



## Reagent preparation before starting experiment

- Dilute the 5x Assay wash buffer to 1x buffer  
40ml 5x Assay wash buffer  
160ml ddH<sub>2</sub>O
- Sample preparation
  - Serum-free cell culture conditioned media can be used directly or dilute 2-fold with 1X Diluent buffer before use. When serum-containing conditioned media is required, be sure to use serum as a control.
  - Cell lysate or tissue lysate can be prepared following the protocol on our website: [http://www.signosisinc.com/pdf/Preparation\\_of\\_Cell\\_Lysates\\_for\\_ELISA.pdf](http://www.signosisinc.com/pdf/Preparation_of_Cell_Lysates_for_ELISA.pdf)
  - Sera or plasma can be used directly or diluted 2-10 fold with 1X Diluent buffer.
- Dilute 100 times of biotin labeled antibody mixture with 1X Diluent buffer.
- Dilute 200 times of streptavidin-HRP with 1X Diluent buffer before use.

## Assay procedure

1. Cut the film over the plate and remove it from the desired columns. Make sure the rest of plate is well sealed.
2. Prepare 3.5ml sample and add 100 µl of sample per well to one section and incubate for 2 hours at room temperature with gentle shaking.
3. Invert the plate over an appropriate container and expel the contents forcibly. Wash the plate by adding 200µl of 1x Assay wash buffer. Repeat the washing process two times for a total of three washes. Complete removal of liquid at each wash by firmly tapping the plate against a pile of clean paper towels.
4. Add 100µl of diluted biotin-labeled antibody mixture to each well and incubate for 1 hour at room temperature with gentle shaking.
5. Repeat the aspiration/wash as in step 3.
6. Add 100 µl of diluted streptavidin-HRP conjugate to each well and incubate for 30 min at room temperature with gentle shaking.
7. Repeat the aspiration/wash as in step 3.

8. Add 100µl substrate to each well and incubate for 30 minutes at least.
9. Add 50µl of Stop solution to each well. The color in the wells should change from blue to yellow.
10. Determine the optical density of each well with a microplate reader at 450 nm within 40 minutes.

## Example of Analysis Data

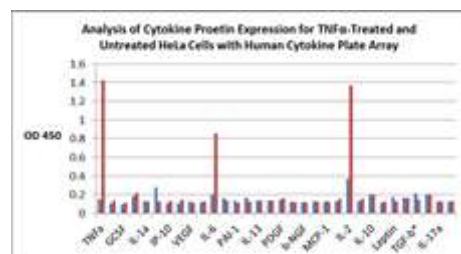


Figure1: Analysis of Cytokine Protein Expression in TNF $\alpha$ -Treated and Untreated HeLa with Human Cytokine ELISA Plate Array

HeLa cells were starved for 24 hours with serum-free medium, subsequently treated the cells with and without 20ng/ul TNF for 16 hours. The serum-free conditioned media were incubated on the plate for 1 hour. After incubating with detection antibody mix and HRP, the plate was detected by a plate reader.

## Diagram of Human Cytokine ELISA Plate Array I

	1	2	3	4	5	6	7	8	9	10	11	12
A	TNF $\alpha$	VEGF	PDGF	IL-10	TNF $\alpha$	VEGF	PDGF	IL-10	TNF $\alpha$	VEGF	PDGF	IL-10
B	IFN $\gamma$	EGF	PLGF	FGFb	IFN $\gamma$	EGF	PLGF	FGFb	IFN $\gamma$	EGF	PLGF	FGFb
C	G-CSF	IL-6	b-NGF	Leptin	G-CSF	IL-6	b-NGF	Leptin	G-CSF	IL-6	b-NGF	Leptin
D	GM-CSF	Resistin	SCF	IGF-1	GM-CSF	Resistin	SCF	IGF-1	GM-CSF	Resistin	SCF	IGF-1
E	IL-1 $\alpha$	PAI-1	MCP-1	TGF- $\beta$	IL-1 $\alpha$	PAI-1	MCP-1	TGF- $\beta$	IL-1 $\alpha$	PAI-1	MCP-1	TGF- $\beta$
F	IL-8	IL-12	MIP-1 $\alpha$	Adipo	IL-8	IL-12	MIP-1 $\alpha$	Adipo	IL-8	IL-12	MIP-1 $\alpha$	Adipo
G	IP-10	IL-13	IL-2	IL-17 $\alpha$	IP-10	IL-13	IL-2	IL-17 $\alpha$	IP-10	IL-13	IL-2	IL-17 $\alpha$
H	Rantes	Eotaxin	IL-4	Blank	Rantes	Eotaxin	IL-4	Blank	Rantes	Eotaxin	IL-4	Blank