



## FewCell ER Stress (UPR) TF Activation Profiling Plate Array

Catalog # FA-1106 (For Research Use Only)

### Introduction

The Unfolded Protein Response (UPR) is a conserved and essential stress response that cells activate to combat Endoplasmic Reticulum (ER) stress, commonly caused by the accumulation of misfolded proteins or failing protein quality control. ER stress is a well-characterized feature of several diseases, such as diabetes, Alzheimer's, Parkinson's, Huntington's, and prion diseases. Early cellular response to ER stress includes the transcriptional upregulation of chaperone proteins, which is mediated by a large number of transcription factors (TFs). **Signosis, Inc.** has developed the *ER Stress/UPR TF Activation Profiling Plate Array*, which can be used to simultaneously monitor 16 ER stress/UPR related TFs including XBP-1, ATF4, ATF6, GADD153/CHOP, CBF/NFY, SREBP1, YY1, PGC-1a, ATF3, AP-1, FOXO1, IRF, p53, NFkB, NRF2/ARE, and HNF4.

### 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 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 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

- PCR machine and PCR tubes
- Microcentrifuge working at 4 °C
- Hybridization incubator at 42°C
- Plate-Shaker
- Plate reader for luminescent detection
- ddH<sub>2</sub>O (DNase-free)
- 8 and 12 Multi-channel pipettes

### Materials Provided with the Kit

Component	Qty	Store at
<b>96-Well Plate (with aluminum adhesive seal)</b>	1	RT
<b>Isolation Columns</b>	6	RT
<b>Elution Buffer</b>	600µL	RT
<b>TF Plate Hybridization Buffer</b>	20mL	RT
<b>5X Plate Hybridization Wash Buffer</b>	30mL	RT
<b>5X Detection Wash Buffer</b>	40mL	RT
<b>Blocking Buffer</b>	30mL	4°C
<b>Filter Wash Buffer</b>	15mL	4°C
<b>Filter Binding Buffer</b>	1.5mL	4°C
<b>Substrate A</b>	1mL	4°C
<b>Substrate B</b>	1mL	4°C
<b>Streptavidin-HRP Conjugate</b>	20µL	4°C
<b>Substrate Dilution Buffer</b>	8mL	4°C
<b>TF Binding Buffer Mix</b>	90µL	-20°C
<b>TF ER Stress Probe Mix</b>	20µL	-20°C
<b>Whole cell lysis buffer</b>	800µL	-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 **200µL** of ddH<sub>2</sub>O in a 1.5mL microcentrifuge tube (per sample; 3 samples would be 600µL ddH<sub>2</sub>O) 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

## The whole cell lysate preparation

1. Wash cells (2000 cells-10,000 cells) with 1X PBS, and add 100ul Whole Cell Lysis Buffer and mix well by pipetting up and down a few times. And incubate for 10 minutes on ice with gently shaking.
2. Centrifuge briefly, then save 90ul of the supernatant. The cleared lysate is now ready for use.

## TF/ DNA Complex Formation

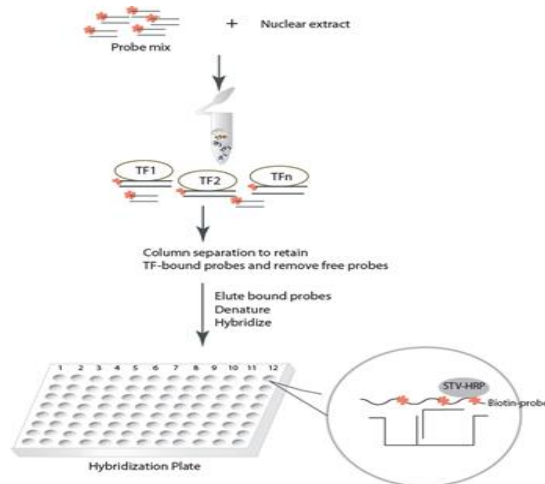
1. Mix the following components for each reaction in a tube  
**15µL TF Binding Buffer Mix**  
**3µL TF Probe mix**  
**XµL Cell Lysate** (10µg-15µg recommended)  
**YµL ddH<sub>2</sub>O** (add up to final volume)  
**30µL Reaction Mix**
2. Incubate the **Reaction Mix** at room temperature (20-23°C) for **30 minutes**.

## Separation of TF DNA Complex from Free Probes

3. Equilibrate an *Isolation Column* by adding **200µL** pre-chilled *Filter Binding Buffer*. Centrifuge the column with the collection tube at **6,000rpm** for **1 minute** in a microcentrifuge at room temperature.
4. Transfer the **30µL 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.
6. Add **500µL** pre-chilled *Filter Wash Buffer* to the *Isolation Column* and incubate for **3 minutes** on ice.
7. 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.
8. Wash the column by adding **500µL** pre-chilled *Filter Wash Buffer* to the *Isolation Column* on ice.
9. Centrifuge the *Isolation Column* with the collection tube for **1 minute** at **6,000rpm** in a microcentrifuge at **4°C**.
10. Repeat steps 8-9 for an additional **3 times** for a total a 4 washes.

## Elution of Bound Probe

11. Add **50µL** 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 1.5mL microcentrifuge tube and centrifuge at **10,000 rpm** for **2 minutes** at room temperature.



13. If you have yet to do so, chill **200µL** ddH<sub>2</sub>O (DNAase free) in a 1.5mL 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 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.**
  - 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 **2mL** pre-warmed *TF Plate Hybridization Buffer* to a dispensing reservoir (DNase free) and then add **200µL** denatured probes. Mix them together by gently shaking the reservoir.
19. Using a 8 multi-channel pipette **100µL** of the mixture from step 18. into the corresponding wells with 8 multi-channel pipette **immediately**.

**Note:** the *96-Well Plate* is divided into 6 sections of two columns each for 6 samples. Two columns are used 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 **100µL** *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 *1X 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.
25. 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*

26. 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.
27. 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.
28. Wash the *96-Well Plate* by adding **200µL** *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 plate-shaker at room temperature. At the last wash, invert plate on clean paper towels for **1 minute** to remove excessive liquid.
28. 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**
29. 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.
30. 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**.

#### Endoplasmic Reticulum Stress TF Activation Profiling Array Diagram

	1	2	3	4	5	6	7	8	9	10	11	12
A	XBP-1	ATF3	XBP-1	ATF3	XBP-1	ATF3	XBP-1	ATF3	XBP-1	ATF3	XBP-1	ATF3
B	ATF4	AP-1	ATF4	AP-1	ATF4	AP-1	ATF4	AP-1	ATF4	AP-1	ATF4	AP-1
C	ATF6	FOXO1	ATF6	FOXO1	ATF6	FOXO1	ATF6	FOXO1	ATF6	FOXO1	ATF6	FOXO1
D	GADD153	IRF	GADD153	IRF	GADD153	IRF	GADD153	IRF	GADD153	IRF	GADD153	IRF
E	CBF/NFY	P53	CBF/NFY	P53	CBF/NFY	P53	CBF/NFY	P53	CBF/NFY	P53	CBF/NFY	P53
F	SREBP1	NFkB	SREBP1	NFkB	SREBP1	NFkB	SREBP1	NFkB	SREBP1	NFkB	SREBP1	NFkB
G	YY1	NRF2/ARE	YY1	NRF2/ARE	YY1	NRF2/ARE	YY1	NRF2/ARE	YY1	NRF2/ARE	YY1	NRF2/ARE
H	ERR	HNF4	ERR	HNF4	ERR	HNF4	ERR	HNF4	ERR	HNF4	ERR	HNF4

#### 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-1004</b>	Cancer Stem Cell TF Activation Profiling Plate Array
<b>FA-1005</b>	Oxidative Stress TF Activation Profiling Plate Array