

TF Activation Profiling Plate Array I

Catalog Number: FA-1001

(For Research Use Only)

Introduction

Transcription factors (TFs) are a group of cellular proteins that play essential roles in regulating gene expression. They act as sensors to monitor cellular change and convert the signals into gene expression. Often a specific cellular signal pathway can activate multiple TFs and the expression of a specific gene is under the control of multiple TFs. Hence, monitoring the activation of multiple TFs simultaneously is critical to understanding the molecular mechanism of cellular regulation underlying cell signaling and gene expression. **Signosis, Inc.'s** *TF Activation Profiling Plate Array I* is used for monitoring 48 different TFs simultaneously in one sample.

Principle of the Assay

Signosis, Inc.'s TF Activation Profiling Plate Array is used for monitoring the activation of multiple TFs simultaneously. In this technology, a series of biotinlabeled probes are made based on the consensus sequences of TF DNA-binding sites. When the probe mix incubates with nuclear extracts, individual probes will find its corresponding TF and form TF/probe complexes, which can be easily separated from free probes through a spin column purification. The bound probes are detached from the complex and analyzed through hybridization with a plate; each well is specifically pre-coated with complementary sequences of the probes. The captured DNA probe is further detected with Streptavidin-HRP Conjugate. Luminescence is reported as relative light units (RLUs) on a microplate luminometer.

Materials Required but Not Provided

- Nuclear Extraction Kit from Signosis (SK-0001)
- PCR machine and PCR tubes
- Microcentrifuge working at 4 °C
- Hybridization incubator at 42°C
- Plate-Shaker
- Plate reader for luminescent detection
- ddH2O (DNAase-free)
- 8 and 12 Multi-channel pipettes

Materials Provided with the Kit

Component	Qty	Store at
96-Well Plate (with aluminum	1	RT
adhesive seal)		
Isolation Columns	2	RT
Elution Buffer	200 µL	RT
TF Plate Hybridization Buffer	12 mL	RT
5X Plate Hybridization Wash	30 mL	RT
Buffer		
5X Detection Wash Buffer	40 mL	RT
Blocking Buffer	30 mL	RT
Filter Wash Buffer	5 mL	4°C
Filter Binding Buffer	1 mL	4°C
Substrate A	1 mL	4°C
Substrate B	1 mL	4°C
Streptavidin-HRP Conjugate	20 µL	4°C
Substrate Dilution Buffer	8 mL	4°C
TF Binding Buffer Mix	30 µL	-20°C
TF Probe Mix I	10 µL	-20°C

Before Starting the Experiment Prepare the Following:

- 1. Place *Filter Binding Buffer* and *Filter Wash Buffer* on **ice** so they are chilled for the assay (for at least **10 minutes**).
- 2. Warm up *TF Plate Hybridization Buffer* and *Hybridization Wash Buffer* **42**°**C** before use.
- 3. Aliquot 500 μ L of ddH₂O in a 1.5 mL microcentrifuge tube **per sample** on ice so that it is chilled for the assay (for at least 10 minutes).
- 4. Dilute **30 mL** of *5X Plate Hybridization Wash Buffer* with **120 mL** of ddH2O before use.
- 5. Dilute **40 mL** of *5X Detection Wash Buffer* with **160 mL** of ddH2O before use.
- 6. Dilute **20 μL** *Streptavidin-HRP* in **10 mL** Blocking Buffer (1:500 dilution).



Please Read the Assay Procedure Before You Begin

www.signosisinc.com Home Page info@signosisinc.com Questions / Comments support@signosisinc.com Technical Support 1(408)-747-0771 Telephone

Assay Procedure

TF/ DNA Complex Formation

- Mix the following components for each reaction in a tube
 15 μL TF Binding Buffer Mix
 3 μL TF Probe Mix I
 X μL Nuclear Extract (5-15 μg recommended)
 Y μL ddH2O (add up to final volume)
 30 μL Reaction Mix [final volume]
- 2. Incubate the **Reaction Mix** at room temperature (20-23°C) for **30 minutes**.

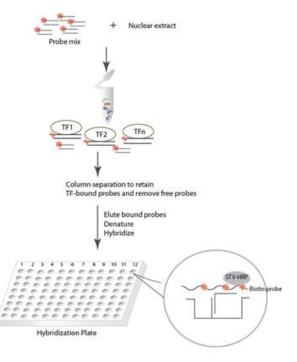
Separation of TF DNA Complex from Free Probes

- Equilibrate an *Isolation Column* by adding 200 μL pre-chilled *Filter Binding Buffer*. Centrifuge the column with the collection tube at 6,000 rpm for 1 minute in a microcentrifuge at room temperature.
- 4. Transfer the **30 μL Reaction Mix** directly onto the filter in the center of the *Isolation Column* (avoiding bubbles).
- 5. Incubate on ice for **30 minutes**. DO NOT incubate longer than 30 minutes; this will result in high background.
- Add 500 µL pre-chilled *Filter Wash Buffer* to the *Isolation Column* and incubate for 3 minutes on ice.
- Centrifuge the *Isolation Column* with the collection tube at 6,000 rpm for 1 minute in a microcentrifuge at 4°C. Discard the flow through from the collection tube.
- Wash the column by adding 500 μL pre-chilled Filter Wash Buffer to the Isolation Column on ice.
- 9. Centrifuge the *Isolation Column* with the collection tube for **1 minute** at **6,000 rpm** in a microcentrifuge at **4°C**. Then discard the flow through.
- 10. Repeat steps 8-9 for an additional **3 times** for a total of 4 washes.

Elution of Bound Probe

- Add 100µL of *Elution Buffer* onto the center of *Isolation Column*, and incubate at room temperature for 5 minutes.
- 12. Place the *Isolation Column* on a new 1.5 mL microcentrifuge tube and centrifuge at **10,000 rpm** for **2 minutes** at room temperature.
- If you have yet to do so, chill 500 µL ddH2O (DNAase free) in a 1.5 mL microcentrifuge tube on ice for at least 10 minutes, and keep on ice.
- 14. Transfer the eluted probe to a PCR tube and denature the eluted probes at **98°C** for **5 minutes**.
- 15. Immediately transfer the denatured probes to the chilled ddH2O from Step 13 and place on ice. The samples are ready for the hybridization phase of the assay. You can store the sample at -20°C for future use. If you decided to store your

www.signosisinc.com Home Page info@signosisinc.com Questions / Comments



sample, go to **step 16**. before proceeding to the hybridization phase.

- 16. <u>Skip this step if you did not freeze your</u> sample for future use.
- A) Thaw your sample back to an aqueous phase at room temperature.
- B) Redistribute the sample into PCR tubes to be reheated at **98°C** for **5 minutes**.
- C) Afterwards, **immediately** place the PCR tubes on ice.
- D) You may now proceed to Step 17.

Hybridization of Eluted Probe with Hybridization Plate

- 17. Remove the clear adhesive film sealing from the provided 96-Well Plate.
- 18. Aliquot **5** mL pre-warmed *TF Plate Hybridization Buffer* to a dispensing reservoir (DNase free) and then add **600** μ L denatured probes. Mix them together by gently shaking the reservoir.
- 19. Using a 8 multi-channel pipette $100 \mu L$ of the mixture from step 18. into the corresponding wells with 8 multi-channel pipette immediately.

Note: the 96-Well Plate is divided into 2 sections of six columns each per sample. If you wish to have a blank to compare your wells against, select one TF you are not interested in and determine its location on the plate by using the diagram on the third page. Add 100 μ L TF Plate Hybridization Buffer only without the eluted probe.

1(408)-747-0771 Telephone 20. Firmly seal the wells with the aluminum adhesive seal to secure well contents. Press the foil over the letters and numbers on the plate to help orient well designations. Hybridize the well contents to the plate by placing the 96-Well Plate in an incubator set at $42^{\circ}C$ overnight.

Detection of Bound Probe

- 21. Remove the aluminum adhesive seal from the experimental wells with a razor blade. Keep the unused wells sealed.
- 22. Invert the *96-Well Plate* over an appropriate container and expel the contents forcibly.
- 23. Wash the plate by adding 200 μL of prewarmed *IX Plate Hybridization Wash Buffer* to each well by row with a 12 multichannel pipette. Incubate the plate for 5 minutes with gentle shaking at room temperature on a plate-shaker. Completely remove at end of 5 minutes by tapping the plate against clean paper towels.
- 24. Repeat step 23 two more times for a total of three washes.
- 25. Add **200 μL** of *Blocking Buffer* to each well by **row** with a **12 multi-channel pipette** and incubate for **5 minutes** at room temperature with gentle shaking on a plate-shaker.
- 26. Invert the plate over an appropriate container to forcibly remove *Blocking Buffer* from the wells.
- If you have yet to do so: add 20 μL of Streptavidin-HRP Conjugate in 10 mL Blocking Buffer (1:500 dilution), enough for the whole plate (6 sections). This is the diluted Streptavidin-HRP Conjugate

- Add 95μ L of *diluted Streptavidin-HRP Conjugate* to each well by row with a 12 multichannel pipette and incubate for 45 minutes at room temperature on a plate-shaker with gentle shaking.
- 29. After the **45 minutes** have elapsed, forcibly remove the *96-Well Plate* contents in an appropriate container. Complete removal of liquid at each wash by firmly tapping the plate against clean paper towels.
- 30. Wash the 96-Well Plate by adding 200 μL 1X Detection Wash Buffer to each well by row with a 12 multi-channel pipette. Incubate the plate for 5 minutes with gentle shaking on a plateshaker at room temperature. Decant the liquid from the wells.
- Repeat step 30 for a total of 3 washes. At the last wash, invert plate on clean paper towels for 1 minute to remove excessive liquid.
- 32. Freshly prepare the *Substrate Solution* in the following ratio:

1 part Substrate A / 1 part Substrate B / 8 parts Substrate Dilution Buffer.

For example, for the entire 96-Well Plate: **1 mL** Substrate A **1 mL** Substrate B **8 mL** Substrate Dilution Buffer

10 mL Substrate Solution

- 33. Add 9 5μL Substrate Solution to each well by row with a 12 multi-channel pipette and incubate the solution in the wells for 1 minute at room temperature.
- 34. Place the plate in the luminometer. Allow plate to sit inside machine for 4 minutes before reading. Set integration time to 1 second with no filter position. For the best results, read the plate within 5-20 minutes.

	1	2	3	4	5	6	7	8	9	10	11	12
A	AP1	CDP	GATA	NF-1	Pit	Stat3	AP1	CDP	GATA	NF-1	Pit	Stat3
в	AP2	CREB	GR/PR	NFAT	PPAR	Stat4	AP2	CREB	GR/PR	NFAT	PPAR	Stat4
С	AR	E2F-1	HIF	NF-E2	PXR	Stat5	AR	E2F-1	HIF	NF-E2	PXR	Stat5
D	ATF2	EGR	HNF4	NFkB	SMAD	Stat6	ATF2	EGR	HNF4	NFkB	SMAD	Stat6
E	Brn-3	ER	IRF	4-Oct	Sp1	TCF/LEF	Brn-3	ER	IRF	4-Oct	Sp1	TCF/LEF
F	C\EBP	Ets	MEF2	p53	SRF	YY1	C\EBP	Ets	MEF2	p53	SRF	YY1
G	CAR	FAST-1	Myb	Pax-5	SATB1	TR	CAR	FAST-1	Myb	Pax-5	SATB1	TR
н	CBF	GAS/ISRE	Myc/Max	Pbx1	Stat1	TFIID	CBF	GAS/ISRE	Myc- Max	Pbx1	Stat1	TFIID

TF Activation Profiling Array I Diagram

info@signosisinc.com Questions / Comments Data Example

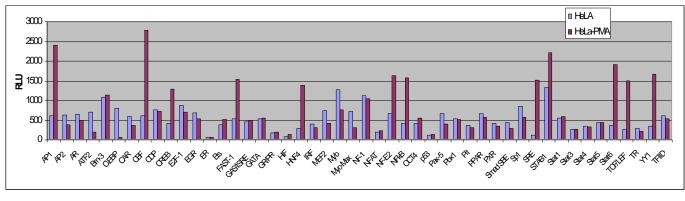


Figure: TF Activation Profiling Plate Array Assay acquired RLUs. HeLA cells were treated with and without PMA. Nuclear Extracts prepared and subjected to the TF Profiling Assay I.

Data analysis notes:

- 1. The TF readings within blank reading $\pm 10\%$ blank reading are considered to be too low for analysis.
- 2. The changes in reading between two samples need to be over 2 fold (increase or decrease) to be significant.
- 3. If you want to choose an internal control for data analysis, select the TF that is not related to your interested pathway and it doesn't show much difference with and without treatment

Gene Description

TF	Gene Description	TF	Gene Description
AP1	Activator protein 1 (JUN/FOS)	NF-1	Nuclear factor 1
AP2	Activator protein 2	NFAT	Nuclear factor of activated T-cells
AR	Androgen receptor	NF-E2	Nuclear factor (erythroid-derived 2)
ATF2	Activating transcription factor 2	NFkB	Nuclear factor of kappa light polypeptide
Brn-3	POU domain, class 4, transcription factor 1	OCT4	POU class 5 homeobox 1
C/EBP	CCAAT/enhancer binding protein (C/EBP),alpha	p53	Tumor protein p53
CAR	Nuclear receptor subfamily 1, group I, member 3	Pax-5	Paired box 5
CBF	CCAAT/enhancer binding protein (C/EBP), zeta	Pbx1	Pre-B cell leukemia transcription factor-1
CDP	Cut-like homeobox 1; CCAAT displacement protein	Pit	Pituitary specific transcription factor 1
CREB	cAMP responsive element binding protein 1	PPAR	Peroxisome proliferator-activated receptor
E2F-1	E2F transcription factor 1	PXR	Pregnane X Receptor
EGR	Early growth response	SMAD	SMAD family
ER	Estrogen receptor	Sp1	SP1 transcription factor
Ets	v-ets erythroblastosis virus E26 oncogene homolog 1	SRF	Serum response factor
FAST-1	Forkhead box H1	SATB1	Special AT-rich sequence binding protein 1
GAS/ISRE	IFN-stimulated response element	Stat1	Signal transducer and activator of transcription 1
GATA	GATA transcription factor	Stat3	Signal transducer and activator of transcription 3
GR/PR	Glucocorticoid receptor/Progesterone receptor	Stat4	Signal transducer and activator of transcription 4
HIF	Hypoxia inducible factor	Stat5	Signal transducer and activator of transcription 5
HNF4	Hepatocyte nuclear factor 4	Stat6	Signal transducer and activator of transcription 6
IRF	Interferon regulatory factor	TCF/LEF	T cell factor / Lymphoid enhancer factor
MEF2	Myocyte enhancer factor 2	YY1	YY1 transcription factor
Myb	v-myb myeloblastosis viral oncogene homolog	TR	Thyroid hormone receptor
Myc-Max	v-myc myelocytomatosis viral oncogene homolog (avian)	TFIID	TATA box binding protein

info@signosisinc.com Questions / Comments support@signosisinc.com Technical Support 1(408)-747-0771 Telephone