

**Description**

CHO-K1 cells stably expressing full-length human recombinant FcGR2B (Fc gamma receptor IIb, FcγRIIB). Cell surface expression of human FcGR2B (Genbank #NM\_004001.4) was confirmed by flow cytometry.

**Background**

FcGR2B (also known as CD32B), is a receptor for the Fc region of immunoglobulin G (IgG) and is known as an immune antibody checkpoint.

The two major forms of FcGR2B, FcGR2B1 and FcGR2B2, arise from either the inclusion or exclusion (respectively) of exon C1 via mRNA splicing, resulting in differing cell type-specific expression and function. Presence of the C1 sequence in isoform 1, which is highly expressed at the surface of B cells, tethers the receptor at the membrane and dramatically increases its half-life at the cell surface. Absence of C1 in isoform 2, expressed in myeloid cells, triggers rapid internalization of the receptor upon ligand binding. FcGR2B induces the phagocytosis of aggregated immunoglobulins. The receptor is also expressed in airways and in liver endothelial cells, where it may act as a “sink” for the removal of IgG immune complexes.

FcGR2B1 operates as a negative regulator of signals induced by antibodies bound to antigens at the surface of cells. Although it acts in concert with dozens of activating receptors, it is the only known negative regulator of B Cell Receptor (BCR)-induced activation of B cells. Thus, the biological function of FcGR2B1 is to tame the antibody-dependent inflammatory response and to clear the circulation of spent immune complexes. Defects in FcGR2B1 signaling lead to overt inflammation and are involved in autoimmune diseases.

In addition, FcGR2B interferes with the efficacy of therapeutic antibodies as it accelerates antibody depletion and decreases B cell responses and antibody production. On the other hand, FcGR2B contributes to the anti-tumor response to antibody checkpoint therapy by boosting CD+ T cells through cross-linking of antibodies directed at stimulatory checkpoints expressed on immune cells such as 4-1BB, OX40, CD40 and GITR.

Therefore, FcGR2B is an important immunotherapy target, both as a direct target for the treatment of B-cell malignancies and in combination with clinically relevant therapeutic monoclonal antibodies to overcome FcGR2B - mediated resistance.

**Application(s)**

- Screen for regulators of antibody-mediated signaling
- Characterize the agonist activity of checkpoint antibodies using FcGR2B receptor-mediated crosslinking

**Materials Provided**

| Components              | Format   |
|-------------------------|--|
| 2 vials of frozen cells | Each vial contains $2 \times 10^6$ cells in 1mL of 10% DMSO in FBS |

**Parental Cell Line**

CHO-K1 cells, Chinese Hamster Ovary, 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 3    | <a href="#">BPS Bioscience #60186</a> |
| Growth Medium 3D | <a href="#">BPS Bioscience #79539</a> |

*Materials Required for Cellular Assay*

| Name   | Ordering Information                  |
|--|---------------------------------------|
| Thaw Medium 1  | <a href="#">BPS Bioscience #60187</a> |
| Assay Medium: Thaw Medium 3                              | <a href="#">BPS Bioscience #60186</a> |
| CD137/NF-κB HEK293 cells                                 | <a href="#">BPS Bioscience #79289</a> |
| Anti-CD137 recombinant human antibody                    | <a href="#">BPS Bioscience #79097</a> |
| ONE-Step™ Luciferase Assay System                        | <a href="#">BPS Bioscience #60690</a> |
| 96 well tissue culture treated white, clear-bottom plate |                                       |
| 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, it is *highly recommended* to use these validated and optimized media from BPS Bioscience. 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 for maintaining the presence of the transfected gene(s) over passages.

Cells should be grown at 37°C with 5% CO<sub>2</sub>. BPS Bioscience's cell lines are stable for at least 15 passages when grown under proper conditions.

*Media Required for Cell Culture*

*Thaw Medium 3 (BPS Bioscience, #60186):*

F-12K medium supplemented with 10% FBS, 1% Penicillin/Streptomycin.

*Growth Medium 3D (BPS Bioscience, #79539):*

F-12K medium supplemented with 10% FBS, 1% Penicillin/Streptomycin plus 1000 µg/ml of Geneticin

### Media Required for Functional Cellular Assay

#### Thaw Medium 1 (BPS Bioscience, #60187):

MEM medium supplemented with 10% FBS, 1% non-essential amino acids, 1 mM Na pyruvate, 1% Penicillin/Streptomycin

#### Thaw Medium 3 (BPS Bioscience, #60186):

F-12K medium supplemented with 10% FBS, 1% Penicillin/Streptomycin

### Cell Culture Protocol

#### Cell Thawing

1. Swirl the vial of frozen cells for approximately 60 seconds in a 37°C water bath. As soon as the 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 3 (**no Geneticin**).

**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 3 (**no Geneticin**).
3. Transfer the resuspended cells to a T25 flask or T75 flask and incubate at 37°C in a 5% CO<sub>2</sub> incubator.
4. After 24 hours of culture, check for cell attachment and viability. Change medium to fresh Thaw Medium 3 (**no Geneticin**), 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 are fully confluent. At first passage and subsequent passages, use Growth Medium 3D (**contains Geneticin**).

#### Cell Passage

1. Aspirate the medium, wash the cells with phosphate buffered saline (PBS), and detach the cells from the culture vessel with 0.25% Trypsin/EDTA.
2. Once the cells have detached, add Growth Medium 3D (**contains Geneticin**) and transfer to a tube. Spin down cells at 300 x g for 5 minutes, remove the medium and resuspend the cells in Growth Medium 3D (**contains Geneticin**). Seed into new culture vessels at the desired sub-cultivation ratio of 1:2 to 1:10 weekly or twice per week.

#### Cell Freezing

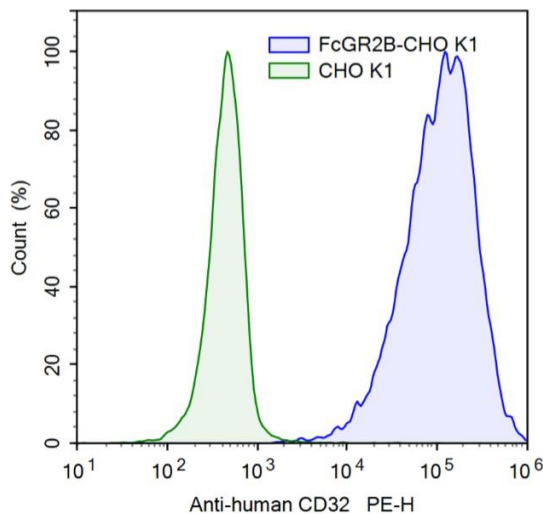
1. Aspirate the medium, wash the cells with phosphate buffered saline (PBS), and detach the cells from the culture vessel with 0.25% Trypsin/EDTA.
2. Once the cells have detached, add Growth Medium 3D and count the cells.
3. Spin down the cells at 300 x g for 5 minutes, remove the medium and resuspend the cells in 4°C Freezing Medium (BPS Bioscience #79796, or 10% DMSO + 90% FBS) at ~2 x 10<sup>6</sup> cells/ml.
4. Dispense 1 ml of cell aliquots into cryogenic vials. Place the vials in an insulated container for slow cooling and store at -80°C overnight.

- Transfer the vials to liquid nitrogen the next day for storage.



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

### A. Validation Data



*Figure 1: Flow cytometry analysis of cell surface expression of FcGR2B in CHO K1 cells. FcGR2B CHO cells (blue) or parental CHO cells (green) were stained with PE-labeled anti-human CD32 antibody (Biolegend, #303206) and analyzed by flow cytometry. Y-axis is the % cell number. X-axis is the intensity of PE.*

### B. Cellular Assay Protocol

Set up each condition at least in triplicate.

- Seed CD137/NF- $\kappa$ B reporter HEK293 cells at a density of  $\sim$ 30,000 cells per well into a white clear-bottom 96-well plate in 100  $\mu$ l of Thaw Medium 1. Leave a couple of wells empty for use as cell-free controls.
- Incubate the plate at 37°C in a CO<sub>2</sub> incubator overnight. Remove 60  $\mu$ l of Thaw Medium 1 from each well.
- 24 hours after seeding, harvest the FcGR2B CHO cells in Thaw Medium 3. Add 50  $\mu$ l of Thaw medium 3 containing  $\sim$ 90,000 FcGR2B-CHO K1 cells to each well of the CD137/NF- $\kappa$ B HEK293 cells.
- Prepare serial dilutions of anti-CD137 antibody in Thaw Medium 1 at concentration 10-fold higher than the desired final concentrations (you will need 10  $\mu$ l/well)
  - Add 10  $\mu$ l of diluted anti-CD137 antibody to the treated wells.
  - Add 10  $\mu$ l of Thaw Medium 1 to control (untreated) wells.
  - Add 50  $\mu$ l of Thaw Medium 1 and 50  $\mu$ l Thaw Medium 3 to cell-free control wells (for determining background luminescence)
- Incubate the plate at 37°C in a CO<sub>2</sub> incubator for  $\sim$ 18 hours.

6. Perform the luciferase assay using the ONE-STEP luciferase assay system:
  - a. Add 100  $\mu$ l of One-Step Luciferase reagent per well and rock at room temperature for ~30 minutes.
  - b. Measure luminescence using a luminometer.
7. Data Analysis: Subtract the average background luminescence (cell-free control wells) from the luminescence reading of all wells. The fold induction of NF- $\kappa$ B luciferase reporter expression is the background-subtracted luminescence of stimulated well divided by the average background-subtracted luminescence of unstimulated control wells.

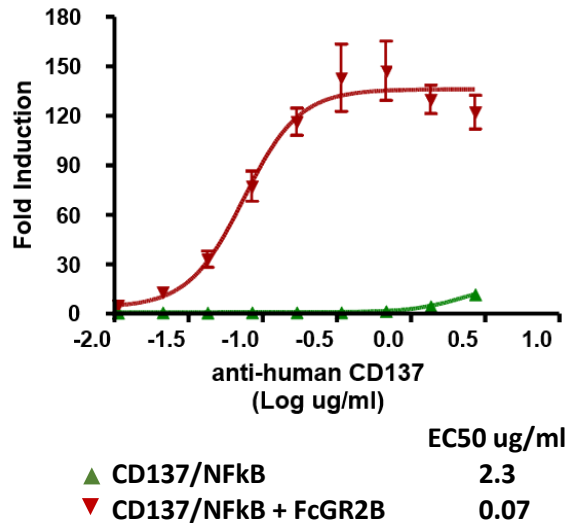


Figure 2: Dose response of anti-CD137 antibody in CD137/NF- $\kappa$ B-reporter HEK293 cells co-cultured with FcGR2B CHO cells. Cross-linking of the anti-CD137 antibody by FcGR2B expressed at the surface of CHO cells allowed activation of NF- $\kappa$ B in CD137-expressing HEK293 cells, indicating that FcGR2B potentiates the efficacy of antibodies directed at checkpoint activator CD137.

### Sequence

Human FcGR2B (GenBank: NM\_004001.4)

MGILSFLPVLATESDWADCKSPQPWGHMLLWTAVLFLAPVAGTPAAPPKAVLKLQPQWINVLQEDSVTLTCRGTHSPESDSIQWF  
 HNGNLIPTHTQPSYRFKANNNDGSEYTCQTGQTSLSDPVHLTVLSEWLVLQTPHLEFQEGETIVLRCHSWKDKPLVKVTFQNGKSK  
 KFSRSDPNFSIPQANHSHSGDYHCTGNIGYTYSSKPVTTITVQAPSSSPMGIIVAVVTGIAVAAIVAAVVALIYCRKKRISALPGYPECRE  
 MGETLPEKPANPTNPDEAD KVGAEANTITYSLLMHPDALEEDDQNR

### References

1. Anania JC, *et al.*, The Human Fc $\gamma$ RII (CD32) Family of Leukocyte FcR in Health and Disease. *Front Immunol.* (2019) **10**: 464.
2. Teige I, *et al.*, Targeting the Antibody Checkpoints to Enhance Cancer Immunotherapy-Focus on Fc $\gamma$ RIIB. *Front Immunol.* (2019) **10**: 481.

### 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> |
|---|------------------|-------------|
| TCR Activator-FcGR2B CHO Cell Line              | 78436            | 2 vials     |
| FcGR2B, Avi-His-Tag Recombinant                 | 100089           | 100 ug      |
| FcGR2B, Avi-His-Tag, Biotin-Labeled Recombinant | 100474           | various     |
| FcGR2B (CD32B) Lentivirus                       | 79877            | 500 µl x2   |