

Description

Recombinant clonal CHO cell line stably expressing full-length human EpCAM (NM_002354). Surface expression of EpCAM was confirmed by flow cytometry. The stable clonal cell line was selected for high levels of EpCAM expression compared to the parental CHO-K1 cell line.

Background

Epithelial cell adhesion molecule (EpCAM, also known as CD326) is a known biomarker of epithelial cells that regulates cell proliferation and differentiation. It is used as an epithelial cell biomarker to detect circulating tumor cells. EpCAM belongs to the small GA₇₃₃ protein family and is a homophilic transmembrane glycoprotein involved in cell-cell adhesion and tissue plasticity. It is a single-chain membrane-spanning protein with three major domains known as EpEX, involved in epidermal growth factor receptor (EGFR) mediated signaling pathways. Due to the versatile function of EpCAM and its crosstalk with other signaling pathways in cancer, EpCAM is considered an attractive target for cancer diagnosis, prognosis, and potential therapies.

Application(s)

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

Materials Provided

Components	Format
2 vials of frozen cells	Each vial contains 2 x 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.

- 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.

Validation Data

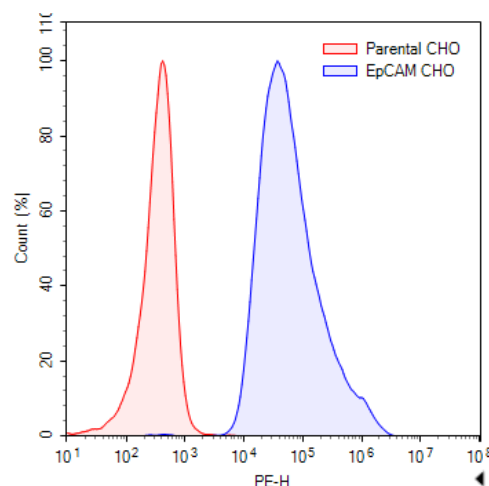


Figure 1: Cell surface expression of EpCAM in EpCAM CHO Cell Line.

EpCAM CHO cell line and parental CHO-K1 were stained with PE-labeled anti-human EpCAM Antibody (Biolegend #324205) and the expression of human EpCAM was analyzed by Flow Cytometry.

Sequence

> NP_002345 *Homo sapiens* epithelial cell adhesion molecule precursor (EpCAM/CD326)
MAPPQVLAFLGLLAAATATFAAAQEEVCENYKLA VNCVFVNNRQCQCTSVGAQNTVICSKLAAKCLVMKAEMNGSKLGRRA
KPEGALQNNNDGLYDPDCDESLFKAKQCNGTSMCWCVNTAGVRRRTDKDTEITCSERVRTYWIIEELKHKAREKPYDSKSLR TALQ
KEITTRYQLDPKFITSILYENNVITIDL VQNSSQKTQNDVDIADVAYYFEKDVKGESLFHSHKMDLTVNGEQLDLDPGQTLIYYVDEK
APEFSMQGLKAGVIAVIVVVVIAVVAGIVVLVISRKKRMAKYEKAEIKEMGEMHRELNA

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.

Reference

Liu, Y., et al. *Exp Hematol Oncol* **11**: 97 (2022).

Related Products

Products	Catalog #	Size
EpCAM, Avi-His-Tag Recombinant	100461	100 µg
EpCAM, Avi-His-Tag, Biotin Labeled Recombinant	100462	various
Claudin-18 Isoform 1 CHO Cell Line	78361	2 vials
Claudin-18 Isoform 2 CHO Cell Line	78533	2 vials