



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 Empty vector control – CHO-K1 Recombinant Cell line Catalog #: 60545

Description

Recombinant CHO-K1 cell transfected with empty expression vector containing the hygromycin resistance gene.

Applications

This cell line serves as a negative control for

TCR Activator - CHO Recombinant Cell line	#60539
DAT CHOK1 Recombinant Cell Line	#60558
CD47 CHO-K1 Cell line	#60602
NET – CHOK1 Recombinant Cell Line	#60557

Format

Two vials containing ~ 2 x 10⁶ cells in 1 ml of 10% DMSO in FBS.

Storage

Store in liquid nitrogen immediately upon receipt. Do not store for long-term at -80°C or on dry ice.

Mycoplasma Testing

This cell line has been screened using the Venor™ GeM Mycoplasma Detection Kit, PCR Based (Sigma, #MP0025) to confirm the absence of Mycoplasma contamination.

Culture Medium:

Thaw Medium 3 (BPS Bioscience, #60186): Ham's F-12 medium (Hyclone # SH30526.01) supplemented with 10% FBS, 1% Penicillin/Streptomycin (Hyclone SV30010.01).

Growth Medium 3B (BPS Bioscience, #79529): Thaw Medium 3, 500 µg/ml Hygromycin B (Thermo Fisher, #10687010).

Recommended Culture Condition:

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 growth medium **without Hygromycin B**, spin down cells, resuspend cells in pre-warmed growth medium **without Hygromycin B**, transfer resuspended cells to T75 flask and culture in 5% CO₂ incubator at 37°C overnight. The next day, replace the medium with fresh growth medium **without Hygromycin B**, and continue growing culture in a 5% 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 **containing Hygromycin**

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

B.

To passage the cells, rinse cells with phosphate buffered saline (PBS), detach cells from culture vessel with 0.05% Trypsin/EDTA, add complete growth medium 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 down the cells, rinse cells with phosphate buffered saline (PBS), detach cells from culture vessel with 0.05% Trypsin/EDTA, add complete growth medium and transfer to a tube, spin down cells, and resuspend in 4°C Freezing Medium (10% DMSO + 90% FBS) at $\sim 2 \times 10^6$ cells/ml. Dispense 1 ml of cell aliquots into cryogenic vials. Place vials in an insulated container for slow cooling and store at -80°C overnight. Transfer to liquid nitrogen the next day for storage.

It is recommended to expand the cells and freeze down more than 10 vials of cells for future use at early passage.

Related Products

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
Thaw Medium 3	60186-1	100 ml
Growth Medium 3B	79529	500 ml
ADCC Bioassay Effector Cell, F variant (Low Affinity)	60540	2 vials
ADCC Bioassay Effector Cell, V variant (High Affinity)	60541	2 vials
TCR Activator - CHO Recombinant Cell line	60539	2 vials
DAT CHOK1 Recombinant Cell Line	60558	2 vials
CD47 CHO-K1 Cell line	60602	2 vials
NET CHOK1 Recombinant Cell Line	60557	2 vials

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