



6042 Cornerstone Court West, Suite B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

RAR γ Reporter (Luc) - HEK293 Cell Line **Catalog #: 60604**

Background

Retinoic acid receptor (RAR) belongs to the family of nuclear receptors and has three subtypes, RAR α , RAR β , and RAR γ . RAR heterodimerizes with retinoic X receptor (RXR) 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 or 9-*cis* retinoic acid, RAR/RXR heterodimer binds to retinoic acid response elements in the promoter region of target genes and recruits coactivator proteins, leading to transcription and expression of the downstream target genes.

Descriptions

The RAR gamma Reporter (Luc)-HEK293 cell line is designed for monitoring the activity of RAR γ . The RAR gamma Reporter (Luc)-HEK293 cell line contains a firefly luciferase gene under the control of retinoic acid response elements stably integrated into HEK293 cells along with full length human RAR γ (accession #P13631-1). This cell line is functionally validated for the response to the stimulation of all-*trans* retinoic acid. The expression of RAR γ is confirmed by western blotting.

Application

- Monitor RAR γ -regulated pathway activity
- Screen agonists or antagonists of RAR γ .

Format

Each vial contains $\sim 2 \times 10^6$ cells in 1 ml of 10% DMSO.

Mycoplasma testing

The cell line has been screened using the PCR-based Venor[®]GeM Mycoplasma Detection kit (Sigma-Aldrich) to confirm the absence of Mycoplasma species.

Storage

Immediately upon receipt, store in liquid nitrogen.

General Culture Conditions

Thaw Medium 6 (BPS Cat. #60183): DMEM medium (Hyclone #SH30243.01) supplemented with 10% FBS (Life technologies #26140-079), 1% Penicillin/Streptomycin (Hyclone SV30010.01).

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court West, Suite B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Complete Growth Medium: Thaw Medium 6 (BPS Cat. #60187) and 400 µg/ml of Geneticin (G418) (Invitrogen #11811031), 1 µg/ml of Puromycin (Hyclone #SV30075.01), and 100 µg/ml Hygromycin (Hyclone #SV30070.01).

Cells should be maintained at 37°C with 5% CO₂ using complete growth medium. If culturing cells in medium from other vendors, it may be necessary to lower the percentage of CO₂ in the incubator depending on the NaHCO₃ level in the basal medium.

To thaw the cells, it is recommended to quickly thaw the frozen cells from liquid nitrogen in a 37°C water-bath, transfer to a tube containing 10 ml of Thaw Medium 6 (**no Geneticin, Puromycin, and Hygromycin**), spin down cells, and resuspend cells in pre-warmed Thaw Medium 6 (**no Geneticin, Puromycin, and Hygromycin**). Transfer resuspended cells to a T25 flask and culture in a 37°C CO₂ incubator. At first passage, switch to complete growth medium (**Thaw Medium 6, Geneticin, Puromycin, and Hygromycin**). Cells should be split before they reach complete confluence.

To passage the cells, rinse cells with phosphate buffered saline (PBS), detach cells from culture vessel with Trypsin/EDTA, and add complete growth medium. Transfer to a tube, spin down cells, resuspend cells and seed appropriate aliquots of cell suspension into new culture vessels. Subcultivation ration: 1:10 to 1:20, twice a week.

To freeze down the cells, rinse cells with phosphate buffered saline (PBS), and detach cells from culture vessel with Trypsin/EDTA. Add complete growth medium and transfer to a tube, spin down cells, and resuspend in freezing medium (10% DMSO + 90% FBS). Place at -80°C overnight and place in liquid nitrogen the next day. Alternatively, vials may be placed directly in liquid nitrogen.

Functional Validation and Assay Performance

The following assays are designed for 96-well format. To perform assay in different tissue culture formats, cell number and reagent volume should be scaled appropriately.

Materials Required but Not Supplied

- all-*trans* retinoic acid (ATRA) (Sigma #R2625): make 1 mM stock solution in DMSO
- Assay medium: phenol red-free DMEM + 10% charcoal stripped FBS (Hyclone #SH3006802) + 1% Pen/Strep
- 96-well tissue culture treated white clear-bottom assay plate (Corning #3610)
- ONE-Step™ Luciferase Assay System (BPS, Cat. #60690)
- Luminometer

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court West, Suite B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

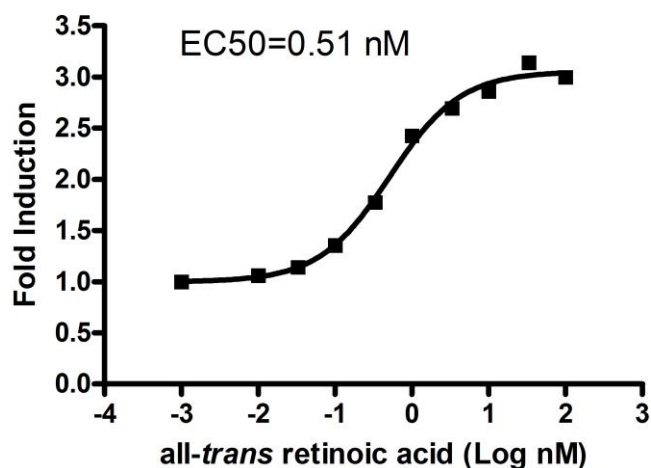
Assay protocol: Dose response of RAR gamma Reporter (Luc) - HEK293 cells to all-trans retinoic acid (ATRA)

1. One day before plating the cells, remove the growth medium from RAR gamma Reporter (Luc)-HEK293 cells and replace with assay medium for 24 hours.
2. Harvest RAR gamma Reporter (Luc)-HEK293 cells and seed cells in 40 μ l of assay medium at a density of \sim 30,000 cells per well into white clear-bottom 96-well microplate.
3. Prepare threefold serial dilution of ATRA in assay medium and add 10 μ l of ATRA solution to each well. The final DMSO concentration is 0.1%.
Add 10 μ l of assay medium with 0.5% DMSO to the unstimulated control wells.
Add 50 μ l of assay medium with 0.1% DMSO to cell-free control wells (for determining background luminescence).
Set up each treatment in at least triplicate.
4. Incubate cells at 37 $^{\circ}$ in a CO₂ incubator for \sim 16 to 24 hours.
5. Perform luciferase assay using ONE-Step™ Luciferase Assay System according to the protocol provided: Add 100 μ l of ONE-Step™ Luciferase reagent per well and rock at room temperature for \sim 10 minutes. Measure luminescence using a luminometer.
If using other luciferase reagents from other vendors, follow the manufacturer's assay protocol.
6. Data Analysis: Subtract average background luminescence (cell-free control wells) from luminescence reading of all wells.
The fold induction of RAR luciferase reporter expression = background-subtracted luminescence of ATRA-stimulated well / average background-subtracted luminescence of unstimulated control wells

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com

Figure 1. Dose response of RAR gamma Reporter (Luc) - HEK293 cells to all-trans retinoic acid (ATRA). Results are shown as fold induction of RAR luciferase reporter expression.



References

1. Petkovich, M, *et al. Nature* (1987) **330(6147)**: 444-450.
2. Allenby, G, *et al. Proc. Natl. Acad. Sci. USA* (1993) **90(1)**: 30-34.

License Disclosure

Purchase of this cell line grants you with a 10-year license to use this cell line in your immediate laboratory, for research use only. This license does not permit you to share, distribute, sell, sublicense, or otherwise make the cell line available for use to other laboratories, departments, research institutions, hospitals, universities, or biotech companies. The license does not permit use of this cell line in humans or for therapeutic or drug use. The license does not permit modification of the cell line in any way. Inappropriate use or distribution of this cell line will result in revocation of the license and result in an immediate cease of sales and distribution of BPS products to your laboratory. BPS does not warrant the suitability of the cell line for any particular use, and does not accept any liability in connection with the handling or use of the cell line. Modifications of this cell line, transfer to another facility, or commercial use of the cells may require a separate license and additional fees; contact sales@bpsbioscience.com for details. Publications using this cell line should reference BPS Bioscience, Inc., San Diego.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com