

Description

Recombinant clonal CHO-K1 stable cell line constitutively expressing full length human CD2 protein. The surface expression of CD2 in this cell line was validated by flow cytometry.

Background

CD2 (Cluster of differentiation 2, also known as T11, SRBC, or LFA-2) is a cell adhesion molecule and surface antigen found on the surface of all peripheral blood T-cells and natural killer (NK) cells. It is used as specific marker of T and NK cells. CD2 interacts with LFA3 (CD58) on antigen-presenting cells to optimize immune recognition. It also functions as a co-stimulator in T and NK cells. It is being investigated as a therapeutic target for the treatment of T cell lymphomas and leukemias.

Application

Screening or performing binding assays of antibodies against human CD2 in a cellular context.

Materials Provided

Components	Format
2 vials of frozen cells	Each vial contains 2×10^6 cells in 1 ml of 10% DMSO

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	BPS Bioscience #60186
Growth Medium 3J	BPS Bioscience #79974

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 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_2 . 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 (Kaighn's Modification of Ham's F-12 Medium) supplemented with 10% FBS, 1% Penicillin/Streptomycin

Growth Medium 3J (BPS Bioscience #79974):

F-12K Medium (Kaighn's Modification of Ham's F-12 Medium) supplemented with 10% FBS, 1% Penicillin/Streptomycin plus 5 µg/ml of Puromycin

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 Puromycin**).
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 Puromycin**).
3. Transfer the resuspended cells to a T25 flask or T75 flask and incubate at 37°C in a 5% CO₂ incubator.
4. After 24 hours of culture, check for cell attachment and viability. Change medium to fresh Thaw Medium 3 (**no Puromycin**), and continue growing in a 5% CO₂ 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 3J (**contains Puromycin**).

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 3J 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 3J (**contains Puromycin**). Seed into new culture vessels at the desired sub-cultivation ratio of 1:10 every 3 days.

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 3J 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 10⁶ 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.
5. 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.

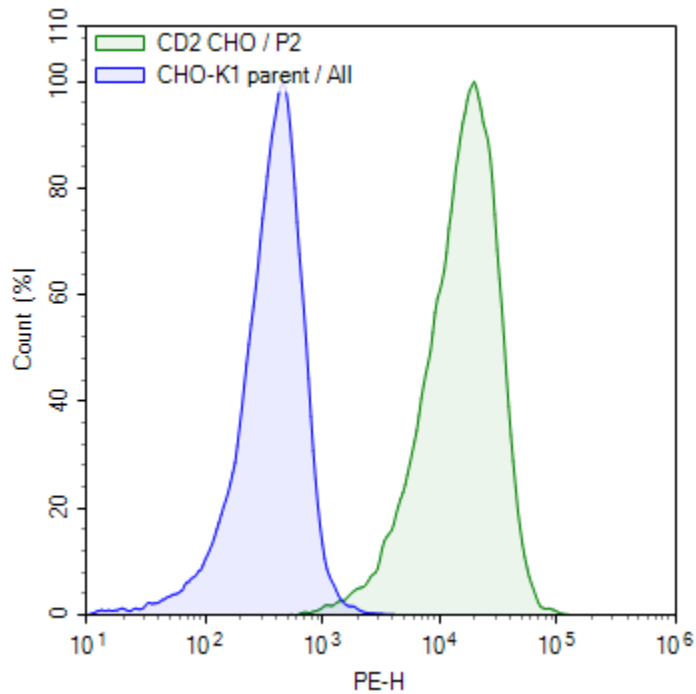
A. Validation Data

Figure 1: Confirmation of CD2 expression on CHO-K1 recombinant cells. Flow cytometry using PE-conjugated anti-human CD2 antibody (BioLegend, #300208) to detect CD2 surface expression on either the CD2 CHO-K1 Recombinant Cell Line (green) or parental CHO-K1 cells (blue).

Sequence

Human CD2 Sequence (Accession Number: NM_001767)

MSFPCKFVASFLIFNVSSKGAVSKEITNALETWGALGQDINLDIPSFQMSDDIDDIKWEKTSKDKKIAQFRKEKETFKEDTYKLFK
 NGTLKIKHLKTDDQDIYKVSIDYDTGKKNVLEKIFDLKIQERVSKPKISWTCINTTLTCEVMNGTDPELNLYQDGKHLKLSQRVITHK
 WTTLSAKFKCTAGNKVSKESSVEPVSCPEKGLDIYLIIGICGGGSLLMVVFALLVFYITKRKKQRSRRNDEELETRAHRVATEERGR
 KPHQIPASTPQNPATSQHPPPPGHRSQAPSHRPPPPGHRVQHQPQKRPPAPSGTQVHQQKGPPLPRPRVQPKPPHGAAENS
 LPSSN

License Disclosure

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

Troubleshooting Guide

Visit bpsbioscience.com/cell-line-faq for detailed troubleshooting instructions. For all further questions, please email support@bpsbioscience.com.

Related Products

<i>Products</i>	<i>Catalog #</i>	<i>Size</i>
CD2, Fc fusion, Avi-Tag, (Human) HiP™	71264	100 µg
CD2, Fc fusion, Biotin-labeled (Human) HiP™	71270	25 µg/50 µg
CD2, Avi-His-Tag	101238	100 µg
CD2, Avi-His-Tag, Biotin-Labeled	101239	25 µg/100 µg
CD58, Fc fusion, Biotin-labeled HiP™	71269	50 µg