

**Description**

MRC2 CHO Cell Line (High Expression) is a CHO-K1 cell line that expresses human MRC2 (mannose receptor C-type 2) protein (accession number NM\_006039.5) under the control of the cytomegalovirus (CMV) promoter. This cell line was generated by lentiviral transduction followed by puromycin selection and limiting dilution. Individual clones were screened based on MRC2 expression levels to obtain this cell line.

This cell line has been validated by flow cytometry.

**Background**

MRC2, also known as mannose receptor C-type 2, uPARAP or Endo180, is a transmembrane receptor characterized by a cysteine-rich N-terminal domain, a fibronectin type II domain, 8 C-type lectin-like domains, and a transmembrane region. MRC2 functions as a recycling endocytic receptor that facilitates collagen uptake and intracellular collagen degradation. It is expressed throughout embryonic development and in adult tissue fibroblasts and macrophages where it plays a role in tissue remodeling and ECM (extracellular matrix) maintenance. MRC2 is upregulated in a variety of cancers including sarcoma and leukemia, as well as in cancer-associated fibroblasts, where it contributes to ECM remodeling and tumor metastasis, making it a promising therapeutic target. Current therapies being developed are focusing primarily on antibody-drug conjugates (ADCs) that leverage MRC2's efficient trafficking as a recycling endocytic receptor to shuttle/internalize cytotoxic drugs into tumors and malignant cells.

**Application**

- Use as target cells to screen antibody-drug conjugates (ADCs) or other biologics targeting MRC2.

**Materials Provided**

Components	Format
2 vials of frozen cells	Each vial contains $\geq 1 \times 10^6$ 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 3L	BPS Bioscience #78104

### 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, the use of validated and optimized media from BPS Bioscience is *highly recommended*. 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 to maintain selective pressure on the cell population expressing the gene of interest.

Cells should be grown at 37°C with 5% CO<sub>2</sub>. BPS Bioscience's cell lines are stable for at least 10 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 3L (BPS Bioscience #78104):*

F-12K medium supplemented with 10% FBS, 1% Penicillin/Streptomycin plus 6 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.

**Note: 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.
3. Transfer the resuspended cells to a T25 flask or T75 flask and incubate at 37°C in a 5% CO<sub>2</sub> incubator.
4. After 24 hours of culture, check for cell attachment and viability. Change medium to fresh Thaw Medium 3 and continue growing in a 5% CO<sub>2</sub> 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 3L.

#### Cell Passage

1. Aspirate the medium, wash the cells with Phosphate Buffered Saline (PBS) without Ca<sup>2+</sup>/Mg<sup>2+</sup>, and detach the cells from the culture vessel with 0.05% Trypsin/EDTA.

## MRC2 CHO Cell Line (High Expression)

2. Once the cells have detached, add Growth Medium 3L and transfer to a tube.
3. Spin down cells at  $300 \times g$  for 5 minutes, remove the medium and resuspend the cells in Growth Medium 3L.
4. Seed into new culture vessels at the recommended sub-cultivation ratio of 1:15 twice per week.

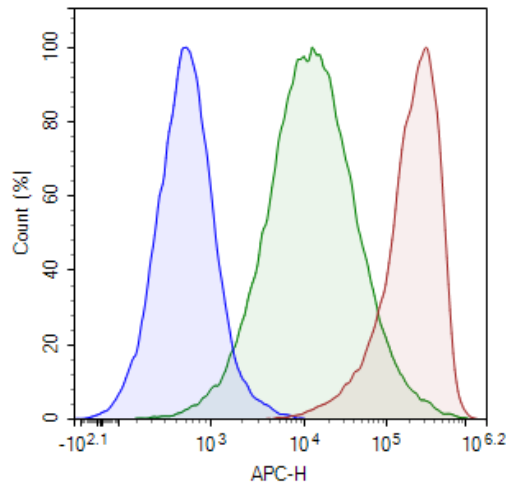
### Cell Freezing

1. Aspirate the medium, wash the cells with PBS without  $\text{Ca}^{2+}/\text{Mg}^{2+}$ , and detach the cells from the culture vessel with 0.05% Trypsin/EDTA.
2. Once the cells have detached, add Growth Medium 3L and count the cells.
3. Spin down the cells at  $300 \times g$  for 5 minutes, remove the medium and resuspend the cells in  $4^{\circ}\text{C}$  Cell Freezing Medium (BPS Bioscience #79796) at  $\sim 2 \times 10^6$  cells/ml.
4. Dispense 1 ml of cell suspension into each cryogenic vial. Place the vials in an insulated container for slow cooling and store at  $-80^{\circ}\text{C}$  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.

### Validation Data



*Figure 1: Cell surface expression of MRC2 in MRC2 CHO Cell Line by flow cytometry.*

High Expressing MRC2 CHO cells (#84205-H) (red), low expressing MRC2 CHO cells (#84205-L) (green), and control parental CHO-K1 cells (blue) were stained with Human MRC2 APC-conjugated Antibody and analyzed by flow cytometry. Y-axis is the % cell number. X-axis is the intensity of APC.

*Data shown is representative.*

## Sequence

Human MRC2 sequence (accession number NM\_006039.5)

MGPGRPAPAPWPRHLLRCVLLLGCLHLGRPGAPGDAALPEPNVFLIFSHGLQGCLQAQGGQVRVTPACNTSLPAQRWKWVSR  
NRLFNLGTMQCLGTGWPGTNTTASLGMYECDREALNLRWHCRTLGDQLSLLLGARTSNISKPGTLERGDQTRSGQWRIYGSEE  
DLCALPYHEVYTIQGNSHGKPCPTIPFKYDNQWFHGCTSTGREDGHLWCATTQDYGKDERWGFCEPIKSNDCETFWDKQDQDLDSC  
YQFNQSTLSWREAWASCEQQGADLLSITEIHEQTYINGLLTGYSSTLWIGLNDLDTSGGWQWSDNSPLKYLNWESDQPDNPS  
EENCGVIRTESSGGWQNRDCSIALPYVCKKPNATAEPTPPDRWANVKVECEPSWQPFQGHYRQLQAEKRSWQESKKAACLRG  
GGDLVSIHSMAELEFITKQIKQEVEELWIGLNDLKLQMNFEWSDGSLVSFTHWHPFEPNNFRDSLEDCVTIWGPEGRWINDSPC  
NQSLPSICKKAGQLSQGAAEEDHGCRKGTWVHSPSCYWLGEDQVTYSEARRLCTDHGSQVLVITNRFEQAFVSSLIYNWEGEYF  
WTALQDLNSTGSFFWLSGDEVMYTHWNRDQPGYSRGGCVLATGSAMGLWEVKNCTSFRRARYICRQSLGTPVTPPELPGPDPT  
PSLTGSCPPQGWASDTKLRYCYKVFSSERLQDKKSWVQAQGACQELGAQLLSLASYEEHFVANMLNKIFGESEPEIHEQHWFWI  
GLNRRDPRGGQSWRWSGQVGFVSYHNFDRSRHDDDDIRGCAVLDLASLQWVAMQCDTQLDWICKIPRGTDVREPDDSPQGR  
REWLRFOEAAYKFFHHSTWAQAQRICWFQAELTSVHSQAELDFLSHNLQKFSRAQEQHWWIGLHTSESDGRFRWTDGSIIN  
FISWAPGKPRPVGKDKKCVYMTASREDWGDQRCLTALPYICKRSNVTKETQPPDLPTTALGGCPSDWIQFLNCKCFVQGGQEPQ  
SRVKWSEAFQSCQEQEAQLVITNPLEQAFITASLPNVTFDLWIGLHASQRDFQWVEQEPLMYANWAPGEPSPGSPAPSGNKP  
TSCAVVLHSPSAHFTGRWDDRSCTEETHGFICQKGTDPSPSPAPALPPAPGTELSYLNGLTFRLLQKPLRWHDALLCESRNASLA  
YVPDPYTAFLTAARGLRTPWIGLAGEEGSRRYSWVSEELNYVGVQDGEPPQPGGCTYVDVDGAWRTTSCDTKLQGAVC  
GVSSGPPPPRRISYHGSCPQGLADSAWIPFREHCYSFHMELLGHKEARQRCRAGGAVLSILDEMENVFVWEHLQSYEGQSRG  
AWLGMNFPNPKGTLVWQDNTAVNYSNWGPPGLGSPMLSHNSCYWIQNSGLWRPGACTNITMGVVCKLPRAEQSSFSPSA  
LPENPAALVVVLMVLLLLLALLTAAILYRRRQSIERGAFFEGARYSRSSSSPTEATEKNILVSDMEMNEQQE

## References

- Evans R., et al., 2023 *Mol Cancer Ther* 22(2):240-253.  
Jungwirth U., et al., 2021 *Nat Commun* 12(1):3516.  
López-Guisa J., et al., 2012 *J Am Soc Nephrol* 23(2):236-251.  
Melander M., et al., 2015 *Int J Oncol* 47(4) :1177-1188.  
Mills T., et al., 2025 *Cancer Letters* 635:218068.  
Yuan Z., et al., 2025 *J Biol Chem* 301(3):108330.

## 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 lot-specific information and all other questions, please visit <https://bpsbioscience.com/contact>.

Version 052726