

Version 3a Last updated 18 January 2024

# ab213760 – Human CD5L / CT-2 ELISA Kit

For the quantitative detection of Human CD5L / CT-2 in cell culture supernatants, cell lysates, serum, and plasma (heparin, EDTA).

This product is for research use only and is not intended for diagnostic use.

# Table of Contents

|                                     |    |
|-------------------------------------|----|
| 1. Overview                         | 1  |
| 2. Protocol Summary                 | 2  |
| 3. Precautions                      | 3  |
| 4. Storage and Stability            | 3  |
| 5. Limitations                      | 4  |
| 6. Materials Supplied               | 4  |
| 7. Materials Required, Not Supplied | 5  |
| 8. Technical Hints                  | 6  |
| 9. Reagent Preparation              | 7  |
| 10. Standard Preparation            | 9  |
| 11. Sample Preparation              | 10 |
| 12. Assay Procedure                 | 12 |
| 13. Calculations                    | 14 |
| 14. Typical data                    | 15 |
| 15. Typical sample values           | 16 |
| 16. Troubleshooting                 | 17 |
| 17. Notes                           | 18 |

# 1. Overview

The Human CD5L / CT-2 Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213760) is designed for the quantitative measurement of Human CD5L / CT-2 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 monoclonal antibody from mouse specific for CD5L / CT-2 has been pre-coated onto 96-well plates. Standards (Expression system for standard: NSO; Immunogen sequence: S20-G347) and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for CD5L / CT-2 is added subsequently and then followed by washing with 1X Wash Buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates are washed away with 1X Wash Buffer. HRP substrate TMB is used to visualize HRP enzymatic reaction. TMB is catalyzed by HRP to produce a blue color product that changed into yellow after adding acidic TMB Stop Solution. The density of yellow is proportional to the Human CD5L / CT-2 amount of sample captured in plate.

CD5 antigen-like, also known as Sp alpha and AIM, is a protein that in humans is encoded by the CD5L gene. It is mapped to 1q21-q23 by fluorescence *in situ* hybridization. It is found that Aim expression is induced in mouse macrophages in response to loading with highly oxidized low density lipoprotein (oxLDL), and that Aim is expressed in foam cells within atherosclerotic lesions. Both the expression of Aim in lesions and its induction by oxLDL require Lxr /Rxr heterodimers. Aim-null macrophages are highly susceptible to oxLDL-induced apoptosis *in vitro* and undergo accelerated apoptosis in atherosclerotic lesions *in vivo*. Double knockout of Aim and Ldlr reduce atherosclerotic lesions. Therefore, it is concluded that AIM expression protects macrophages from apoptosis within atherosclerotic lesions, promoting early lesion development.

## 2. Protocol Summary

Prepare all reagents, samples, and standards as instructed



Add 100  $\mu$ L standard or sample to appropriate wells

Incubate at 37°C for 90 minutes



Discard plate content. Do not wash.

Add 100  $\mu$ L biotinylated Antibody in to all wells

Incubate at 37°C for 60 minutes



Wash each well three times with 300  $\mu$ L of 1X Wash Buffer



Add 100  $\mu$ L ABC working solution

Incubate at 37°C for 30 minutes



Wash each well five times with 300  $\mu$ L of 1X Wash Buffer

Add 90  $\mu$ L of prepared TMB Color Developing Agent

Incubate at 37°C in dark for 15-25 minutes



Add 100  $\mu$ L TMB Stop Solution and read OD at 450 nm within 30 minutes

### 3. Precautions

**Please read these instructions carefully prior to beginning the ELISA assay.**

- All kit components have been formulated and quality control tested to function successfully as a kit.
- We understand that, occasionally, experimental protocols might need to be modified to meet unique experimental circumstances. However, we cannot guarantee the performance of the product outside the conditions detailed in this protocol booklet.
- Reagents should be treated as possible mutagens and should be handled with care and disposed of properly. Please review the Safety Datasheet (SDS) provided with the product for information on the specific components.
- Observe good laboratory practices. Gloves, lab coat, and protective eyewear should always be worn. Never pipette by mouth. Do not eat, drink or smoke in the laboratory areas.
- All biological materials should be treated as potentially hazardous and handled as such. They should be disposed of in accordance with established safety procedures.

### 4. Storage and Stability

**Store ELISA kit at -20°C immediately upon receipt.**

Refer to list of materials supplied for storage conditions of individual components. Observe the storage conditions for individual prepared components in the Materials Supplied section.

Aliquot components in working volumes before storing at the recommended temperature.

## 5. Limitations

- ELISA kit intended for research use only. Not for use in diagnostic procedures.
- Do not mix or substitute reagents or materials from other kit lots or vendors. Kits are QC tested as a set of components and performance cannot be guaranteed if utilized separately or substituted.

## 6. Materials Supplied

| Item                                                       | Quantity          | Storage Condition<br>(Before prep) | Storage Condition<br>(After prep) |
|------------------------------------------------------------|-------------------|------------------------------------|-----------------------------------|
| Anti-Human CD5L / CT-2 Coated Microplate<br>(12 x 8 wells) | 1 x 96 well plate | -20°C                              | -20°C                             |
| Lyophilized recombinant Human CD5L / CT-2 Standard         | 2 x 1 vial        | -20°C                              | -20°C                             |
| Biotinylated anti- Human CD5L / CT-2 Antibody              | 100 µL            | -20°C                              | -20°C                             |
| Avidin-Biotin-Peroxidase Complex (ABC)                     | 100 µL            | -20°C                              | -20°C                             |
| Sample Diluent Buffer                                      | 30 mL             | -20°C                              | -20°C                             |
| Antibody Diluent Buffer                                    | 12 mL             | -20°C                              | -20°C                             |
| ABC Diluent Buffer                                         | 12 mL             | -20°C                              | -20°C                             |
| TMB Color Developing Agent                                 | 10 mL             | -20°C                              | -20°C                             |
| TMB Stop Solution                                          | 10 mL             | -20°C                              | -20°C                             |
| Adhesive Plate Seal                                        | 4                 | -20°C                              | -20°C                             |
| Wash Buffer (25X)                                          | 20 mL             | -20°C                              | -20°C                             |

## 7. Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully perform this assay:

- Microplate reader capable of measuring absorbance at 450 nm.
- Automated plate washer.
- Multi- and single-channel pipettes.
- Clean tubes and Eppendorf tubes.

## 8. Technical Hints

- Samples generating values higher than the highest standard should be further diluted in the appropriate sample dilution buffers.
- Avoid foaming or bubbles when mixing or reconstituting components.
- Avoid cross contamination of samples or reagents by changing tips between sample, standard and reagent additions.
- Ensure plates are properly sealed or covered during incubation steps.
- Don't let the 96-well plate dry, for a dry plate will inactivate active components on plate.
- Complete removal of all solutions and buffers during wash steps is necessary to minimize background.
- All samples should be mixed thoroughly and gently.
- Avoid multiple freeze/thaw of samples.
- When generating positive control samples, it is advisable to change pipette tips after each step.
- Before using the kit, spin tubes and bring down all components to the bottom of tubes.
- 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 Color Developing Agent solution will be pre-warmed in 37°C for 30 minutes before using.
- To avoid high background always add samples or standards to the well before the addition of the antibody cocktail.
- This kit is sold based on number of tests. A 'test' simply refers to a single assay well. The number of wells that contain sample, control or standard will vary by product. Review the protocol completely to confirm this kit meets your requirements. Please contact our Technical Support staff with any questions.



## 9. Reagent Preparation

- Equilibrate all reagents to room temperature (18-25°C) prior to use. The kit contains enough reagents for 96 wells.
- Prepare only as much reagent as is needed on the day of the experiment.

### 9.1 Anti-human CD5L / CT-2 coated Microplate (12 x 8 wells)

One plate of 96 wells. Ready to use. Store at -20°C.

### 9.2 Lyophilized recombinant Human CD5L / CT-2 standard (2 x 10 ng)

- 9.2.1 CD5L / CT-2 standard solution should be prepared no more than 2 hours prior to the experiment. Two tubes of CD5L / CT-2 standard (2 x 10 ng) are included in each kit. Use one tube for each experiment.
- 9.2.2 Add 1 mL sample diluent buffer into one tube to create 10,000 pg/mL of Human CD5L / CT-2 stock solution. Keep the tube at room temperature for 10 minutes and mix thoroughly.

### 9.3 Biotinylated anti- Human CD5L / CT-2 antibody

The solution should be prepared no more than 2 hours prior to the experiment.

- 9.3.1 The total volume should be: 100 µL/well x (the number of wells). (Allowing 100 µL – 200 µL more than total volume)
- 9.3.2 Biotinylated anti-Human CD5L / CT-2 antibody should be diluted in 1:100 with the antibody diluent buffer and mixed thoroughly. (i.e. Add 1 µL Biotinylated Anti-Human CD5L / CT-2 antibody to 99 µL antibody diluent buffer.)

### 9.4 Avidin-Biotin-Peroxidase Complex (ABC)

The solution should be prepared no more than 1 hour prior to the experiment.

- 9.4.1 The total volume should be: 100 µL/well x (the number of wells). (Allowing 100 µL - 200 µL more than total volume)
- 9.4.2 Avidin- Biotin-Peroxidase Complex (ABC) should be diluted in 1:100 with the ABC dilution buffer and mixed thoroughly. (i.e. Add 1 µL ABC to 99 µL ABC diluent buffer.)

**9.5 Sample diluent buffer**

30 mL. Ready to use. Store at -20°C.

**9.6 Antibody diluent buffer**

12 mL. Ready to use. Store at -20°C.

**9.7 ABC diluent buffer**

12 mL. Ready to use. Store at -20°C.

**9.8 TMB Color Developing Agent**

10 mL. Ready to use. Store at -20°C.

**9.9 TMB Stop Solution**

10 mL. Ready to use. Store at -20°C.

**9.10 1X Wash Buffer**

Prepare 500 mL of working 1X Wash Buffer by diluting 20 ml of the supplied Wash Buffer (25X) with 480 ml of deionized or distilled water. If crystals have formed in the concentrate, warm to room temperature and mix it gently until crystals have completely dissolved.

## 10. Standard Preparation

- 10.1** To prepare standards, label 6 Eppendorf tubes with 5,000 pg/mL, 2,500 pg/mL, 1,250 pg/mL, 625 pg/mL, 312.5 pg/mL and 156.25 pg/mL respectively.
- 10.2** Aliquot 300 µL of the sample diluent buffer into each tube.
- 10.3** Add 300 µL of the above 10,000 pg/mL CD5L / CT-2 solution into 1<sup>st</sup> tube and mix.
- 10.4** Transfer 300 µL from 1<sup>st</sup> tube to 2<sup>nd</sup> tube and mix. Transfer 300 µL from 2<sup>nd</sup> tube to 3<sup>rd</sup> tube and mix, and so on.

| Tube # | Volume to dilute                      | Volume of diluent | Concentration (pg/mL) |
|--------|---------------------------------------|-------------------|-----------------------|
| 1      | 300 µL of 10,000 pg/mL stock solution | 300 µL            | 5,000                 |
| 2      | 300 µL of tube #1                     | 300 µL            | 2,500                 |
| 3      | 300 µL of tube #2                     | 300 µL            | 1,250                 |
| 4      | 300 µL of tube #3                     | 300 µL            | 625                   |
| 5      | 300 µL of tube #4                     | 300 µL            | 312.5                 |
| 6      | 300 µL of tube #5                     | 300 µL            | 156.25                |

**Δ Note:** The standard solutions are best used within 2 hours. The 10,000 pg/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.

## 11. Sample Preparation

Store samples to be assayed within 24 hours at 4°C. For long-term storage, aliquot and freeze samples at -20°C. Avoid repeated freeze-thaw cycles.

- Cell lysates: After sufficient splitting, there should be no obvious cell sediment. Centrifuge cell lysates at approximately 10,000 x g for 5 minutes. Collect the cell lysate supernatants to go ahead.
- Serum: Allow the serum to clot in a serum separator tube (about 4 hours) at room temperature. Centrifuge at approximately 1,000 x g for 15 minutes. Analyze the serum immediately or aliquot and store samples at -20°C.
- Cell culture supernatant: Remove particulates by centrifugation, assay immediately or aliquot and store samples at -20°C.
- Plasma: Collect plasma using heparin or EDTA as an anticoagulant. Centrifuge for 15 minutes at 1,500 x g within 30 minutes of collection. Assay immediately or aliquot and store samples at -20°C.

### Activate the sample:

CD5 combines with its ligand (CD5L / CT-2) to form complex in samples. Free CD5L / CT-2 needs to be released by activation.

### Activating Reagent:

**Solution A:** 1N HCl: add 8.33 mL of 12N HCl into 91.67 mL of H<sub>2</sub>O.

**Solution B:** 1.2N NaOH/0.5M HEPES: add 12 mL of 10N NaOH and 11.9 g HEPES into 75 mL of H<sub>2</sub>O, add H<sub>2</sub>O to adjust volume to 100 mL.

- Cell culture supernatant: add activating reagent pro rate, i.e. add 20 µL of Solution A into 40 µL of sample, 10 minutes later, add 20 µL of Solution B. pH7.0-7.6.
- Serum: add activating reagent pro rate, i.e. add 20 µL of Solution A into 40 µL of sample, 10 minutes later, add 20 µL of Solution B. pH7.0-7.6. Prior to the assay, the activated sample requires a 2000-fold dilution. Add 10 µL of activated sample into 390 µL wash buffer (1:40). And then add 4 µL of above diluted reagent into 196 µL sample diluent buffer (1:50).

- Plasma (heparin, EDTA): add activating reagent pro rate, i.e. add 20  $\mu\text{L}$  of Solution A into 40  $\mu\text{L}$  of sample, 10 minutes later, add 20  $\mu\text{L}$  of Solution B. pH7.0-7.6. Prior to the assay, the activated sample requires a 4000-fold dilution. Add 10  $\mu\text{L}$  of activated sample into 790  $\mu\text{L}$  wash buffer (1:80). And then add 4  $\mu\text{L}$  of above diluted reagent into 196  $\mu\text{L}$  sample diluent buffer (1:50). Optimal dilutions should be determined by end users.
  - It is unnecessary to activate the recombinant CD5L / CT-2.
- Δ Note:** Samples are diluted partly by addition of activating reagent, so please pay attention to this when calculating target protein concentration.

It is recommended 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 (100,000 pg/mL-1,000,000 pg/mL). The working dilution is 1:100. i.e. Add 1  $\mu\text{L}$  sample into 99  $\mu\text{L}$  sample diluent buffer.
- Medium target protein concentration (10,000 pg/mL-100,000 pg/mL). The working dilution is 1:10. i.e. Add 10  $\mu\text{L}$  sample into 90  $\mu\text{L}$  sample diluent buffer.
- Low target protein concentration (156 pg/mL-10,000 pg/mL). The working dilution is 1:2. i.e. Add 50  $\mu\text{L}$  sample to 50  $\mu\text{L}$  sample diluent buffer.
- Very Low target protein concentration (0 pg/mL-156 pg/mL). No dilution necessary, or the working dilution is 1:2.

## 12. Assay Procedure

- It is recommended to assay all standards, controls and samples in duplicate.
  - The ABC working solution and TMB color developing agent must be kept warm at 37°C for 30 minutes before use. When diluting samples and reagents, they must be mixed completely and evenly. Standard CD5L / CT-2 detection curve should be prepared for each experiment. The user will decide sample dilution fold by crude estimation of CD5L / CT-2 amount in samples.
- 12.1** Aliquot 100 µL per well of the 10,000 pg/mL, 5,000 pg/mL, 2,500 pg/mL, 1,250 pg/mL, 625 pg/mL, 312.5 pg/mL, 156.25 pg/mL Human CD5L / CT-2 standard solutions into the pre-coated 96-well plate.
  - 12.2** Add 100 µL of the sample diluent buffer into the control well (Zero well).
  - 12.3** Add 100 µL of each properly diluted sample of Human cell culture supernatants, cell lysates, tissue homogenates, serum or plasma (heparin, EDTA) to each empty well. See "Sample Preparation" above for details. It is recommended that each Human CD5L / CT-2 standard solution and each sample be measured in duplicate.
  - 12.4** Seal the plate with a new adhesive cover provided and incubate at 37°C for 90 minutes.
  - 12.5** 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.
  - 12.6** Add 100 µL of biotinylated anti-Human CD5L / CT-2 antibody working solution into each well, seal the plate with a new adhesive cover provided and incubate at 37°C for 60 minutes.
  - 12.7** Wash plate 3 times with 1X Wash Buffer, and each time let washing buffer stay in the wells for 1 minute. 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 300 µL of 1X Wash 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 1X Wash Buffer, overfilling wells with 1X Wash Buffer. Blot the plate onto paper towels or other absorbent material.)

- 12.8** Add 100  $\mu$ L of prepared ABC working solution into each well, seal the plate with a new adhesive cover provided and incubate at 37°C for 30 minutes.
- 12.9** Wash plate 5 times with 1X Wash Buffer, and each time let washing buffer stay in the wells for 1-2 minutes. Discard the washing buffer and blot the plate onto paper towels or other absorbent material. (See Step 12.7 for plate washing method.)
- 12.10** Add 90  $\mu$ L of prepared TMB color developing agent into each well, seal the plate with a new adhesive cover and incubate at 37°C in dark for 15-25 minutes.

**Δ Note:** 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 Human CD5L / CT-2 standard solutions; the other wells show no obvious color.

- 12.11** Add 100  $\mu$ L of prepared TMB Stop Solution into each well. The color changes into yellow immediately.
- 12.12** Read the O.D. absorbance at 450 nm in a microplate reader within 30 minutes after adding the TMB Stop Solution.

## 13. Calculations

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 Human CD5L / CT-2 concentration of the samples can be interpolated from the standard curve.

(the relative O.D.450) = (the O.D.450 of each well) – (the O.D.450 of Zero well).

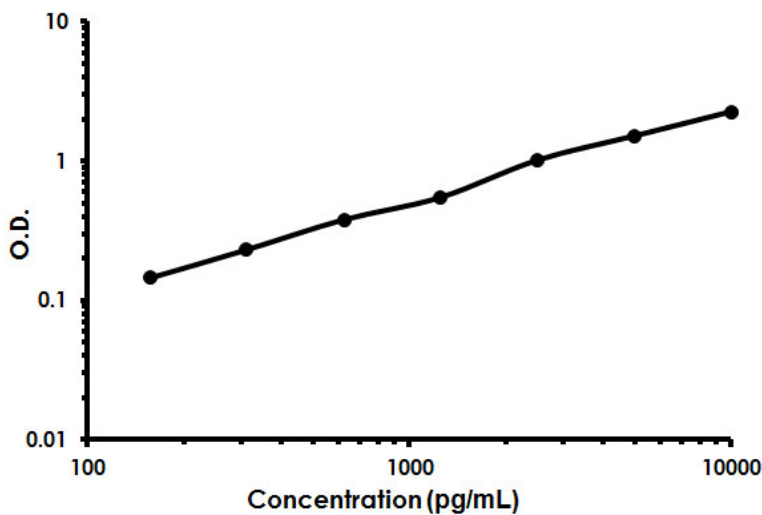
**Δ Note:** if the samples measured were diluted, multiply the dilution factor to the concentrations from interpolation to obtain the concentration before dilution.



# 14. Typical data

**Typical standard** curve – Data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.

| Sample | Human CD5L / CT-2 (pg/mL) | O.D.  |
|--------|---------------------------|-------|
| 1      | 0                         | 0.091 |
| 2      | 156                       | 0.145 |
| 3      | 312                       | 0.231 |
| 4      | 625                       | 0.379 |
| 5      | 1,250                     | 0.547 |
| 6      | 2,500                     | 1.009 |
| 7      | 5,000                     | 1.503 |
| 8      | 10,000                    | 2.241 |



**Figure 1.** Human CD5L / CT-2 ELISA Kit (ab213760) Standard Curve.

## 15. Typical sample values

### Sensitivity –

The biological sensitivity of the assay is <10 pg/mL.

The range is 156 pg/mL – 10,000 pg/mL.

### Precision –

**Intra-assay precision:** (Precision within an assay) Three samples of known concentration were tested on one plate to assess intra-assay precision.

| Sample | Number of measures | Mean (pg/mL) | Standard Deviation | CV% |
|--------|--------------------|--------------|--------------------|-----|
| 1      | 16                 | 245          | 10.04              | 4.1 |
| 2      | 16                 | 1,172        | 79.69              | 6.8 |
| 3      | 16                 | 5,450        | 414.2              | 7.6 |

**Inter-assay precision:** (Precision between assays) Three samples of known concentration were tested in separate assays to assess inter-assay precision.

| Sample | Number of assays | Mean (pg/mL) | Standard Deviation | CV% |
|--------|------------------|--------------|--------------------|-----|
| 1      | 24               | 224          | 13.21              | 5.9 |
| 2      | 24               | 1,288        | 91.44              | 7.1 |
| 3      | 24               | 5,192        | 415.36             | 8.0 |

### Specificity:

Natural and recombinant Human CD5L / CT-2.

### Cross-reactivity:

There is no detectable cross-reactivity with other relevant proteins.

## 16.Troubleshooting

| Problem             | Cause                                                      | Solution                                                                                                   |
|---------------------|------------------------------------------------------------|------------------------------------------------------------------------------------------------------------|
| Poor standard curve | Inaccurate Pipetting                                       | Check Pipettes                                                                                             |
|                     | Improper standard dilution                                 | Prior to opening, briefly spin the stock standard tube and dissolve the powder thoroughly by gentle mixing |
| Low Signal          | Incubation times too brief                                 | Ensure sufficient incubation times standard/sample incubation                                              |
|                     | Inadequate reagent volumes or improper dilution            | Check Pipettes and ensure correct preparation                                                              |
|                     | Incubation times with TMB Color Developing Agent too brief | Ensure sufficient incubation time until blue color develops prior addition of TMB Stop Solution            |
| Large CV            | Plate is insufficiently washed                             | Review manual for proper wash technique. If using a plate washer, check all ports for obstructions.        |
|                     | Contaminated wash buffer                                   | Prepare fresh wash buffer                                                                                  |
| Low sensitivity     | Improper storage of the ELISA kit                          | All components -20°C.<br>Keep TMB substrate solution protected from light.                                 |

## 17. Notes



## Technical Support

Copyright © 2024 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

[www.abcam.com/contactus](http://www.abcam.com/contactus)

[www.abcam.cn/contactus](http://www.abcam.cn/contactus) (China)

[www.abcam.co.jp/contactus](http://www.abcam.co.jp/contactus) (Japan)