

ab119523 – E-selectin Monkey ELISA Kit

Instructions for Use

For the quantitative measurement of Monkey E-selectin concentrations in cell culture supernatant and serum.

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

Table of Contents

INTRODUCTION

- 1. BACKGROUND** 2
- 2. ASSAY SUMMARY** 4

GENERAL INFORMATION

- 3. PRECAUTIONS** 5
- 4. STORAGE AND STABILITY** 5
- 5. MATERIALS SUPPLIED** 5
- 6. MATERIALS REQUIRED, NOT SUPPLIED** 6
- 7. LIMITATIONS** 6
- 8. TECHNICAL HINTS** 7

ASSAY PREPARATION

- 9. REAGENT PREPARATION** 8
- 10. STANDARD PREPARATIONS** 10
- 11. SAMPLE COLLECTION AND STORAGE** 12
- 12. PLATE PREPARATION** 13

ASSAY PROCEDURE

- 13. ASSAY PROCEDURE** 14

DATA ANALYSIS

- 14. CALCULATIONS** 16
- 15. TYPICAL DATA** 17
- 16. TYPICAL SAMPLE VALUES** 18
- 17. ASSAY SPECIFICITY** 19

RESOURCES

- 18. TROUBLESHOOTING** 20
- 19. NOTES** 21

1. BACKGROUND

Abcam's E-selectin monkey *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for accurate quantitative measurement of monkey E-selectin concentrations in cell culture supernatant and serum.

E-selectin specific antibodies have been precoated onto 96-well plates. Standards and test samples are added to the wells along with an HRP-conjugated E-selectin detection antibody. The microplate is then incubated at room temperature. After the removal of unbound proteins by washing TMB is added and catalyzed by HRP to produce a blue color product that changes to yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the amount of E-selectin captured on the plate.

E-selectin (ELAM-1, Endothelial Leukocyte Adhesion Molecule-1) belongs to the selectin family of adhesion molecules. Together with LECAM-1 (L-selectin) and GMP-140 (P-selectin), E-selectin mediates the initial interactions of leukocytes and platelets with endothelial cells.

Molecular structure: The extracellular part of all selectins consists of an amino terminal c-type lectin domain which specifically binds to carbohydrate ligands. This is followed by an EGF-like domain, and, in the case of E-selectin, by 6 short consensus repeats. The transmembrane portion of the molecule is followed by a short cytoplasmic tail.

Selectins guide non-activated polymorphonuclear cells to the areas of inflammation in creating first, loose contacts with the endothelial layer. The potential binding partner of E-selectin contains sialyl LewisX oligosaccharide. Other suitable ligands for the lectin domain of E-selectin are sialylated, fucosylated lactosaminoglycans. Together with GMP-140, E-selectin is expressed on cytokine-activated endothelial cells, and contributes to the adhesion of still resting leukocytes to the endothelium. This initial binding event is an essential prerequisite for

the activation of the immune cells via different inflammatory mediators. In contrast to GMP-140, E-selectin is maximally expressed 2-4 hours after cell activation. Within the next 24-48 hours E-selectin is again eliminated from the cytoplasmic membrane by shedding into the circulation. The circulating form or soluble (E-selectin) of this selectin exerts chemotactical signals on neutrophils and additionally activates the 2-integrins - E-selectin assists in preparing the migration capacity of these cells.

Determination of monkey E-selectin could provide more detailed insights into the pathological modifications during various diseases.

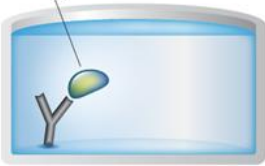
2. ASSAY SUMMARY

Primary capture antibody



Equilibrate all reagents to room temperature. Prepare all the reagents, samples, and standards as instructed.

Sample



Add standard or sample to each well. Incubate the plate.

HRP conjugated antibody



Wash and add prepared HRP labeled secondary detector antibody. Incubate at room temperature

Substrate **Colored product**



Wash and add TMB Substrate Solution to each well. Add Stop Solution to each well. Read immediately.

3. PRECAUTIONS

Please read these instructions carefully prior to beginning the assay.

All kit components have been formulated and quality control tested to function successfully as a kit. Modifications to the kit components or procedures may result in loss of performance.

4. STORAGE AND STABILITY

Upon receipt, store kit immediately at 2-8°C.

Refer to list of materials supplied for storage conditions of individual components. Observe the storage conditions for individual prepared components in section 9 Reagent Preparation.

5. MATERIALS SUPPLIED

Item	Amount	Storage Condition (Before Preparation)
Microplate coated with monoclonal antibody to E-selectin (12 x 8 wells)	96 wells	2-8 °C
HRP Conjugated anti-E-selectin monoclonal antibody	150 µL	2-8 °C
E-selectin Standard lyophilized	2 Vials	2-8 °C
Sample Diluent	12 mL	2-8 °C
20X Assay Buffer Concentrate	5 mL	2-8 °C
20X Wash Buffer Concentrate	50 mL	2-8 °C
TMB Substrate Solution	15 mL	2-8 °C
Stop Solution (1M Phosphoric acid)	15 mL	2-8 °C

6. MATERIALS REQUIRED, NOT SUPPLIED

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

- 5 mL and 10 mL graduated pipettes
- 5 μ L to 1000 μ L adjustable single channel micropipettes with disposable tips
- 50 μ L to 300 μ L adjustable multichannel micropipette with disposable tips
- Multichannel micropipette reservoir
- Beakers, flasks, cylinders necessary for preparation of reagents
- Device for delivery of wash solution (multichannel wash bottle or automatic wash system)
- Microplate strip reader capable of reading at 450 nm (620 nm as optional reference wave length)
- Glass-distilled or deionized water
- Statistical calculator with program to perform regression analysis

7. LIMITATIONS

- Assay kit intended for research use only. Not for use in diagnostic procedures
- Do not use kit or components if it has exceeded the expiration date on the kit labels
- 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

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
- Complete removal of all solutions and buffers during wash steps.
- As exact conditions may vary from assay to assay, a standard curve must be established for every run.
- Disposable pipette tips, flasks or glassware are preferred, reusable glassware must be washed and thoroughly rinsed of all detergents before use.
- Improper or insufficient washing at any stage of the procedure will result in either false positive or false negative results. Empty wells completely before dispensing fresh wash solution, fill with Wash Buffer as indicated for each wash cycle and do not allow wells to sit uncovered or dry for extended periods.
- **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 and samples to room temperature (18-25°C) prior to use.

9.1 1X Wash Buffer

Prepare 1X Wash Buffer by diluting the 20X Wash Buffer Concentrate with distilled or deionized water. To make 500 mL 1X Wash Buffer, combine 25 mL 20X Wash Buffer Concentrate with 475 mL distilled or deionized water. Mix thoroughly and gently to avoid foaming.

Note: The 1X Wash Buffer should be stored at 2-8 °C and is stable for 30 days.

9.2 1X Assay Buffer

Prepare 1X Assay Buffer by diluting the 20X Assay Buffer Concentrate with distilled or deionized water. To make 50 mL 1X Assay Buffer, combine 2.5 mL 20X Assay Buffer Concentrate with 47.5 mL distilled or deionized water. Mix thoroughly and gently to avoid foaming.

Note: The 1X Assay Buffer should be stored at 2-8 °C and is stable for 30 days.

9.3 1X HRP Conjugated Antibody

To prepare the HRP Conjugated Antibody, dilute the anti-HRP Conjugated Antibody 100-fold with 1X Assay Buffer. Use the following table as a guide to prepare as much 1X HRP Conjugated Antibody as needed by adding the required volume (μL) of the HRP Conjugated Antibody to the required volume (mL) of 1X Assay Buffer. Mix gently and thoroughly.

Number of strips	Volume of HRP-Conjugate anti-Monkey E-selectin antibody (μL)	Volume of 1X Assay Buffer (mL)
1 - 6	30	2.97
7 - 12	60	5.94

Note: The 1X Biotin-Conjugated Antibody should be used within 30 minutes after dilution.

10. STANDARD PREPARATIONS

Prepare serially diluted standards immediately prior to use. Always prepare a fresh set of standards for every use.

- 10.1 Prepare a 120 U/mL **Stock Standard** by reconstituting one vial of the E-selectin standard with the volume of distilled water stated on the label. Hold at room temperature for 10-30 minutes. The 120 U/mL **Stock Standard** cannot be stored for later use.
- 10.2 Label eight tubes with numbers 1 - 7.
- 10.3 Add 225 μ L Sample Diluent to all tubes.
- 10.4 Prepare a 60 U/mL **Standard 1** by transferring 225 μ L of the 120 U/mL Stock Standard to tube 1. Mix thoroughly and gently.
- 10.5 Prepare **Standard 2** by transferring 225 μ L from Standard 1 to tube 2. Mix thoroughly and gently.
- 10.6 Prepare **Standard 3** by transferring 225 μ L from Standard 2 to tube 3. Mix thoroughly and gently.
- 10.7 Using the table below as a guide, repeat for tubes 4 through to 6.
- 10.8 **Standard 7** contains no protein and is the Blank control

ASSAY PREPARATION

Standard	Sample to Dilute	Volume to Dilute (μL)	Volume of Diluent (μL)	Starting Conc. (U/mL)	Final Conc. (U/mL)
1	Stock	225	225	120	60
2	Standard 1	225	225	60	30
3	Standard 2	225	225	30	15
4	Standard 3	225	225	15	7.5
5	Standard 4	225	225	7.5	3.8
6	Standard 5	225	225	3.8	1.9
7	None	-	225	-	0



11. SAMPLE COLLECTION AND STORAGE

- Cell culture supernatant and serum were tested with this assay. Other biological samples might be suitable for use in the assay. Remove serum from the clot as soon as possible after clotting.
- Samples containing a visible precipitate must be clarified prior to use in the assay. Do not use grossly hemolyzed or lipemic specimens.
- Samples should be aliquoted and must be stored frozen at -20°C to avoid loss of bioactive monkey E-selectin. If samples are to be run within 24 hours, they may be stored at 2° to 8°C.
- Avoid repeated freeze-thaw cycles. Prior to assay, the frozen sample should be brought to room temperature slowly and mixed gently.
- Aliquots of serum samples (spiked or unspiked) were stored at -20°C and thawed 5 times, and the monkey E-selectin levels determined. There was no significant loss of monkey E-selectin immunoreactivity detected by freezing and thawing.
- Aliquots of serum samples (spiked or unspiked) were stored at -20°C, 2-8°C, room temperature (RT) and at 37°C, and the monkey E-selectin level determined after 24 h. There was no significant loss of monkey E-selectin immunoreactivity detected during storage under above conditions.

12. PLATE PREPARATION

- The 96 well plate strips included with this kit are supplied ready to use
- Unused well strips should be returned to the plate packet and stored at 2-8°C
- For statistical reasons, we recommend each sample should be assayed with a minimum of two replicates (duplicates)
- Well effects have not been observed with this assay

13. ASSAY PROCEDURE

- **Equilibrate all materials and prepared reagents to room temperature prior to use.**
 - **It is recommended to assay all standards, controls and samples in duplicate.**
- 13.1. Prepare all reagents, working standards, and samples as directed in the previous sections. Determine the number of microplate strips required to test the desired number of samples plus appropriate number of wells needed for running blanks and standards.
 - 13.2. Wash the microplate twice with approximately 400 μ L 1X Wash Buffer per well with thorough aspiration of microplate contents between washes. Allow the 1X Wash Buffer to remain in the wells for about 10 - 15 seconds before aspiration. Take care not to scratch the surface of the microplate.
 - 13.3. After the last wash step, empty wells and tap microplate on absorbent pad or paper towel to remove excess 1X Wash Buffer. Use the microplate strips immediately after washing. Alternatively the microplate strips can be placed upside down on a wet absorbent paper for not longer than 15 minutes. Do not allow wells to dry.
 - 13.4. Add 100 μ L of prepared standards (including the no standard blank control) to the appropriate wells.
 - 13.5. Add 80 μ L of Sample Diluent to the sample wells.
 - 13.6. Add 20 μ L of samples to appropriate wells.
 - 13.7. Add 50 μ L of 1X HRP Conjugated Antibody to all wells.
 - 13.8. Cover with adhesive film and incubate at room temperature (18° to 25°C) for 2 hours (microplate can be incubated on a shaker set at 400 rpm).
 - 13.9. Remove adhesive film and empty wells. Wash microplate strips 5 times according to step 13.2. Proceed immediately to step 13.10.

13.10. Pipette 100 μ L of TMB Substrate Solution to all wells.

13.11. Incubate the microplate strips at room temperature (18 to 25°C) for 10 minutes. Avoid direct exposure to intense light.

Note: The color development on the plate should be monitored and the substrate reaction stopped (see step 13.12) before the signal in the positive wells becomes saturated. Determination of the ideal time period for color development should be done individually for each assay. It is recommended to add the stop solution when the highest standard has developed a dark blue color. Alternatively the color development can be monitored by the ELISA reader at 620 nm. The substrate reaction should be stopped as soon as Standard 1 has reached an OD of 0.9 - 0.95.

13.12. Stop the enzyme reaction by adding 100 μ L of Stop Solution into each well.

Note: It is important that the Stop Solution is mixed quickly and uniformly throughout the microplate to completely inactivate the enzyme. Results must be read immediately after the Stop Solution is added or within one hour if the microplate strips are stored at 2 - 8°C in the dark.

13.13. Read absorbance of each microplate on a spectrophotometer using 450 nm as the primary wave length (optionally 620 nm as the reference wave length; 610 nm to 650 nm is acceptable). Blank the plate reader according to the manufacturer's instructions by using the blank wells. Determine the absorbance of both the samples and the standards.

Note: In case of incubation without shaking the obtained O.D. values may be lower than indicated below. Nevertheless the results are still valid.

14. CALCULATIONS

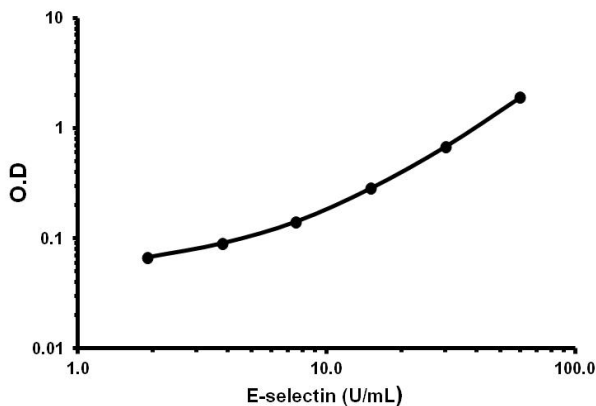
Average the duplicate readings for each standard, sample and control blank. Subtract the no protein control blank from all mean readings. Plot the mean standard readings against their concentrations and draw the best smooth curve through these points to construct a standard curve. Most plate reader software or graphing software can plot these values and curve fit. A five parameter algorithm (5PL) usually provides the best fit, though other equations can be examined to see which provides the most accurate (e.g. linear, semi-log, log/log, 5-parameter logistic). Extrapolate protein concentrations for unknown and control samples from the standard curve plotted. Samples producing signals greater than that of the highest standard should be further diluted in appropriate buffer and reanalyzed, then multiplying the concentration found by the appropriate dilution factor.

If samples have been diluted 1:5, as stated in step 13.6, the concentration obtained from the standard curve must be multiplied by the dilution factor ($\times 5$) to obtain an accurate value, in addition to any initial sample dilution factor.

Calculation of samples with a concentration exceeding standard 1 may result in incorrect, low monkey E-selectin levels. Such samples require further external predilution according to expected monkey E-selectin values with Sample Diluent in order to precisely quantitate the actual monkey E-selectin level.

15. TYPICAL DATA

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



Standard Curve Measurements			
Conc. (U/mL)	O.D. 450 nm		Mean
	1	2	O.D.
0	0.049	0.051	0.050
1.9	0.070	0.065	0.067
3.8	0.095	0.095	0.090
7.5	0.150	0.131	0.141
15.0	0.288	0.279	0.283
30	0.681	0.668	0.676
60	1.955	1.823	1.889

Figure 1. Example of Monkey E-selectin protein standard curve.

16. TYPICAL SAMPLE VALUES

SENSITIVITY -

The limit of detection of monkey E-selectin defined as the analyte concentration resulting in an absorbance significantly higher than that of the dilution medium (mean plus 2 standard deviations) was determined to be 0.4 U/mL (mean of 6 independent assays).

RECOVERY –

The spike recovery was evaluated by spiking 3 levels of monkey E-selectin into pooled normal monkey serum samples. Recoveries were determined in 3 independent experiments with 4 replicates each. The amount of endogenous monkey E-selectin in unspiked serum was subtracted from the spike values. The overall mean recovery was 85%.

LINEARITY OF DILUTION –

Serum samples with different levels of Monkey E-selectin were analyzed at serial 2 fold dilutions with 4 replicates each. The recovery ranged from 97.4% to 119.2% with an overall recovery of 105.2%

PRECISION –

Intra- and Inter-assay reproducibility was determined by measuring samples containing different concentrations of Monkey E-selectin.

	Intra-Assay	Inter-Assay
n=	4	4
%CV	< 5	< 10

17. ASSAY SPECIFICITY

The interference of circulating factors of the immune system was evaluated by spiking these proteins at physiologically relevant concentrations into a monkey E-selectin positive serum. There was no cross reactivity detected.

18. TROUBLESHOOTING

Problem	Cause	Solution
Poor standard curve	Inaccurate pipetting	Check pipettes
	Improper standards 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; change to overnight standard/sample incubation
	Inadequate reagent volumes or improper dilution	Check pipettes and ensure correct preparation
Samples give higher value than the highest standard	Starting sample concentration is too high.	Dilute the specimens and repeat the assay
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 kit	Store the all components as directed.

19. NOTES

UK, EU and ROW

Email: technical@abcam.com | Tel: +44-(0)1223-696000

Austria

Email: wissenschaftlicherdienst@abcam.com | Tel: 019-288-259

France

Email: supportscientifique@abcam.com | Tel: 01-46-94-62-96

Germany

Email: wissenschaftlicherdienst@abcam.com | Tel: 030-896-779-154

Spain

Email: soportecientifico@abcam.com | Tel: 911-146-554

Switzerland

Email: technical@abcam.com

Tel (Deutsch): 0435-016-424 | Tel (Français): 0615-000-530

US and Latin America

Email: us.technical@abcam.com | Tel: 888-77-ABCAM (22226)

Canada

Email: ca.technical@abcam.com | Tel: 877-749-8807

China and Asia Pacific

Email: hk.technical@abcam.com | Tel: 400 921 0189 / +86 21 2070 0500

Japan

Email: technical@abcam.co.jp | Tel: +81-(0)3-6231-0940

www.abcam.com | www.abcam.cn | www.abcam.co.jp