

## Description

Recombinant clonal stable CHO cell line constitutively expressing full length human CD38 protein (also known as ADPRC1, Genbank accession #NM\_001775) and human CD19 protein (also known as B4 or CVID3, Genbank accession #NM\_001770). This cell line was derived from our CHO-K1 Luciferase cells (BPS Bioscience, #79725), therefore it also constitutively expresses the firefly luciferase reporter. Surface expression of CD38 and CD19 were confirmed by flow cytometry.

## Background

The CD38 protein is a dimeric, non-lineage-restricted, type II transmembrane glycoprotein that synthesizes and hydrolyzes the second messengers cyclic ADP-ribose and NADP. CD38 is highly expressed by lymphoid and myeloid cells, particularly plasma cells. Increased CD38 expression on chronic lymphocytic leukemia (CLL) cells is linked to aggressive disease features and poor clinical outcome. CD38 is used as a prognostic marker for patients with CLL and multiple myeloma (MM), and is an ideal target for immunotherapy in CLL and MM.

B-lymphocyte antigen CD19 (Cluster of Differentiation 19), also known as B-Lymphocyte Surface Antigen B4 and CVID3, is a transmembrane protein expressed in follicular dendritic cells and all B lineage cells except plasma cells. CD19 plays two major roles in human B cells. It acts as an adaptor protein to recruit cytoplasmic signaling proteins to the membrane and it works within the CD19/CD21 complex to decrease the threshold for B cell receptor signaling pathways. Due to its presence on all B cells, it is a biomarker for B lymphocyte development and lymphoma diagnosis and can be used as a target for leukemia immunotherapies. CD19-targeted therapies based on T cells that express CD19-specific chimeric antigen receptors (CARs) have been utilized for their antitumor abilities in patients with CD19+ lymphoma and leukemia, such as Non-Hodgkin's Lymphoma (NHL), CLL and ALL.

## Application

1. Useful for validation of anti-CD38 and anti-CD19 bispecific antibody.
2. Useful as CD38- and/or CD19-expressing target cells in co-culture assay with CD38- and/or CD19-CAR-T cells, for both CD38/CD19-specific cell killing assay and cytokine production assay.
3. Useful for screening and validating antibodies against CD38 or CD19 and anti CD38 or anti-CD19 CAR-T for immunotherapy research and drug discovery.

## Materials Provided

Components	Format
2 vials of frozen cells	2 x 10 <sup>6</sup> cells in 1 ml of 10% DMSO

## Host Cell

CHO K1 cell line, Chinese Hamster Ovary, epithelial-like cells, adherent

## Mycoplasma Testing

The cell line has been screened using the MycoAlert™ Mycoplasma Detection kit (Lonza, #LT07-218) to confirm the absence of Mycoplasma species.

## Materials Required but Not Supplied



These materials are not supplied with this cell line but are necessary for cell culture and cellular assays. BPS Bioscience reagents systems 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.

*Materials Required for Cell Culture*

Name	Ordering Information
Thaw Medium 3	<a href="#">BPS Bioscience #60186</a>
Growth Medium 3K	<a href="#">BPS Bioscience #78041</a>

*Materials Required but Not Supplied*

Name	Ordering Information
ONE-Step™ Luciferase Assay System	<a href="#">BPS Bioscience #60690</a>
96-well tissue culture-treated white clear-bottom assay plate	
Luminometer	

**Storage Conditions**

Cells will arrive upon 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. To formulate a comparable but not BPS validated media, formulation components can be found below.



Note: Thaw Media does *not* contain selective antibiotics. However, Growth Media *does* contain selective antibiotics, which are used for maintaining cell lines over many passages. Cells should be grown at 37°C with 5% CO<sub>2</sub> using Growth Medium 3K.

*Media Required for Cell Culture*

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

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

*Growth Medium 3K (BPS Bioscience #78041):*

F-12K Medium supplemented with 10% FBS, 1% Penicillin/Streptomycin plus 1000 µg/ml Geneticin, 5 µg/ml Puromycin and 500 µg/ml of Hygromycin B to ensure recombinant expression.

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

**Cell Culture Protocol***Cell Thawing*

1. To thaw the cells, it is recommended to quickly thaw the frozen cells from liquid nitrogen in a 37°C water-bath, then transfer the entire contents of the vial to a tube containing 10 ml of Thaw Medium 3 (**no Geneticin, Puromycin or Hygromycin B**).
2. Spin down the cells, remove supernatant and resuspend cells in 5 ml of pre-warmed Thaw Medium 3 (**no Geneticin, Puromycin or Hygromycin B**).
3. Transfer the resuspended cells to a T25 flask and incubate at 37°C in a 5% CO<sub>2</sub> incubator.
4. After 24 hours of culture, add an additional ~3 ml of Thaw Medium 3 (**no Geneticin, Puromycin or Hygromycin B**) and continue growing culture in a CO<sub>2</sub> incubator at 37°C until the cells are ready to be split.

- Cells should be split before they are fully confluent. At first passage, switch to Growth Medium 3K (**contains Geneticin, Puromycin and Hygromycin B**).

#### Cell Passage

- To passage the cells, remove the medium, rinse cells with phosphate buffered saline (PBS), and detach cells from culture vessel with 0.25% Trypsin/EDTA.
- After detachment, add Growth Medium 3K (**contains Geneticin, Puromycin and Hygromycin B**) and transfer to a tube, spin down cells, resuspend cells in Growth Medium 3K and seed appropriate aliquots of cell suspension into new culture vessels. Sub cultivation ration: about 1:20 every 5 days.

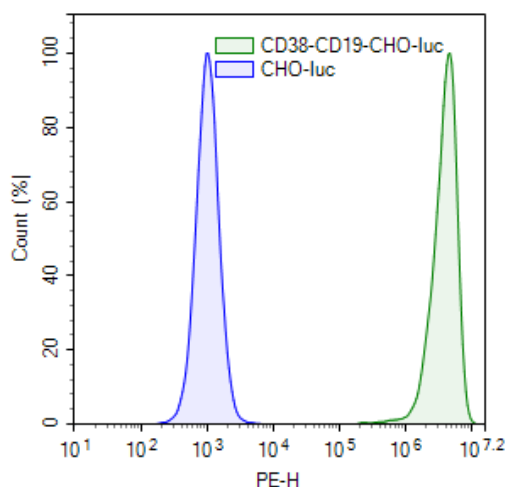
#### Cell Freezing

- To freeze down the cells, remove the medium, rinse cells with phosphate buffered saline (PBS), and detach cells from culture vessel with 0.25% Trypsin/EDTA.
- After detachment, add Thaw Medium 3 (**no Geneticin, Puromycin or Hygromycin B**) and count the cells, then transfer to a tube, spin down cells, and resuspend in 4°C Freezing Medium (BPS Bioscience, #79796) 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.



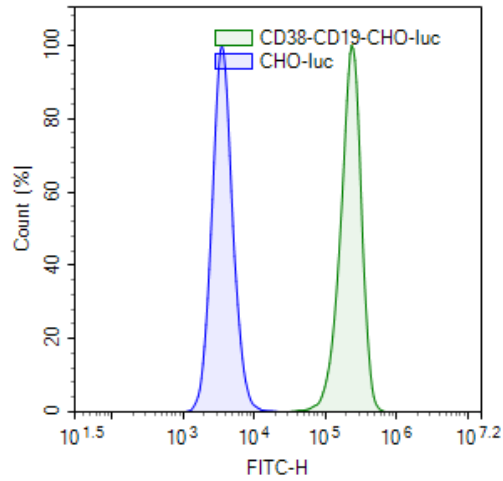
Note: It is recommended to expand the cells and freeze down at least 10 vials of cells at an early passage for future use.

#### Validation Data



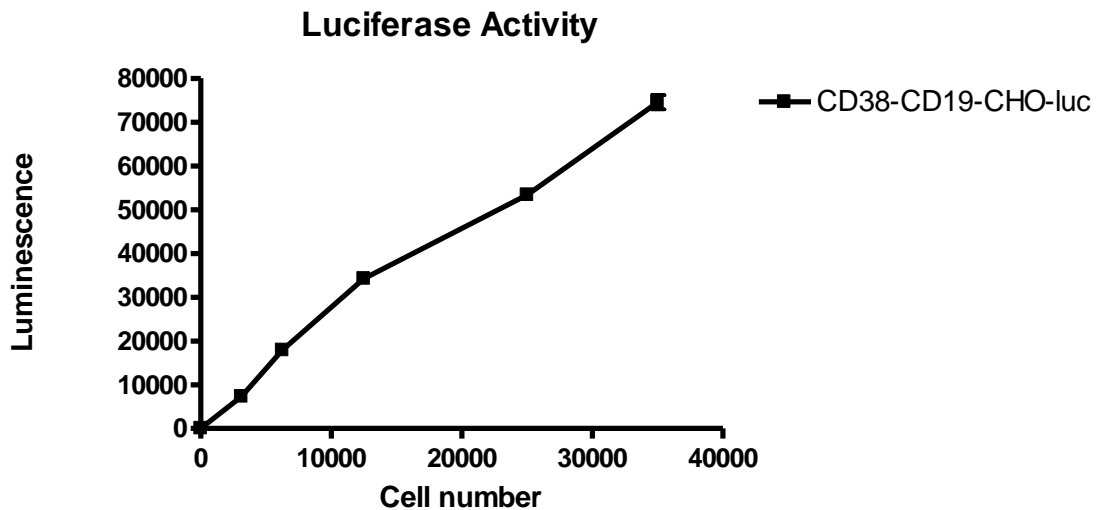
#### Figure 1. Expression of CD38 validated by flow cytometry.

Flow cytometry using PE-conjugated anti-human CD38 antibody (BioLegend, #303506) to detect CD38 surface expression on either the CD38 / CD19 / Firefly Luciferase - CHO Recombinant Cell Line (green) or parental CHO-luc cells (blue).



**Figure 2. Expression of CD19 validated by flow cytometry.**

Flow cytometry using FITC-conjugated anti-human CD19 antibody (BioLegend, #392508) to detect CD19 surface expression on either the CD38 / CD19 / Firefly Luciferase - CHO Recombinant Cell Line (green) or parental CHO-luc cells (blue).



**Figure 3. Luciferase activity of CD38 / CD19 / Firefly Luciferase - CHO Recombinant Cells.**

CD38 / CD19 / Firefly Luciferase - CHO Recombinant Cells were seeded in a 96-well plate at various densities. After four hours, luciferase activity under CMV promoter was measured using the ONE-Step luciferase assay system (BPS Bioscience, #60690).

**Sequence**

Human CD38 sequence (accession number NM\_001775)

MANCEFSPVSGDKPCCRLSRAQLCLGVSILVLILVVVLAVVVPRWRQQWSGPGTTKRFPETVLARCVMKYTEIHPMRHVDCQS  
VWDAFKGAFISKHPCNITEEDYQPLMKLGTQTVPCNKILLWSRIKDLAHQFTQVQRDMFTLEDTLGLYLADDLTWCGEFNTSKIN  
YQSCPDWRKDCSNNPVSVFWKTVSRRFAEAACDVVHVMLNGSRSKIFDKNSTFGSVEVHNLQPEKVQTLAWVIHGGREDSR  
DLCQDPTIKELESIISKRNQIFSCKNIRPDKFLQCVKNPEDSSCTSEI

Human CD19 sequence (accession number NM\_001770)

MPPPRLFFLLFLTPMEVRPEEPLVVKVEEGDNAVLQCLKGTS DGPTQQLTWSRESPLKPFLLKLSGLPGLGIHMRPLAIWLFIFNV  
SQQMGGFYLCQPGPPSEKAWQPGWTVNVEGSGELFRWNVSDLGGLGCGLNRSSEGPSSPSGKLMSPKLYVWAKDRPEIWE  
GEPCLPPRDSLNSQLSQDLTMAPGSTLWLSGCVPPDSVSRGPLSWTHVHPKPGPKSLLSLELKD RPARDMWVVMETGLLLPRA  
TAQDAGKYYCHRGNTMSFHLAITARPVLWHWLLRTGGWKVSAVTLAYLIFCLCSLVGILHLQRALVLRKRKRMTDPTRRRFFKV  
TPPPGSGPQNQYGNVLSLPTPTSGLGRAQRWAAGLGGTAPSYGNPSSDVQADGALGSRSPPGVGPEEEEEGEGYEEPDSEEDSE  
FYENDSNLGDQSLQDGSQGYENPEDEPLGPEDEDSFSNAESYENEDELTQPVARTMDFLSPHGS AWDPSREATSLGSQSYED  
MRGILYAAPQ LRSIRGQPGPNHEEDADSYENMDNPDGPDPAWGGGGRMGTWSTR

**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>
CD19 CHO Recombinant Cell Line (Various Expression)	79561	2 vials
CD19/Luciferase-CHO Recombinant Cell line	79714	2 vials
CD38-CHO Recombinant Cell Line (Various Expression)	79615	2 vials
Human CD19, Fc-Fusion, Avi-Tag HiP™	79472	50 µg
Human CD19, Fc-fusion (IgG1), Avi-Tag, Biotin-Labeled HiP™	79475	50 µg
Human CD19, Fc-Avi-Tag, PE-labeled	100732	50 µg
Human CD38, Avi-His-Tag	100346	100 µg
Human CD38, His-Tag (Human), HiP™	71277	100ug
Human CD38, His-Tag, APC-labeled	71883	100 µg
Human CD38, His-Tag, PE-labeled	71882	50 µg
CD38 Inhibitor Screening Assay Kit (Hydrolase Activity)	79287	96 reactions
CD38 Inhibitor Screening Assay Kit (Cyclase Activity)	71275	96 reactions