

Description

Recombinant clonal CHO cell line stably expressing full-length human CEACAM5 (NM_004363). Surface expression of CEACAM5 was confirmed by flow cytometry. The stable clonal cell line was selected for different levels of CEACAM5 expression (High, Medium, and Low) compared to the parental CHO-K1 cell line.

Background

Carcinoembryonic antigen-related cell adhesion molecule 5 (CEACAM5, also known as CD66e) is a cell surface glycoprotein that serves as a cell adhesion protein. It has been used as a clinical biomarker to detect liver metastasis from gastrointestinal cancers and to predict gastrointestinal cancer relapse. CEACAM5 was recently identified as a potential target antigen for CAR T-cell therapy. Additionally, CEACAM5 may be involved in the inhibition of cell differentiation, apoptosis, and cell polarity.

Application(s)

- Screen and validate antibodies against CEACAM5 for drug discovery and research.
- Screen for compounds that regulate or inhibit CEACAM5 signaling in a cellular context.
- Perform binding assays to screen for potential CEACAM5 ligands.

Materials Provided

Components	Format
2 vials of frozen cells	Each vial contains 2×10^6 cells in 1 ml of cell freezing medium (BPS Bioscience #79796)

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% 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 3J (BPS Bioscience #79974):

F-12K 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 1:6 to 1:8 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 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

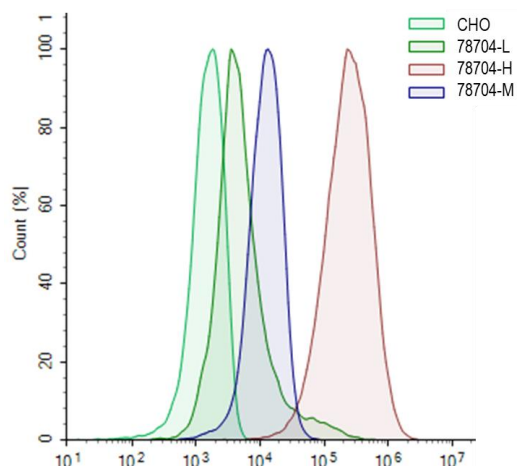


Figure 1: Cell surface expression of CEACAM5 in CEACAM5 CHO Cell Line.

CEACAM5 CHO cell line along with control parental CHO-K1 were stained with Alexa Fluor® 647 anti-human CD66d/e Antibody (Biolegend #392805) and the expression of human CEACAM5 was analyzed by flow cytometry.

Sequence

> NP_004354. *Homo sapiens* CEA cell adhesion molecule 5 (CEACAM5)

MESPSAPPHRWCI PWQRLLLTASLLTFWNPTTAKLTIESTPFNV AEGKEVLLVHNL PQHLFGYSWYKGERVDGNRQIIGYVIGT
 QQATPGPAYSGREIIPNASLLIQNIIQNDTGFYTLHVIKSDLVNEEATGQFRVYPELPKPSISSNNSKPVEDKDAVAFTCEPETQDA
 TYLWWWVNNQSLPVPSPRLQLSNGNRTLTLFNVTRNDTASYKCETQNPVSARRSDSVILNVLYGPDAPTISPLNTSYRSGENLNLSC
 AASNPPAQYSWVFNQSTQELFIPNITVNNSGSYTCQAHNSDTGLNRTT VTTITVYAEPKPFITSNNSNPVEDEDAVALTCE
 PEIQNTTYLWWWVNNQSLPVPSPRLQLSNDNRTLLSVTRNDVGPYECGIQNELSVDHSDPVLNVLYGPDPTISPSYTYRPGVN
 LLSLSCHAASNPPAQYSWLIDGNIQQHTQELFISNITEKNSGLYTCQANNSASGHSRTTVKTITVSAELPKPSISSNNSKPVEDKDAV
 AFTCEPEAQNTTYLWWWVNGQSLPVPSPRLQLSNGNRTLTLFNVTRNDARAYVCGIQNSVSANRSDPVTLDVLYGPDTPHSPDSS
 YLSGANLNLSCHSANPSPQYSWRINGIPQHTQVLFIAKITPNNGTYACFVSNLATGRNNSIVKSITVSASGTSPGLSAGATVGI
 MIGVLVGVALI

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

Products	Catalog #	Size
CEACAM5, Avi-His-Tag Recombinant	100509	100 µg
CEACAM, Avi-His-Tag Recombinant	70206	100 µg
CEACAM, Avi-His-Tag, Biotin Labeled Recombinant	70201	25 µg/50 µg
Claudin-18 Isoform 1 CHO Cell Line	78361	2 vials
Claudin-18 Isoform 2 CHO Cell Line	78533	2 vials
EpCAM CHO Cell Line	78683	2 vials