

(Medium Expression)

Description

Recombinant CD5-CHO K1 cell line stably expressing a full-length human CD5 receptor (accession number: NM_014207.3). Surface expression of hCD5 was confirmed by flow cytometry. Each stable clonal cell line was selected for different levels of hCD5 expression (High, Medium and Low) to mimic different stages of cancer target cells with various CD5 expression levels.

Background

CD5 is a type-1 transmembrane glycoprotein with an extracellular region composed of three scavenger receptor, cysteine-rich domains. This lymphocytic-specific scavenger receptor functions at the interphase of the innate and adaptive immune responses. It is involved in: (i) microbial-associated pattern recognition; and (ii) modulation of intracellular signals mediated by antigen-specific receptor present in T and B cells. CD5 is expressed by all T cells and some B cell subsets (B1a and Breg or B10). T cells usually express higher levels of CD5 than B cells, and CD5 is upregulated in T cells upon strong activation. CD5 is a signaling co-receptor, associated with the antigen-specific receptor complex of T and B cells, and down-modulates the activation/differentiation signals delivered upon specific antigen recognition. The signaling pathway used by CD5 is only partially known.

Application

- Screen for activators or inhibitors of antibody-mediated signaling for immunotherapy research and drug discovery.
- Characterize CD5 antibodies and ligands.

Materials Provided

Components	Format
2 vials of frozen cells	~2 x 10 ⁶ cells in 1 ml of 90% FBS, 10% DMSO

Host Cell

CHO K1

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 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	BPS Bioscience #60186
Growth Medium 3B	BPS Bioscience #79529

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 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% CO₂. 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 3B (BPS Bioscience #79529):

F-12K Medium supplemented with 10% FBS, 1% Penicillin/Streptomycin plus 500 µg/ml of Hygromycin B.

Validation

Cell surface expression of human CD5 in CHO K1 cells was confirmed by flow cytometry.

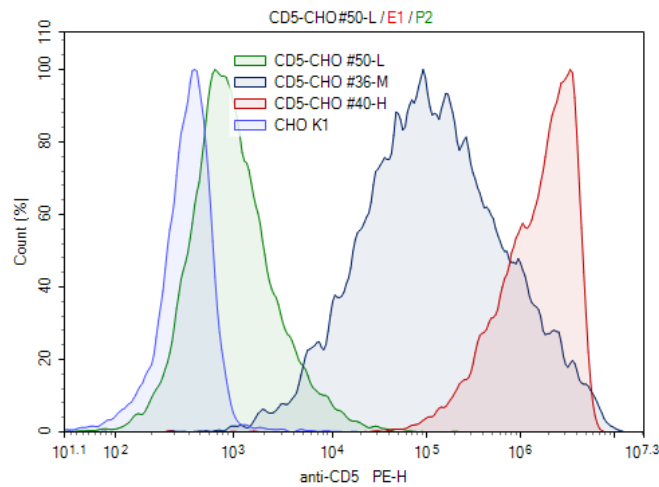


Figure 1. Flow cytometry analysis of cell surface expression of hCD5 in CHO K1 cells. CD5-CHO K1 cells or control CHO K1 cells were stained with PE-labeled anti-human CD5 antibody (Biolegend, #980358) and analyzed by flow cytometry. Y-axis is the % cell number. X-axis is the intensity of PE.

Green: low expression cell line; black: medium expression cell lines; red: high expression cell lines; blue: parental control CHO K1 cells.

Sequence

Human CD5 sequence (accession number: NM_014207.3)

MPMGSLQPLATLYLLGMLVASCLGRLSWYDPDFQARLTRNSKCGQLEVVYKDGWHMVCSSQSWGRSSKQWEDPSQASKVC
 QRLNCGVPLSLGPFVLYTTPQSSIICYGQLGFSFNCSHSRNDMCHSLGLTCLEPQKTTPTTRPPPTTTPEPTAPPRLQLVAQSGGQ
 HCAGVVEFYSGSLGGTISYEAQDKTQDLENFLCNLQCGSFLKHLPETEAGRAQDPGEPREHQPLPIQWKIQNSSCTSLEHCFRKI
 KPQKSGRVLALLCSGFQPKVQSRLVGGSSICEGTVEVRQGAQWAALCDSSSARSSLRWEEVCREQQCGSVNSYRVLDAGDPTSR
 GLFCPHQKLSQCHELWERNYSYCKKVFVTCQDPNPAGLAAGTVASIILALVLLVLLVCGPLAYKKLKFRQKKQRQWIGPTGM
 NQNMSFHRNHTATVRSNENPTASHVDNEYSQPPRNSHLSAYPALEGALHRSSMQPDNSSSDSYDLHGAQRL

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>
CD5 (Human) CRISPR/Cas9 Lentivirus (Non-Integrating)	78198	500 µl x 2
CD5 (Human) CRISPR/Cas9 Lentivirus (Integrating)	78119	500 µl x 2
CD5 – CHO K1 Recombinant Cell Line (Low Expression)	78290-L	2 vials
CD5 – CHO K1 Recombinant Cell Line (High Expression)	78290-H	2 vials
CD5, Fc Fusion, Avi-Tag, Biotin-labeled, HiP™	101006	20 µg/50 µg
CD5L, Avi-His-Tag	101007	100 µg/1 mg