



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

Data Sheet

Empty vector control - HEK293 Recombinant Cell line Catalog #: 90334

Product description

Recombinant HEK293 cell transfected with empty expression vector containing the neomycin resistance gene. This cell line serves as a negative control for

- TMEM16A-HEK293 Cell Line Cat. #90230
- TRPC3-HEK293 Cell Line Cat. #90130
- TRPC7-HEK293 Cell Line Cat. #90030
- IKCA1 (KCNN4) - HEK293 Cell Line Cat. #90330
- ERG (Kv11.1) - HEK293 Cell Line Cat. #60619

Format

Each vial contains 1.5×10^6 cells in 1 ml of 10% DMSO.

Storage

Immediately upon receipt, store in liquid nitrogen.

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.

Culture conditions

Thaw Medium 1 (BPS Cat. #60187): MEM medium (Hyclone #SH30024.01) supplemented with 10% FBS (Invitrogen #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 1B (BPS Cat. #79531): Thaw Medium 1 (BPS Cat. #60187) plus 400 µg/ml of Geneticin (life Technologies #11811031) to ensure the recombinant expression is maintained.

Cells should be grown using at 37°C with 5% CO₂ using Growth Medium 1B. hERG-HEK293 cells should exhibit a typical cell division time of ~24 hours.

To thaw the cells, it is recommended to quickly thaw the frozen cells from liquid nitrogen in a 37°C water-bath and transfer to a tube containing 10 ml of Thaw Medium 1 (**no Geneticin**). Spin down cells, resuspend cells in pre-warmed Thaw Medium 1 (**no Geneticin**), and transfer resuspended cells to T25 flask. Culture at 37°C in a CO₂ incubator overnight. The next day, replace the medium with fresh Thaw Medium 1 (**no Geneticin**) and continue growing culture in

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 W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

a CO₂ incubator at 37°C until the cells are ready to be split. Cells should be split before they reach complete confluence. At first passage, switch to Growth Medium 1B (**contains Geneticin**).

To passage the cells, rinse cells with phosphate buffered saline (PBS) and detach cells from culture vessel with 0.05% Trypsin/EDTA. Add Growth Medium 1B and 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 weekly or twice a week.

To freeze cells, rinse cells with phosphate buffered saline (PBS) and detach cells from culture vessel with 0.05% Trypsin/EDTA. Add Growth Medium 1B and transfer to a tube, spin down cells and resuspend in freezing medium (90% FBS, 10% DMSO). Freeze cells at -80°C overnight and transfer to liquid nitrogen the next day.

License Disclosure

Visit bpsbioscience.com/license for the label license and other key information about this product.

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**](mailto:info@bpsbioscience.com)

Please visit our website at: [**www.bpsbioscience.com**](http://www.bpsbioscience.com)