



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

Nrf2 Antioxidant Pathway **ARE Reporter – Hep G2 Cell line** **Catalog #: 60513**

Product Description

The Nrf2 antioxidant response pathway plays an important role in cellular antioxidant defense. Nrf2, a basic leucine zipper transcription factor, induces the expression of antioxidant and phase II enzymes by binding to the ARE (antioxidant response element) region of the gene promoter. Under basal conditions, Nrf2 is retained in the cytosol by binding to the cytoskeletal protein Keap1. Upon exposure to oxidative stress or other ARE activators, Nrf2 is released from Keap1 and translocates to the nucleus, where it can bind to the ARE, leading to the expression of antioxidant and phase II enzymes that protect the cell from oxidative damage.

The ARE Reporter – Hep G2 cell line contains a firefly luciferase gene under the control of ARE stably integrated into Hep G2 cells. This cell line is validated for the response to the stimulation of tert-butylhydroquinone and sulforaphane.

Application

- Monitor Nrf2 antioxidant response pathway activity.
- Screen for activators or inhibitors of Nrf2 antioxidant response pathway.

Format

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

Storage

Immediately upon receipt, store in liquid nitrogen.

General Culture Conditions

Thaw Medium 1 (BPS Cat. #60187): MEM medium (Hyclone #SH30024.01) supplemented with 10% FBS (Life Technologies #26140-079), 1% non-essential amino acids (Hyclone #SH30238.01), 1 mM Na-pyruvate (Hyclone #SH30239.01), 1% Penicillin/Streptomycin (Hyclone SV30010.01).

Growth Medium 1K (BPS Cat. #79533): Thaw Medium 1 (BPS Cat. #60187) and 600 $\mu\text{g/ml}$ of Geneticin (Life Technologies #11811031).

Cells should be grown at 37°C with 5% CO₂ using Growth Medium 1K (Thaw Medium 1 plus Geneticin). It may be necessary to adjust the percentage of CO₂ in the incubator depending on the NaHCO₃ level in the basal medium.

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

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 1 (**no geneticin**), spin down cells, and resuspend cells in pre-warmed Thaw Medium 1 (**no geneticin**). Transfer the resuspended cells to a T25 flask and incubate at 37°C in a 5% CO₂ incubator. At first passage, switch to Growth Medium 1K (**contains geneticin**). 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 0.05% Trypsin/EDTA, add Growth Medium 1K and transfer to a tube. Spin down cells, resuspend cells, and seed appropriate aliquots of cell suspension into new culture vessels. Subcultivation ration: 1:5 to 1:10 weekly.

To freeze down the cells, rinse cells with phosphate buffered saline (PBS), and detach cells from culture vessel with Trypsin/EDTA. Add Growth Medium 1K 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 the assay in different tissue culture formats, the cell number and reagent volumes should be scaled appropriately.

Materials Required but Not Supplied

- Tert-butylhydroquinone (Sigma # 112941). Prepare stock solution in DMSO.
- DL-Sulforaphane (Sigma # S4441). Prepare stock solution in DMSO.
- Assay medium: Thaw Medium 1 (BPS Cat. #60187)
- Growth Medium 1K (BPS Cat. #79533)
- 96-well tissue culture plate or 96-well tissue culture-treated white clear-bottom assay plate
- ONE-Step™ Luciferase Assay System (BPS Cat. #60690).
- Luminometer

Mycoplasma testing

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

A. Response of ARE Reporter - Hep G2 cells to antioxidant inducers.

1. Harvest ARE Reporter – Hep G2 cells from culture in Growth Medium 1K and seed cells at a density of ~ 40,000 cells per well into white clear-bottom 96-well microplate in 45 µl of assay medium.

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

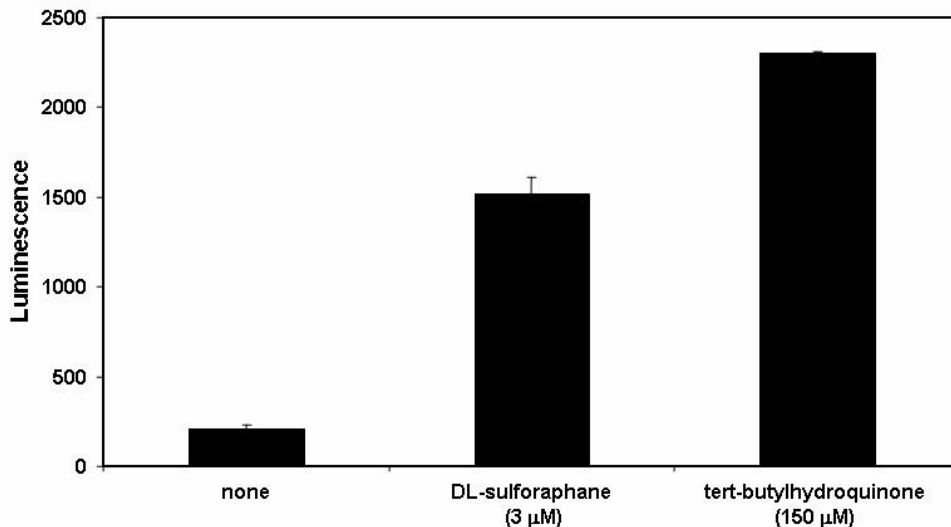
Please visit our website at: www.bpsbioscience.com



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

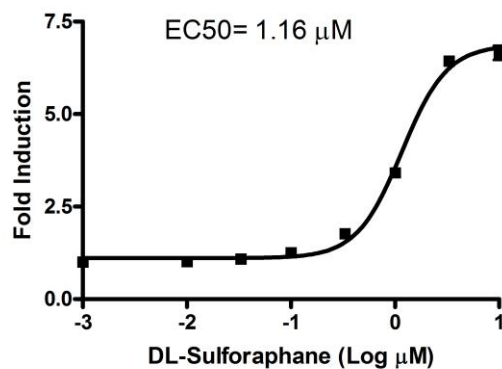
- Dilute antioxidant inducer (tert-butylhydroquinone or DL-Sulforaphane) into assay medium and add 5 μ l of dilution to each well. The final DMSO concentration can be up to 0.5%.
Add 5 μ l of assay medium with same concentration of DMSO but without the antioxidant inducer to the unstimulated control wells.
Add 50 μ l of assay medium with DMSO to cell-free control wells (for determining background luminescence).
Set up each treatment in at least triplicate.
- Incubate cells at 37°C in a CO₂ incubator for 15 to 18 hours.
- The next day, perform luciferase assay using the ONE-Step™ Luciferase Assay System: Add 100 μ l of ONE-Step™ Luciferase Assay reagent per well and rock at room temperature for ~15 minutes. Measure luminescence using a luminometer.
If using luciferase reagents from other vendors, follow the manufacturer's assay protocol.
- Data Analysis: Subtract the average background luminescence (cell-free control wells) from the luminescence reading of all wells.
The fold induction of ARE luciferase reporter expression = background-subtracted luminescence of stimulated well / average background-subtracted luminescence of unstimulated control wells.

Figure 1. ARE Reporter – Hep G2 cell response to antioxidant inducer.



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 2. Dose response of ARE Reporter – Hep G2 cells to DL-Sulforaphane. The results are shown as fold induction of ARE luciferase reporter expression.



References

1. Lee, JM *et. al.* (2004) An important role of Nrf2-ARE pathway in the cellular defense mechanism. *J Biochem Mol Biol.* **37(2)**: 139-143.
2. Dinkova-Kostova, AT *et.al.* (2002) Direct evidence that sulfhydryl groups of Keap1 are the sensors regulating induction of phase 2 enzymes that protect against carcinogens and oxidants. *Proc Natl Acad Sci U S A.* **99(18)**: 11908-11913.

Related Products

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
ERK Signaling Pathway SRE Reporter – HEK293 Cell Line	60406	1 vial
Hedgehog Pathway Gli Reporter – NIH3T3 Cell Line	60409	1 vial
JAK/STAT Signaling Pathway ISRE Reporter – HEK293 Cell Line	60510	1 vial
JNK Signaling Pathway AP1 Reporter – HEK293 Cell Line	60405	1 vial
NK-kB Reporter (Luc) – HEK293 Cell Line	60650	1 vial
RARalpha Reporter (Luc) – HEK293 Cell Line	60503	1 vial
Wnt Signaling Pathway TCF/LEF Reporter – HEK293 Cell Line	60501	1 vial
ONE-Step™ Luciferase Assay System	60690-1	10 ml
ONE-Step™ Luciferase Assay System	60690-2	100 ml

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

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