

Human sICAM-1 Coated ELISA Kit

PRODUCT INFORMATION & MANUAL

Enzyme-linked Immunosorbent Assay for quantitative detection of human sICAM-1.

Catalog Number **REF** *BMS201CE, BMS201TENCE*



For in-vitro diagnostic use. Not for therapeutic procedures.

Information in this document is subject to change without notice.

DISCLAIMER

TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information

These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

Corporate entity

Life Technologies | Carlsbad, CA 92008 USA | Toll free in USA 1.800.955.6288

Trademarks

All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. All other trademarks are the property of their respective owners.

© 2017 Thermo Fisher Scientific Inc. All rights reserved.



Bender MedSystems GmbH Campus Vienna Biocenter 2 1030 Vienna, Austria

TABLE OF CONTENTS

1. Intended Use	4
2. Summary	4
3. Principles of the Test	6
4. Reagents Provided.....	7
5. Storage Instructions – ELISA Kit.....	8
6. Specimen Collection and Storage Instructions.....	8
7. Materials Required But Not Provided	8
8. Precautions for Use	9
9. Preparation of Reagents	10
10. Test Protocol	12
11. Calculation of Results	14
12. Limitations	15
13. Performance Characteristics.....	16
14. Reagent Preparation Summary	19
15. Test Protocol Summary	20
PRODUKTINFORMATION UND HANDBUCH (Deutsch)	21
1. Mitgelieferte Reagenzien	21
2. Lagerhinweise	22
3. Sicherheitsvorkehrungen für den Gebrauch	22
4. Vorbereitung der Reagenzien.....	23
5. Testprotokoll	25
INFORMACIÓN Y MANUAL DEL PRODUCTO (Espanol).....	28
1. Reactivos Suministrados	28
2. Instrucciones de Conservación.....	29
3. Precauciones de uso.....	29
4. Preparación de los Reactivos	30
5. Protocolo de Ensayo	32
INFORMATIONS SUR LE PRODUIT ET MANUEL (Français)	34
1. Réactifs Fournis	34
2. Instruction de Stockage	35
3. Préventions de Sécurité pour l'Usage.....	35
4. Préparation des Réactifs	36
5. Protocole de Test	38
INFORMAZIONI SUL PRODOTTO E MANUALE (Italiano)	40

1. Reagenti Forniti	40
2. Istruzioni di Conservazione	41
3. Precauzioni per l'Uso	41
4. Preparazione dei Reagenti	42
5. Procedura del Test.....	44
Documentation and support.....	46

1. Intended Use

The human sICAM-1 ELISA is an enzyme-linked immunosorbent assay for the quantitative detection of human sICAM-1. **The human sICAM-1 ELISA is for in vitro diagnostic use. Not for use in therapeutic procedures.**

2. Summary

Intercellular Adhesion Molecule-1 (ICAM-1) is a member of the immunoglobulin supergene family and functions as a ligand for the Lymphocyte Function-Associated Antigen-1 (LFA-1), an alpha-beta- complex that is a member of the leukocyte integrin family of cell-cell and cell-matrix receptors. This family consists of the leukocyte adhesion glycoproteins LFA-1 which mediates lymphocyte adhesion, Mac-1 which mediates granulocyte adhesion and p150,95.

ICAM-1 is a single-chain glycoprotein with a polypeptide core of 55kD that can be expressed on non-hematopoietic cells of many lineages such as vascular endothelial cells, thymic epithelial cells, other epithelial cells and fibroblasts and on hematopoietic cells such as tissue macrophages, mitogen-stimulated T-lymphoblasts, germinal center B- cells and dendritic cells in tonsils, lymph nodes and Peyer's patches. ICAM-1 is inducible on fibroblasts and endothelial cells by inflammatory mediators such as IL-1, TNF and IFN gamma within few hours and is correlated to the infiltration of lymphocytes into inflammatory lesions. ICAM-1 seems to be the initial marker of inflammatory reactions and is expressed prior to, and to a greater extent than is HLA-DR.

The role of ICAM-1 as a disease marker has been demonstrated for a number of different indications and pathological situations.

ICAM-1 upregulation in **allergic airway inflammation** is responsible for the recruitment of activated leukocytes and the pathogenesis of allergic rhinitis.

In **allergic contact dermatitis** ICAM-1 on keratinocytes was induced already 4 hours after application of the allergic patch test.

In bladder cancer there is a direct correlation between constitutive ICAM-1 expression and the histopathologic grade of the tumor. Sera of **GI-cancer** patients with liver metastasis showed significant higher sICAM-1 levels than those of patients without metastasis.

ICAM-1 is expressed on malignant cells in myeloid as well as **B lymphoid malignancies**. In **lymphoproliferative disorders** ICAM-1 is related to the degree of malignancy. In HTLV-1 associated myelopathy, and adult T-cell leukemia sICAM-1 serum levels are elevated.

Patients with malignant **melanoma** have significantly increased serum levels of sICAM-1, which is of prognostic importance.

Significantly elevated concentrations of sICAM-1 are detected in **HIV-1** infected persons.

In **malaria tropica** ICAM-1 serves in the adhesion of infected erythrocytes to the capillary endothelium which event is important in the pathogenesis of cerebral malaria.

sICAM-1 is a good prognostic parameter of responsiveness of **hepatitis B** infection to IFN beta therapy.

ICAM-1 seems to provide the mechanism crucial to **allograft rejection of the cornea**.

Expression of ICAM-1 is also increased during rejection on the capillary endothelium, the myocardial membrane, and the endocardium of the **transplanted heart**.

sICAM-1 serum levels significantly increased with acute renal graft rejection. Measuring sICAM-1 is helpful in discriminating rejection from Cyclosporine-A intoxication of the **transplanted kidney**.

Strong expression of ICAM-1 is also seen in patients with acute rejection versus stable **liver transplants**, or patients with non-rejection complications.

Serum levels of circulating ICAM-1 and L-selectin were found elevated in insulin-dependent **diabetes mellitus** (IDDM) and in subjects at risk of IDDM.

Significant elevation of serum ICAM-1 has been demonstrated in anterior **uveitis** in intermediate uveitis and in patients with sarcoidosis. In the first 12-24 hours of monitoring acute **myocardial infarction** a decrease of sICAM-1 is measurable. This can provide prognostic significance to sICAM-1 also for myocardial **ischemia** and **reperfusion**.

Increased glomerular ICAM-1 expression is seen in early cases of different forms of **glomerulonephritis** and tubular *de novo* expression of ICAM-1 shows a strong correlation with disease activity.

In **asthma** ICAM-1 is upregulated on inflamed airway epithelium and bronchial endothelium, thereby mediating eosinophil adhesion. sICAM-1 is significantly elevated in the sera of patients with **idiopathic pulmonary fibrosis** or **sarcoidosis**.

sICAM-1 is a reliable marker for an inflammatory process within the **central nervous system** which is associated with blood: CSF barrier disturbance.

Soluble ICAM-1 is not detectable in most midtrimester amniotic fluid samples but when present is significantly related to **intrauterine growth retardation** and elevated midtrimester levels of maternal serum alpha fetoprotein.

Elevated levels of sICAM-1 correlate with the activity of the **rheumatoid arthritis**.

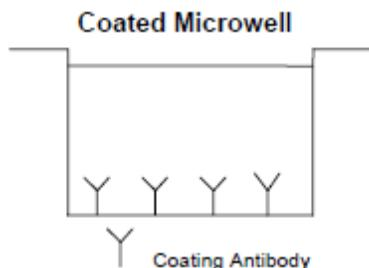
In **psoriasis** ICAM-1 on keratinocytes shows strong correlation with severity of disease and decreases under successful therapy. Before treatment sICAM-1 levels are significantly elevated compared to healthy controls.

For literature update refer to [www.Thermo Fisher.com](http://www.ThermoFisher.com).

3. Principles of the Test

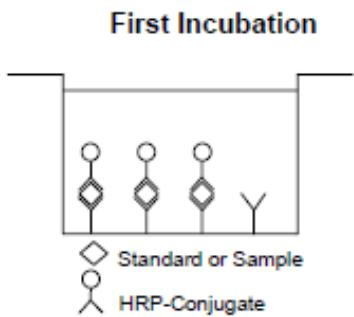
An anti-human sICAM-1 coating antibody is adsorbed onto microwells.

Figure 1



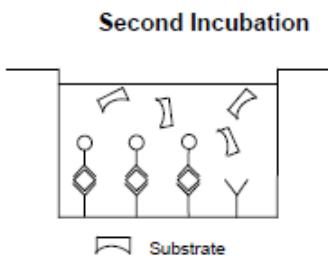
Human sICAM-1 present in the sample or standard binds to antibodies adsorbed to the microwells and the HRP-conjugated anti-human sICAM-1 antibody is added and binds to human sICAM-1 captured by the first antibody.

Figure 2



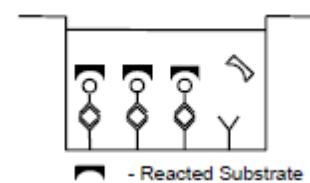
Following incubation unbound HRP-conjugated anti-human sICAM-1 is removed during a wash step, and substrate solution reactive with HRP is added to the wells.

Figure 3



A colored product is formed in proportion to the amount of human sICAM-1 present in the sample or standard. The reaction is terminated by addition of acid and absorbance is measured at 450 nm. A standard curve is prepared from 5 human sICAM-1 standard dilutions and human sICAM-1 concentration determined.

Figure 4



4. Reagents Provided

4.1 Reagents for human sICAM-1 ELISA BMS201CE (96 tests)

- 1 aluminum pouch with a **Microwell Plate coated** with monoclonal antibody to human sICAM-1
- 1 vial (100 µl) **HRP-Conjugate** anti-human sICAM-1 monoclonal antibody
- 2 vials (500 µl) human sICAM-1 **Standard**, 100 ng/ml
- 1 vial **Control high**, lyophilized
- 1 vial **Control low**, lyophilized
- 1 vial (12 ml) **Sample Diluent**
- 1 vial (5 ml) **Assay Buffer Concentrate** 20x (PBS with 1% Tween 20, 10% BSA)
- 1 bottle (50 ml) **Wash Buffer Concentrate** 20x (PBS with 1% Tween 20)
- 1 vial (15 ml) **Substrate Solution** (tetramethyl-benzidine)
- 1 vial (15 ml) **Stop Solution** (1M Phosphoric acid)
- 2 **Adhesive Films**

4.2 Reagents for human sICAM-1 ELISA BMS201TENCE (10x96 tests)

- 10 aluminum pouches with a **Microwell Plate coated** with monoclonal antibody to human sICAM-1
- 10 vials (100 µl) **HRP-Conjugate** anti-human sICAM-1 monoclonal antibody
- 10 vials (500 µl) human sICAM-1 **Standard**, 100 ng/ml 10 vials **Control high**, lyophilized
- 10 vials **Control low**, lyophilized 10 vials (12 ml) **Sample Diluent**
- 1 vial (5 ml) **Assay Buffer Concentrate** 20x (PBS with 1% Tween 20, 10% BSA)
- 3 bottles (50 ml) **Wash Buffer Concentrate** 20x (PBS with 1% Tween 20)
- 10 vials (15 ml) **Substrate Solution** (tetramethyl-benzidine)
- 1 vial (100 ml) **Stop Solution** (1M Phosphoric acid)
- 10 **Adhesive Films**

5. Storage Instructions – ELISA Kit

Store kit reagents between 2° and 8°C except controls. Store lyophilized controls at -20°C. Immediately after use remaining reagents should be returned to cold storage (2° to 8°C), or to -20°C, respectively. Expiry of the kit and reagents is stated on labels.

Expiry of the kit components can only be guaranteed if the components are stored properly, and if, in case of repeated use of one component, this reagent is not contaminated by the first handling.

6. Specimen Collection and Storage Instructions

Cell culture supernatant, serum, plasma (EDTA, citrate, heparin), amniotic fluid, spontaneous urine and bile were tested with this assay. Other biological samples might be suitable for use in the assay. Remove serum or plasma from the clot or cells as soon as possible after clotting and separation.

Pay attention to a possible “**Hook Effect**” due to high sample concentrations (see chapter 11).

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 human sICAM-1. If samples are to be run within 24 hours, they may be stored at 2° to 8°C (for sample stability refer to 13.5).

Avoid repeated freeze-thaw cycles. Prior to assay, the frozen sample should be brought to room temperature slowly and mixed gently.

To measure human sICAM-1 in spontaneous urine use undiluted samples.

Cell culture media without serum components are not suitable for human sICAM-1 determination with the ELISA.

7. Materials Required But Not Provided

- 5 ml and 10 ml graduated pipettes
- 5 µl to 1000 µl adjustable single channel micropipettes with disposable tips
- 50 µl to 300 µl 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)
- Microwell 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

8. Precautions for Use

- All reagents should be considered as potentially hazardous. We therefore recommend that this product is handled only by those persons who have been trained in laboratory techniques and that it is used in accordance with the principles of good laboratory practice. Wear suitable protective clothing such as laboratory overalls, safety glasses and gloves. Care should be taken to avoid contact with skin or eyes. In the case of contact with skin or eyes wash immediately with water. See material safety data sheet(s) and/or safety statement(s) for specific advice.
- Reagents are intended for in vitro diagnostic use and are not for use in therapeutic procedures.
- Do not mix or substitute reagents with those from other lots or other sources.
- Do not use kit reagents beyond expiration date on label.
- Do not expose kit reagents to strong light during storage or incubation.
- Do not pipette by mouth.
- Do not eat or smoke in areas where kit reagents or samples are handled.
- Avoid contact of skin or mucous membranes with kit reagents or specimens.
- Rubber or disposable latex gloves should be worn while handling kit reagents or specimens.
- Avoid contact of substrate solution with oxidizing agents and metal.
- Avoid splashing or generation of aerosols.
- In order to avoid microbial contamination or cross-contamination of reagents or specimens which may invalidate the test use disposable pipette tips and/or pipettes.
- Use clean, dedicated reagent trays for dispensing the conjugate and substrate reagent.
- Exposure to acid inactivates the conjugate.
- Glass-distilled water or deionized water must be used for reagent preparation.
- Substrate solution must be at room temperature prior to use.
- Decontaminate and dispose specimens and all potentially contaminated materials as they could contain infectious agents. The preferred method of decontamination is autoclaving for a minimum of 1 hour at 121.5°C.
- Liquid wastes not containing acid and neutralized waste may be mixed with sodium hypochlorite in volumes such that the final mixture contains 1.0% sodium hypochlorite. Allow 30 minutes for effective decontamination. Liquid waste containing acid must be neutralized prior to the addition of sodium hypochlorite.

9. Preparation of Reagents

Buffer Concentrates should be brought to room temperature and should be diluted before starting the test procedure. If crystals have formed in the **Buffer Concentrates**, warm them gently until they have completely dissolved.

9.1. Wash Buffer (1x)

Pour entire contents (50 ml) of the **Wash Buffer Concentrate (20x)** into a clean 1000 ml graduated cylinder. Bring to final volume of 1000 ml with glass-distilled or deionized water. Mix gently to avoid foaming.

Transfer to a clean wash bottle and store at 2° to 25°C. Please note that Wash Buffer (1x) is stable for 30 days.

Wash Buffer (1x) may also be prepared as needed according to the following table:

Number of Strips	Wash Buffer Concentrate (20x) (ml)	Distilled Water (ml)
1 - 6	25	475
1 - 12	50	950

9.2. Assay Buffer (1x)

Pour the entire contents (5 ml) of the **Assay Buffer Concentrate (20x)** into a clean 100 ml graduated cylinder. Bring to final volume of 100 ml with distilled water. Mix gently to avoid foaming.

Store at 2° to 8°C. Please note that the Assay Buffer (1x) is stable for 30 days.

Assay Buffer (1x) may also be prepared as needed according to the following table:

Number of Strips	Assay Buffer Concentrate (20x) (ml)	Distilled Water (ml)
1 - 6	2.5	47.5
1 - 12	5.0	95.0

9.3. HRP-Conjugate

Please note that the HRP-Conjugate should be used within 30 minutes after dilution.

Make a 1:100 dilution of the concentrated **HRP-Conjugate** solution with Assay Buffer (1x) in a clean plastic tube as needed according to the following table:

Number of Strips	HRP-Conjugate (ml)	Assay Buffer (1x) (ml)
1 - 6	0.03	2.97
1 - 12	0.06	5.94

9.4. Human sICAM-1 Standard

Standard dilutions can be prepared directly on the microwell plate (see 10.c) or alternatively in tubes (see 9.4.1).

9.4.1. External Standard Dilution

Label 4 tubes, one for each standard point.

S2, S3, S4, S5

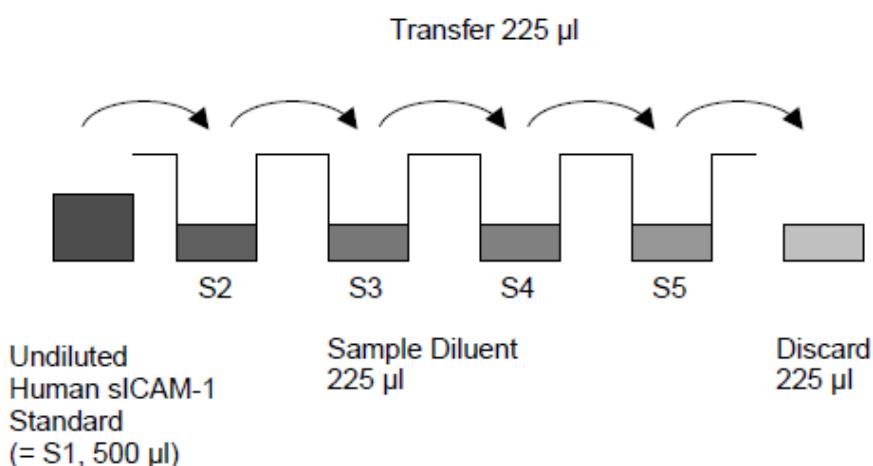
Then prepare 1:2 serial dilutions for the standard curve as follows: Pipette 225 µl of Sample Diluent into tubes S2 – S5.

Pipette 225 µl of undiluted standard (serves as the highest standard S1, concentration of standard 1 = 100 ng/ml) into the first tube, labeled S2, and mix thoroughly before the next transfer (concentration of standard 2 = 50 ng/ml).

Repeat serial dilutions 3 more times thus creating the points of the standard curve (see Figure 5).

Sample Diluent serves as blank.

Figure 5



9.5. Controls

Reconstitute by adding 100 µl distilled water to lyophilized controls (10- 30 minutes). Further treat the controls like your samples in the assay. For control range please refer to certificate of analysis or vial label. Store reconstituted controls aliquoted at -20°C. Avoid repeated freeze and thaw cycles.

10. Test Protocol

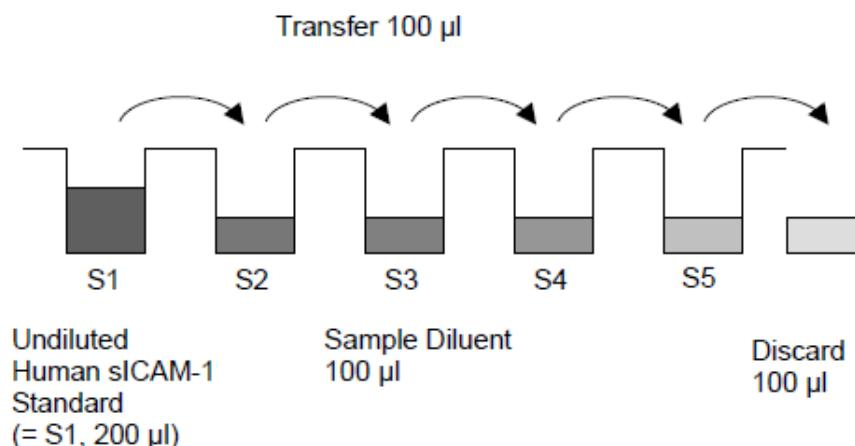
1. Determine the number of microwell strips required to test the desired number of samples plus appropriate number of wells needed for running blanks and standards. Each sample, standard, blank and optional control sample should be assayed in duplicate. Remove extra microwell strips from holder and store in foil bag with the desiccant provided at 2°-8°C sealed tightly.
2. Wash the microwell strips twice with approximately 400 µl **Wash Buffer** per well with thorough aspiration of microwell contents between washes. Allow the Wash Buffer to sit in the wells for about **10 – 15 seconds** before aspiration. Take care not to scratch the surface of the microwells.

After the last wash step, empty wells and tap microwell strips on absorbent pad or paper towel to remove excess Wash Buffer. Use the microwell strips immediately after washing. Alternatively microwell strips can be placed upside down on a wet absorbent paper for not longer than 15 minutes. **Do not allow wells to dry.**

3. **Standard dilution on the microwell plate** (Alternatively the standard dilution can be prepared in tubes - see 9.4.1):

Add 100 µl of Sample Diluent in duplicate to **standard wells B1/2- E1/2**, leaving A1/A2 empty. Pipette 200 µl of undiluted **standard** (see Preparation of Standard 9.4, concentration of standard 1, S1 = 100.0 ng/ml) in duplicate into well A1 and A2 (see Table 1). Transfer 100 µl to wells B1 and B2. Mix the contents of wells B1 and B2 by repeated aspiration and ejection (concentration of standard 2, S2 = 50.0 ng/ml), and transfer 100 µl to wells C1 and C2, respectively (see Figure 6). Take care not to scratch the inner surface of the microwells. Continue this procedure 2 times, creating two rows of human sICAM-1 standard dilutions ranging from 100.0 to 6.3 ng/ml. Discard 100 µl of the contents from the last microwells (E1, E2) used.

Figure 6



In case of an **external standard dilution** (see 9.4.1), pipette 100 µl of these standard dilutions (S1 – S5) in the standard wells according to Table 1.

Table 1

Table depicting an example of the arrangement of blanks, standards and samples in the microwell strips:

	1	2	3	4
A	Standard 1 (100.0 ng/ml)	Standard 1 (100.0 ng/ml)	Sample 3	Sample 3
B	Standard 2 (50.0 ng/ml)	Standard 2 (50.0 ng/ml)	Sample 4	Sample 4
C	Standard 3 (25.0 ng/ml)	Standard 3 (25.0 ng/ml)	Sample 5	Sample 5
D	Standard 4 (12.5 ng/ml)	Standard 4 (12.5 ng/ml)	Sample 6	Sample 6
E	Standard 5 (6.3 ng/ml)	Standard 5 (6.3 ng/ml)	Sample 7	Sample 7
F	Blank	Blank	Sample 8	Sample 8
G	Sample 1	Sample 1	Sample 9	Sample 9
H	Sample 2	Sample 2	Sample 10	Sample 10

4. Add 100 µl of Sample Diluent in duplicate to the blank wells.
5. Add 90 µl of Sample Diluent to the sample wells.
6. Add 10 µl of each sample in duplicate to the sample wells.
7. Prepare HRP-Conjugate (see Preparation of HRP-Conjugate 9.3).
8. Add 50 µl of HRP-Conjugate to all wells.
9. Cover with an adhesive film and incubate at room temperature (18 to 25°C) for 1 hour, if available on a microplate shaker set at 400 rpm.
10. Remove adhesive film and empty wells. Wash microwell strips 3 times according to point b. of the test protocol. Proceed immediately to the next step.
11. Pipette 100 µl of TMB Substrate Solution to all wells.
12. Incubate the microwell strips at room temperature (18° to 25°C) for about 10 min. Avoid direct exposure to intense light.

The development on the plate should be monitored and the substrate reaction stopped (see next point of this protocol) before positive wells are no longer properly recordable. Determination of the ideal time period for color development has to 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. Stop the enzyme reaction by quickly pipetting 100 µl of Stop Solution into each well. It is important that the Stop Solution is spread quickly and uniformly throughout the microwells to completely inactivate the enzyme. Results must be read immediately after the Stop Solution is added or within one hour if the microwell strips are stored at 2 - 8°C in the dark.
14. Read absorbance of each microwell on a spectro-photometer 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.

11. Calculation of Results

- Calculate the average absorbance values for each set of duplicate standards and samples. Duplicates should be within 20 per cent of the mean value.
- Create a standard curve by plotting the mean absorbance for each standard concentration on the ordinate against the human sICAM-1 concentration on the abscissa. Draw a best fit curve through the points of the graph (a 5-parameter curve fit is recommended).
- To determine the concentration of circulating human sICAM-1 for each sample, first find the mean absorbance value on the ordinate and extend a horizontal line to the standard curve. At the point of intersection, extend a vertical line to the abscissa and read the corresponding human sICAM-1 concentration.
- **If instructions in this protocol have been followed, samples have been diluted 1:10 (10 µl sample + 90 µl Sample Diluent) and the concentration read from the standard curve must be multiplied by the dilution factor (x 10).**
- **Calculation of samples with a concentration exceeding standard 1 will result in incorrect, low human sICAM-1 levels (Hook Effect). Such samples require further external predilution according to expected human sICAM-1 values with Sample Diluent in order to precisely quantitate the actual humans ICAM-1 level.**
- It is suggested that each testing facility establishes a control sample of known human sICAM-1 concentration and runs this additional control with each assay. If the values obtained are not within the expected range of the control, the assay results may be invalid.
- A representative standard curve is shown in Figure 7. This curve cannot be used to derive test results. Each laboratory must prepare a standard curve for each group of microwell strips assayed.

Figure 7

Representative standard curve for human sICAM-1 ELISA. Human sICAM-1 was diluted in serial 2-fold steps in Sample Diluent. Do not use this standard curve to derive test results. A standard curve must be run for each group of microwell strips assayed.

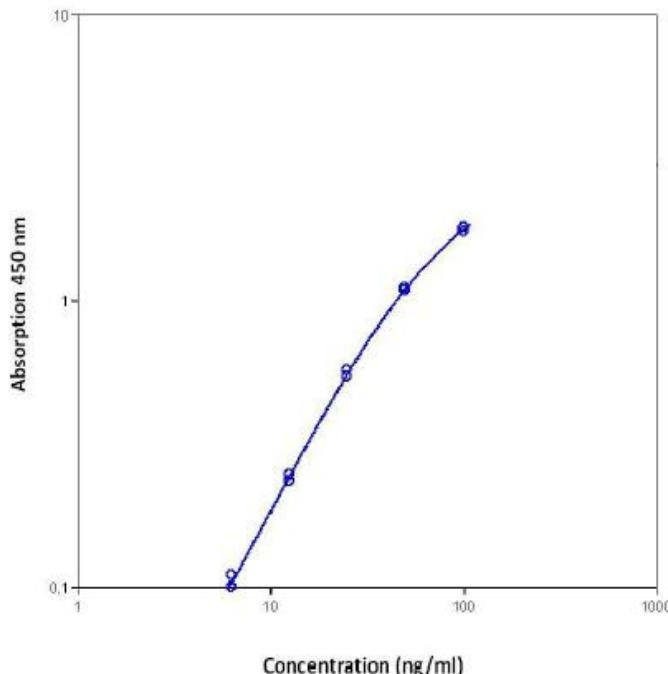


Table 2

Typical data using the human sICAM-1 ELISA

Measuring wavelength: 450 nm

Reference wavelength: 620 nm

Standard	Human sICAM-1 Concentration (ng/ml)	O.D. at 450 nm	Mean O.D. at 450 nm	C.V. (%)	
1	100.0	1.737	1.772	2.8	
		1.806			
2	50.0	1.114	1.101	1.7	
		1.087			
3	25.0	0.570	0.558	3.0	
		0.546			
4	12.5	0.247	0.241	3.5	
		0.235			
5	6.3	0.110	0.105	6.7	
		0.100			
Blank		0	0.005	20.0	
		0.006			

The OD values of the standard curve may vary according to the conditions of assay performance (e.g. operator, pipetting technique, washing technique or temperature effects). Furthermore shelf life of the kit may affect enzymatic activity and thus color intensity. Values measured are still valid.

12. Limitations

- Since exact conditions may vary from assay to assay, a standard curve must be established for every run.
- Bacterial or fungal contamination of either screen samples or reagents or cross-contamination between reagents may cause erroneous results.
- 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.
- The use of radioimmunotherapy has significantly increased the number of patients with human anti-mouse IgG antibodies (HAMA). HAMA may interfere with assays utilizing murine monoclonal antibodies leading to both false positive and false negative results. Serum samples containing antibodies to murine immunoglobulins can still be analysed in such assays when murine immunoglobulins (serum, ascitic fluid, or monoclonal antibodies of irrelevant specificity) are added to the sample.

13. Performance Characteristics

13.1. Sensitivity

The limit of detection of human sICAM-1 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 2.2 ng/ml (mean of 6 independent assays).

13.2. Reproducibility

13.2.1. Intra-assay

Reproducibility within the assay was evaluated in 2 independent experiments. Each assay was carried out with 6 replicates of 8 serum samples containing different concentrations of human sICAM-1. 2 standard curves were run on each plate. Data below show the mean human sICAM-1 concentration and the coefficient of variation for each sample (see Table 3). The calculated overall intra-assay coefficient of variation was 4.1%.

Table 3

The mean human sICAM-1 concentration and the coefficient of variation for each sample

Sample	Experiment	Mean Human sICAM-1 Concentration (ng/ml)	Coefficient of Variation (%)
1	1	330.1	7.8
	2	385.3	2.4
2	1	169.4	4.1
	2	187.3	3.4
3	1	518.9	7.2
	2	560.1	6.8
4	1	939.7	7.1
	2	1054.0	2.2
5	1	378.2	2.5
	2	404.1	4.2
6	1	318.3	2.1
	2	349.4	2.3
7	1	213.5	2.7
	2	231.3	3.3
8	1	149.6	3.4
	2	164.8	4.4

13.2.2. Inter-assay

Assay to assay reproducibility within one laboratory was evaluated in 2 independent experiments. Each assay was carried out with 6 replicates of 8 serum samples containing different concentrations of human sICAM-1. 2 standard curves were run on each plate. Data below show the mean human sICAM-1 concentration and the coefficient of variation calculated on 18 determinations of each sample (see Table 4). The calculated overall inter-assay coefficient of variation was 7.7%.

Table 4

The mean human sICAM-1 concentration and the coefficient of variation of each sample

Sample	Mean Human sICAM-1 Concentration (ng/ml)	Coefficient of Variation (%)
1	376.3	11.1
2	184.6	7.2
3	529.1	5.1
4	934.7	8.2
5	363.2	9.6
6	323.4	7.5
7	217.7	5.8
8	154.2	6.8

13.3. Spike Recovery

The spike recovery was evaluated by spiking 4 levels of human sICAM-1 into Sample Diluent (serum matrix). Recoveries were determined in 3 independent experiments with 4 replicates each.

The recovery ranged from 82% to 109% with an overall mean recovery of 99%.

13.4. Dilution Parallelism

Four serum samples with different levels of human sICAM-1 were analysed at serial two-fold dilutions with four replicates each.

The recovery ranged from 80% to 102% with an overall recovery of 93% (see Table 5).

Table 5

Sample	Dilution	Expected Human sICAM-1 Concentration (ng/ml)	Observed Human sICAM-1 Concentration (ng/ml)	Recovery of Expected Human sICAM-1 Concentration (%)
1	1:10	--	368	--
	1:20	184	148	80
	1:40	92	82	89
	1:80	46	46	100
2	1:10	--	163	--
	1:20	82	76	93
	1:40	41	40	98
	1:80	21	17	85
3	1:10	--	530	--
	1:20	265	238	90
	1:40	133	125	94
	1:80	66	67	102
4	1:10	--	602	--
	1:20	301	296	98
	1:40	151	135	89
	1:80	75	74	99

13.5. Sample Stability

13.5.1. Freeze-Thaw Stability

Aliquots of serum samples (spiked or unspiked) were stored at -20°C and thawed 5 times, and the human sICAM-1 levels determined. There was no significant loss of human sICAM-1 immunoreactivity detected by freezing and thawing.

13.5.2. Storage Stability

Aliquots of serum samples (spiked or unspiked) were stored at -20°C, 2- 8°C, room temperature (RT) and at 37°C, and the human sICAM-1 level determined after 24, 48 and 96 h. There was no significant loss of human sICAM-1 immunoreactivity detected during storage under above conditions.

13.6. Comparison of Serum and Plasma

From 3 individuals, serum as well as EDTA, citrate and heparin plasma obtained at the same time point were evaluated. Human sICAM-1 levels were not significantly different and therefore all these blood preparations are suitable for human sICAM-1 determinations.

13.7. Specificity

The interference of circulating factors of the immune system was evaluated by spiking these proteins at physiologically relevant concentrations into a human sICAM-1 positive serum.

There was no cross reactivity detected with TNF-R (60 kDa), TNF-R (80 kDa), IL-8/NAP-1, TNF alpha, TNF beta, IFN gamma, IFN alpha 2C, IFN omega, IL-6, IL-2R, ELAM-1 and L-selectin.

13.8. Expected Values

Panels of 40 serum as well as EDTA, citrate and heparin plasma samples from randomly selected apparently healthy donors (males and females) were tested for human sICAM-1.

For detected human sICAM-1 levels see Table 6.

Table 6

Sample Matrix	Number of Samples Evaluated	Range (ng/ml)	Mean (ng/ml)	Standard Deviation (ng/ml)
Serum	40	302 - 1115	504	171
Plasma (EDTA)	40	249 - 966	543	168
Plasma (Citrate)	40	59 - 529	305	103
Plasma (Heparin)	40	133 - 623	385	122

14. Reagent Preparation Summary

14.1. Wash Buffer (1x)

Add **Wash Buffer Concentrate 20x** (50 ml) to 950 ml distilled water.

Number of Strips	Wash Buffer Concentrate (ml)	Distilled Water (ml)
1 - 6	25	475
1 - 12	50	950

14.2. Assay Buffer (1x)

Add **Assay Buffer Concentrate 20x** (5 ml) to 95 ml distilled water.

Number of Strips	Assay Buffer Concentrate (ml)	Distilled Water (ml)
1 - 6	2.5	47.5
1 - 12	5.0	95.0

14.3. HRP-Conjugate

Make a 1:100 dilution of **HRP-Conjugate** in Assay Buffer (1x):

Number of Strips	HRP-Conjugate (ml)	Assay Buffer (1x) (ml)
1 - 6	0.03	2.97
1 - 12	0.06	5.94

14.4. Controls

Add 100 µl distilled water to lyophilized **controls**.

15. Test Protocol Summary

1. Determine the number of microwell strips required.
2. Wash microwell strips twice with Wash Buffer.
2. Standard dilution on the microwell plate: Add 100 µl Sample Diluent, in duplicate, to all standard wells leaving the first wells empty. Pipette 200 µl undiluted standard into the first wells and create standard dilutions by transferring 100 µl from well to well. Discard 100 µl from the last wells.
Alternatively external standard dilution in tubes (see 9.4.1): Pipette 100 µl of these standard dilutions in the microwell strips.
4. Add 100 µl Sample Diluent, in duplicate, to the blank wells.
5. Add 90 µl Sample Diluent to sample wells.
6. Add 10 µl sample in duplicate, to designated sample wells.
7. Prepare HRP-Conjugate.
8. Add 50 µl HRP-Conjugate to all wells.
9. Cover microwell strips and incubate 1 hour at room temperature (18° to 25°C).
10. Empty and wash microwell strips 3 times with Wash Buffer.
11. Add 100 µl of TMB Substrate Solution to all wells.
12. Incubate the microwell strips for about 10 minutes at room temperature (18° to 25°C).
13. Add 100 µl Stop Solution to all wells.
14. Blank microwell reader and measure color intensity at 450 nm.

Note: If instructions in this protocol have been followed, samples have been diluted 1:10 (10 µl sample + 90 µl Sample Diluent) and the concentration read from the standard curve must be multiplied by the dilution factor (x 10).

PRODUKTINFORMATION UND HANDBUCH (Deutsch)

1. Mitgelieferte Reagenzien

1.1. Mitgelieferte Reagenzien für human sICAM-1 ELISA BMS201CE (96 Tests)

- 1 Aluminiumbeutel mit **Mikrotiterplatte, beschichtet** mit Antikörper (monoklonal) gegen human sICAM-1
- 1 Fläschchen (100 µl) **HRP-Konjugat**, monoklonaler anti-human sICAM-1 Antikörper
- 2 Fläschchen (500 µl) human sICAM-1-**Standard**, 100 ng/ml
- 1 Fläschchen lyophilisierte **Kontrolle, hoch konzentriert**
- 1 Fläschchen lyophilisierte **Kontrolle, niedrig konzentriert**
- 1 Fläschchen (12 ml) **Verdünnungslösung**
- 1 Fläschchen (5 ml) **Probenpufferkonzentrat** 20x (PBS mit 1% Tween 20, 10% BSA)
- 1 Flasche (50 ml) **Waschpufferkonzentrat** 20x (PBS mit 1% Tween 20)
- 1 Fläschchen (15 ml) **Substratlösung** (Tetramethylbenzidin)
- 1 Fläschchen (15 ml) **Stopplösung** (1 M Phosphorsäure)
- 2 Klebefolien

1.2. Mitgelieferte Reagenzien für human sICAM-1 ELISA BMS201TENCE (10x96 Tests)

- 10 Aluminiumbeutel mit **Mikrotiterplatte, beschichtet** mit Antikörper (monoklonal) gegen human sICAM-1
- 10 Fläschchen (100 µl) **HRP-Konjugat**, monoklonaler anti-human sICAM-1 Antikörper
- 10 Fläschchen (500 µl) human sICAM-1-**Standard**, 100 ng/ml 10 Fläschchen lyophilisierte **Kontrolle, hoch konzentriert**
- 10 Fläschchen lyophilisierte **Kontrolle, niedrig konzentriert**
- 10 Fläschchen (12 ml) **Verdünnungslösung**
- 1 Fläschchen (5 ml) **Probenpufferkonzentrat** 20x (PBS mit 1% Tween 20, 10% BSA)
- 3 Flaschen (50 ml) **Waschpufferkonzentrat** 20x (PBS mit 1% Tween 20)
- 10 Fläschchen (15 ml) **Substratlösung** (Tetramethylbenzidin)
- 1 Fläschchen (100 ml) **Stopplösung** (1 M Phosphorsäure)
- 10 Klebefolien

2. Lagerhinweise

Lagern Sie den Inhalt des Kits mit Ausnahme der Kontrollen bei 2°-8°C. Lagerung der lyophilisierten Kontrollen bei -20°C. Verbliebene Reagenzien nach Verwendung sofort wieder auf 2°-8°C, bzw. auf -20°C kühlen.

Das Ablaufdatum des Kits und der Reagenzien ist auf den Etiketten angegeben.

Die Haltbarkeit des Kits und der Komponenten kann nur bei fachgerechter Lagerung garantiert werden, sowie bei mehrfacher

Verwendung nur dann, wenn die Reagenzien bei der ersten

Verwendung nicht kontaminiert wurden.

3. Sicherheitsvorkehrungen für den Gebrauch

- Alle enthaltenen Reagenzien sollten als potenziell gefährlich betrachtet werden. Daher wird empfohlen, dass dieses Produkt nur von Personen mit labortechnischer Erfahrung und in Übereinstimmung mit GLP Richtlinien verwendet wird. Passende Schutzbekleidung, wie Labormäntel, Sicherheitsbrillen und Laborhandschuhe müssen getragen werden. Vermeiden Sie jeden Kontakt der Reagenzien mit Haut oder Augen. Im Falle des Kontaktes von Reagenzien mit Haut oder Augen, sofort mit Wasser spülen. Bitte entnehmen Sie weitere spezifische Hinweise den Sicherheitsdatenblättern und/oder den Sicherheitsbestimmungen.
- Die Reagenzien sind ausschließlich für Diagnosezwecke bestimmt und nicht für den Einsatz bei Therapien.
- Reagenzien aus verschiedenen Chargen oder anderer Herkunft nicht mischen oder untereinander austauschen.
- Verwenden Sie die Kitreagenzien nicht nach dem Ablaufdatum (siehe Etikett).
- Setzen Sie die Kitreagenzien während der Lagerung oder Inkubation keiner starken Lichteinstrahlung aus.
- Nicht mit dem Mund pipettieren.
- In Bereichen, in denen mit Kitreagenzien oder Proben hantiert wird, nicht essen, trinken oder rauchen.
- Vermeiden Sie den Kontakt der Haut/Schleimhäute mit Kitreagenzien/Proben.
- Tragen Sie während des Hantierens mit Kitreagenzien oder Proben geeignete Gummi- oder Einweghandschuhe.
- Vermeiden Sie den Kontakt zwischen Substratlösung und Oxidationsmitteln/Metallen.
- Vermeiden Sie Verspritzen von Flüssigkeit oder Bildung von Aerosolen.
- Zur Vermeidung von Kontamination mit Mikroben oder Kreuzkontamination der Reagenzien oder Proben, die den Test ungültig machen könnten, verwenden Sie Einwegpipettenspitzen und/oder Einwegpipetten.
- Verwenden Sie saubere, geeignete Reagenzgefäße für das Dispensieren von Konjugat und Substratreagenzien.
- Vermeiden Sie Kontakt mit Säuren, da dadurch Konjugate inaktiviert werden.
- Für die Reagensherstellung muss destilliertes oder entionisiertes Wasser verwendet werden.
- Die Substratlösung muss vor der Verwendung auf Raumtemperatur gebracht werden.
- Dekontaminieren und entsorgen Sie Proben sowie alle möglicherweise kontaminierten Materialien so, als ob sie Infektionserreger enthalten könnten. Die bevorzugte Dekontaminationsmethode ist Autoklavieren für mind. eine Stunde bei 121.5°C.
- Flüssige Abfälle, die kein Säure enthalten, sowie neutralisierte Abfälle werden zur Dekontamination mit Natrium Hypochlorit versetzt (Endkonzentration von Natrium Hypochlorit 1.0%). Nach 30 min ist eine effektive Dekontamination erreicht. Flüssige Abfälle, die Säure enthalten, müssen vor der Dekontamination neutralisiert werden.

4. Vorbereitung der Reagenzien

Bringen Sie die **Pufferkonzentrate** auf Raumtemperatur und stellen Sie die Verdünnung vor Beginn des Tests her. Sollten sich in den **Pufferkonzentraten** Kristalle gebildet haben, erwärmen Sie diese vorsichtig bis zur vollständigen Auflösung der Kristalle.

4.1. Waschpuffer (1x)

Leeren Sie den gesamten Inhalt (50 ml) des **Waschpufferkonzentrats** (20x) in einen sauberen 1000-ml-Messzylinder. Füllen Sie mit destilliertem oder entionisiertem Wasser auf, bis ein Endvolumen von 1000 ml erreicht ist. Mischen Sie vorsichtig um Schäumen zu vermeiden.

Füllen Sie in eine saubere Waschflasche um und lagern Sie den Waschpuffer (1x) bei 2° bis 25°C lagern. Bitte beachten Sie, dass dieser 30 Tage haltbar ist.

Der benötigte Waschpuffer (1x) kann auch entsprechend der untenstehenden Tabelle hergestellt werden:

Anzahl der Streifen	Waschpufferkonzentrat (20x) (ml)	Destilliertes Wasser (ml)
1 - 6	25	475
1 - 12	50	950

4.2. Probenpuffer (1x)

Leeren Sie den gesamten Inhalt (5 ml) des **Probenpufferkonzentrates** (20x) in einen sauberen 100-ml-Messzylinder. Füllen Sie mit destilliertem oder entionisiertem Wasser auf, bis ein Endvolumen von 100 ml erreicht ist. Mischen Sie vorsichtig um Schäumen zu vermeiden.

Probenpuffer (1x) bei 2° bis 8°C lagern. Bitte beachten Sie, dass der Probenpuffer (1x) 30 Tage haltbar ist.

Der benötigte Probenpuffer (1x) kann auch entsprechend der untenstehenden Tabelle hergestellt werden:

Anzahl der Streifen	Probenpufferkonzentrat (20x) (ml)	Destilliertes Wasser (ml)
1 - 6	2.5	47.5
1 - 12	5.0	95.0

4.3. HRP-Konjugat

Bitte beachten Sie, dass die HRP-Konjugatlösung nach der Verdünnung nur 30 Minuten haltbar ist.

Stellen Sie eine 1:100 Verdünnung der konzentrierten HRP- Konjugatlösung in Probenpuffer (1x) in einem sauberen Gefäß entsprechend der untenstehenden Tabelle her.

Anzahl der Streifen	HRP-Konjugat (ml)	Probenpuffer (1x) (ml)
1 - 6	0.03	2.97
1 - 12	0.06	5.94

4.4. Human sICAM-1-Standard

Die **Standardverdünnungen** können direkt auf den Mikrotiterplatten (siehe 5.c) oder in Reaktionsgefäß en (siehe 4.4.1) hergestellt werden.

4.4.1. Externe Standardverdünnung

Beschriften Sie 4 Gefäße, jedes für einen Standardpunkt:

S2, S3, S4, S5

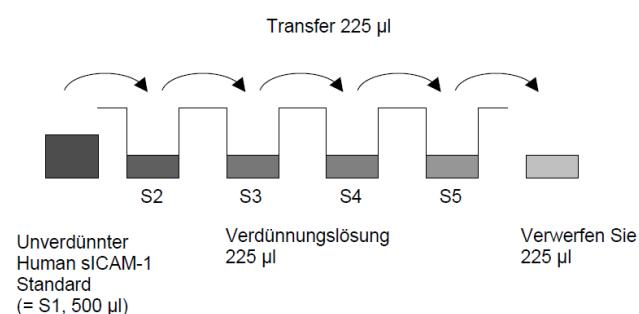
Danach stellen Sie eine 1:2 Verdünnungsreihe für die Standardkurve her: Pipettieren Sie in die Gefäße S2 – S5 225 µl der Verdünnungslösung.

Pipettieren Sie 225 µl des unverdünnten Standards (ist Standard mit der höchsten Konzentration S1 = 100 ng/ml) in das erste Gefäß mit der Beschriftung S2 und mischen Sie (Konzentration des Standard 2 = 50 ng/ml).

Wiederholen Sie diese Verdünnungsschritte 3x. Die so hergestellte Verdünnungsreihe dient zur Erstellung der Standardkurve (siehe Abbildung 1).

Verdünnungslösung dient als Blindwert.

Abbildung 1



4.5. Kontrollen

Lösen Sie die **Kontrollen** durch Zugabe von 100 µl destilliertem Wasser auf. Für die Kontrollen 10-30 min Rekonstitutionszeit einhalten. Mixen oder schütteln Sie die Fläschchen vorsichtig um eine vollständige Lösung zu erreichen. Verfahren Sie in der Folge mit den Kontrollen analog zu den Proben. Der Konzentrationsbereich der Kontrollen ist am Analysenzertifikat oder am Flaschenetikett angegeben.

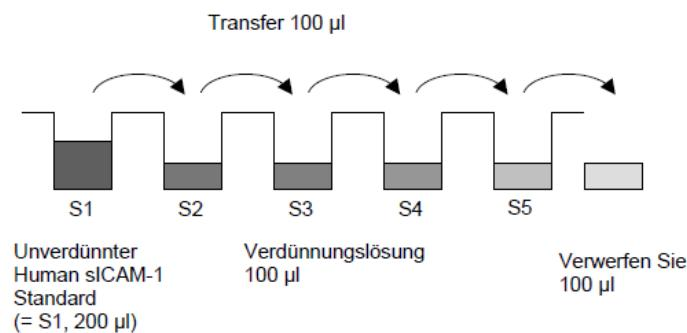
Lagern sie die rekonstituierten Kontrollen aliquotiert bei -20°C. Vermeiden Sie wiederholtes Frieren und Tauen.

5. Testprotokoll

- 1 Bestimmen Sie die Anzahl der Mikrowellstreifen die für das Testen der gewünschten Anzahl von Proben benötigt werden, sowie die Mikrowellstreifen für Blindwert und Standards. Probe, Standard, Blindwert und optionale Kontrollproben immer jeweils doppelt testen. Entfernen Sie die zusätzlichen Mikrowellstreifen von der Halterung und bewahren Sie diese mit dem mitgelieferten Trockenmittel in dem Folienbeutel fest verschlossen bei 2°-8°C auf.
- 2 Waschen Sie die Mikrowellstreifen 2 mal mit ca. 400 µl **Waschpuffer** pro Vertiefung; zwischen den Waschgängen den Inhalt der Vertiefungen gründlich absaugen. Vor dem Absaugen Waschpuffer **10-15 Sekunden** einwirken lassen. Achten Sie darauf, die Oberfläche der Vertiefungen nicht zu zerkratzen.
Leeren Sie die Vertiefungen nach dem letzten Waschschritt und klopfen Sie die Mikrowellstreifen auf einem Saug- oder Papiertuch aus um überschüssigen Waschpuffer zu entfernen. Verwenden Sie die Mikrowellstreifen sofort nach dem Waschen, oder legen Sie diese für maximal 15 min umgedreht auf ein nasses Saugtuch. **Lassen Sie die Vertiefungen nicht austrocknen.**
- 3 **Standardverdünnung auf der Mikrotiterplatte** (Wahlweise können die Standardverdünnungen auch in Reaktionsgefäßern hergestellt werden – siehe 4.4.1)

Pipettieren Sie 100 µl Verdünnungslösung in die **Standardvertiefungen B1/2-E1/2**, wobei A1/2 leer gelassen wird. Pipettieren Sie 200 µl des unverdünnten **Standards** (siehe Herstellung des Standards 4.4.1, Konzentration des Standards, S1 = 100.0 ng/ml) in die Vertiefungen A1 und A2 (Doppelbestimmung, siehe Tabelle 1). Transferieren Sie 100 µl in die Vertiefungen B1 und B2. Mischen Sie den Inhalt der Vertiefungen B1 und B2 durch wiederholtes Aufsaugen und Zugeben (Konzentration des Standards S2 = 50.0 ng/ml) und transferieren Sie 100 µl in die Probenvertiefungen C1 und C2 (siehe Abbildung 2). Achten Sie darauf, die Oberfläche der Vertiefungen nicht zu zerkratzen. Wiederholen Sie diese Verdünnungsschritte 2x, wodurch zwei human sICAM-1 Verdünnungsreihen mit den Konzentrationen von 100.0 bis 6.3 ng/ml hergestellt werden. Verwerfen Sie 100 µl aus den letzten Standardvertiefungen (E1/2). Die so hergestellten Verdünnungsreihen dienen zur Erstellung der Standardkurve.

Abbildung 2



Falls sie eine **externe Standardverdünnungsreihe** erstellen (siehe 4.4.1), pipettieren Sie 100 µl der Standardverdünnungen (S1 –S5) in die Standardvertiefungen (entsprechend Tabelle 1).

Tabelle 1

Diagramm mit Beispiel für die Anordnung von Blindwert, Standards und Proben in den Mikrowellstreifen:

	1	2	3	4
A	Standard 1 (100.0 ng/ml)	Standard 1 (100.0 ng/ml)	Probe 3	Probe 3
B	Standard 2 (50.0 ng/ml)	Standard 2 (50.0 ng/ml)	Probe 4	Probe 4
C	Standard 3 (25.0 ng/ml)	Standard 3 (25.0 ng/ml)	Probe 5	Probe 5
D	Standard 4 (12.5 ng/ml)	Standard 4 (12.5 ng/ml)	Probe 6	Probe 6
E	Standard 5 (6.3 ng/ml)	Standard 5 (6.3 ng/ml)	Probe 7	Probe 7
F	Blindwert	Blindwert	Probe 8	Probe 8
G	Probe 1	Probe 1	Probe 9	Probe 9
H	Probe 2	Probe 2	Probe 10	Probe 10

- 4 Pipettieren Sie in alle Blindwertvertiefungen (Doppelbestimmung), 100 µl Verdünnungslösung.
- 5 Pipettieren Sie in alle Probenvertiefungen 90 µl Verdünnungslösung.
- 6 Pipettieren Sie je 10 µl von jeder Probe (Doppelbestimmung) in die Probenvertiefungen und mischen Sie den Inhalt durch.
- 7 Stellen Sie das HRP-Konjugat (siehe: Vorbereitung der Reagenzien HRP-Konjugat 4.3) her.
- 8 Pipettieren Sie in alle Vertiefungen, einschließlich der Blindwertvertiefungen 50 µl HRP-Konjugat.
- 9 Mit einer Klebefolie abdecken und bei Raumtemperatur (18° bis 25°C) für 1 Stunde inkubieren, wenn möglich auf einem Schüttler bei 400 U/min.
- 10 Entfernen Sie die Klebefolie und entleeren Sie die Vertiefungen. **Waschen** Sie die Mikrowellstreifen 3 mal wie in Punkt b des Testprotokolls beschrieben. Verwenden Sie die Mikrowellstreifen sofort nach dem Waschen.
- 11 Pipettieren Sie in alle Vertiefungen, einschließlich der Blindwertvertiefungen, 100 µl TMB-Substratlösung.
- 12 Inkubieren Sie die Mikrowellstreifen bei Raumtemperatur (18° bis 25°C) für ca. 10 Minuten. Vermeiden Sie direkte, starke Lichteinstrahlung.

Die Farbentwicklung innerhalb der einzelnen Vertiefungen muss beobachtet und die Substratreaktion gestoppt werden (siehe nächster Protokollpunkt), bevor die gefärbten Vertiefungen nicht mehr richtig gemessen werden können. Die optimale Inkubationszeit für die Farbentwicklung muß bei jedem Versuch neu bestimmt werden.

Es wird empfohlen, die Stopplösung zuzugeben, wenn der höchste Standardpunkt eine dunkelblaue Farbe angenommen hat. Alternativ kann die Farbentwicklung auch mit einem Photometer bei 620 nm verfolgt werden. Die Substratreaktion sollte gestoppt werden, wenn der höchste Standardpunkt eine OD von 0.9 -0.95 erreicht.

- 13 Stoppen Sie die Enzymreaktion durch rasche Zugabe von 100 µl Stopplösung in jede Vertiefung, einschließlich der Blindwertvertiefungen. Für eine vollständige Inaktivierung der Enzyme ist es wichtig, die Stopplösung rasch und gleichmäßig in den Vertiefungen zu verteilen. Die OD Werte müssen sofort nach Beigabe der Stopplösung oder innerhalb einer Stunde nach Lagerung der Mikrowellstreifen in Dunkelheit bei 2-8°C gemessen werden.
- 14 Messen Sie die Absorption jeder Vertiefung mit einem Spektrophotometer. Verwenden Sie dabei 450 nm als primäre Wellenlänge (optional 620 nm als Referenzwellenlänge; 610 nm bis 650 nm sind möglich). Stellen Sie das Plattenmessgerät nach Anleitung des Herstellers und unter Verwendung der Blindwertvertiefungen auf den Leerwert ein. Bestimmen Sie die Absorption der Proben wie auch der human sICAM-1-Standards.

Die Proben wurden im Zuge der Testdurchführung 1:10 verdünnt. Daher muß der aus der Standardkurve berechnete Wert mit dem Verdünnungsfaktor multipliziert werden (x 10).

Anmerkung: Falls die Platte während der Inkubation nicht geschüttelt wurde, können die erreichten OD Werte niedriger als die unten angeführten sein. Die Ergebnisse sind trotzdem gültig.

INFORMACIÓN Y MANUAL DEL PRODUCTO (Espanol)

1. Reactivos Suministrados

1.1. Reactivos Suministrados para human sICAM-1 ELISA BMS201CE (96 tests)

- 1 bolsa de aluminio con **una placa de micropocillos recubiertos** con anticuerpos monoclonales anti-human sICAM-1
- 1 vial (100 µl) con **conjugado de HRP** (anticuerpos monoclonales anti- human sICAM-1)
- 2 viales (500 µl) con **Estándar** human sICAM-1, 100 ng/ml
- 1 vial del **control alto**, liofilizado
- 1 vial del **control bajo**, liofilizado
- 1 vial (12 ml) de **diluyente de muestra**
- 1 vial (5 ml) de **concentrado de tampón de ensayo** 20x (PBS con Tween 20 al 1% y BSA al 10%)
- 1 frasco (50 ml) de **concentrado de tampón de lavado** 20x (PBS con Tween 20 al 1%)
- 1 vial (15 ml) de **solución de sustrato** (tetrametil-bencidina)
- 1 vial (15 ml) de **solución de parada** (ácido fosfórico 1M)
- 2 **tapas para placas**, adhesives

1.2. Reactivos Suministrados para human sICAM-1 ELISA BMS201TENCE (10x96 tests)

- 10 bolsas de aluminio con **una placa de micropocillos recubiertos** con anticuerpos monoclonales anti-human sICAM-1
- 10 viales (100 µl) con **conjugado de HRP** (anticuerpos monoclonales anti-human sICAM-1)
- 10 viales (500 µl) con **Estándar** human sICAM-1, 100 ng/ml 10 viales del **control alto**, liofilizado
- 10 viales del **control bajo**, liofilizado
- 10 viales (12 ml) de **diluyente de muestra**
- 1 vial (5 ml) de **concentrado de tampón de ensayo** 20x (PBS con Tween 20 al 1% y BSA al 10%)
- 3 frascos (50 ml) de **concentrado de tampón de lavado** 20x (PBS con Tween 20 al 1%)
- 10 viales (15 ml) de **solución de sustrato** (tetrametil-bencidina)
- 1 vial (100 ml) de **solución de parada** (ácido fosfórico 1M)
- 10 **tapas para placas**, adhesives

2. Instrucciones de Conservación

Conservar los reactivos del kit a una temperatura comprendida entre 2 y 8°C excepto los controles. Conservar los controles liofilizados a -20°C. Inmediatamente después de utilizarlos deberá volver a conservar los reactivos a dicha temperatura (2° to 8°C), controles a -20°C. En las etiquetas figuran las fechas de caducidad del kit y de los reactivos.

Sólo se podrá garantizar la fecha de caducidad de los componentes del kit si se conservan adecuadamente y, en caso de uso reiterado de un mismo componente, si el reactivo no queda contaminado en la primera manipulación.

3. Precauciones de uso

- Todos los productos químicos deben considerarse potencialmente peligrosos. Por tanto, recomendamos que este producto sea manipulado únicamente por aquellas personas que hayan sido entrenadas en técnicas de laboratorio que sea usado de acuerdo con los principios de buenas prácticas de laboratorio. Se debe llevar ropa de protección apropiada como puedan ser las batas de laboratorio, gafas de seguridad y guantes. Se debe trabajar con cuidado para evitar cualquier contacto con piel y ojos. En el caso de que tenga lugar un contacto con piel u ojos, proceder de forma inmediata a lavar la parte afectada con abundante agua. Véase la(s) hoja(s) de seguridad y/o declaraciones de seguridad para recomendaciones específicas.
- Los reactivos están destinados para un uso en diagnóstico in vitro y no se deben usar en procedimientos terapéuticos.
- No mezclar o sustituir los reactivos por los equivalentes de otros lotes u otras fuentes.
- No usar reactivos caducados.
- No exponer los reactivos del kit a una luz intensa durante su almacenamiento o incubación.
- No pipetear con la boca.
- No se recomienda comer o fumar en las zonas donde se manipulen muestras o reactivos.
- Evitar el contacto de los reactivos del kit o de las muestras con piel o mucosas.
- Se recomienda el uso de guantes desechables de goma o látex durante la manipulación de las muestras y reactivos.
- Evitar el contacto de la solución de sustrato con agentes oxidantes y metales.
- Evitar salpicaduras y la generación de aerosoles.
- Con el propósito de evitar una contaminación microbiológica o contaminaciones cruzadas de reactivos y muestras que puedan invalidar el test se recomienda el uso de pipetas y/o puntas de pipetas de un solo uso.
- Usar recipientes limpios y específicos de reactivos para la dispensación de reactivos de sustrato.
- La exposición a los ácidos inactiva el conjugado.
- Se debe usar agua destilada o desionizada en la preparación de los reactivos.
- La solución de sustrato debe de estar a temperatura ambiente antes de su uso.
- Descontaminar y disponer las muestras y todos los materiales potencialmente contaminados como si pudieran contener agentes infecciosos. El método preferente de descontaminación es un autoclavado durante un mínimo de 1 hora a 121.5°C.
- Los residuos líquidos que no contengan ácido y los residuos neutralizados pueden ser mezclados con hipoclorito sódico en volúmenes tales que la mezcla final contenga 1.0% de hipoclorito sódico. Dejar actuar durante 30 minutos para una efectiva descontaminación. Los residuos líquidos que contengan ácido deben ser neutralizados previamente a la adición de hipoclorito sódico.

4. Preparación de los Reactivos

Los **tampones concentrados** debe de alcanzar la temperatura ambiente y ser diluidos antes de iniciar el procedimiento del test. Si en el concentrado de **tampones concentrados** se han formado cristales, caliente suavemente hasta su completa disolución.

4.1. Tampón de Lavado (1x)

Vierta todo el contenido (50 ml) del **concentrado de tampón de lavado (20x)** en un matraz aforado de 1000 ml limpio. Enrase en matraz con agua destilada o desionizada. Mezcle suavemente para evitar la formación de espuma.

Transfiera la solución a un frasco de lavado limpio y consérvela a una temperatura entre 2°C y 25°C. El tampón de lavado permanece estable durante 30 días.

En función de la cantidad que vaya a necesitar, prepare el tampón de lavado de acuerdo a la siguiente tabla:

Número de tiras	Tampón de lavado (20x) (ml)	Agua destilada (ml)
1 - 6	25	475
1 - 12	50	950

4.2. Tampón de Ensayo (1x)

Vierta todo el contenido (5 ml) del **concentrado de tampón de ensayo (20x)** en un matraz aforado de 100 ml limpio. Enrase en matraz con agua destilada o desionizada. Mezcle suavemente para evitar la formación de espuma.

Conserve la solución a una temperatura de entre 2°C y 8°C. El tampón de trabajo permanece estable durante 30 días.

En función de la cantidad que vaya a necesitar, prepare el tampón de ensayo de acuerdo a la siguiente tabla:

Número de tiras	Tampón de Ensayo (20x) (ml)	Aqua destilada (ml)
1 - 6	2.5	47.5
1 - 12	5.0	95.0

4.3. Conjugado de HRP

Se utilizará el conjugado de HRP antes de transcurridos 30 minutos desde su dilución.

Justo antes de utilizar el conjugado de HRP, se debe diluirlo con Tampón de Ensayo (1x) en un tubo de ensayo de plástico limpio, en una proporción de 1:100.

En función de la cantidad que vaya a necesitar, prepare el conjugado de HRP de acuerdo a la siguiente tabla:

Número de tiras	Conjugado de biotina (ml)	Tampón de Ensayo (1x) (ml)
1 - 6	0.03	2.97
1 - 12	0.06	5.94

4.4. Dilución estándar human sICAM-1

Las diluciones estándar pueden ser preparadas directamente en la placa multipocillo (véase 5.b) o alternativamente en tubos (véase 4.4.1).

4.4.1. Dilución Estándar Externa

Rotular 4 tubos, uno para cada punto de la curva estándar.

S2, S3, S4, S5.

Acto seguido, preparar diluciones seriadas 1:2 para la curva estándar como se indica a continuación:

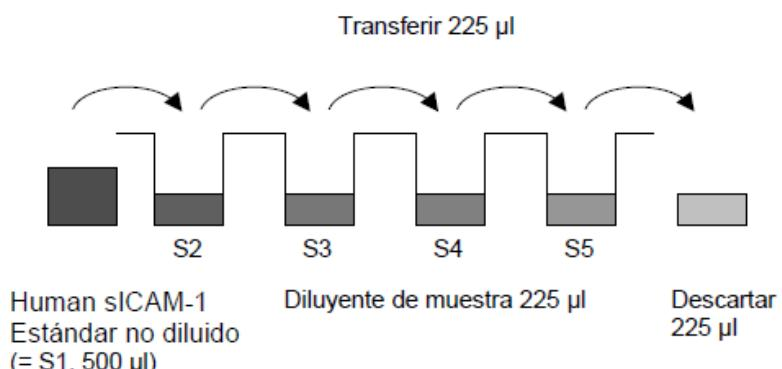
Pipetear 225 µl de Diluyente de muestra a tubos S2-S5.

Pipetear 225 µl de estándar rno diluido (sirve como el estándar más alto S1, concentración del estándar 1 = 100 ng/ml) en el primer tubo, etiquetado como S2, y mezclar (concentración del estándar 2 = 50 ng/ml).

Repetir la serie de diluciones 3 veces más de manera que se obtengan los diferentes puntos de la curva estándar (véase Figura 1).

Diluyente de muestra sirve como blanco.

Figura 1



4.5. Controles

Solubilizar añadiendo 100 µl de agua destilada al **controles** liofilizados. Permitir que el controles liofilizados se asiente durante 10-30 minutos. Mezclar bien previamente a realizar las diluciones. Vortear concienzudamente para asegurar una solubilización homogénea y completa. A partir de aquí, tratar el controles de la misma forma que las muestras. El rango del control viene indicado en el certificado de calidad y en la etiqueta del vial. Almacenar el controles reconstituidos y alicuotado a -20°C. Evitar ciclos repetidos de congelación y descongelación.

5. Protocolo de Ensayo

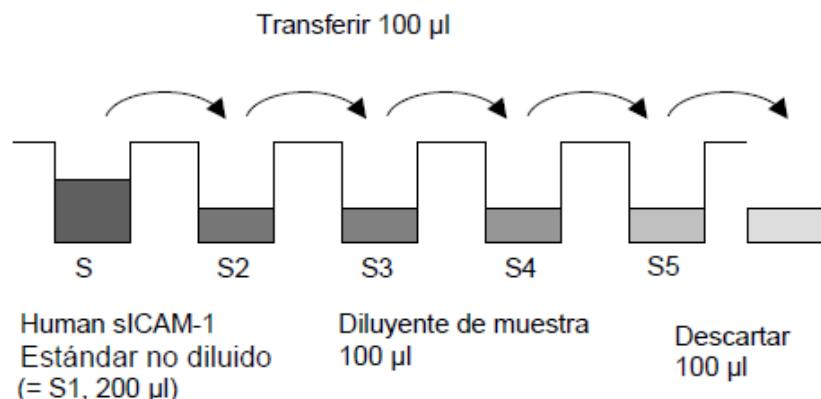
1. Determine el número de tiras necesarias para analizar el número deseado de muestras y además añada las tiras para blancos y patrones (de color). Todas las muestras, estándares, blancos y las posibles muestras de control deben ser analizadas por duplicado. Retire del soporte las tiras sobrantes y consérvelas, junto con el desecante suministrado en una bolsa metalizada y cerrada herméticamente, a una temperatura de 2°-8° C. Coloque las tiras que contienen la curva de valoración en las posiciones A1/A2 a H1/H2 (véase la Figura 2).
2. Lave 2 veces las tiras con aproximadamente 400 µl de **tampón de lavado** por cada pocillo, aspirando completamente el contenido de los pocillos entre cada lavado. Permitir que el tampón de lavado permanezca en los pocillos durante **10-15 segundos** antes de su aspiración. Evite rayar la superficie de los pocillos.

Tras el último lavado, golpee suavemente las tiras contra un papel absorbente o una toallita de papel para eliminar el exceso de tampón de lavado. Utilice las tiras inmediatamente después de lavadas o bien colóquelas boca abajo sobre un papel absorbente húmedo durante como máximo 15 minutos. **No deje secar los pocillos.**

3. **Dilución de los Estándares en la placa multipocillo** (Alternativamente, la dilución de los estándares puede ser preparada en tubos – véase 4.4.1)

Añadir 100 µl de a **pocillos estándar B1/2-E1/2**, dejando A1/A2 vacíos. Pipetear 200 µl de **estándar no diluido** (véase Preparación del Estándar 4.4, concentración del estándar 1, S1 = 100.0 ng/ml) por duplicado en los pocillos A1 y A2 (véase Tabla 1). Transferir 100 µl a los pocillos B1 y B2. Mezclar el contenido de los pocillos B1 y B2 por repetidas aspiraciones y expulsiones del contenido con la pipeta (concentración del estándar 2, S2 = 50.0 ng/ml), y transferir 100 µl a los pocillos C1 y C2, respectivamente (véase Figura 2). Levar cuidado de no rascar la superficie interior de los micropocillos con la punta de la pipeta. Continuar este procedimiento 2 veces, formando dos filas de diluciones estándar del human sICAM-1 ordenadas des de 100.0 a 6.3 ng/ml. Descartar 100 µl de los contenidos de los últimos micropocillos (E1, E2) usados.

Figura 2



En caso de **una dilución estándar externa** (véase 4.4.1), pipetear 100 µl de estas diluciones estándar (S1 – S5) en los pocillos correspondientes al estándar de acuerdo con la Tabla 1.

4. Añada 100 µl **Diluyente de muestra** a los pocillos del blanco, por duplicado.
5. Añada 90 µl de Diluyente de muestra a los pocillos con muestras.

6. Por duplicado, añada 10 µl de cada muestra a los pocillos designados.

Tabla 1

Tabla que describe un ejemplo de la disposición de los blancos, estándares y muestras en los micropocillos de las tiras:

	1	2	3	4
A	Estándar 1 (100.0 ng/ml)	Estándar 1 (100.0 ng/ml)	Muestra 3	Muestra 3
B	Estándar 2 (50.0 ng/ml)	Estándar 2 (50.0 ng/ml)	Muestra 4	Muestra 4
C	Estándar 3 (25.0 ng/ml)	Estándar 3 (25.0 ng/ml)	Muestra 5	Muestra 5
D	Estándar 4 (12.5 ng/ml)	Estándar 4 (12.5 ng/ml)	Muestra 6	Muestra 6
E	Estándar 5 (6.3 ng/ml)	Estándar 5 (6.3 ng/ml)	Muestra 7	Muestra 7
F	Blanco	Blanco	Muestra 8	Muestra 8
G	Muestra 1	Muestra 1	Muestra 9	Muestra 9
H	Muestra 2	Muestra 2	Muestra 10	Muestra 10

7. Prepare conjugado de HRP (véase la preparación de conjugado de HRP 4.3.)
8. Añada 50 µl **conjugado de HRP** a todos los pocillos.
9. Cubra la placa con una tapa e incúbela a temperatura ambiente (18°C - 25°C) durante 1 hora (en un agitador mecánico a 400 rpm, si es posible).
10. Retire la tapa y vacíe los pocillos. Lavar los micropocillos de las tiras 3 veces de acuerdo al punto b del protocolo del test. Proseguir inmediatamente después al próximo paso.
11. Pipetee 100 µl de solución de sustrato TMB y viértalos en todos los pocillos, incluidos los del blanco.
12. Incube las tiras a temperatura ambiente (18°C - 25°C) durante aproximadamente 10 minutos. Evite la exposición directa a la luz intensa.

Deben monitorizarse los valores DO de la placa para detener la reacción del sustrato (véase el siguiente punto de este protocolo) antes de que deje de ser posible registrar correctamente los pocillos positivos.

La determinación del tiempo adecuado para el desarrollo del color, debe realizarse de forma individual para cada ensayo.

Se recomienda añadir la solución de parada cuando el estándar más alto presente un color azul oscuro. Alternativamente el desarrollo de color puede ser monitorizado con un lector de placas de ELISA a 620 nm. La reacción del substrato debería ser parada cuando este estándard alcance una DO entre 0.9 y 0.9 5.

13. Detenga la reacción enzimática pipeteando rápidamente 100 µl de solución de parada en cada pocillo, incluidos los del blanco. Es importante dispensar la solución de parada de forma rápida y uniforme en todos los pocillos para inactivar totalmente la enzima. Los resultados deben leerse inmediatamente después de añadir la solución de parada o, como máximo, en el plazo de 1 hora si las tiras se conservan a una temperatura entre 2 - 8°C en un lugar oscuro.
14. Lea la absorbancia de cada pocillo en un espectrofotómetro utilizando 450 nm como longitud de onda principal (opcionalmente 620 nm como longitud de onda de referencia; los valores comprendidos entre 610 nm y 650 nm son aceptables). Utilizando los pocillos de blanco, haga el blanco del lector de placas de acuerdo con las instrucciones del fabricante. Determine la absorbancia de las muestras y de los human sICAM-1.

Las muestras han sido diluidas 1:10, por tanto la concentración leída a partir de la curva estándar debe ser multiplicada por el factor de dilución (x 10).

Nota: En caso de incubar sin agitar, los valores de D.O. pueden ser inferiores a los indicados más abajo. De todas formas los resultados siguen siendo válidos.

INFORMATIONS SUR LE PRODUIT ET MANUEL (Français)

1. Réactifs Fournis

1.1. Reactifs pour ELISA de human sICAM-1 BMS201CE (96 essais)

- 1 pochette en aluminium contenant **une plaque de microtitration** recouverte d'anticorps monoclonaux anti-human sICAM-1
- 1 flacon (70 µl) de **conjugué HRP** anti-human sICAM-1 (anticorps monoclonaux anti-human sICAM-1)
- 2 flacons (500 µl) d'**étalon** human sICAM-1, 100 ng/ml
- 1 flacon de **contrôle élevé**, lyophilisé
- 1 flacon de **contrôle bas**, lyophilisé
- 1 flacon (12 ml) **diluant d'échantillon**
- 1 flacon (5 ml) **tampon d'essai concentré** 20 x
(PBS avec Tween 20 1% et de la sérumalbumine bovine 10%)
- 1 flacon (50 ml) de **tampon de lavage concentré** 20 x (PBS avec du Tween 20 1 %)
- 1 flacon (15 ml) de **solution de substrat** (tétraméthyle-benzidine)
- 1 flacon (15 ml) de **solution d'arrêt** (acide phosphorique 1 M)
- 2 **couvercle-plaques** adhésifs

1.2. Reactifs pour ELISA de human sICAM-1 BMS201TENCE (10x96 essais)

- 10 pochettes en aluminium contenant **une plaque de microtitration** recouverte d'anticorps monoclonaux anti-human sICAM-1
- 10 flacons (70 µl) de **conjugué HRP** anti-human sICAM-1 (anticorps monoclonaux anti-human sICAM-1)
- 10 flacons (500 µl) d'**étalon** human sICAM-1, 100 ng/ml 10 flacons de **contrôle élevé**, lyophilisé
- 10 flacons de **contrôle bas**, lyophilisé 10 flacons (12 ml) **diluant d'échantillon**
- 1 flacon (5 ml) **tampon d'essai concentré** 20 x
(PBS avec Tween 20 1% et de la sérumalbumine bovine 10%)
- 3 flacons (50 ml) de **tampon de lavage concentré** 20 x (PBS avec du Tween 20 1 %)
- 10 flacons (15 ml) de **solution de substrat** (tétraméthyle-benzidine)
- 1 flacon (100 ml) de **solution d'arrêt** (acide phosphorique 1 M) 10 **couvercle-plaques** adhésifs

2. Instruction de Stockage

Conserver les réactifs du kit entre 2° et 8°C et les contrôles à -20°C. Immédiatement après l'utilisation, les réactifs doivent être rangés au frais (2° à 8°C), les contrôles à -20°C. La date de péremption du kit est spécifiée sur les étiquettes.

Le délai de péremption du kit ne peut être garanti que si les composants sont conservés correctement et si, en cas d'utilisation répétée d'un composant, le réactif n'a pas été contaminé lors d'une première utilisation.

3. Préventions de Sécurité pour l'Usage

- Tout réactifs doivent être considérés comme potentiellement dangereux. Pour cela il est recommandé que ce produit est utilisé que par des personnes ayant une qualification de laboratoire et qu'il soit utilisé à l'avenant au code GLP. Une tenue correspondante comme une blouse de travail, des lunettes protectrices et des gants de travail doivent-être portés. Evitez tous contacts de réactifs avec la peau ou les yeux. En cas de contact avec les yeux ou la peau rincez immédiatement avec de l'eau. Veuillez consulter tous conseils spécifiques dans les fiches de données de sécurité et/ou les règles de sécurité.
- Les réactifs sont réservés exclusivement au diagnostique et non pas au thérapeutique.
- Evitez de mélanger et d'échanger les réactifs de lots différents et de provenance différents.
- Evitez l'utilisation des réactifs perimés (voyez étiquette).
- N'exposez pas les réactifs à la lumière pendant le stockage ou l'incubation.
- Ne pas pipeter avec la bouche
- Ne pas manger, boire ou fumer dans les zones de manipulation de réactifs et d'échantillons.
- Evitez le contact de la peau et des muqueuses avec les réactifs.
- Pendant le travail avec les réactifs, utilisez des gants appropriés.
- Evitez le contact de substrats avec des métaux/oxydant.
- Evitez de gicler des liquides et la formation d'Aérosoles.
- A fin d'éviter des contaminations avec microbes ou contaminations de reactifs et d'échantillons qui pourraient rendre le test sans valeur, veuillez utiliser des pointes de pipettes jetables.
- Utilisez des tubes appropriés pour dispenser le conjugué et le substrat.
- Toute exposition aux acides inactive le conjugué.
- Pour la préparation des réactifs de l'eau déstilée ou déionisé doit être utilisée.
- La solution de substrat doit être rendue à température ambiante avant usage.
- Décontaminez et éliminez les échantillons et tous matériaux contaminés de manière comme si ils contenaient des germes de maladies infectieuses. La méthode préférée de décontamination est par l'autoclave pour au moins une heure à 121.5 °C.
- Traitez les déchets liquides non-acidiques tel que des déchets neutralisés par l'hypochlorite de sodium (concentration finale d'hypochlorite: 1.0%). Après 30 minutes la décontamination effective est atteinte. Les déchets liquides contenant de l'acide doivent être neutralisés avant la décontamination.

4. Préparation des Réactifs

Placer **les concentrés** de tampon à une température ambiante et diluer avant de commencer le test. Si des cristaux se sont formés dans les **concentrés de tampon**, chauffer doucement ces derniers jusqu'à fin de les dissoudre la dissolution des cristaux totale.

4.1. Tampon de Lavage (1x)

Verser tout le contenu (50 ml) du concentré de **tampon de lavage** (20x) dans un cylindre gradué propre de 1000 ml. Porter le volume final à 1000 ml avec de l'eau distillée ou déionisée dans un alambic en verre. Mélanger doucement pour éviter la formation de mousse.

Transférer tout dans une bouteille de lavage et conserver à une température comprise entre 2° et 25°C. Noter que le tampon de lavage reste stable pendant 30 jours.

Le tampon de lavage peut être préparé selon le tableau suivant:

Nombre de bandes	Tampon de lavage (20x) (ml)	Eau distillée (ml)
1 - 6	25	475
1 - 12	50	950

4.2. Tampon d'essai (1x)

Bien mélanger le contenu de la bouteille. Ajouter le contenu du **tampon d'essai concentré** (20x) (5.0 ml) aux 95 ml d'eau distillée ou déionisée et mélanger doucement pour éviter la formation de mousse. Stocker le tout entre 2° et 8°C. Noter que le tampon d'essai reste stable pendant 30 jours.

Le tampon d'essai peut être préparé selon le tableau suivant :

Nombre de bandes	Tampon d'essai (20x) (ml)	Eau distillée (ml)
1 - 6	2.5	47.5
1 - 12	5.0	95.0

4.3. Préparation du conjugué HRP

Noter que le conjugué HRP doit être utilisé dans les 30 minutes qui suivent la dilution.

Le conjugué HRP doit être dilué au 1:100 avec le Tampon d'essai (1x) juste avant l'utilisation dans un tube à essais en plastique propre.

Le conjugué HRP peut être préparé selon le tableau suivant :

Nombre de bandes	Conjugué HRP (ml)	Tampon d'essai (1x) (ml)
1 - 6	0.03	2.97
1 - 12	0.06	5.94

4.4. Étalon human sICAM-1

Des **dilutions d' étalon** peuvent être préparées directement sur la plaque de microtitration (voir 5.c) ou comme alternative dans des tubes (voir 4.4.1).

4.4.1. Dilution d' étalon externe

Etiquetter les tubes 4, une pour chaque point d' étalon.

S2, S3, S4, S5

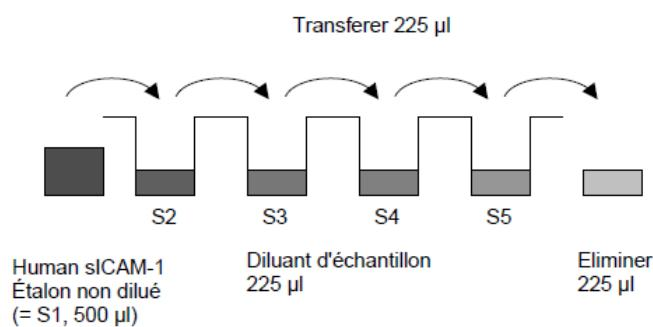
Puis préparer séries de dilutions 1:2 pour la courbe d' étalonnage de manière suivante: Pipeter 225 µl de Diluant d'échantillon dans les tubes S2 – S5.

Pipeter 225 µl d' étalon non dilué (sert comme étalon le plus élevé S1, concentration d' étalon 1 = 100 ng/ml) dans un premier tube marqué S2 et agiter (concentration d' étalon 2 = 50 ng/ml).

Répéter des séries de dilutions 3 fois pour créer les dilutions d' étalon pour la courbe d' étalonnage (voir Figure 1).

Diluant d'échantillon sert comme contrôle vide.

Figure 1



4.5. Contrôles

Solubiliser en ajoutant 100 µl d'eau distillée aux **contrôles** lyophilisés. Laisser reconstituer les contrôles pendant 10-30 min. Agiter doucement jusqu'à la dissolution complète et homogène. Traiter ensuite les contrôles comme les échantillons dans le test. Pour la gamme étalon référez au certificat d'analyse ou l'étiquette de l'ampoule. Conserver les contrôles réconstitués aliquotés à -20°C. Eviter la congélation et le dégel répété.

5. Protocole de Test

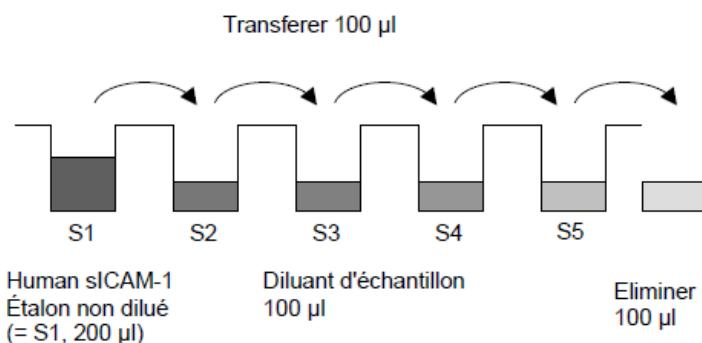
1. Déterminer le nombre de barrettes de puits de microtitration nécessaires pour tester le nombre souhaité d'échantillons plus les barrettes nécessaires aux contrôles vides et aux étalons. Chaque échantillon, étalon, contrôle vide et contrôle doit être testé en double. Retirer les barrettes de microtitration inutiles du support et les stocker à 2°-8°C dans une pochette hermétiquement refermée, avec le dessiccatif fourni.
2. Laver deux fois les barrettes de puits avec environ 400 µl de **tampon de lavage** pour chaque puits et en aspirant à fond le contenu des puits entre les lavages. Laisser le Tampon de lavage dans les puits pendant **10 - 15 secondes** avant l'aspiration. Veiller à ne pas rayer la surface des puits de microtitration.

Après le dernier lavage, vider les barrettes de puits et les tapoter sur un tampon absorbant ou une serviette en papier pour éliminer l'excès de tampon de lavage. Utiliser les barrettes de micropuits immédiatement après le lavage ou les placer renversées sur un papier absorbant pendant 15 minutes au maximum. **Ne pas laisser sécher les puits.**

3. **Dilution d'étalon sur la plaque de microtitration** (Comme alternative des dilutions d'étalon peuvent être préparées dans des tubes –voir 4.4.1)

Ajouter en double 100 µl de Diluant d'échantillon dans les **puits d'étalon** B1/2- E1/2, en laissant les puits A1/2 vide. Pipeter en double 200 µl d'**étalon** non dilué (voir Préparation d'étalon 4.4, concentration d'étalon 1, S1 = 100.0 ng/ml) dans les puits A1 et A2 (voir Tableau 1). Transférer 100 µl dans les puits B1 et B2 (concentration d'étalon 2, S2 = 50.0 ng/ml). Mélanger bien le contenu des puits B1 et B2 par aspiration et ejection répétée et transférer 100 µl dans les puits C1 et C2, respectivement. (voir Figure 2) Veiller à ne pas rayer la surface des puits de microtitration. Continuer la procédure 2 fois en préparant deux séries de dilutions d'étalon human sICAM-1, de 100.0 à 6.3 ng/ml. Eliminer 100 µl du contenu des derniers puits (E1, E2).

Figure 2



Dans le cas d'**une dilution d'étalon externe** (voir 4.4.1), pipeter 100 µl de ces dilutions d'étalon (S1 –S5) dans les puits de façon montrée dans Tableau 1.

- Ajouter 100 µl de **Diluant d'échantillon** dans tous les **puits de contrôle vide**.

Tableau 1

Exemple d'arrangement d'échantillons, d'étalons et de contrôles vides dans les barrettes de puits de microtitration.

	1	2	3	4
A	Étalon 1 (100.0 ng/ml)	Étalon 1 (100.0 ng/ml)	Échantillon 3	Échantillon 3
B	Étalon 2 (50.0 ng/ml)	Étalon 2 (50.0 ng/ml)	Échantillon 4	Échantillon 4
C	Étalon 3 (25.0 ng/ml)	Étalon 3 (25.0 ng/ml)	Échantillon 5	Échantillon 5
D	Étalon 4 (12.5 ng/ml)	Étalon 4 (12.5 ng/ml)	Échantillon 6	Échantillon 6
E	Étalon 5 (6.3 ng/ml)	Étalon 5 (6.3 ng/ml)	Échantillon 7	Échantillon 7
F	Contrôle vide	Contrôle vide	Échantillon 8	Échantillon 8
G	Échantillon 1	Échantillon 1	Échantillon 9	Échantillon 9
H	Échantillon 2	Échantillon 2	Échantillon 10	Échantillon 10

- Ajouter 90 µl de **Diluant d'échantillon** dans tous les puits d'échantillon.
- Ajouter 10 µl de chaque échantillon, en double, dans les puits d'échantillon.
- Préparer du Conjugué HRP (se reporter à la préparation des réactifs Conjugué HRP 4.3).
- Ajouter 50 µl de Conjugué HRP dans tous les puits.
- Recouvrir avec un couvre-plaque et incuber à température ambiante (entre 18° et 25°C) pendant 1 heure, si possible sur un agitateur rotateur réglé à 400 tr/min.
- Retirer le couvre-plaque et vider les puits. **Laver** 3 fois les barrettes de puits de microtitration comme indiqué à point b de ce protocol. Utiliser les barrettes de micropuits immédiatement après le lavage.
- Pipeter 100 µl de solution de substrat TMB dans chaque puits, y compris les puits de contrôle vide.
- Incuber les puits de microtitration à température ambiante (entre 18 et 25 C) pendant environ 10 minutes. Éviter toute exposition directe à une source de lumière intense.

Les valeurs de densité optique au niveau de la plaque doivent être surveillées et la réaction du substrat stoppée (voir le point prochain) avant que les puits positifs ne soient plus correctement mesurables.

La duree de l'incubation pour le developpement de couleur doit être determiné pour chaque essai individuellement.

Il est recommandé d'ajouter la solution stop quand une couleur bleu sombre se développe à la concentration la plus haute de la gamme étalon. Une autre alternative consiste à suivre le développement de la couleur par lecteur ELISA à 620 nm. La réaction du substrat doit être arrêtée dès que la DO atteint 0.9 à 0.95.

- Arrêter la réaction enzymatique en pipetant rapidement 100 µl de **solution d'arrêt** dans chaque puits, y compris les puits de contrôle vide. Il est important que la solution d'arrêt soit répandue rapidement et uniformément dans les puits pour inactiver complètement l'enzyme. Les résultats doivent être lus immédiatement après l'ajout de la solution d'arrêt ou dans l'heure qui suit si les barrettes de microtitration sont conservées à l'obscurité entre 2 et 8 °C.
- Lire l'absorbance de chaque puits sur un spectrophotomètre avec 450 nm comme longueur d'onde primaire (éventuellement 620 nm comme longueur d'onde de référence; 610 à 650 nm sont acceptables). Mesurer le contrôle vide du lecteur de plaque conformément aux instructions du fabricant, en utilisant les puits de contrôle vide. Déterminer l'absorbance des échantillons et des human sICAM-1.

Les échantillons ont été dilués 1:10 en cours de test. Pour cette raison, la valeur de concentration déterminée par la gamme étalon doit être multipliée par le facteur de dilution (x 10).

Remarque: Si la plaque n'est pas agitée pendant l'incubation, les valeurs de densité optique peuvent être inférieur aux valeurs indiqués plus haut. Néanmoins ces valeurs sont valables.

INFORMAZIONI SUL PRODOTTO E MANUALE (Italiano)

1. Reagenti Forniti

1.1. Reagenti Forniti per human sICAM-1 ELISA BMS201CE (96 tests)

- 1 busta d'alluminio con **Piastra Micropozzetti rivestita** con anticorpo monoclonale anti human sICAM-1
- 1 flaconcino (100 µl) di anticorpo **HRP-Coniugato** (anticorpo policlonale human sICAM-1)
- 2 flaconcini (500 µl) human sICAM-1 **Standard**, 100 ng/ml
- 1 flaconcino di **Controllo basso**, liofilizzato
- 1 flaconcino di **Controllo alto**, liofilizzato
- 1 flaconcino (12 ml) con **Diluente del Campione**
- 1 flaconcino (5 ml) con **Tampone del Saggio concentrata** 20x (PBS con 1% Tween 20 e 10% BSA)
- 1 bottiglia (50 ml) con **Tampone di Lavaggio concentrato** 20x (PBS con 1% Tween 20)
- 1 flaconcino (15 ml) di **Soluzione Substrato** (tetrametilbenzidina)
- 1 flaconcino (15 ml) di **Soluzione bloccante** (acido fosforico 1M)
- 2 **Copripiasta** adesivi

1.2. Reagenti Forniti per human sICAM-1 ELISA BMS201TENCE (10x96 tests)

- 10 buste d'alluminio con **Piastra Micropozzetti rivestita** con anticorpo monoclonale anti human sICAM-1
- 10 flaconcini (100 µl) di anticorpo **HRP-Coniugato** (anticorpo monoclonale verso human sICAM-1)
- 10 flaconcini (500 µl) human sICAM-1 **Standard**, 100 ng/ml 10 flaconcini di **Controllo basso**, liofilizzato
- 10 flaconcini di **Controllo alto**, liofilizzato
- 10 flaconcini (12 ml) con **Diluente del Campione**
- 1 flaconcino (5 ml) con **Tampone del Saggio concentrata** 20x (PBS con 1% Tween 20 e 10% BSA)
- 3 bottiglie (50 ml) con **Tampone di Lavaggio concentrato** 20x (PBS con 1% Tween 20)
- 10 flaconcini (15 ml) di **Soluzione Substrato** (tetrametilbenzidina)
- 1 flaconcino (100 ml) di **Soluzione bloccante** (acido fosforico 1M) 10 **Copripiasta** adesivi

2. Istruzioni di Conservazione

Conservare i reagenti del kit a 2°-8° C e i controlli a -20° C. Subito dopo l'uso riporre i reagenti nel luogo di conservazione a 2°-8° C e i controlli a -20° C. La scadenza del kit e dei reagenti è indicata sulle etichette.

La data di scadenza dei componenti del kit può essere garantita solo se questi sono conservati correttamente e, in caso di uso ripetuto di un componente, il reagente non è stato contaminato durante la prima manipolazione.

3. Precauzioni per l'Uso

- Tutti i prodotti chimici vanno considerati come potenzialmente pericolosi. Raccomandiamo, perciò, l'utilizzo di questo prodotto solo da personale addestrato alle tecniche di laboratorio e che siano avvezze alle comuni pratiche di laboratorio. Indossare abbigliamento idoneo come camici, guanti ed occhiali. Attenzione ad evitare contatto con la pelle e gli occhi. Nel caso di contatto con pelle o occhi, immediatamente lavare con acqua. Consultare la scheda di sicurezza del prodotto per specifici consigli.
- I reagenti sono per uso in vitro diagnostico e non sono per uso terapeutico.
- Non mischiare tra loro reagenti di diversi lotti o provenienza.
- Non usare i kit dopo la data di scadenza.
- Non esporre i reagenti del kit, durante la conservazione e incubazione a forti fonti di luce.
- Non pipettare utilizzando la bocca.
- Non mangiare o fumare nell'area dove sono utilizzati i reagenti del kit o i campioni.
- Evitare il contatto dei reagenti o campioni con la pelle o le mucose.
- Guanti di gomma o lattice dovrebbero essere sempre indossati quando si usano reagenti e campioni.
- Evitare il contatto tra il substrato del kit e agenti ossidanti e metallo.
- Evitare schizzi o produzione di aerosol.
- Per evitare contaminazione microbica o cross-contaminazione dei reagenti o dei campioni che invaliderebbero il test, usare sempre pipette e puntali mono-uso.
- Usare vaschette pulite e dedicate per la dispensare il reagente substrato.
- L'esposizione agli acidi inattiva il coniugato.
- Acqua distillata o de-ionizzata deve essere utilizzata per la preparazione dei reagenti.
- La soluzione di substrato deve essere portata a temperatura ambiente prima dell'utilizzo.
- Decontaminare ed eliminare i campioni e tutto il materiale potenzialmente contaminante perché potrebbero contenere agenti infettanti. Il metodo preferito per la decontaminazione è l'autoclavaggio per minimo 1 ora a 121.5°C.
- Gli scarti liquidi, non contenenti acido e gli scarti neutralizzati possono essere mischiati con sodio ipoclorito in un volume finale di 1.0%. Lasciare minimo 30 minuti per l'effettiva decontaminazione. Scarti liquidi contenenti acido devono essere neutralizzati prima dell'aggiunta di sodio ipoclorito.

4. Preparazione dei Reagenti

Prima di cominciare con le procedure del test i **concentrati** dei tamponi devono essere portati a temperatura ambientale e diluiti alle concentrazioni adeguate. Se i **concentrati dei tamponi** presenta cristalli in sospensione, riscaldare lievemente i tamponi fino a ottenere la completa dissoluzione dei cristalli.

4.1. Tampone di Lavaggio (1x)

Versare l'intero contenuto (50 ml) del **tampone di lavaggio concentrato** (20x) in un cilindro graduato pulito da 1000 ml. Portare il volume finale a 1000 ml utilizzando acqua distillata o acqua deionizzata. Mescolare delicatamente per evitare la formazione di schiuma.

Trasferire il prodotto in una bottiglia pulita e conservare a temperature comprese fra 2°C e 25°C. Il tampone di lavaggio è stabile per 30 giorni.

Se necessario, è possibile preparare il tampone di lavaggio secondo la tabella seguente:

Numero di strip	Tampone di lavaggio (20x) (ml)	Acqua distillata (ml)
1 - 6	25	475
1 - 12	50	950

4.2. Tampone del Saggio (1x)

Versare l'intero contenuto (5 ml) del **tampone del saggio concentrato** (20x) in un cilindro graduato pulito da 100 ml. Portare il volume finale a 100 ml utilizzando acqua distillata o acqua deionizzata. Mescolare delicatamente per evitare la formazione di schiuma.

Conservare a temperatura compresa fra 2°C e 8°C. La soluzione tampone diluita è stabile per 30 giorni.

Se necessario, è possibile preparare la soluzione tampone secondo la tabella seguente:

Numero di strip	Tampone del Saggio (20x) (ml)	Acqua distillata (ml)
1 - 6	2.5	47.5
1 - 12	5.0	95.0

4.3. HRP-Coniugato

Il HRP-Coniugato deve essere utilizzato entro 30 minuti dalla diluizione.

Il HRP-Coniugato deve essere diluito 1:100 con Tampone del Saggio (1x) in una provetta di plastica pulita secondo la tabella seguente:

Numero di strip	HRP-Coniugato (ml)	Tampone del Saggio (1x) (ml)
1 - 6	0.03	2.97
1 - 12	0.06	5.94

4.4. Human sICAM-1 Standard

La diluizione dello standard può essere fatto direttamente nella piastra (vedi 5.c.) oppure nei tubi (vedi 4.4.1).

4.4.1. Diluizione degli Standard esterni

Etichettare 4 tubi, uno per ogni punto dello standard.

S2, S3, S4, S5

Preparare diluizione seriali 1:2 per lo standard nel seguente modo:

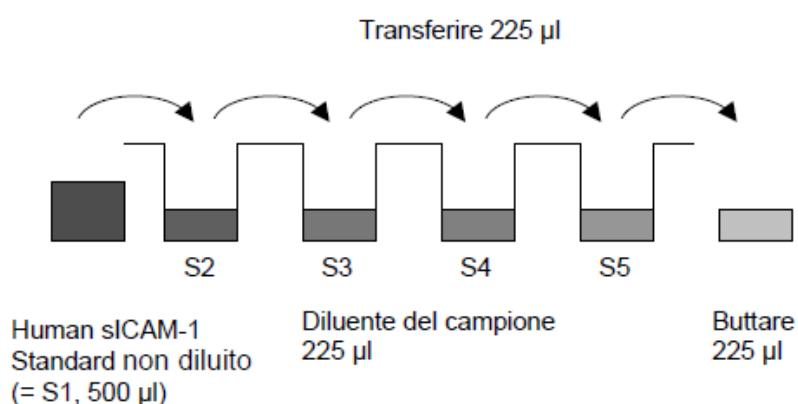
Pipettare 225 ul di Diluente del campione nei tubi S2 – S5.

Pipettare 225 ul di ricostituito non diluito (serve come standard più alto S1, concentrazione dello standard 1=100 ng/ml) nel primo tubo, etichettato S2, e mescolare (concentrazione dello standard 2=50.0 ng/ml).

Ripetere le 3 diluizioni seriali in modo da creare i punti della curva di calibrazione (vedere Figura 1)

Diluente del campione serve come bianco.

Figura 1



4.5. Controlli

Solubilizzare aggiungendo 100 µl di acqua distillata ai controlli liofilizzati. Permettere ai controlli di riposare per 10-30 minuti. Agitare o mescolare delicatamente per assicurare una solubilizzazione completa ed omogenea. In seguito considerare i controlli allo stesso modo dei campioni del dosaggio. Per il range dei valori del controllo si rimanda al certificato di analysis o all'etichetta presente sulla fiala. Conservare i controlli ricostituiti in aliquote a -20°C. Evitare ripetuti cicli di scongelamento.

5. Procedura del Test

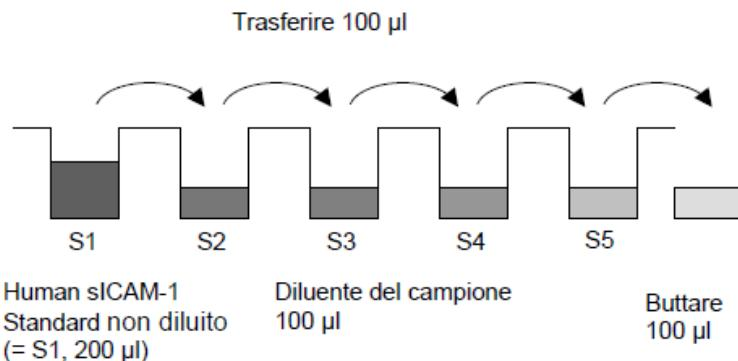
1. Stabilire il numero di strip dei micropozzetti necessarie per analizzare la quantità desiderata di campioni più le strip per i bianchi e gli standard. Tutti i campioni, gli standardi, il bianco e i campioni di controllo opzionali devono essere processati in duplicato. Rimuovere dal supporto le strip micropozzetti non utilizzate e conservarle nella bustina metallica contenente la polvere essiccatrice, mantenendole a 2°-8°C e perfettamente sigillate.
2. Lavare due volte le strip micropozzetti utilizzando circa 400 µl di **tampone di lavaggio** per pozzetto, aspirando accuratamente il contenuto dei micropozzetti tra un lavaggio e l'altro. Permettere al tampone di lavaggio di rimanere, nei pozzetti, circa **10-15 secondi** prima dell'aspirazione. Evitare di scalfire la superficie dei micropozzetti.

Dopo l'ultimo lavaggio, asciugare le strip micropozzetti con un tampone o carta assorbente per rimuovere il tampone di lavaggio in eccesso. Utilizzare le strip subito dopo il lavaggio o sistemarle capovolte su carta assorbente umida per non più di 15 min. **Non lasciar asciugare i pozzetti.**

3. **Diluizione dello standard in micropozzetti** (alternativamente la diluizione dello standard può avvenire in tubi – vedi 4.4.1) Aggiungere 100 ul di Diluente del campione in duplicato a **pozzetti standard B1/2- E1/**, lasciando A1/A2 vuoti. Pipettare 200 ul **standard** non diluito (vedi preparazione dello standard 4.4.1, concentrazione dello standard 1, S1 = 100 ng/ml) in duplicato nei pozzetti A1 e A2 (vedi Tabla 1). Trasferire 100 ul ai pozzetti B1 e B2 (concentrazione dello standard 2, S2 = 50 ng/ml). Mescolare il contenuto dei pozzetti B1 e B2 attraverso ripetute aspirazione ed iniezioni e trasferire 100 ul, rispettivamente, nei pozzetti C1 e C2 (vedere Figura 2). Fare attenzione a non graffiare la parte interna dei pozzetti. Continuare questa procedura per 2 volte, creando due colonne di standard in diluizione con concentrazione da 100.0 a 6.3 ng/ml. Buttare 100 µl del contenuto degli ultimi pozzetti (E1 e E2).

In caso di **diluizione esterna dello standard** (vedi 4.4.1) pipettare 100 ul di queste diluizioni standard (S1 –S5) nei pozzetti degli standard come da Tabla 1.

Figura 2



4. Dispensare 100 µl di **Diluente del campione** in duplicato ai **pozzetti de bianco**.
5. Dispensare 90 µl di **Diluente del campione** in duplicato ai **pozzetti dei campioni**.
6. Dispensare 10 µl di **campione** in duplicato ai **pozzetti dei campioni**.

Tabla 1

Tabella rappresenta un esempio dell'organizzazione dei bianchi, standardi e campioni nei pozzetti:

	1	2	3	4
A	Standard 1 (100.0 ng/ml)	Standard 1 (100.0 ng/ml)	Campione 3	Campione 3
B	Standard 2 (50.0 ng/ml)	Standard 2 (50.0 ng/ml)	Campione 4	Campione 4
C	Standard 3 (25.0 ng/ml)	Standard 3 (25.0 ng/ml)	Campione 5	Campione 5
D	Standard 4 (12.5 ng/ml)	Standard 4 (12.5 ng/ml)	Campione 6	Campione 6
E	Standard 5 (6.3 ng/ml)	Standard 5 (6.3 ng/ml)	Campione 7	Campione 7
F	Bianco	Bianco	Campione 8	Campione 8
G	Campione 1	Campione 1	Campione 9	Campione 9
H	Campione 2	Campione 2	Campione 10	Campione 10

7. Preparare la **HRP-coniugato** (consultare la sezione HRP-coniugato 4.3 sulla preparazione dei reagenti).
8. Dispensare 50 µl di **HRP-coniugato** a ciascun pozzetto.
9. Coprire con un copripiasta e incubare a temperatura ambiente (18°- 25°C) per 1 ora utilizzando, se disponibile, un vortex a 400 rpm.
10. Rimuovere il copripiasta e svuotare i pozzetti. **Lavare** le strip della pozzetti 3 volte come descritto in punto 5.b. del protocollo. Procedere immediatamente al punto successivo.
11. Pipettare 100 µl di **soluzione substrato TMB** in tutti i pozzetti, inclusi quelli del blank.
12. Incubare le strip a temperatura ambiente (18°-25° C) per circa 10 minuti. Evitare l'esposizione diretta a luci intense.

È necessario monitorare i valori O.D. a livello della piastra e interrompere la reazione del substrato (vedi il punto prossimo del protocollo) prima che i pozzetti positivi cessino di essere appropriatamente registrabili.

La determinazione del tempo necessario per lo sviluppo del colore dev'essere fatto per ogni singolo parametro.

Si raccomanda di aggiungere la soluzione di stop quando lo standard più elevato ha sviluppato un colore blu scuro.

Alternativamente lo sviluppo del colore può essere monitorato con un lettore ELISA a 620 nm. La reazione del substrato deve essere bloccata non appena viene misurato un valore delle OD di 0.9 - 0.95.

13. Interrompere la reazione enzimatica pipettando rapidamente 100 µl di **soluzione bloccante** in ciascun pozzetto, inclusi i pozzetti del bianco. È importante che la soluzione bloccante si diffonda rapidamente e uniformemente attraverso i micropozzetti per inattivare completamente l'enzima. I risultati devono essere letti immediatamente dopo l'aggiunta della soluzione bloccante o entro 1 ora se le strip sono conservate in un luogo buio a 2°-8° C.
14. Leggere l'assorbanza di ciascun micropozzetto su uno spettrofotometro che utilizza 450 nm come lunghezza d'onda primaria (620 nm come lunghezza d'onda di riferimento alternativa; valori da 610 nm a 650 nm sono accettabili). Azzerare il lettore della piastra secondo le istruzioni del produttore e utilizzando i pozzetti del bianco. Determinare l'assorbanza sia dei campioni, sia degli standard di human sICAM-1.

I campioni sono stati diluiti 1: 10, quindi la concentrazione dalla curva standard risultante deve essere moltiplicata per il fattore di diluizione (x 10).

Annotazione: In caso di incubazione senza agitazione i valori di densità ottica (O.D.) potranno essere più bassi di quanto indicato sotto. Tuttavia i risultati saranno da ritenersi validi.

Documentation and support

Obtaining support

Technical support	For the latest services and support information for all locations, visit www.thermofisher.com . At the website, you can: <ul style="list-style-type: none">• Access worldwide telephone and fax numbers to contact Technical Support and Sales facilities• Search through frequently asked questions (FAQs)• Submit a question directly to Technical Support (thermofisher.com/support)• Search for user documents, SDSs, vector maps and sequences, application notes, formulations, handbooks, certificates of analysis, citations, and other product support documents• Obtain information about customer training• Download software updates and patches
Safety Data Sheets (SDS)	Safety Data Sheets (SDSs) are available at thermofisher.com/support .
Limited product warranty	Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.thermofisher.com/us/en/home/global/terms-and-conditions.html . If you have any questions, please contact Life Technologies at www.thermofisher.com/support .

For support visit thermofisher.com/support or email techsupport@lifetech.com
thermofisher.com

23 January 2017

ThermoFisher
SCIENTIFIC