



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

CD123 (IL3Ra)-CHO Recombinant Cell Line (High Expression) Catalog # 79640-H

Description

Stable, recombinant clonal CHO cell line constitutively expressing full length human CD123 protein (Genbank #NM_002183.2). Surface expression of CD123 was confirmed by flow cytometry. This clonal cell line was selected for high level expression of CD123. Clones exhibiting lower levels of CD123 expression are also available (#79640-M and 79640-L). Each stable clonal cell line was selected for different levels of CD123 expression to mimic different stages of cancer target cells with various CD123 expression levels.

Background

CD123, also known as interleukin 3 receptor alpha (IL3Ra), is a transmembrane protein normally expressed on basophils and plasmacytoid dendritic cells and overexpressed on multiple blood cancers and leukemic hematopoietic stem cells. It normally plays a role in cell proliferation. CAR-Ts and other therapies targeting CD123 are currently being evaluated in clinical trials. CD123 may be especially valuable in treating B cell acute lymphoblastic leukemia patients who have relapsed after CD19 CAR-T treatment.

Application

Useful for screening and validating antibodies against CD123 and CD123 CAR-T for immunotherapy research and drug discovery. Also useful for CD123 binding assays to screen for CD123 ligands.

Host Cell

CHO K1 cell line, Chinese Hamster Ovary

Format

Each vial contains ~ 2 x 10⁶ cells in 1 ml of FBS + 10% DMSO.

Storage

Store in liquid nitrogen immediately upon receipt.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Cell Culture

Thaw Medium 3 (BPS Bioscience, #60186): Ham's F-12K medium (Hyclone, #SH30526.01) supplemented with 10% FBS (Life Technologies, #26140-079), 1% Penicillin/Streptomycin (Hyclone, #SV30010.01).

Growth Medium 3B (BPS Bioscience, #79529): Thaw Medium 3 (BPS Bioscience, #60186) plus 500 µg/ml Hygromycin (Thermo Fisher/Life Technologies, #10687010).

Recommended Culture Condition

Thawing cells: Prepare a 15 ml conical tube with 10 ml of pre-warmed Thaw Medium 3 (**no hygromycin**). Quickly thaw cells in a 37°C water bath with constant and slow agitation. Clean the outside of the vial with 70% ethanol and immediately transfer the entire content to Thaw Medium 3 (**no hygromycin**). Avoid pipetting up and down, and gently rock the conical tube.

Spin the cells down at 150 x g for 5 minutes. Discard the medium and re-suspend the cell pellet in fresh Thaw Medium 3 (**no hygromycin**). Transfer the entire content to a T75 flask to distribute the cells. Incubate the cells in a humidified 37°C incubator with 5% CO₂. After 48-72 hours of incubation, change to fresh Thaw Medium 3 (**no hygromycin**), without disturbing the attached cells. Switch to Growth Medium 3B at the first passage.

Subculture: When cells reach 90% confluency, remove the medium and wash twice with PBS (without magnesium or calcium). Treat cells with 1 ml of 0.05% trypsin/EDTA and incubate for 2-3 minutes at 37°C. After confirming cell detachment by light microscopy, add 14 ml pre-warmed medium and gently pipette up and down to dissociate cell clumps. Dispense 0.5 ml of the cell suspension into a new T75 flask containing 9.5 ml pre-warmed media. Incubate cells in a humidified 37°C incubator with 5% CO₂. Cells should be split twice per week at a 1:30 split ratio. Freeze cells in Thaw Medium 3 + 10% DMSO. Cells have been demonstrated to be stable for at least 20 passages; BPS recommends preparing frozen stocks at an early passage.

Mycoplasma Testing

This cell line has been screened using the Quick Test Mycoplasma Detection Kit (Biotool.com, #B39032) to confirm the absence of mycoplasma contamination.

Application References

1. Pizzitola, I., *et al.* Chimeric antigen receptors against CD33/CD123 antigens efficiently target primary acute myeloid leukemia cells in vivo. *Leukemia*. 2014 Aug, **28(8)**: 1596-1605.
2. Mardiros, A., *et al.* T cells expressing CD123 chimeric antigen receptors for treatment of acute myeloid leukemia. *Curr Opin Hematol*. 2015 Nov, **22(6)**: 484-488.
3. Ruella, M., *et al.* Dual CD19 and CD123 targeting prevents antigen-loss relapses after CD19-directed immunotherapies. *J. Clin. Invest.* 2016 Oct, **126(10)**: 3814-3826.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Vector and Sequence

Human CD123 (NM_002183.2) was cloned into pCMV3.

```
ATGGTCCTCCTTTGGCTCACGCTGCTCCTGATCGCCCTGCCCTGTCTCCTGCAAACGAAG
GAAGATCCAAACCCACCAATCACGAACCTAAGGATGAAAGCAAAGGCTCAGCAGTTGACCT
GGGACCTTAACAGAAATGTGACCGATATCGAGTGTGTTAAAGACGCCGACTATTCTATGCC
GGCAGTGAACAATAGCTATTGCCAGTTTGGAGCAATTTCTTATGTGAAGTGACCAACTAC
ACCGTCCGAGTGGCCAACCCACCATTCTCCACGTGGATCCTCTTCCCTGAGAACAGTGGG
AAGCCTTGGGCAGGTGCGGAGAATCTGACCTGCTGGATTTCATGACGTGGATTTCTTGAGC
TGCAGCTGGGCGGTAGGCCCGGGGGCCCCCGCGGACGTCCAGTACGACCTGTACTTGAA
CGTTGCCAACAGGCGTCAACAGTACGAGTGTCTTCACTACAAAACGGATGCTCAGGGAAC
ACGTATCGGGTGTCTGTTTCGATGACATCTCTCGACTCTCCAGCGGTTCTCAAAGTTCCCAC
ATCCTGGTGCGGGGCAGGAGCGCAGCCTTCGGTATCCCCTGCACAGATAAGTTTGTCTGTC
TTTTACAGATTGAGATATTAACCTCCACCCAACATGACTGCAAAGTGTAATAAGACACATTC
CTTTATGCACTGGAAAATGAGAAAGTCATTTCAATCGCAAATTTTCGCTATGAGCTTCAGATAC
AAAAGAGAATGCAGCCTGTAATCACAGAACAGGTCAGAGACAGAACCTCCTTCCAGCTACT
CAATCCTGGAACGTACACAGTACAAATAAGAGCCCGGGAAAGAGTGTATGAATTCTTGAGC
GCCTGGAGCACCCCCCAGCGCTTCGAGTGCGACCAGGAGGAGGGCGCAAACACACGTGC
CTGGCGGACGTGCTGCTGATCGCGCTGGGGACGCTGCTGGCCCTGGTCTGTGTCTTCG
TGATCTGCAGAAGGTATCTGGTGATGCAGAGACTCTTTCCCCGCATCCCTCACATGAAAGA
CCCCATCGGTGACAGCTTCCAAAACGACAAGCTGGTGGTCTGGGAGGCGGGCAAAGCCG
GCCTGGAGGAGTGTCTGGTGACTGAAGTACAGGTCGTGCAGAAAACCTTAA
MVLWLTLALLIALPCLLQTKEDPNPPITNLRMKAKAQLTWDLNRNVTDIECVKDADYSMPAVN
NSYCQFGAISLCEVTNYTVRVANPPFSTWILFPENSGKPWAGAENLTCWIHDVDFLSCSWAVG
PGAPADVQYDLYLNVANRRQQYECLHYKTDAQGTRIGCRFDDISRLSSGSQSSHILVRGRSAA
FGIPCTDKFVFSQIEILTPPNMTAKCNKTHSFMHWKMRSHFNRFYELQIQKRMQPVITEQV
RDRTSFQLLNPGTYTVQIRARERVYEFLSAWSTPQRFECDEEGANTRAWRTSLLIALLGTLALL
VCVFVICRRYLVMQRLFPRIHPMKDPIGDSFQNDKLVVWEAGKAGLEECLVTEVQVVQKT
```

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com

Quality Assurance

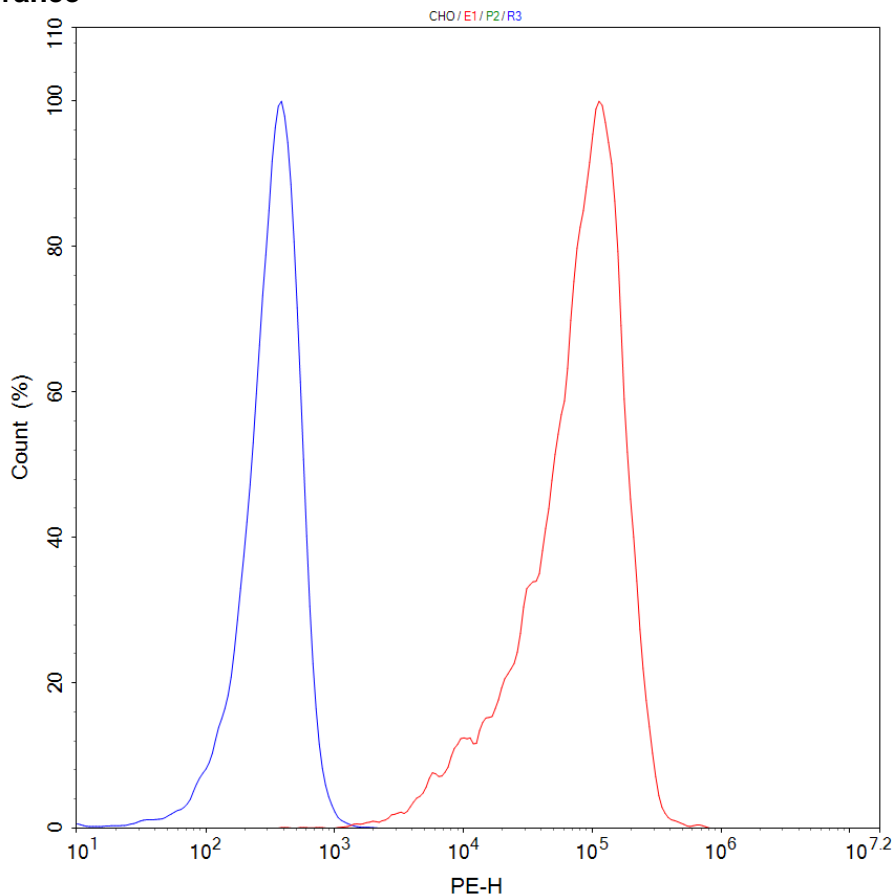


Figure 1. Expression of CD123 validated by flow cytometry. Flow cytometry using PE-conjugated anti-human CD123 antibody (Biolegend, #306006) detects CD123 on the surface of CD123-CHO Recombinant Cell Line, #79640-H. (CD123-CHO, red; CHO parental, blue).

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

License Disclosure: Purchase of this cell line grants you with a 10-year license to use this cell line in your immediate laboratory, for research use only. This license does not permit you to share, distribute, sell, sublicense, or otherwise make the cell line available for use to other laboratories, departments, research institutions, hospitals, universities, or biotech companies. The license does not permit the use of this cell line in humans or for therapeutic or drug use. The license does not permit modification of the cell line in any way. Inappropriate use or distribution of this cell line will result in revocation of the license and result in an immediate cease of sales and distribution of BPS products to your laboratory. BPS does not warrant the suitability of the cell line for any particular use and does not accept any liability in connection with the handling or use of the cell line. Modifications of this cell line, transfer to another facility, or commercial use of the cells may require a separate license and additional fees; contact sales@bpsbioscience.com for details. Publications using this cell line should reference BPS Bioscience, Inc., San Diego.

Related Products

Product	Cat. #	Size
CD123 (IL3Ra)-CHO Recombinant Cell Line (Medium Expression)	79640-M	2 vials
CD123 (IL3Ra)-CHO Recombinant Cell Line (Low Expression)	79640-L	2 vials
Growth Medium 3B	79529	500 ml
Thaw Medium 3	60186	100, 500ml
FOLH1 (PSMA)-CHO Recombinant Cell Line (High Expression)	79641-H	2 vials
FOLH1 (PSMA)-CHO Recombinant Cell Line (Medium Expression)	79641-M	2 vials
FOLH1 (PSMA)-CHO Recombinant Cell Line (Low Expression)	79641-L	2 vials
CD19-CHO Recombinant Cell Line (High Expression)	79561-H	2 vials
CD19-CHO Recombinant Cell Line (Medium Expression)	79561-M	2 vials
CD19-CHO Recombinant Cell Line (Low Expression)	79561-L	2 vials
CD22-CHO Recombinant Cell Line (High Expression)	79557-H	2 vials
CD22-CHO Recombinant Cell Line (Medium Expression)	79557-M	2 vials
BCMA-CHO Recombinant Cell Line (Medium Expression)	79500-M	2 vials
BCMA-CHO Recombinant Cell Line (High Expression)	79500-H	2 vials
BCMA-CHO Recombinant Cell Line (Low Expression)	79500-L	2 vials

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: [**info@bpsbioscience.com**](mailto:info@bpsbioscience.com)

Please visit our website at: [**www.bpsbioscience.com**](http://www.bpsbioscience.com)