity
No detectable cross-reactivity with other relevant proteins.

Resources

References

1. Haaparanta, T.; Uitto, J.; Ruoslahti, E.; Engvall, E. Molecular cloning of the cDNA encoding human laminin A chain. *Matrix* 11: 151-160, 1991.
2. Sasaki, M.; Kato, S.; Kohno, K.; Martin, G. R.; Yamada, Y. Sequence of the cDNA encoding the laminin B1 chain reveals a multidomain protein containing cysteine-rich repeats. *Proc. Nat. Acad. Sci.* 84: 935-939, 1987.
3. Nagayoshi, T.; Mattei, M.-G.; Passage, E.; Knowlton, R.; Chu, M.-L.; Uitto, J. Human laminin A chain (LAMA) gene: chromosomal mapping to locus 18p11.3. *Genomics* 5: 932-935, 1989.
4. Jaye, M.; Modi, W. S.; Ricca, G. A.; Mudd, R.; Chiu, I.-M.; O'Brien, S. J.; Drohan, W. N. Isolation of a cDNA clone for the human laminin-B1 chain and its gene localization. *Am. J. Hum. Genet.* 41: 605-615, 1987.
5. Sasaki, M.; Yamada, Y. The laminin B2 chain has a multidomain structure homologous to the B1 chain. *J. Biol. Chem.* 262: 17111-17117, 1987.
6. Mattei, M.-G.; Weil, D.; Pribula-Conway, D.; Bernard, M. P.; Passage, E.; Van Cong, N.; Timpl, R.; Chu, M.-L. cDNA cloning, expression and mapping of human laminin B2 gene to chromosome 1q31. *Hum. Genet.* 79: 235-241, 1988.

Plate Layout

12								
11								
10								
9								
8								
7								
6								
5								
4								
3								
2								
1								
	A	B	C	D	E	F	G	H