

**Description**

RAR $\alpha$  Luciferase Reporter HEK293 Cell Line are engineered HEK293 cells expressing a conditional firefly luciferase reporter under the control of retinoid acid response elements (RARE), and constitutively expressing full length human Retinoic Acid Receptor alpha (RAR $\alpha$ , NM\_0.00964). This cell line is functionally validated to respond to *all-trans* retinoic acid (ATRA) stimulation.

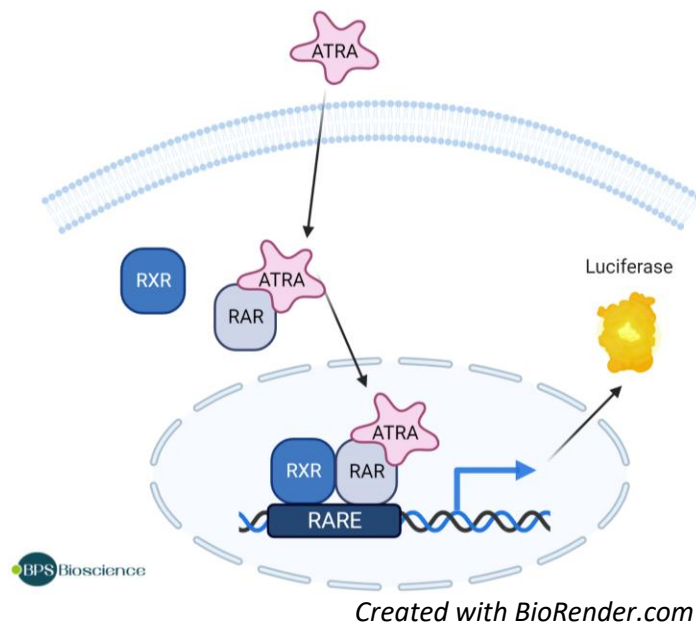


Figure 1: Activation of retinoid acid-dependent luciferase reporter in the RAR $\alpha$ /RAR $\alpha$  Luciferase Reporter HEK293 Cell Line.

**Background**

RAR (Retinoic Acid Receptor) belongs to a family of nuclear receptors and has three subtypes: RAR $\alpha$ , RAR $\beta$ , and RAR $\gamma$ . RAR heterodimerizes with RXR (Retinoid X Receptor) and acts as a transcription factor that regulates the growth and differentiation of both normal and malignant cells. When RAR binds to its ligands, *all-trans* retinoic acid (ATRA) or *9-cis* retinoic acid, the RAR/RXR heterodimer binds to the retinoic acid response elements in the promoter region of target genes. This recruits coactivator proteins, leading to transcription and expression of the downstream target genes. Retinoic acid (a derivative of Vitamin A) signaling plays critical roles in embryonic and immune system development. Mutations in RAR $\alpha$  can lead to Acute Promyelocytic Leukemia (APL), characterized by the accumulation of progenitor cells from the blood lineage (promyelocytes), and other cancer types. Use of RA may prove beneficial in cancer therapy.

**Application**

- Monitor RAR $\alpha$ -regulated pathway activity.
- Screen RAR $\alpha$  agonists and antagonists.

**Materials Provided**

Components	Format
2 vials of frozen cells	Each vial contains >1 x 10 <sup>6</sup> cells in 1 ml of Cell Freezing Medium (BPS Bioscience #79796)

**Parental Cell Line**

HEK293, Human Embryonic Kidney, epithelial-like cells, adherent.

**Mycoplasma Testing**

The cell line has been screened to confirm the absence of Mycoplasma species.

**Materials Required but Not Supplied**

These materials are not supplied with the cell line but are necessary for cell culture and cellular assays. BPS Bioscience's reagents are validated and optimized for use with this cell line and are highly recommended for best results. Media components are provided in the Media Formulations section below.

*Media Required for Cell Culture*

Name	Ordering Information
Thaw Medium 6	<a href="#">BPS Bioscience #60183</a>
Growth Medium 6A	<a href="#">BPS Bioscience #79542</a>

*Materials Required for Cellular Assay*

Name	Ordering Information
ATRA	Sigma #R2625
Assay Medium 6A	<a href="#">BPS Bioscience #82211</a>
96-well tissue culture treated white clear-bottom assay plate	Corning #3610
ONE-Step™ Luciferase Assay System	<a href="#">BPS Bioscience #60690</a>
Luminometer	

**Storage Conditions**

Cells are shipped in dry ice and should immediately be thawed or stored in liquid nitrogen upon receipt. Do not use a -80°C freezer for long term storage. Contact technical support at [support@bpsbioscience.com](mailto:support@bpsbioscience.com) if the cells are not frozen in dry ice upon arrival.

**Media Formulations**

For best results, the use of validated and optimized media from BPS Bioscience is *highly recommended*. Other preparations or formulations of media may result in suboptimal performance.



Note: Thaw Media do *not* contain selective antibiotics. However, Growth Media *do* contain selective antibiotics, which are used to maintain selective pressure on the cell population expressing the gene of interest.

Cells should be grown at 37°C with 5% CO<sub>2</sub>. BPS Bioscience's cell lines are stable for at least 10 passages when grown under proper conditions.

*Media Required for Cell Culture*

*Thaw Medium 6 (BPS Bioscience #60183):*

DMEM supplemented with 10% FBS and 1% Penicillin/Streptomycin.

*Growth Medium 6A (BPS Bioscience #79542):*

DMEM supplemented with 10% FBS, and 1% Penicillin/Streptomycin plus 400  $\mu\text{g/ml}$  of Geneticin (G418), 1  $\mu\text{g/ml}$  Puromycin and 100  $\mu\text{g/ml}$  Hygromycin.

*Media Required for Functional Cellular Assay*

*Assay Medium 6A (BPS Bioscience #82211):*

Phenol red-free DMEM, 10% Charcoal Stripped FBS, and 1% Penicillin/Streptomycin.

**Cell Culture Protocol**

**Note: HEK293 cells are derived from human material and thus the use of adequate safety precautions is recommended.**

*Cell Thawing*

1. Swirl the vial of frozen cells for approximately 60 seconds in a 37°C water bath. Once cells are thawed (it may be slightly faster or slower than 60 seconds), quickly transfer the entire contents of the vial to a tube containing 10 ml of pre-warmed Thaw Medium 6.  
**Leaving the cells in the water bath at 37°C for too long will result in rapid loss of viability.**
2. Immediately spin down the cells at 300 x *g* for 5 minutes, remove the medium and resuspend the cells in 5 ml of pre-warmed Thaw Medium 6.
3. Transfer the resuspended cells in Thaw Medium 6 to a T25 flask and incubate at 37°C in a 5% CO<sub>2</sub> incubator.
4. After 48-72 hours in culture, check for cell viability, change to fresh Thaw Medium 6, and continue growing in a 5% CO<sub>2</sub> incubator at 37°C until the cells are ready to passage.
5. Cells should be passaged before they reach 100% confluency. Switch to Growth Medium 6A for passage.

*Cell Passage*

1. Aspirate the medium, wash the cells with phosphate buffered saline (PBS) without Ca<sup>2+</sup>/Mg<sup>2+</sup>, and detach the cells from the culture vessel with 0.25% Trypsin/EDTA following volumes recommended for the cell vessel being used.
2. Once the cells have detached, add Growth Medium 6A and transfer to a tube.
3. Spin down cells at 300 x *g* for 5 minutes, remove the medium and resuspend the cells in Growth Medium 6A.
4. Seed into new culture vessels at the recommended sub-cultivation ratio of 1:10 to 1:20 twice a week.

*Cell Freezing*

1. After detachment, spin down the cells at 300 x *g* for 5 minutes.
2. Remove the medium and resuspend the cell pellet in 4°C Cell Freezing Medium (BPS Bioscience #79796) at a density of  $\sim 2 \times 10^6$  cells/ml.

3. Dispense 1 ml of cell suspension into each cryogenic vial. Place the vials in an insulated container for slow cooling and store at -80°C overnight.
4. Transfer the vials to liquid nitrogen the next day for long term storage.



Note: It is recommended to expand the cells and freeze at least 10 vials at an early passage for future use.

### Validation Data

#### Dose response of RAR $\alpha$ Reporter – HEK293 Recombinant Cell Line to *all-trans* retinoic acid (ATRA)

##### Assay Protocol

- The following assay is designed for 96-well format. To perform the assay in different tissue culture formats, cell number and reagent volume should be scaled appropriately.
- All samples and controls should be performed in triplicate.
- This assay should include “Cell-Free Control”, “Unstimulated Control” and “Test Compound” conditions.

##### Assay Medium:

##### Assay Medium 6A.

1. 24 hours before plating the cells remove Growth Medium 6A and replace with Assay Medium.
2. Dissociate the cells and seed ~30,000 cells/well in 90  $\mu$ l of Assay Medium in a 96-well white clear-bottom plate.
3. Add 90  $\mu$ l of Assay Medium to “Cell Free Control” wells (for determining background luminescence).
4. Prepare ATRA (10  $\mu$ l/ well): start by preparing a stock solution of ATRA in 100% DMSO at 2000-fold the highest desired concentration. The final volume is 100  $\mu$ l.

Dilute 200-fold in Assay Medium to prepare an intermediate 10x stock solution with 0.5% DMSO.

Prepare a serial dilution of ATRA in 0.5% DMSO in Assay Medium at concentrations 10-fold higher than the desired final concentrations. Prepare Assay Medium with 0.5% DMSO (Diluent Solution) to use in controls (Diluent Solution).

*Note: The final concentration of DMSO in the assay should not exceed 0.1%.*

5. Add 10  $\mu$ l of diluted ATRA to each well.
6. Add 10  $\mu$ l of Diluent Solution to the “Unstimulated Control” wells “Cell-Free Control” wells (background).

Component	Cell Free Control	Unstimulated Control	Test Compound
Assay Medium	90 $\mu$ l	-	-
Cell suspension	-	90 $\mu$ l	90 $\mu$ l
Diluted ATRA	-	-	10 $\mu$ l
Diluent Solution	10 $\mu$ l	10 $\mu$ l	-
<b>Total</b>	<b>100 <math>\mu</math>l</b>	<b>100 <math>\mu</math>l</b>	<b>100 <math>\mu</math>l</b>

- Incubate cells at 37°C in a CO<sub>2</sub> incubator for ~16 to 24 hours.
- Add 100  $\mu$ l of ONE-Step™ Luciferase reagent per well and gently agitate at Room Temperature for ~15 minutes.
- Measure the luminescence using a luminometer.
- Subtract average background luminescence (cell-free control wells) from the luminescence reading of all wells. Fold induction of RAR $\alpha$  luciferase reporter expression can be calculated by dividing the ATRA-stimulated Luminescence with the Unstimulated Control Luminescence.

$$\text{Fold induction} = \frac{\text{ATRA stimulated Lumin.} - \text{Background}}{\text{Unstimulated Lumin.} - \text{Background}}$$

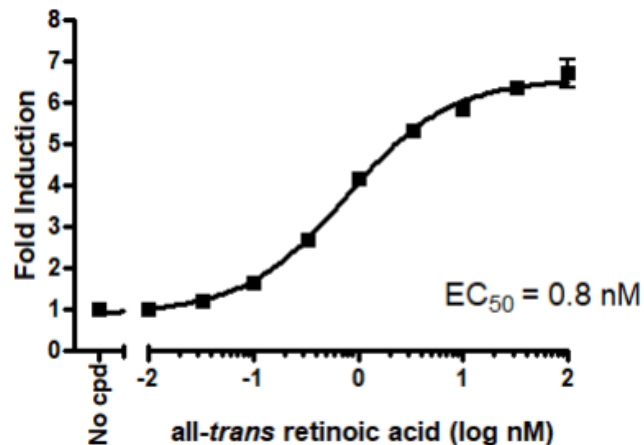


Figure 2: Dose response of RAR $\alpha$  Luciferase Reporter HEK293 Cell Line to all-trans retinoic acid (ATRA).

Cells were incubated with increasing amounts of ATRA. Luciferase activity was measured using the ONE-Step™ Luciferase Assay System. Results are expressed as Fold Induction (calculated against the “Unstimulated Cells Control”, which was set to 1).

Data shown is representative. For lot-specific information, please contact BPS Bioscience, Inc. at [support@bpsbioscience.com](mailto:support@bpsbioscience.com).

**References**

Petkovich M., *et al.*, 1987 *Nature*. 330 (6147):440-50.  
 Allenby J., *et al.*, 1993 *Proc Natl Acad Sci USA* 90 (1):30-4.  
 Peng Z., *et al.*, 2023 *J Immunotherapy Cancer* 11 (3): e006325.

**License Disclosure**

Visit [bpsbioscience.com/license](https://bpsbioscience.com/license) for the label license and other key information about this product.

**Troubleshooting Guide**

Visit [bpsbioscience.com/cell-line-faq](https://bpsbioscience.com/cell-line-faq) for detailed troubleshooting instructions. For all further questions, please email [support@bpsbioscience.com](mailto:support@bpsbioscience.com).

**Related Products**

<i>Products</i>	<i>Catalog #</i>	<i>Size</i>
RAR $\gamma$ Reporter - HEK293 Recombinant Cell Line	60604	2 vials
RAR $\beta$ Reporter – HEK293 Recombinant Cell Line	60603	2 vials
RAR $\alpha$ Reporter Cellular Assay Pack	79322	2 vials
RAR $\beta$ Reporter Cellular Assay Pack	79323	2 vials
RAR $\gamma$ Reporter Cellular Assay Pack	79324	2 vials

*Version 112723*