



## Cancer Stem Cell TF Activation Profiling Plate Array

Catalog Number FA-1004

(For Research Use Only)

### Introduction

Transcription factors (TF) have been considered as master switches for cell fate determination. The cancer stem cell paradigm postulates that decontrolled tissue-specific stem cells or progenitor cells are precursors for cancer biogenesis. Cancer stem cells are a subset of cancer cells within a tumor that have stem cell-like characteristics. During tumorigenesis, deregulated transcription factor expression or activation can promote abnormal self-renewal, proliferation, and differentiation of the cells. For example, OCT4, Sox2, KLF4, Myc and Nanog have been shown to possess the capability to reprogram somatic cells into pluripotent cells by reactivating pluripotent and silencing somatic genes. Understanding the activation patterns of the related transcription factors in a tumor can help to understand the transcriptional mechanism of cancer stem cell and provide better approaches for targeting and eliminating cancer stem cells.

### Principle of the Assay

Signosis, Inc.'s *Cancer Stem Cell TF activation Profiling Plate Array* is used for monitoring the activation of multiple TFs simultaneously. With our array, a series of biotin-labeled 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 spin column purification. The bound probes are detached from the complex and analyzed through hybridization to the 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 set to 42°C
- Plate-shaker
- Plate reader for luminescent detection
- ddH<sub>2</sub>O (DNAase-free)
- 8 and 12 Multi-channel pipettes

### Materials Provided with the Kit

Component	Qty	Store at
96-Well Plate (with aluminum adhesive seal)	1	RT
Isolation Columns	4	RT
Elution Buffer	600µL	RT
TF Plate Hybridization Buffer	20mL	RT
5X Plate Hybridization Wash Buffer	30mL	RT
5X Detection Wash Buffer	40mL	RT
Blocking Buffer	30mL	4°C
Filter Wash Buffer	15mL	4°C
Filter Binding Buffer	1.5mL	4°C
Substrate A	1mL	4°C
Substrate B	1mL	4°C
Streptavidin-HRP Conjugate	20µL	4°C
Substrate Dilution Buffer	8mL	4°C
TF Binding Buffer Mix	90µL	-20°C
TF Cancer Stem Cell Probe Mix I	20µL	-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*, *Blocking Buffer*, and *Hybridization Wash Buffer* **42°C** before use.
3. Aliquot **300µL** of ddH<sub>2</sub>O in a 1.5mL microcentrifuge tube **per sample** on ice so that it is chilled for the assay (for at least **10 minutes**).
4. Dilute **30mL** of *5X Plate Hybridization Wash Buffer* with **120mL** of ddH<sub>2</sub>O before use.
5. Dilute **40mL** of *5X Detection Wash Buffer* with **160mL** of ddH<sub>2</sub>O before use.
6. Dilute **20µL** *Streptavidin-HRP* in **10mL** *Blocking Buffer* (1:500 dilution).



**Please Read the  
Assay Procedure  
Before You Begin**

# Assay Procedure

## TF/ DNA Complex Formation

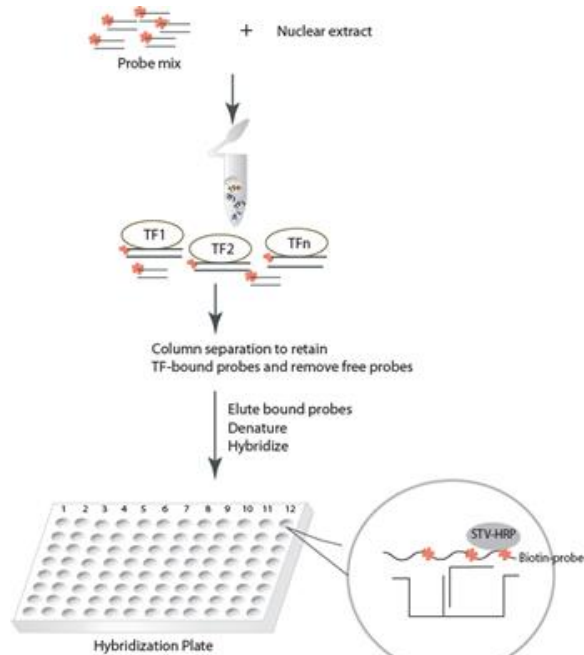
- Mix the following components for each reaction in a tube  
**15 $\mu$ L** *TF Binding Buffer Mix*  
**3 $\mu$ L** *TF Cancer Stem Cell Probe Mix I*  
**X $\mu$ L** Nuclear Extract (5 $\mu$ g-15 $\mu$ g recommended)  
**Y $\mu$ L** ddH<sub>2</sub>O (add up to final volume)  
**30 $\mu$ L** **Reaction Mix** [final volume]
- 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 $\mu$ L** pre-chilled *Filter Binding Buffer*. Centrifuge the column with the collection tube at **6,000rpm** for **1 minute** in a microcentrifuge at room temperature.
- Transfer the **30 $\mu$ L** **Reaction Mix** directly onto the filter in the center of the *Isolation Column* (avoiding bubbles).
- Incubate on ice for **30 minutes**. **DO NOT** incubate longer than 30 minutes; this will result in high background.
- Add **500 $\mu$ L** 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 $\mu$ L** pre-chilled *Filter Wash Buffer* to the *Isolation Column* on ice.
- Centrifuge the *Isolation Column* with the collection tube for **1 minute** at **6,000rpm** in a microcentrifuge at **4°C**. Then discard the flow through.
- Repeat steps 8-9 for an additional **3 times** for a total a 4 washes.

## Elution of Bound Probe

- Add **75 $\mu$ L** of *Elution Buffer* onto the center of *Isolation Column*, and incubate at room temperature for **5 minutes**.
- Place the *Isolation Column* on a new 1.5mL microcentrifuge tube and centrifuge at **10,000 rpm** for **2 minutes** at room temperature.
- If you have yet to do so, chill **300 $\mu$ L** ddH<sub>2</sub>O (DNAase free) in a 1.5mL microcentrifuge tube on ice for at least **10 minutes**, and **keep on ice**.
- Transfer the eluted probe to a PCR tube and denature the eluted probes at **98°C** for **5 minutes**.
- Immediately** transfer the denatured probes to the chilled ddH<sub>2</sub>O 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



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

### 16. **Skip this step if you did not freeze your sample for future use.**

- Thaw your sample back to an aqueous phase at room temperature.
- Redistribute the sample into PCR tubes to be reheated at **98°C** for **5 minutes**.
- Afterwards, **immediately** place the PCR tubes on ice.
- You may now proceed to Step 17.

## Hybridization of Eluted Probe with Hybridization Plate

- Remove the clear adhesive film sealing from the provided *96-Well Plate*.
- Aliquot **3mL** pre-warmed *TF Plate Hybridization Buffer* to a dispensing reservoir (DNase free) and then add **300 $\mu$ L** denatured probes. Mix them together by gently shaking the reservoir.
- 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 4 sections of three columns. 1 sample is used per section (Columns 1→3 is one 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 $\mu$ L** *TF Plate Hybridization Buffer* only **without** the eluted probe.

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°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 **200µL** of pre-warmed *IX Plate Hybridization Wash Buffer* to each well by **row** with a **12 multi-channel 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 **200µL** 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.
27. If you have yet to do so: add **20µL** of *Streptavidin-HRP Conjugate* in **10mL Blocking Buffer** (1:500 dilution), enough for the whole plate (6 sections). This is the *diluted Streptavidin-HRP Conjugate*
28. Add **95µL** of *diluted Streptavidin-HRP Conjugate* to each well by **row** with a **12**

- multi-channel 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 **200µL IX 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 plate-shaker at room temperature. Decant the liquid from the wells.
31. 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:  
**1mL Substrate A**  
**1mL Substrate B**  
**8mL Substrate Dilution Buffer**  
**10mL Substrate Solution**
33. Add **95µL 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**.

#### Cancer Stem Cell TF Activation Profiling Array Diagram

	1	2	3	4	5	6	7	8	9	10	11	12
A	Androgen R	KLF4	Snail	Androgen	KLF4	Snail	Androgen	KLF4	Snail	Androgen	KLF4	Snail
B	AP1(Jun)	Myc	SOX2	AP1(Jun)	Myc	SOX2	AP1(Jun)	Myc	SOX2	AP1(Jun)	Myc	SOX2
C	AP2	Nanog	SOX9	AP2	Nanog	SOX9	AP2	Nanog	SOX9	AP2	Nanog	SOX9
D	CREB	NFkB1	STAT3	CREB	NFkB1	STAT3	CREB	NFkB1	STAT3	CREB	NFkB1	STAT3
E	ER	NKX3.1	TBX3	ER	NKX3.1	TBX3	ER	NKX3.1	TBX3	ER	NKX3.1	TBX3
F	FoxO3	Oct-3/4	Twist-1	FoxO3	Oct-3/4	Twist-1	FoxO3	Oct-3/4	Twist-1	FoxO3	Oct-3/4	Twist-1
G	GLI	p53	WT1	GLI	p53	WT1	GLI	p53	WT1	GLI	p53	WT1
H	HIF-1	PRDM14	Blank	HIF-1	PRDM14	Blank	HIF-1	PRDM14	Blank	HIF-1	PRDM14	Blank

#### Related Products

Catalog #	Product Description
<b>FA-1001</b>	TF Activation Profiling Plate Array I
<b>FA-1002</b>	TF Activation Profiling Plate Array II
<b>FA-1003</b>	Stem Cell TF Activation Profiling Plate Array
<b>FA-1005</b>	Oxidative Stress TF Activation Profiling Plate Array