



6405 Mira Mesa Blvd Ste 100  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

## **Data Sheet**

### **CD22 CHO Recombinant Cell Line (High Expression)**

### **Catalog #79557-H**

#### **Description**

Recombinant clonal stable CHO cell line constitutively expressing full length human CD22 protein (Genbank #NM\_001771). Surface expression of CD22 was confirmed by flow cytometry. This clonal cell line was selected for high level expression of CD22. A clone exhibiting lower levels of CD22 expression is also available (#79557-M). Please contact BPS Bioscience for additional information.

#### **Background**

CD22, also known as Siglec-2, is expressed on the membrane of B-cells. It is reported to act as an inhibitory co-receptor of the B-cell receptor to control the body's B-cell response. In 2017 the FDA approved inotuzumab ozogamicin (Besponsa), an antibody-drug conjugate targeting CD22, for patients with B-cell acute lymphoblastic leukemia (ALL). Additional therapies targeting CD22 are under evaluation.

#### **Application**

Useful for screening and validating antibodies against CD22 and different CD22 CAR-T for immunotherapy research and drug discovery. Also useful for CD22 binding assays to screen for CD22 ligands.

#### **Host Cell**

CHO K1 cell line, Chinese Hamster Ovary

#### **Format**

Each vial contains ~ 2.5 x 10<sup>6</sup> cells in 1 ml of Thaw Media + 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 **1.858.202.1401** Fax **1.858.481.8694**

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

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



6405 Mira Mesa Blvd Ste 100

San Diego, CA 92121

Tel: 1.858.202.1401

Fax: 1.858.481.8694

Email: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

### Cell Culture

**Thaw Medium 3 (BPS Bioscience, #60186):** F-12K Medium supplemented with 10% FBS, 1% Penicillin/Streptomycin

**Growth Medium 3D (BPS Bioscience, #79539):** F-12K Medium supplemented with 10% FBS, 1% Penicillin/Streptomycin plus 1 mg/ml G418.

### Recommended Culture Condition

Thawing cells: Prepare a 15 ml conical tube with 10 ml of pre-warmed Thaw Medium 3 (**no G418**). 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 G418**). 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 G418**). Transfer the entire content to a T75 flask to distribute the cells. Incubate the cells in a humidified 37°C incubator with 5% CO<sub>2</sub>. After 48-72 hours of incubation, change to fresh Thaw Medium 3 (**no G418**), without disturbing the attached cells. Switch to Growth Medium 3D 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 0.25% trypsin/EDTA and incubate for 2-3 minutes at 37°C. After confirming cell detachment by light microscopy, add 10 ml pre-warmed medium and gently pipette up and down to dissociate cell clumps. Dispense 1 ml of the cell suspension into a new T75 flask containing 9 ml pre-warmed media. Incubate cells in a humidified 37°C incubator with 5% CO<sub>2</sub>. Cells should be split twice per week at a 1:10 split ratio. Freeze cells in Thaw Medium 3 + 10% DMSO. Cells have been demonstrated to be stable for at least 15 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. Poe J, *et al.* CD22 and Siglec-G in B cell function and tolerance. *Trends Immunol.* 2012 Aug; **33(8)**:413-420.
2. Dorner T, *et al.* The mechanistic impact of CD22 engagement with epratuzumab on B cell function: Implications for the treatment of systemic lupus erythematosus. *Autoimmunity Reviews.* 2015 **14**:1079-1086.
3. Wei G., *et al.* Novel immunotherapies for adult patients with B-lineage acute lymphoblastic leukemia. *Journal of Hematology and Oncology.* 2017 **10**:1-13.

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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

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

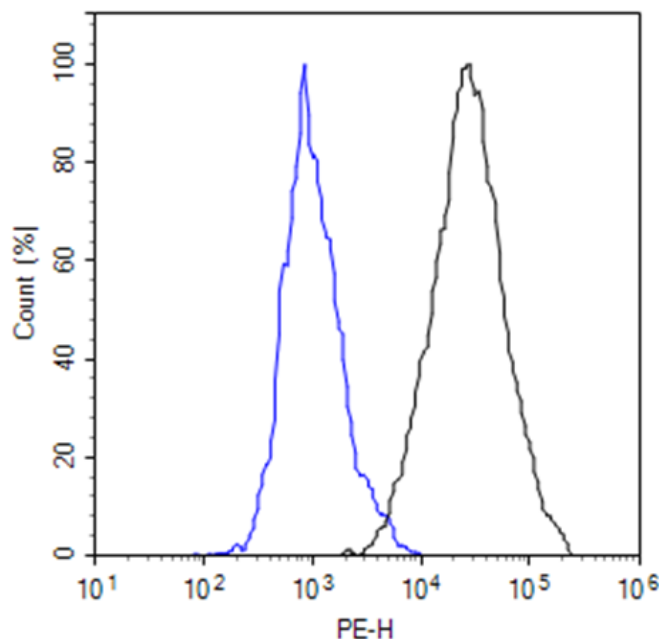
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)

### Vector and Sequence

Human CD22 (NM\_001771) was cloned into pIRESneo3.

TYRALDGDLESFILFHNPEYNKNTSKFDGTRLYESTKDGKVPSEQKRVQFLGDKNKNCTLSIHP  
 VHLNDSGQLGLRMESKTEKWMERIHNLVNSERPFPPHIQLPPEIQESQEVTLTCLLNFSYGYPI  
 QLQWLLEGVPMRQAAVTSTSLTIKSVFTRSELKFSQWSSHGKIVTCQLQDADGKFLSNDTVQ  
 LNVKHTPKLEIKVTPSDAIVREGDSVTMTCEVSSSNPEYTTVSWLKDGTSLKKQNTFTLNLREV  
 TKDQSGKYCCQVSNDVGPGRSEEVFLQVQYAPEPSTVQILHSPAVEGSQVEFLCMSLANPLP  
 TNYTWYHNGKEMQGRTEEKVHIPKILPWHAGTYSCVAENILGTGQRGPGAELDVQYPPKKVTT  
 VIQNPMPPIREGDTVTLSCNYNSSNPSVTRYEWKPHGAWEEPSLGLVKIQNVGWDNTTIACAAC  
 NSWCSWASPVALNVQYAPRDVVRKIKPLSEIHSNGNSVSLQCDFSSSHPKEVQFFWEKNGRL  
 LGKESQLNFDISPEDAGSYSCWVNNSIGQTASKAWTLEVLYAPRRLRVSMSPGDQVMGKS  
 ATLTCESDANPPVSHYTWFDWNNQSLPYHSQKLRLEPVKVQHSGAYWCQGTNSVGKGRSPL  
 STLTVYYSPETIGRRVAVGLGSCLAAILAICGLKLQRRWKRTQSQQGLQENSSGQSFFVRNKK  
 VRRAPLSEGPLSLGCYNPMMEDGISYTTLRFPEMNIPRTGDAESSEMQRPPPDCDDTDTYSA  
 LHKRQVGDYENVIPDFPEDEGIHYSELIQFGVGERPQAQENVVDYVILKH

### Quality Assurance



**Figure 1. Expression of CD22 validated by flow cytometry.** Flow cytometry using PE-conjugated anti-human CD22 antibody (Biolegend, #302506) detects CD22 on the surface of CD22 CHO Recombinant Cell Line, #79557-H. (CD22 CHO, black; 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 **1.858.202.1401** Fax **1.858.481.8694**

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

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



6405 Mira Mesa Blvd Ste 100

San Diego, CA 92121

**Tel:** 1.858.202.1401

**Fax:** 1.858.481.8694

**Email:** [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

**License Disclosure:**

Visit [bpsbioscience.com/license](http://bpsbioscience.com/license) for the label license and other key information about this product.

**Related Products**

**Product**

Growth Medium 3D

Thaw Medium 3

CD22, Fc-fusion (IgG1), Avi-Tag HiP™

CD22, Fc-fusion (IgG1), Avi-Tag HiP™

CD22 CHO Recombinant Cell Line (Medium Expression)

BCMA-CHO Recombinant Cell Line (Medium Expression)

BCMA-CHO Recombinant Cell Line (High Expression)

BCMA-CHO Recombinant Cell Line (Low Expression)

**Cat. #**

79539

60186

79464

79466

79557-M

79500-M

79500-H

79500-L

**Size**

500 ml

100, 500ml

100 µg

50 µg

2 vials

2 vials

2 vials

2 vials

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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

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

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