

ab193660 – Mouse Cytokine Antibody Array - Membrane (144 Targets)

Instructions for Use

For the simultaneous detection of 144 mouse Cytokine proteins in serum, plasma, cell culture media and other liquid samples 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 6
- 6. MATERIALS REQUIRED, NOT SUPPLIED 7
- 7. LIMITATIONS 7
- 8. TECHNICAL HINTS 8

ASSAY PREPARATION

- 9. REAGENT PREPARATION 10
- 10. SAMPLE PREPARATION AND STORAGE 11
- 11. ARRAY MAP 12

ASSAY PROCEDURE

- 12. ASSAY PROCEDURE 14

DATA ANALYSIS

- 13. CALCULATIONS 17
- 14. TYPICAL DATA 19

RESOURCES

- 16. TROUBLESHOOTING 20
- 17. NOTES 22

1. BACKGROUND

Abcam's Mouse Cytokine Antibody Array - Membrane (144 Targets) ab193660 can be used for the simultaneous detection of 144 mouse proteins in serum, plasma, cell culture media and other liquid samples types.

Targets: 4-1BB (TNFRSF9/CD137), 6CKine (CCL21), ACE, ALK-1, Amphiregulin, Axl, bFGF, BLC (CXCL13), Cardiotrophin-1 (CT-1), CD27 (TNFRSF7), CD27 Ligand (TNFSF7), CD30 Ligand (TNFSF8), CD30 (TNFRSF8), CD36 (SR-B3), CD40 (TNFRSF5), CD40 Ligand (TNFSF5), Chordin, CRG-2, CTACK (CCL27), CTLA-4 (CD152), CXCL16, Decorin, DKK-1, CD26 (DPPIV), Dtk, E-Cadherin, EGF, Endoglin (CD105), Eotaxin-1 (CCL11), Eotaxin-2 (MPIF-2/CCL24), Epigen, Epiregulin, E-Selectin, Fas Ligand (TNFSF6), Fc gamma RIIB (CD32b), Flt-3 Ligand, Fractalkine (CX3CL1), Galectin-1, GCSF, GITR (TNFRSF18), GITR Ligand (TNFSF18), GM-CSF, Granzyme B, Gas 1, Gas 6, HAI-1, HGF, HGFR, ICAM-1 (CD54), IFN-gamma, IGFBP-2, IGFBP-3, IGFBP-5, IGFBP-6, IGF-1, IGF-2, IL-1 beta (IL-1 F2), IL-1 R4 (ST2), IL-10, IL-11, IL-12 p40/p70, IL-12 p70, IL-13, IL-15, IL-17A, IL-17B, IL-17 RB, IL-17E (IL-25), IL-17F, IL-1 alpha (IL-1 F1), IL-1 RA (IL-1 F3), IL-2, IL-2 R alpha, IL-20, IL-21, IL-28A (IFN-LAMBDA 2), IL-3, IL-3 R beta, IL-4, IL-5, IL-6, IL-6 R, IL-7, IL-9, I-TAC (CXCL11), JAM-A (CD321/F11R), KC (CXCL1), Leptin, Leptin R, LIX, L-Selectin (CD62L), Lungkine (CXCL15), Lymphotactin (XCL1), MAdCAM-1, MCP-1, MCP-5, M-CSF, MDC, MFG-E8, MIG (CXCL9), MIP-1 alpha (CCL3), MIP-1 gamma, MIP-2, MIP-3 beta (CCL19), MIP-3 alpha (CCL20), MMP-2, MMP-3, Neprilysin, Osteopontin (SPP1), Osteoprotegerin (TNFRSF11B), Pentraxin-3 (TSG-14), Platelet Factor 4 (CXCL4), Prolactin, Pro-MMP-9, P-Selectin, RAGE, RANTES (CCL5), Resistin, SCF, SDF-1 alpha (CXCL12 alpha), Sonic Hedgehog N-Terminal (Shh-N), TNF RI (TNFRSF1A), TNF RII (TNFRSF1B), TACI (TNFRSF13B), TARC (CCL17), I-309 (TCA-3/CCL1), TECK (CCL25), TCK-1 (CXCL7), TIMP-1, TIMP-2, TNF alpha, Thrombopoietin (TPO), TRANCE (TNFSF11), TREM-1, TROY (TNFRSF19), TSLP, TWEAK (TNFSF12), TWEAK R

(TNFRSF12), VCAM-1 (CD106), VEGF-A, VEGFR1, VEGFR2, VEGFR3, VEGF-D.

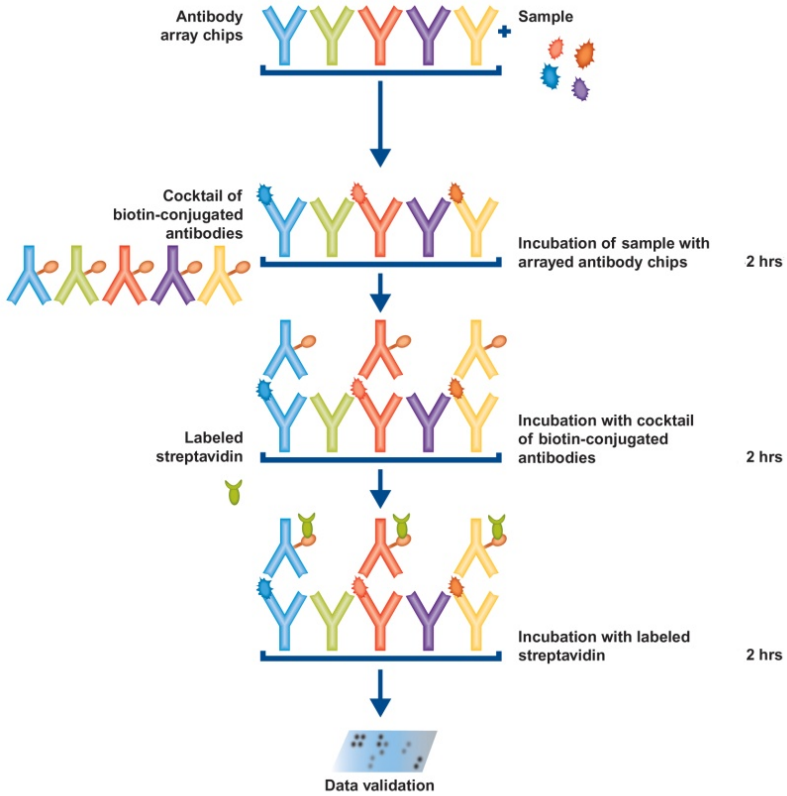
New techniques such as cDNA microarrays have enabled us to analyze global gene expression. However, almost all cell functions are executed by proteins, which cannot be studied simply through DNA and RNA techniques. Experimental analysis clearly shows disparity can exist between the relative expression levels of mRNA and their corresponding proteins⁴. Therefore, analysis of the proteomic 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 distinct from classic hormones or neurotransmitters, play important roles in inflammation, 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 cell signaling pathways. 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.

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, for short-term storage, store array membranes and 1X Blocking Buffer at $\leq -20^{\circ}\text{C}$, and all other component at 2-8°C.

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

5. MATERIALS SUPPLIED

Item	Quantity			Storage Condition Before Preparation
	2X Membranes	4X Membranes	8X Membranes	
Mouse Cytokine Antibody Arrays Membranes (C3)	2X C3 Membranes	4X C3 Membranes	8X C3 Membranes	-20°C
Mouse Cytokine Antibody Arrays Membranes (C4)	2X C4 Membranes	4X C4 Membranes	8X C4 Membranes	-20°C
Mouse Cytokine Antibody Arrays Membranes (C5)	2X C5 Membranes	4X C5 Membranes	8X C5 Membranes	-20°C
Biotinylated Antibody Cocktail (C3)	1X C3 Vials	2X C3 Vials	4X C3 Vials	-20°C
Biotinylated Antibody Cocktail (C4)	1X C4 Vials	2X C4 Vials	4X C4 Vials	-20°C
Biotinylated Antibody Cocktail (C5)	1X C5 Vials	2X C5 Vials	4X C5 Vials	-20°C
1,000X HRP-Conjugated Streptavidin	1X 50 µL	2X 50 µL	4X 50 µL	-20°C
1X Blocking Buffer	2X 25 mL	3X 25 mL	8X 25 mL	-20°C
20X Wash Buffer I	1X 20 mL	2X 20 mL	4X 20 mL	-20°C
20X Wash Buffer II	1X 20 mL	2X 20 mL	4X 20 mL	-20°C
2X Cell Lysis Buffer	1X 16 mL	2X 16 mL	4X 16 mL	-20°C
Detection Buffer C	1X 2.5 mL	2X 2.5 mL	4X 2.5 mL	-20°C
Detection Buffer D	1X 2.5 mL	2X 2.5 mL	4X 2.5 mL	-20°C
8-Well Incubation Tray with Lid	1 Unit	2 Units	4 Units	-20°C
Plastic sheets	1 Unit	2 Units	2 Units	-20°C

The kit also includes plastic sheets, a booklet, an array template and a packing list.

6. MATERIALS REQUIRED, NOT SUPPLIED

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

- Pipettors, pipette tips and other common lab consumables
- Orbital shaker or oscillating rocker
- Tissue paper, blotting paper or chromatography paper
- Adhesive tape or plastic wrap
- Distilled or de-ionized water
- Chemiluminescent blot documentation system (CCD Camera, X-Ray Film and a suitable film processor, Gel documentation system, or other chemiluminescent detection system capable of imaging a Western blot)

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 $\leq -20^{\circ}\text{C}$.
- Completed membranes may be stored up to 5 days in 1X Wash Buffer II or 1X Blocking Buffer at $2-8^{\circ}\text{C}$. 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 .

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 following incubations may be done overnight at $2-8^{\circ}\text{C}$: all wash steps, Step 12.4. (sample incubation), Step 12.8. (biotin-Ab incubation) and Step 12.11. (HRP-Conjugated Streptavidin incubation).

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

Thaw all reagents to room temperature immediately before use. If wash buffers contain visible crystals, warm to room temperature and mix gently until dissolved.

The Biotinylated Antibody Cocktail and the HRP-Streptavidin Concentrate vials should be briefly centrifuged (~1,000 x g) before opening to ensure maximum recovery and mixed well as precipitates may form during storage.

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 Biotinylated Antibody Cocktail

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 1,000X HRP stock vial well before use. Dilute 1,000X HRP-Conjugated Streptavidin 1,000-fold with 1X Blocking Buffer to prepare the 1X working concentration.

9.5 1X Cell Lysis Buffer

Dilute 2X Cell Lysis Buffer Concentrate 2-fold with distilled or deionized water. Only use for preparing cell or tissue lysates.

- 1X Wash Buffers can be stored at 2-8°C for up to 1 month
- The Blocking buffer and 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
- **Other Body Fluids and Liquids:** Neat or 2-fold to 5-fold dilution
- **Cell and Tissue Lysates:** load 50 to 500 μg of total protein (after a 5-fold to 10-fold dilution to minimize the effect of any detergent(s). Therefore the original lysate concentration should be 1 to 5 mg/mL.

11. ARRAY MAP

POS – Positive Control, NEG – Negative Control, BLANK – No Antibody

Array Map for Mouse Cytokine Antibody Array – Membrane C3 (144

Mouse Cytokine Antibody Array C3

	A	B	C	D	E	F	G	H	I	J	K	L	M	N
1	POS	POS	NEG	NEG	BLANK	Axl	BLC	CD30 L	CD30	CD40	CRG-2	CTACK	CXCL16	Eotaxin1
2	POS	POS	NEG	NEG	BLANK	Axl	BLC	CD30 L	CD30	CD40	CRG-2	CTACK	CXCL16	Eotaxin1
3	Eotaxin 2	Fas Ligand	CX3CL1	GCSF	GM-CSF	IFN gamma	IGFBP 3	IGFBP 5	IGFBP 6	IL-1 alpha	IL-1 beta	IL-2	IL-3	IL-3 Rb
4	Eotaxin 2	Fas Ligand	CX3CL1	GCSF	GM-CSF	IFN gamma	IGFBP 3	IGFBP 5	IGFBP 6	IL-1 alpha	IL-1 beta	IL-2	IL-3	IL-3 Rb
5	IL-4	IL-5	IL-6	IL-9	IL-10	IL12 p40/70	IL-12 p70	IL-13	IL-17	KC	Leptin R	Leptin	LUX	L Selectin
6	IL-4	IL-5	IL-6	IL-9	IL-10	IL12 p40/70	IL-12 p70	IL-13	IL-17	KC	Leptin R	Leptin	LUX	L Selectin
7	Ltn/ XCL1	MCP 1	MCP-5	M-CSF	MIG	MIP-1 alpha	MIP-1 gamma	MIP-2	MIP-3 beta	MIP-3 alpha	PF-4	P Selectin	RANTES	SCF
8	Ltn/ XCL1	MCP 1	MCP-5	M-CSF	MIG	MIP-1 alpha	MIP-1 gamma	MIP-2	MIP-3 beta	MIP-3 alpha	PF-4	P Selectin	RANTES	SCF
9	SDF-1 alpha	TARC	TCA-3	TECK	TIMP-1	TNF alpha	sTNFRI	sTNFRII	TPO	VCAM-1	VEGF	BLANK	BLANK	POS
10	SDF-1 alpha	TARC	TCA-3	TECK	TIMP-1	TNF alpha	sTNFRI	sTNFRII	TPO	VCAM-1	VEGF	BLANK	BLANK	POS

Targets) ab193660

Array Map for Mouse Cytokine Antibody Array – Membrane C4 (144

Mouse Cytokine Antibody Array C4

	A	B	C	D	E	F	G	H	I	J	K	L
1	POS	POS	NEG	NEG	BLANK	bFGF	CD26	Dtk	E Selectin	Fc gamma RIIB	Fit-3 Ligand	GITR
2	POS	POS	NEG	NEG	BLANK	bFGF	CD26	Dtk	E Selectin	Fc gamma RIIB	Fit-3 Ligand	GITR
3	HGF R	ICAM-1	IGFBP 2	IGF-1	IGF-2	IL-15	IL-17B R	IL-7	I-TAC	Lungkine	MDC	MMP-2
4	HGF R	ICAM-1	IGFBP 2	IGF-1	IGF-2	IL-15	IL-17B R	IL-7	I-TAC	Lungkine	MDC	MMP-2
5	MMP-3	OPN	OPG	Pro MMP-9	Resistin	Shh-N	TCK-1	TIMP-2	TRANCE	TROY	TSLP	VEGF R1
6	MMP-3	OPN	OPG	Pro MMP-9	Resistin	Shh-N	TCK-1	TIMP-2	TRANCE	TROY	TSLP	VEGF R1
7	VEGF R2	VEGF R3	VEGF D	BLANK	BLANK	BLANK	BLANK	BLANK	BLANK	BLANK	BLANK	POS
8	VEGF R2	VEGF R3	VEGF D	BLANK	BLANK	BLANK	BLANK	BLANK	BLANK	BLANK	BLANK	POS

Targets) ab193660

ASSAY PREPARATION

Array Map for Mouse Cytokine Antibody Array – Membrane C5 (144 Targets) ab193660

Mouse Cytokine Antibody Array C5

	A	B	C	D	E	F	G	H	I	J	K	L	M	N
1	POS	POS	NEG	NEG	4-1BB	6Ckine	ACE	ALK-1	AREG	CT-1	CD27	CD27 Ligand	CD36	CD40 Ligand
2	POS	POS	NEG	NEG	4-1BB	6Ckine	ACE	ALK-1	AREG	CT-1	CD27	CD27 Ligand	CD36	CD40 Ligand
3	Chordin	CTLA 4	Decorin	DKK-1	E Catherin	EGF	Endoglin	Epigen	EPR	Galectin 1	GAS 1	GAS 6	GITR Ligand	Granzyme B
4	Chordin	CTLA 4	Decorin	DKK-1	E Catherin	EGF	Endoglin	Epigen	EPR	Galectin 1	GAS 1	GAS 6	GITR Ligand	Granzyme B
5	HAI-1	HGF	IL-1 R4	IL-11	IL-17 B	IL-17 E	IL-17 F	IL-1ra	IL-2R alpha	IL-20	IL-21	IL-28	IL-6 R	JAM-A
6	HAI-1	HGF	IL-1 R4	IL-11	IL-17 B	IL-17 E	IL-17 F	IL-1ra	IL-2R alpha	IL-20	IL-21	IL-28	IL-6 R	JAM-A
7	MAdCAM-1	MFG E8	NEP	PTX3	PRL	RAGE	TAC1	TREM 1	TWEAK	TWEAK R	NEG	NEG	NEG	POS
8	MAdCAM-1	MFG E8	NEP	PTX3	PRL	RAGE	TAC1	TREM 1	TWEAK	TWEAK R	NEG	NEG	NEG	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.

NOTE: The antibody printed side is marked by a dash (-) or number (#) in the upper left corner.

- 12.2. Block membranes by incubating with 2 mL 1X Blocking Buffer at room temperature (RT) for 30 minutes.
- 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 1): Longer incubations can help maximize the spot signal intensities. However, doing so can also increase the background response so complete liquid removal and washing is critical.

NOTE 2): If sample volume is limited, one C3, one C4, and one C5 membrane can be incubated together in a single well. For 3 membranes per well, use 1.2 mL of sample per well. Rotate bottom membrane to the top every 30 minutes and make sure sample is pipetted in between membranes to ensure even coverage

- 12.5. Aspirate samples from each well.
- 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 Biotinylated Antibody Cocktail into each well and incubate for 1.5-2 hours at RT or overnight at 2-8°C.

NOTE: Ensure only C3 antibody vials are used with C3 membranes, C4 antibody vials are used with C4 membranes, and C5 antibody vials are used with C5 membranes

- 12.9. Aspirate 1X Biotinylated Antibody Cocktail 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 2-8°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.

- 12.16. Transfer and place the membranes, printed side up, onto a plastic sheet (provided) lying on a flat surface.

NOTE: Multiple membranes can be placed next to each other and fit onto a single plastic sheet. Use additional plastics sheets if necessary.

- 12.17. Into a single, clean tube, pipette equal volumes (1:1) of Detection Buffer C and Detection Buffer D. For 1 membrane add 250 µL of Detection Buffer C and 250 µL of Detection Buffer D into tube. Mix well.

- 12.18. Gently pipette 500 µL of the Detection Buffers mixture onto each membrane (e.g., 500 µL = 250 µL of Detection Buffer C and 250 µL of Detection Buffer D) and incubate for

2 minutes at room temperature (**Do not rock or shake**). Immediately afterwards, proceed to Step 12.19.

NOTE: Exposure should ideally start within 5 minutes after finishing Step 12.18 and completed within 20 minutes as chemiluminescence signals will fade over time.

- 12.19. 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.

- 12.20. 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.

- 12.21. 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.

- 12.22. 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's Cytokine Antibody Membrane 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	Chemiluminescent imager is not working properly	Contact image manufacturer
	Too short exposure	Expose the membranes longer
	Degradation of components due to improper storage	Store entire kit at $\leq -20^{\circ}\text{C}$. Do not use kit after expiration date. See storage guidelines.
	Improper preparation or dilution of the HRP-Streptavidin	Centrifuge vial briefly before use, mix well, and do not dilute more than 1,000-fold
	Waiting too long before exposing	The entire detection process should be completed in 10-15 minutes
Positive controls spots signals visible but no other spots	Low sample protein levels	Decrease sample dilution, concentrate samples, or load more protein initially
	Skipped sample incubation step	Samples must be loaded after the blocking step
	Too short incubations	Ensure the incubations are performed for the appropriate time or try the optional overnight incubation(s)
Uneven signal or background	Bubbles present on or below membrane	Don't rock/rotate the tray too vigorously or pipette the sample or reagent with excessive force
	Insufficient sample or reagent volume	Load enough sample and reagent to completely cover the membrane
	Insufficient mixing of reagents	Gently mix all reagents before loading onto the membrane, especially the HRP-Streptavidin and Biotin Antibody Cocktail
	Rocking/rotating on an uneven surface while incubating	Rock/rotate on a flat surface or the sample or reagent can "pool" to one side

RESOURCES

High background signal	Too much HRP-Streptavidin or Biotinylated Antibody Cocktail	Prepare these signal enhancing components precisely as instructed
	Membranes dried out	Do not let the membranes dry out during the experiment. Cover the incubation tray with the lid to minimize evaporation
	Too high sample protein concentration	Increase dilution of the sample or load less protein
	Exposed too long	Decrease exposure time
	Insufficient washing	Ensure all the wash steps are carried out and the wash buffer is removed completely after each wash step
	Non-specific binding	Ensure the blocking buffer is stored and used properly

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)