

ab169812 – Human Chemokine Antibody Array – Membrane

Instructions for Use

For the simultaneous detection of 38 Human Chemokine levels in all sample types

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

Table of Contents

INTRODUCTION

- 1. BACKGROUND 2
- 2. ASSAY SUMMARY 4

GENERAL INFORMATION

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

ASSAY PREPARATION

- 9. REAGENT PREPARATION 9
- 10. SAMPLE PREPARATION AND STORAGE 10
- 11. ARRAY MAP 15

ASSAY PROCEDURE

- 12. ASSAY PROCEDURE 17

DATA ANALYSIS

- 13. CALCULATIONS 20
- 14. TYPICAL DATA 22

RESOURCES

- 16. TROUBLESHOOTING 23
- 17. NOTES 25

1. BACKGROUND

New techniques such as cDNA microarrays have enabled us to analyze global gene expression. However, almost all cell functions are executed by proteins, which cannot be studied simply through DNA and RNA techniques. Experimental analysis clearly shows a disparity between the relative expression levels of mRNA and their corresponding proteins. Therefore, analysis of the protein profile is critical.

The conventional approach to analyzing multiple protein expression levels has been to use 2-D SDS-PAGE coupled with mass spectrometry. However, these methods are slow, expensive, labor-intensive and require specialized equipment. Thus, effective study of multiple protein expression levels can be complicated, costly and time-consuming. Moreover, these traditional methods of proteomics are not sensitive enough to detect most cytokines (typically at pg/mL concentrations).

Cytokines, broadly defined as secreted cell–cell signaling proteins and peptides distinct from classic hormones or neurotransmitters, play an important role in innate immunity, apoptosis, angiogenesis, cell growth and differentiation. They are involved in most disease processes, including cancer, obesity and inflammatory and cardiac diseases.

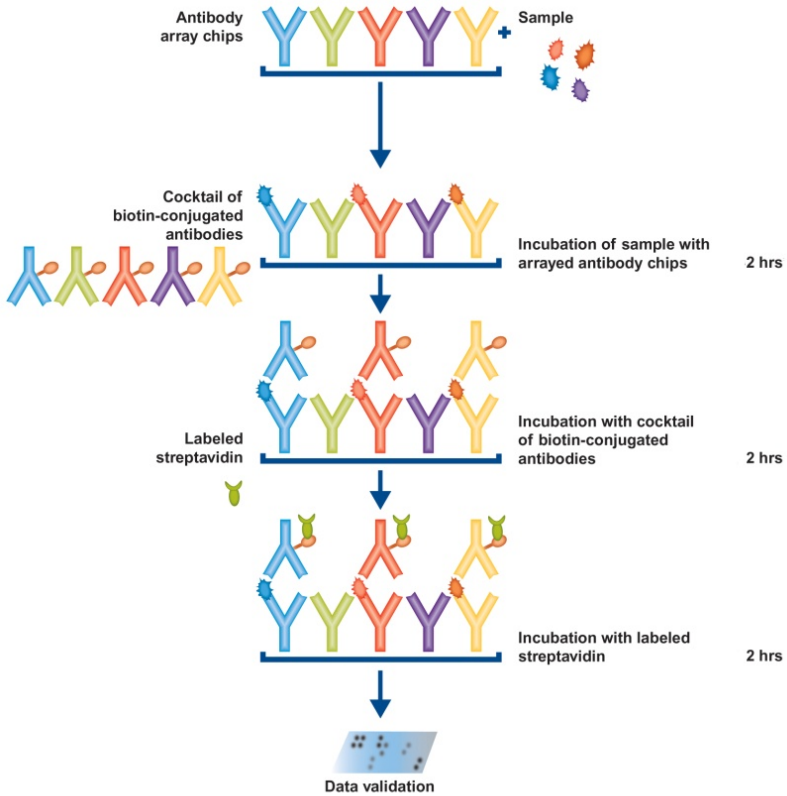
Simultaneous detection of multiple cytokines undoubtedly provides a powerful tool to study cytokines. Regulation of cellular processes by cytokines is a complex, dynamic process, often involving multiple proteins. Positive and negative feedback loops, pleiotrophic effects and redundant functions, spatial and temporal expression of or synergistic interactions between multiple cytokines, even regulation via release of soluble forms of membrane-bound receptors, all are common mechanisms modulating the effects of cytokine signaling. As such, unraveling the role of individual cytokines in physiologic or pathologic

processes generally requires consideration and detection of multiple cytokines rather than of a single cytokine.

Abcam's Cytokine Antibody Arrays have several advantages over detection of cytokines using single-target ELISA kits:

1. More Data, Less Sample: Antibody arrays provide high-content screening using the same sample volume as for ELISA.
2. Global View of Cytokine Expression: Antibody array screening improves the chances for discovering key factors, disease mechanisms or biomarkers related to cytokine signaling.
3. Greater Sensitivity: As little as 4 pg/ml of MCP-1 can be detected using the Cytokine array format. In contrast, our similar MCP-1 ELISA assay has a sensitivity of 40 pg/ml of MCP-1.
4. Increased Range of Detection: ELISA assays typically detect a concentration range of 100- to 1000-fold, however, Abcam arrays can detect a range of 10,000-fold.
5. Better Precision: As determined by densitometry, the inter-array spot-to-spot Coefficient of Variation (CV) ranges from 5 to 10%, which compare favorably with the typical CV in ELISA testing of 10-15%.

2. ASSAY SUMMARY



3. PRECAUTIONS

Please read these instructions carefully prior to beginning the assay.

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

4. STORAGE AND STABILITY

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

Once thawed, store array membranes and 1X Blocking Buffer at $\leq -20^{\circ}\text{C}$, and all other component at 4°C .

Array kits are robust and will retain full activity even if stored for up to 24 hours at room temperature.

5. MATERIALS SUPPLIED

Item	Quantity		Storage Condition (Before Preparation)
Chemokine Antibody Array Membranes	4 Membranes	8 Membranes	-20°C
2000X Biotin-Conjugated Anti-Cytokines	2 Vials	4 Vials	-20°C
1000X HRP-Conjugated Streptavidin	1 Vial	1 Vial	-20°C
1X Blocking Buffer	25 mL	50 mL	-20°C
20X Wash Buffer I	10 mL	20 mL	-20°C
20X Wash Buffer II	10 mL	20 mL	-20°C
2X Cell Lysis Buffer	10 mL	16 mL	-20°C
Detection Buffer C	1.5 mL	2.5 mL	-20°C
Detection Buffer D	1.5 mL	2.5 mL	-20°C
8-Well Plastic Tray	1 Unit	1 Unit	-20°C
Plastic sheets	1 Unit	1 Unit	-20°C

6. MATERIALS REQUIRED, NOT SUPPLIED

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

- Small plastic boxes or containers.
- Pipettors, pipet tips and other common lab consumables.
- Distilled or De-ionized Water.
- Tissue paper, blotting paper or chromatography paper.
- Orbital shaker or oscillating rocker.
- Saran Wrap or similar plastic film.
- A chemiluminescent blot documentation system.
- X-ray Film and a suitable film processor.

7. LIMITATIONS

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

8. TECHNICAL HINTS

Handling Array Membranes

- Array membranes are fragile when dry. Handle with care.
- Wet or dry, grasp membranes by the edges using forceps.
- Do not allow membranes to dry out during experiments.
- The printed side of each membrane is denoted by a dash mark (-) or array number in the upper left corner.
- Unused membranes should be stored at -20°C or -80°C.
- Completed membranes may be stored up to 5 days in 1X Wash Buffer II or 1X Blocking Buffer at 4°C. Be sure to cover container to prevent evaporation during storage.
- For longer term storage of completed membranes, place wetted membranes between 2 plastic sheets (provided in kit), wrap in plastic wrap and store at -20°C or -80°C.

Incubation and Washes

- Perform all incubation and wash steps under gentle rotation or rocking motion (~0.5 to 1 cycle/sec) to ensure complete and even solution coverage as well as to avoid foaming or bubbles from appearing on the membrane surface.
- All washes and incubations (except for detection buffers incubation step) should be performed using the 8-Well Incubation Tray provided in the kit.
- Cover the 8-Well Incubation Tray with the lid for all incubation steps to avoid evaporation.
- Completely cover the membranes with sample or reagent during each incubation. Avoid forceful pipetting directly onto the membrane, instead gently pipette in a corner of each well.
- The sample, biotin-conjugate antibody, and HRP-Streptavidin incubation steps may be performed overnight at 4°C. Overnight incubations are the most effective method of increasing spot intensities but may also increase background noise.

- If you perform overnight sample incubations, we recommend adding the optional “Large Volume Wash” described in Step 12.7 to minimize background signals.

Chemiluminescence Detection

- Trying multiple exposure times is recommended to obtain optimum results. Anywhere from a few seconds to 10 minutes is common with 30 seconds to 1 minute being suitable for most samples.
- If the signals are too weak, increase exposure time (e.g. 2-10 minutes). If the signals are too strong, reduce exposure time (e.g. 3-30 seconds).
- Blot documentation systems that use CCD cameras to detect chemiluminescence are ideal for imaging Abcam array membranes. They can easily be programmed to take multiple exposures, and the dynamic range of these detectors tends to be 2-3 orders of magnitude greater than that of X-ray film or and much more sensitive to chemiluminescence than phosphoimaging systems.

9. REAGENT PREPARATION

Keep all reagents on ice during preparation. Reagents should only be used in their 1X working concentration.

9.1. 1X Wash Buffer I

Dilute 20X Wash Buffer I 20-fold with distilled or deionized water to prepare the 1X Wash Buffer I.

9.2. 1X Wash Buffer II

Dilute 20X Wash Buffer II 20-fold with distilled or deionized water to prepare the 1X Wash Buffer II.

9.3. 1X Biotin-Conjugated Anti-Cytokines

Briefly centrifuge each vial (1 vial is enough to test 2 membranes) and reconstitute by adding 2 mL of 1X Blocking Buffer. Mix thoroughly and gently.

9.4. 1X HRP-Conjugated Streptavidin

Mix the 1000X HRP stock vial well before use. Dilute 1000X HRP-Conjugated Streptavidin 1000-fold with 1X Blocking Buffer to prepare the 1X working concentration.

- 1X Wash Buffers can be stored at 4°C for up to 1 month
- Detection Buffers C and D are supplied at working concentrations.

10. SAMPLE PREPARATION AND STORAGE

10.1. **General Considerations**

- Freeze samples as soon as possible after collection.
- Avoid multiple freeze-thaw cycles. If possible, sub-aliquot your samples prior to initial storage.
- Spin samples hard (5-10 minutes at 10,000 to 15,000 RPM) immediately prior to incubation of samples with array.
- Optimal sample concentrations may need to be determined empirically based on the signal intensities of spots and background signals obtained with each sample.
- If spot intensities are weak, increase sample concentration in subsequent experiments.
- If background or spot intensities are too strong, decrease sample concentration in subsequent experiments.
- Most samples will not need to be concentrated. If concentration is required, we recommend using a spin-column concentrator with a chilled centrifuge.
- Unless otherwise noted, dilute all samples using the same dilution factor in 1X Blocking Buffer.

10.2. **Recommended Sample Volumes and Dilution Factors**

NOTE: All sample dilutions should be made using the 1X Blocking Buffer provided in this kit. For all sample types, final sample volume = 1.0 mL per membrane

- **Cell Cultured Media:** Neat (no dilution needed)
- **Serum & Plasma:** 2-fold to 10-fold dilution
- **Most other Body Fluids:** Neat or 2-fold to 5-fold dilution
- **Cell and Tissue Lysates:** Minimum 5-fold through 10 fold to equal concentrations of total protein in lysate. Optimal sample concentration of cell and tissue lysates should be determined empirically. For more details, please see below (Cell and Tissue Lysates/Homogenates) below.

- Other Liquid Sample Types: Most often Neat or 2-fold to 5-fold. However, optimal dilutions should be determined empirically. For more details, please see Other Liquid Samples section below.

10.3. Preparing Serum and Plasma Samples

- Prepare samples according to established protocols or collection tube manufacturer's instructions. Sub-aliquot into plastic tubes. Store at -20°C or -80°C .
- We do not recommend comparing results between serum and plasma samples or between plasma prepared using different anticoagulants.
- For most applications, you may test plasma samples prepared using any anticoagulant (ie, Heparin, EDTA or Citrate). However, EDTA-prepared plasma may interfere with optimal detection of MMPs and other metal-binding proteins.
- If possible, avoid testing hemolyzed Serum or Plasma samples, as these samples may generate anomalous cytokine expression patterns and/or high background signals.

10.4. Preparing Cell-Cultured Media

- Expression of proteins in cell culture may depend on many variables, including cell type, starting cell number, media composition and growth conditions.
- To start, we recommend seeding $\geq 1 \times 10^6$ cells in a 100 mm culture dish. However, you should consult the scientific literature for tips on how to cultivate your particular cell type.
- Cell type and experimental protocols can profoundly influence cytokine expression. Please consult the scientific literature for details on the effectiveness of various treatments to elicit a desired response, optimal timeframes for growing cells prior to treatment, optimal concentrations and exposure times for treatments and the timing of sample collection

- If possible, use media that is free of recombinant or purified growth factors. If you must add them, we strongly recommend testing an uncultured media aliquot as a sample “blank” to assess baseline signal response for comparison with cultured media samples.
- Serum-containing media rarely produce a baseline signal response with this product. Nevertheless, an ideal experimental design would be to test uncultured media as a sample “blank” to assess baseline signal responses.

10.5. Preparing Cell and Tissue Lysates/Homogenates

IMPORTANT: Lysate sample volumes required must be determined empirically and will depend upon the total protein concentration of each lysate and the intensity of background signals for each sample.

- You must determine the total protein concentration of each lysate/homogenate. We recommend using the BCA method, as it is insensitive to detergents commonly found in lysis buffers.
- Minimum Recommended Total Lysate Protein Concentration (prior to sample dilution) = 1.0 $\mu\text{g}/\mu\text{L}$
- Minimum Recommended Dilution of Lysates (prior to sample incubation): 5 to 10 fold with 1X Blocking Buffer. Dilute all lysate samples to the same final concentration of total lysate protein in 1X Blocking Buffer to 1 mL final volume.
- For your first experiment, we recommend using 200-250 μg of total protein in 1 mL of 1X Blocking Buffer (final volume) for each array membrane.
- Optimal amounts of total lysate protein may range from 50-1000 μg per array membrane. Based upon the signal intensities of background and spots obtained with each sample, you may need to increase or decrease the volume of lysate used in subsequent experiments.

- We recommend preparing cell or tissue lysates using 2X Cell Lysis Buffer which is provided. Be sure to properly dilute Cell Lysis Buffer (1:1 with deionized H₂O) prior to use.
- Other lysis buffers can be used if they are non-denaturing, non-reducing, total salt concentration ≤ 700 mM), contain $\leq 2\%$ total detergent (v/v) and are free of sodium azide (NaN₃).
- We strongly recommend adding protease inhibitors to your cell lysis buffer. Any broad-spectrum protease inhibitor cocktail intended for preparation of mammalian cell/tissue lysates should be sufficient, but please consult the scientific literature before deciding upon the exact composition of cocktail to use.
- Optimal protocols for mechanical disruption vary for different cell and tissue types. Please consult the scientific literature for examples of successful detection of proteins using ELISA or multiplex ELISA techniques in lysates made from cell or tissue samples similar to yours.
- CAUTION. Sonication can quickly heat volumes of 1 mL or less and denature proteins in your samples!
- After extraction, clarify your lysates by centrifugation and save the supernatant for your experiment. Preparing sub-aliquots is strongly recommended.

10.6. Other Liquid Samples

- Abcam Cytokine Antibody Arrays are compatible with most liquid samples, including extracts, perfusates and lavages, as well as body fluids, such as CSF, Sputum, Saliva, Tears, and Urine.
- Be sure to measure the total protein concentrations of these samples prior to sample incubation.
- For samples that have fairly consistent concentrations of total protein between samples (CV $\leq 20\%$), dilute samples using equal volumes for each sample (ie, use the same dilution factor, v/v).

ASSAY PREPARATION

- For samples exhibiting wider ranges of total protein content, dilute to equal concentrations of total protein (eg, 200 µg of total protein) in 1X Blocking Buffer to a final volume of 1 mL.

11. ARRAY MAP

POS – Positive Control

NEG – Negative Control

BLANK – No Antibody

	A	B	C	D	E	F
1	POS	POS	NEG	NEG	BLC	CCL28
2	POS	POS	NEG	NEG	BLC	CCL28
3	Eotaxin-3	Fractalkine	GCP-2	GRO	GRO α	HCC-4
4	Eotaxin-3	Fractalkine	GCP-2	GRO	GRO α	HCC-4
5	MCP-2	MCP-3	MCP-4	MDC	MIG	MIP-1 α
6	MCP-2	MCP-3	MCP-4	MDC	MIG	MIP-1 α
7	PARC	RANTES	SDF-1 α	SDF-1 β	TARC	TECK
8	PARC	RANTES	SDF-1 α	SDF-1 β	TARC	TECK

ASSAY PREPARATION

	G	H	I	J	K	L
1	Ck β 8-1	CTACK	CXCL16	ENA-78	Eotaxin	Eotaxin-2
2	Ck β 8-1	CTACK	CXCL16	ENA-78	Eotaxin	Eotaxin-2
3	I-309	I-TAC	IL-8	IP-10	Lymphotactin	MCP-1
4	I-309	I-TAC	IL-8	IP-10	Lymphotactin	MCP-1
5	MIP-1 β	MIP-1 δ	MIP-3 α	MIP-3 β	MPIF-1	NAP 2
6	MIP-1 β	MIP-1 δ	MIP-3 α	MIP-3 β	MPIF-1	NAP 2
7	BLANK	BLANK	BLANK	BLANK	BLANK	POS
8	BLANK	BLANK	BLANK	BLANK	BLANK	POS

12. ASSAY PROCEDURE

Please prepare all reagents immediately prior to use. All incubations and washes must be performed under gentle rotation/rocking.

- 12.1. Place each membrane printed side up into the 8-well tray provided in the kit.
- 12.2. Block membranes by incubating with 2 mL 1X Blocking Buffer at room temperature (RT) for 30 min.
- 12.3. Aspirate 1X Blocking Buffer from each well.
- 12.4. Pipette 1 mL of diluted or undiluted sample into each well and incubate for 1.5 - 2 hours at room temperature.

NOTE: When trying the protocol for the first time or when the signal is low, we strongly recommend incubating samples overnight at 4°C.

- 12.5. Aspirate samples from each well.

NOTE: After the overnight incubation, we strongly recommend an additional step with Large Volume Wash: After step 12.4, place membranes into clean container(s). Add 20-30 mL of Wash Buffer I per membrane, and wash at RT with gentle shaking or rocking for 30-45 minutes. Return membranes to the 8-well tray.

- 12.6. Pipette 2 mL of 1X Wash Buffer I into each well and incubate for 5 minutes at RT. Repeat this 2 more times for a total of 3 washes using fresh buffer and aspirating each time.
- 12.7. Pipette 2 mL of 1X Wash Buffer II into each well and incubate for 5 minutes at RT. Repeat this 1 more time for a total of 2 washes using fresh buffer and aspirating each time.
- 12.8. Pipette 1 mL of 1X Biotin-Conjugated Anti-Cytokines into each well and incubate for 1.5 - 2 hours at RT or overnight at 4°C.

- 12.9. Aspirate 1X Biotin-Conjugated Anti-Cytokines from each well.
- 12.10. Wash membranes as directed in Steps 12.6 and 12.7.
- 12.11. Pipette 2 mL of 1X HRP-Conjugated Streptavidin into each well and incubate for 2 hours at room temperature or overnight at 4°C.
- 12.12. Aspirate 1X HRP-Conjugated Streptavidin from each well.
- 12.13. Wash membranes as directed in Steps 12.6 and 12.7.
- 12.14. Transfer and place each membrane printed side up onto a sheet of chromatography paper, tissue paper, or blotting paper lying on a flat surface, such as a benchtop.
- 12.15. Remove any excess wash buffer by blotting the membrane edges with another piece of chromatography, blotting, or tissue paper. Transfer and place the membranes, printed side up, onto a plastic sheet (provided) lying on a flat surface.
- 12.16. Into a single, clean tube, pipette equal volumes (1:1) of Detection Buffer C and Detection Buffer D. For 1 membrane add 250 μ L of Detection Buffer C and 250 μ L of Detection Buffer D into tube. Mix well.
- 12.17. Gently pipette 500 μ L of the Detection Buffers mixture onto each membrane (e.g., 500 μ L = 250 μ L of Detection Buffer C and 250 μ L of Detection Buffer D) and incubate for 2 minutes at room temperature (**Do not rock or shake**). Immediately afterwards, proceed to Step 12.17.

NOTE: Exposure should ideally start within 5 minutes after finishing Step 12.17 and completed within 20 minutes as chemiluminescence signals will fade over time.
- 12.18. Place another plastic sheet on top by starting at one end and gently “rolling” the flexible plastic across the surface to the opposite end to smooth out any air bubbles. The membranes should now be “sandwiched” between two plastic sheets.

NOTE: Avoid sliding the plastic sheet along the membranes' printed surface.

- 12.19. If using a CCD camera (recommended), transfer the sandwiched membranes to the imaging system and expose. (See tips for obtaining array images in Section C).

NOTE: If using X-ray film remove the top plastic sheet covering the printed side so that the membranes can be directly exposed to the film.

- 12.20. Try multiple exposures to obtain an image with low background and strong positive control signal spots that do not bleed into one another. Typical exposure times are between few seconds to 2 minutes.
- 12.21. To store, without direct pressure, gently sandwich the membranes between 2 plastic sheets (if not already), tape the sheets together or wrap in plastic wrap to secure them, and store at $\leq 20^{\circ}\text{C}$.

13. CALCULATIONS

Obtaining Densitometry Data:

Visual comparison of array images may be sufficient to see differences in relative protein expression. However, most researchers will want to perform numerical comparisons of the signal intensities (or more precisely, signal *densities*), using 2-D densitometry. Gel/Blot documentation systems and other chemiluminescent or phosphorescent detection systems are usually sold as a package with compatible densitometry software.

To obtain densitometry data from an X-ray film, one must first scan the film to obtain a digitized image using an ordinary office scanner with resolution of 300 dpi or greater. Any densitometry software should be sufficient to obtain spot signal densities from your scanned images. One such software program, ImageJ, is available for free from the NIH (for more info, visit <http://rsbweb.nih.gov/ij/>).

We suggest using the following guidelines when extracting densitometry data from our array images:

- For each array membrane, identify a single exposure that the exhibits low background signal intensity and strong Positive Control signals that do not “bleed” into one another. Exposure times do not need to be identical for each array, but Positive Control signals on each image should have similar intensities.
- Measure the density of each spot using a circle that is roughly the size of one of the largest spots. Be sure to use the same circle (area and shape) for measuring the signal densities on every array for which you wish to compare the results.
- For each spot, use the summed signal density across the entire circle (i.e., total signal density per unit area)

Once you have obtained the raw densitometry data, you should subtract the background and normalize to the Positive Control signals before proceeding to analysis.

Background Subtraction:

On each array, several "Negative Control" and/or "Blank" spots will be included. Blank spots are literally blank; nothing has been printed there. Negative Control spots are printed with the same buffer used to dilute antibodies printed on the array. Thus, the signal intensities of the Negative Controls represent the background plus non-specific binding to the printed spots. We recommend subtracting the mean of 4 or more Negative Control spots for background correction.

Normalization of Array Data:

The amount of biotin-conjugated IgG protein printed for each Positive Control spot is consistent from array to array. As such the intensity of these Positive Control signals can be used to normalize signal responses for comparison of results across multiple arrays, much like housekeeping genes and proteins are used to normalize results of PCR gels and Western Blots, respectively.

To normalize array data, one array is defined as "reference" to which the other arrays are normalized. This choice can be arbitrary. You can calculate the normalized values as follows:

$$X(Ny) = X(y) * P1/P(y)$$

Where:

P1 = mean signal density of Positive Control spots on reference array

P(y) = mean signal density of Positive Control spots on Array "y"

X(y) = mean signal density for spot "X" on Array for sample "y"

X(Ny) = normalized signal intensity for spot "X" on Array "y"

After normalization to Positive Control signal intensities, you can compare the relative expression levels, analyte-by-analyte, among or between your samples or groups. By comparing these signal intensities, one can determine relative differences in cytokine expression in each sample.

14. TYPICAL DATA

Typical results obtained with Abcam Antibody Arrays:



The preceding figure presents typical images obtained with Abcam Human Cytokine Antibody Array. These membranes were probed with conditioned media from two different cell lines. Membranes were exposed to film at room temperature for 1 minute.

Note the strong signals of the Positive Control spots, provided by biotin-conjugated IgG printed directly onto the array membrane in the upper-left and lower-right corners. These Positive Control spots are useful for proper orientation of the array image.

The signal intensity for each antigen-specific antibody spot is proportional to the relative concentration of the antigen in that sample. Comparison of signal intensities for individual antigen-specific antibody spots between and among array images can be used to determine relative differences in expression levels of each analyte sample-to-sample or group-to-group.

15. TROUBLESHOOTING

Problem	Cause	Recommendation
No signal for any spots, including Positive Controls	Complete detection failure	Repeat incubation with HRP-Streptavidin and Detection Buffers.
Weak or no signals antigen-specific spots	Sample is too dilute	Repeat experiment using higher sample concentration.
	Improper dilution of HRP-Streptavidin	Tube may contain precipitants. Repeat detection, mix 1000X HRP-Streptavidin well before diluting reagent.
	Waiting too long to detect chemiluminescent signals.	Repeat detection, making sure to complete this process within 20 minutes.
	Other Tips	Incubate with sample O/N at 4°C.
		Increase concentration of HRP-Streptavidin.
Increase concentration of Biotin-conjugated Anti-Cytokine.		
Extend exposure time (may go overnight).		
Uneven signal or background	Bubbles present on membrane during incubations	Be sure to completely remove all bubbles from membrane surface.
	Membranes were not evenly covered during washes/incubations or allowed to dry out	Completely cover membranes with solution, use a rocker or shaker during washes and incubations.

RESOURCES

Problem	Cause	Recommendation
High background signals	Overexposure	Decrease exposure time.
	Sample is too concentrated	Repeat experiment using more dilute sample.
	NOTE: To reduce background on completed membrane, wash overnight at 4°C in 1X Wash Buffer II, then re-incubate with HRP-Streptavidin and repeat detection.	

16. NOTES

Technical Support

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

For all technical or commercial enquiries please go to:

www.abcam.com/contactus

www.abcam.cn/contactus (China)

www.abcam.co.jp/contactus (Japan)