

CAR-T CELL TOOLS

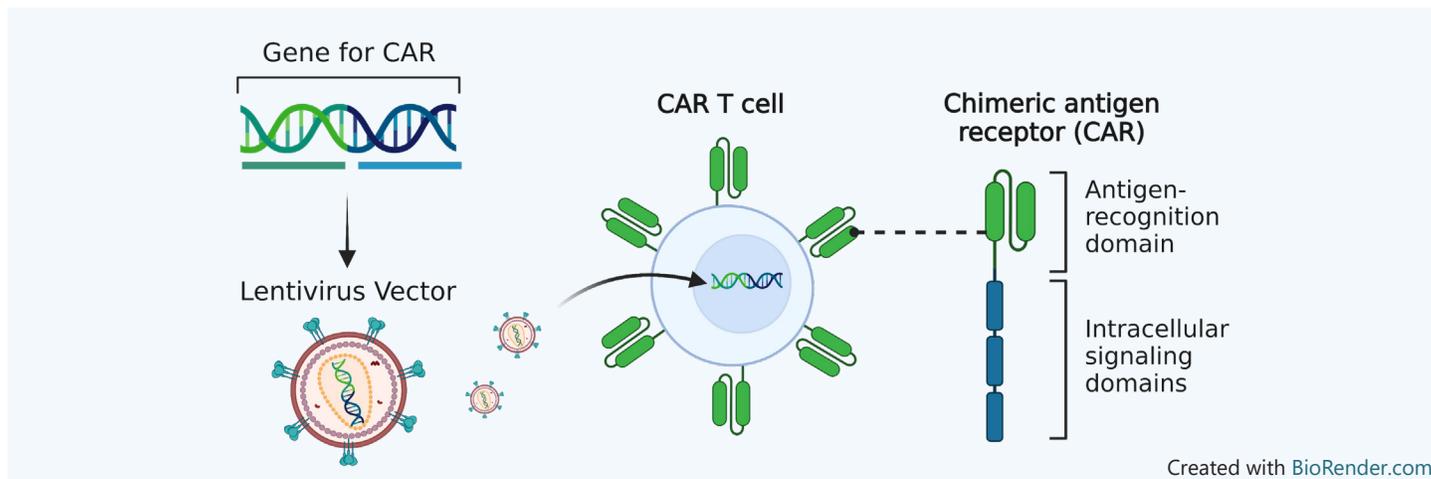
FOR ENGINEERING THE IMMUNE SYSTEM

Cell Lines | Primary Cells | Proteins | Antibodies | Viruses | Services



Engineering Effective CAR-T Cells

The development of Chimeric Antigen Receptor (CAR)-expressing T cells has made significant progress in the treatment of some types of cancer, with potential for applications in transplantation and chronic inflammation as well. More work is required to improve and expand therapies, and to limit harmful side-effects. The future is bright for CAR-T research, and BPS Bioscience continues to develop unique cell lines and other tools to help researchers create, evaluate, and enhance CAR-T cells for the improvement of human health.



Our Advantages



Produced In-House

- Made in the USA at our San Diego, CA laboratory
- Experience customized, personal support directly from our scientists



Committed to Excellence

- ISO 9001:2015-certified Quality Management System
- Lot-specific quality control testing



Expansive Portfolio

- Choose from ready-to-use proteins, cell lines, primary cells, antibodies, BiTE® (Bispecific T-cell Engager) molecules, lentivirus, and AAV
- Consistently launching new and novel products to meet research needs



CAR-T Cell Service Platform

- Lentivirus vector design and construction
- Preparation and validation of functional CAR-T cells
- Cytotoxic (cell killing) assays
- Generation of BiTE® constructs and production of BiTE® molecules
- Antibody screening
- Custom cell line development

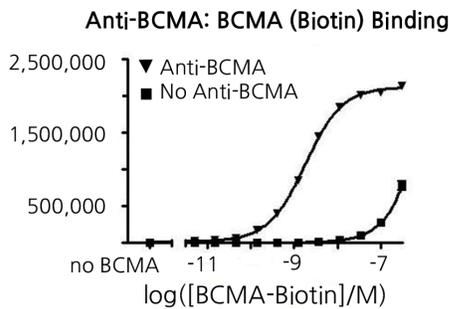
Proteins and Antibodies

● Recombinant Proteins

We specialize in the development and manufacture of bioactive enzymes and proteins, including a large selection related to CAR-T targets and T cell function. Our High-Purity (HiP™), low-aggregation protein products are optimal for generating clear and consistent research data. High purity means lower amounts of byproducts and contaminants from the manufacturing process and higher amounts of the full, expected length protein, which enables more accurate, better-quality results. Low aggregation means improved, more precise measurements for binding studies. BPS Bioscience maintains the highest standards for protein aggregation in drug discovery research.



- Functional proteins for binding and blocking assays
- CAR-T target proteins
- Immune checkpoint inhibitor proteins
- Cell activating and co-stimulatory molecules
- Epitope tagged, biotin-labeled, or fluorophore-labeled



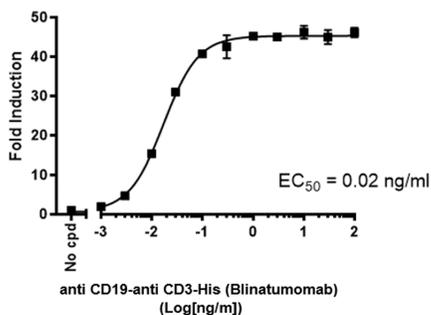
Anti-BCMA antibody (single-chain variable fragment) (#100173) was used to coat wells at 100 ng/well. Biotin-labeled BCMA recombinant protein (#79467) was titrated from 0-300 nM with a 3-fold dilution series and added to wells. Binding was detected with Streptavidin-HRP (#79742) and developed with a chemiluminescence readout.

● Antibodies



- Bispecific T cell Engager (BiTE®) molecules and trispecific antibodies, including anti-CD19-anti-CD3
- CAR-T targets, such as anti-BCMA, anti-CD19, and more
- T cell agonist antibodies, such as anti-CD3 and anti-CD28
- Recombinant production ensures consistent performance
- Human Ig isotype controls

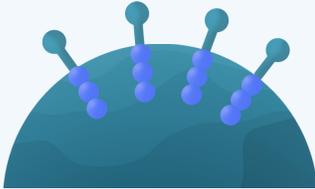
Activation of Jurkat Reporter Cells by Anti-CD19-Anti-CD3 BiTE® in Presence of CD19-positive Raji Cells



Anti-CD19-Anti-CD3 BiTE (equivalent to Blinatumomab) (#100441) was added at increasing concentrations to NFAT Reporter Luciferase Jurkat cells (#60641) in the presence of CD19-positive Raji cells. Luciferase induction was measured using the ONE-Step™ Luciferase Assay System (#60690).

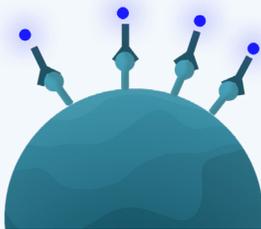
Cell Lines and Primary Cells

● CAR-T Cells



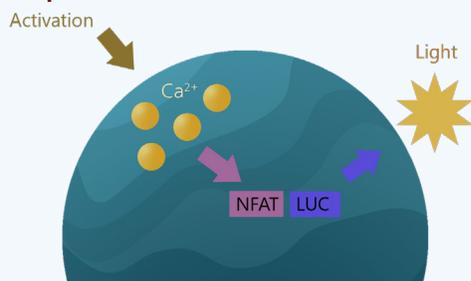
- For design of co-culture studies
- Positive controls for developing new CARs
- Cell lines and primary cells

● Overexpression/Knockout Cell Lines



- Overexpressing CAR-target cell lines, including BCMA, CD19, and more
- Antibody screening and binding studies
- TCR, B2M, and CIITA knockout cell lines to model universal CAR-T cells

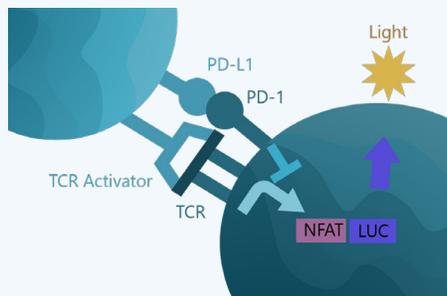
● Reporter Cell Lines



Measure:

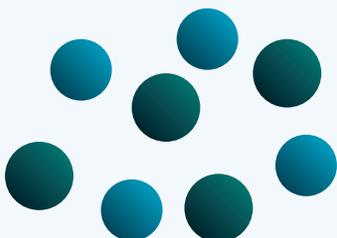
- Signaling activity
- Promoter activity
- Transcription factor activity

● Co-Stimulatory Cell Lines



- Inhibitor screening
- Antibody affinity
- Co-culture studies

● Primary Cells

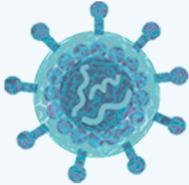


- Human $CD4^+$ and $CD8^+$ T cells, negatively selected
- Anti-BCMA and anti-CD19 CAR $CD4^+$ / $CD8^+$ T cells
- PBMCs
- Isolated from peripheral blood of healthy donors

Lentivirus and AAV Vectors

Virus-based tools such as lentiviruses and AAV are critical reagents for cell-engineering, particularly in CAR-T, gene therapy, and other personalized medicine. We have designed a suite of ready-to-use lentivirus and AAV vectors for CAR-T research and development.

Lentiviruses

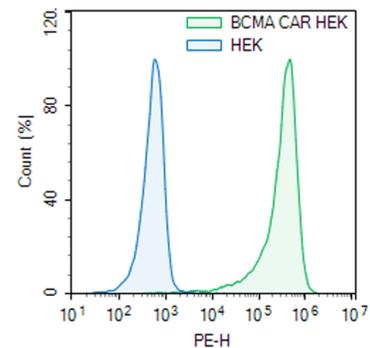
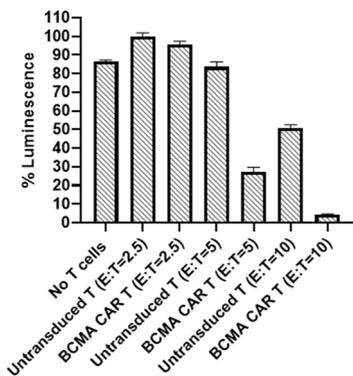


Lentiviruses are a popular tool for transducing CAR genes into primary T cells. Our replication-incompetent lentiviruses have been VSV-G pseudotyped, making these virus particles safe, stable and especially useful to target a wide range of cell types, particularly T cells in culture. Lentiviruses confer a number of advantages over other transduction methods.

Advantages

- Can infect actively dividing and non-dividing cells
- Can infect a wide range of cell stages
- Size of inserted DNA can be up to 10 kb
- Long term stable expression of a transgene
- Low cellular toxicity
- High transduction efficiency

Anti-BCMA CAR Lentivirus (Clone C11D5.3 ScFv-CD8-4-1BB-CD3ζ) (#78655)



Anti-BCMA CAR Lentivirus-transduced CD4⁺ and CD8⁺ T cells induce killing of Firefly Luciferase-RPMI8226 target cells.

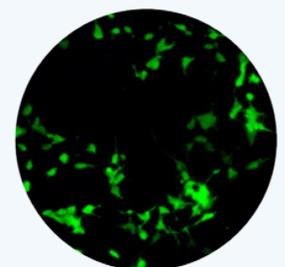
Anti-BCMA CAR Lentivirus-transduced HEK293 cells express CAR molecules that bind to biotinylated recombinant BCMA and PE-streptavidin.

AAV Reporter Vectors



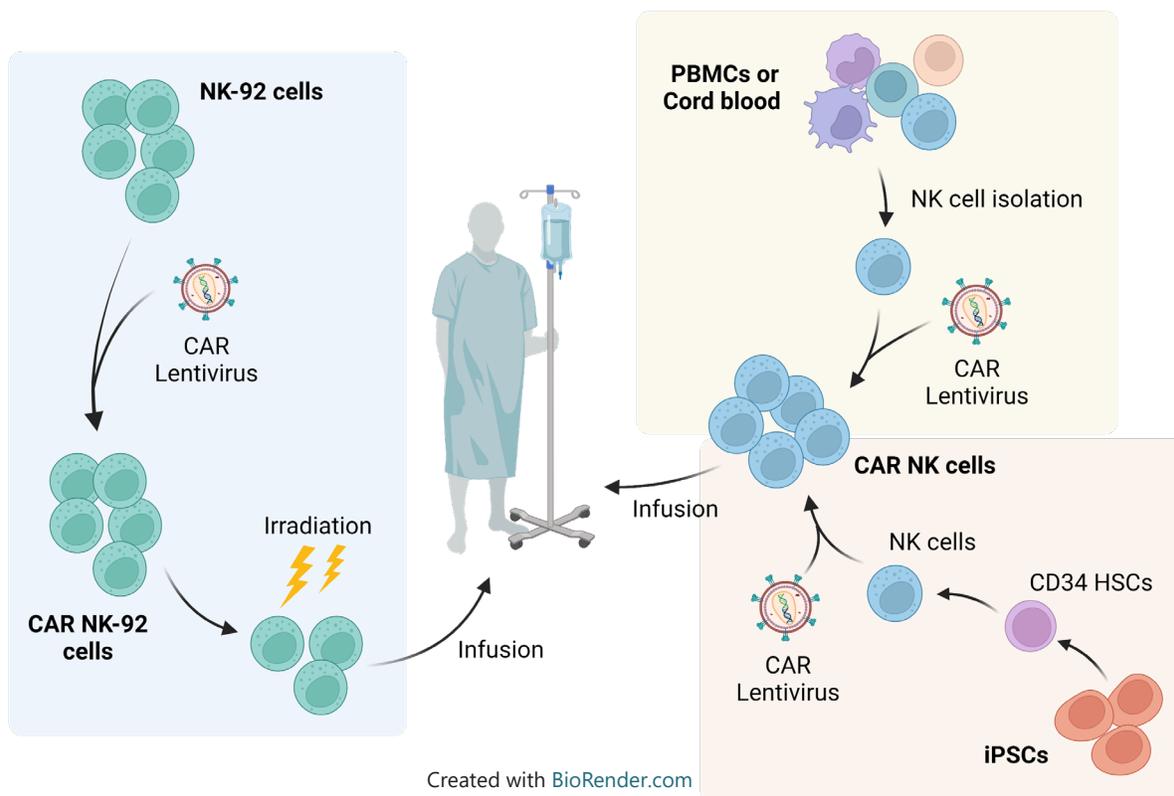
Adeno-Associated Virus (AAV) is a small non-enveloped dependovirus consisting of an icosahedral capsid containing a short, single-stranded DNA genome. Demonstrated as being safe for use in humans, AAV has been used for gene therapy to engineer cells using the viral genome to deliver the gene of interest.

We offer a growing line of AAV reporter vectors, such as luciferase or fluorescent markers for use as transduction controls, to track transgene expression over time, or for optimization of transduction and experimental conditions.



CAR-NK Cell-Related Products

Engineered CAR-NK cells are the next generation in CAR-expressing cell therapies. CAR-NK cells deliver a number of advantages over the existing CAR-T cell therapy, including fewer, less harmful side-effects, high feasibility for off-the-shelf manufacturing, which improves on delivery times, multiple mechanisms for activating cytotoxicity, and potential to be derived from multiple cell sources.



● NK-92 Cells

- Firefly Luciferase (#78400) or eGFP (#78399) constitutive expression
- Useful for NK cell killing assay controls or as a platform for CAR-NK cells

● Recombinant Proteins

- CD16A
- CD38
- KIR2DL1
- KIR2DL2
- KIR3DS1
- NKG2A
- NKG2D
- Nkp46
- Functional proteins for binding, blocking, and enzymatic assays
- Epitope tagged, Fc-fusion, or biotin-labeled
- Bulk production and customization

● Recombinant Cell Lines

- IL-15 Responsive Luciferase Reporter Cell Line (#78402)
- KIR3DL3/IL-2 Luciferase Reporter Jurkat Cell Line (#78322)
- FcGR3A (CD16A) CHO Cell Line (#78332)

● Lentiviruses

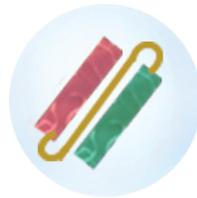
- Ideal for introducing transgenes into primary cells
- Stable integration for long term expression
- Custom production available

Custom CAR-T Cells

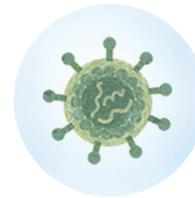
BPS Bioscience provides full service production of CAR-T cells to your desired specifications. With our milestone-measured process, you can monitor your steps to successful custom CAR-T cell generation.



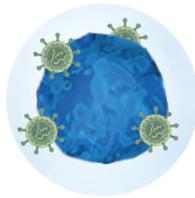
1
Researcher provides Ab sequence against antigen



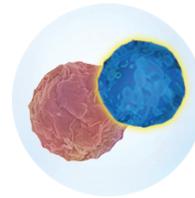
2
Engineering & validation of ScFv for specificity and affinity



3
CAR Lentivirus production and initial validation



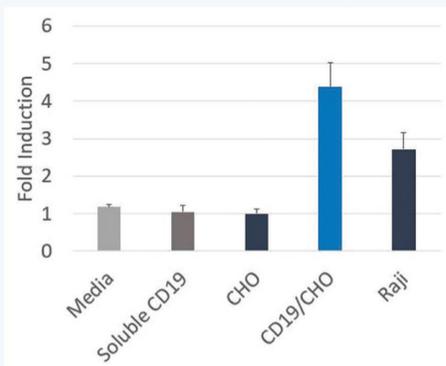
4
T cell preparation & transduction



5
Functional validation of CAR-T cells

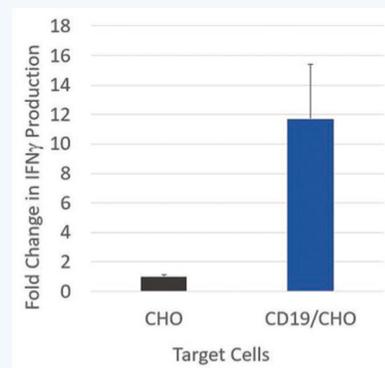
● Functional Validation

Primary Screening & Verification of CAR Activity Using a Reporter Cell Line



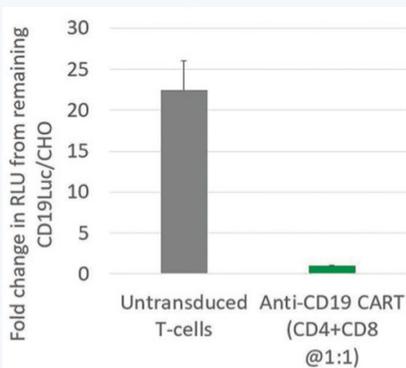
Luciferase activity in a stable cell pool of anti-CD19 CAR-expressing NFAT luciferase Jurkat cells co-cultured with the indicated targets and controls.

IFN γ Cytokine Detection from Activated CAR-T Cells



IFN γ production from Anti-CD19 CAR-T cells induced by CD19-expressing CHO cells (effector:target = 10:1). IFN γ was measured by ELISA (#79777).

Target Cell Killing Assays



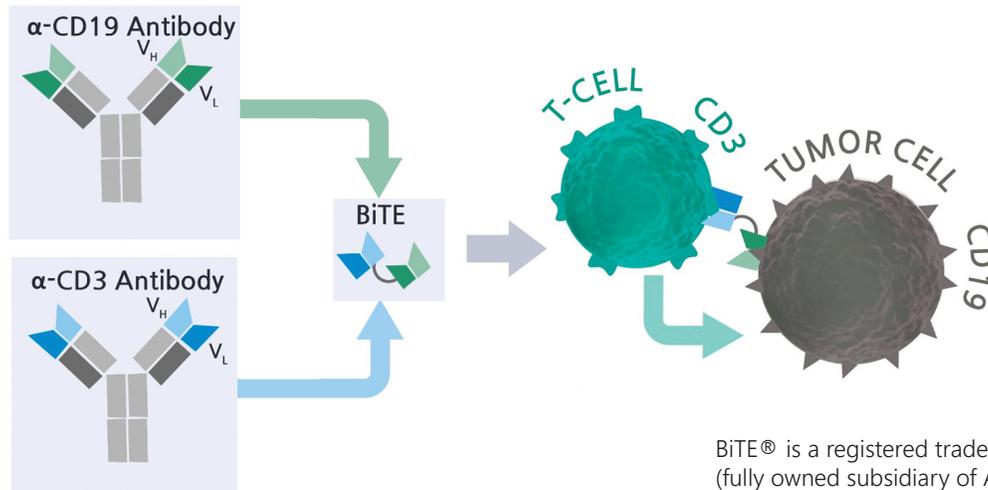
CD19 Luciferase CHO cells (79714) were targeted and killed by anti-CD19 CAR-T cells (effector:target = 10:1).

Additional Assays

- Flow cytometry to confirm CAR expression
- Mycoplasma testing

BiTE Molecules & Services

Bispecific T cell Engager (BiTE®) molecules are bivalent structures derived from two distinct antibodies designed as an immune-based therapeutic, by simultaneously engaging CD3 on T cells and a tumor-associated antigen expressed on cancer cells. This cell-to-cell ligation enables T cell targeting and killing of the tumor cells.



Our Capabilities

- Designing and generating BiTE constructs and producing BiTE molecules
- Measuring the affinity of BiTE binding to antigen targets using interferometry (Gator™, Probe Life) or ELISA-based assay
- Assessing T cell activation using reporter cell-based assays and measuring of K_d values using in vitro assay kits
- Evaluating bispecific constructs in reporter cell-based assays using our selection of over 100 antigen-expressing cancer cell lines

Comparing CAR-T vs BiTE Therapy



CAR-T

- Ex vivo engineered T cells derived from patients, complex production
- Indicated for hematologic cancers
- MHC/TCR-independent, independent of endogenous T cell populations, long lasting
- Lacks efficacy against solid tumors, antigen expression-dependent



BiTE

- In vitro engineered protein, not patient derived, relatively easy production
- Useful for hematologic cancers with potential for solid tumors
- MHC/TCR-independent, dependent on endogenous T cell populations
- Antigen-dependent, requires continuous/repeated administration due to short half-life

Antibodies	Catalog#
Anti-BCMA Antibody	101219
Anti-BCMA Antibody (Single-Chain Variable Fragment), His-Tag	100173
Anti-BCMA-Anti-CD19-Anti-CD3 Trispecific Molecule	100761
Anti-BCMA-Anti-CD3 Bispecific Molecule	100689
Anti-BCMA-Anti-CD3 IgG Bispecific Antibody	101968
Anti-CD19 Antibody, FITC-Labeled	101863
Anti-CD19 IgG Antibody	100981
Anti-CD19 IgG Antibody, Biotin-Labeled	101093
Anti-CD19-Anti-CD3 Bispecific Molecule	100441
Anti-CD19-Anti-CD3 IgG format Bispecific Antibody	101076
Anti-CD20 Antibody, FITC-Labeled	101864
Anti-CD20 Antibody, PE-Labeled	101672
Anti-CD20 Functional Antibody	71209
Anti-CD20 IgG Antibody, Biotin-labeled	101207
Anti-CD20-Anti-CD3 Bispecific Antibody	100836
Anti-CD20-Anti-CD3 IgM format Bispecific Antibody	100860
Anti-Claudin-18 Isoform 2 Antibody, FITC-Labeled	101866
Anti-Claudin-18 Isoform 2 Antibody, PE-Labeled	101676
Anti-Claudin-18 Isoform 2 IgG Antibody	101564
Anti-Claudin-18 Isoform 2 IgG Antibody, Biotin-Labeled	101565
Anti-Claudin-18 Isoform2-Anti-CD3 IgG Bispecific Antibody	101541
Anti-IL-2RA (CD25) Neutralizing Antibody	101593
Anti-PSMA Antibody	101695
Anti-PSMA, Biotin Label Antibody	101757
Anti-PSMA-Anti-CD3 IgG format Bispecific Antibody	101242

Biochemical Assay Kits	Catalog#
BAFF:BCMA[Biotinylated] Inhibitor Screening Assay Kit	79667
BCMA:APRIL[Biotinylated] Inhibitor Screening Assay Kit	79722
CD38 (Dog) Inhibitor Screening Assay Kit (Hydrolase Activity)	78108
CD38 (Mouse) Fluorogenic Assay Kit (Cyclase Activity)	78285
CD38 (Mouse) Inhibitor Screening Assay Kit (Hydrolase Activity)	79682
CD38 (Pig) Fluorogenic Assay Kit (Hydrolase Activity)	78178
CD38 Inhibitor Screening Assay Kit (Cyclase Activity)	71275
CD38 Inhibitor Screening Assay Kit (Hydrolase Activity)	79287

Biochemical Assay Kits	Catalog#
CD38 Inhibitor Screening Assay Kit (Hydrolase Activity)	79672
CD47:SIRP- α [Biotinylated] Inhibitor Screening Assay Kit	72044
CD47:SIRP- γ [Biotinylated] Inhibitor Screening Assay Kit	72059
Cytotoxicity Dye Kit (CFSE, 7-ADD)	82296

Cell Isolation Kits/Components	Catalog#
CD19 Positive Cell Isolation Kit	78564

Cell Lines	Catalog#
Anti-BCMA CAR /NFAT (Luciferase) Reporter Jurkat Cell Line	79694
Anti-CD19 CAR / NFAT (Luciferase) Reporter Jurkat Cell Line (CD19 SCFV-CD28-4-1BB-CD3 ζ)	79853
Anti-CD19 CAR Negative Control/NFAT (Luciferase) Reporter Jurkat Cell Line (CD19 SCFV-CD28 Transmembrane Motif)	79854
B2M Knockout Jurkat Cell Line	78342
B2M Knockout NFAT Luciferase Reporter Jurkat Cell Line	78363
B2M Knockout THP-1 Cell Line	78389
B2M/CIITA Double Knockout THP-1 Cell Line	78391
BCMA / CD20 / Firefly Luciferase CHO Cell Line	78185
BCMA / GLuc - CHO Recombinant Cell Line	79830
BCMA / Luciferase - CHO Recombinant Cell Line	79724
BCMA / NF- κ B - Reporter HEK293 Recombinant Cell Line	79755
BCMA CHO Recombinant Cell Line (High or Low Expression)	79500
CD19 / BCMA / Firefly Luciferase - CHO Recombinant Cell Line	78030
CD19 / CD20 / Firefly Luciferase CHO Cell Line	78186
CD19 / Firefly Luciferase - CHO Recombinant Cell Line	79714
CD19 CHO Recombinant Cell Line (Low, Medium and High Expression)	79561
CD20 CHO Recombinant Cell Line (High or Medium Expression)	79624
CD20/Firefly Luciferase CHO Cell Line	78620
CD22 / Luciferase - CHO Recombinant Cell Line	79715
CD22 CHO Recombinant Cell Line (Medium and High Expression)	79557
CD38 / BCMA / Firefly Luciferase CHO Recombinant Cell Line	78148
CD38 / CD19 / Firefly Luciferase CHO Recombinant Cell Line	78149
CD38 CHO Recombinant Cell Line (High, Medium or Low Expression)	79615
CD47 - HEK293 Cell Line	71249
CD7 CHO Cell Line (Medium or High Expression)	78324
CD8+ TCR Knockout NFAT-Luciferase Reporter Jurkat Cell Line	78757

Cell Lines	Catalog#	Lentiviruses	Catalog#
CIITA Knockout THP-1 Cell Line	78390	Anti-BCMA CAR Lentivirus (VHH1/VHH2 ScFv-CD8-4-1BB-CD3ζ)	78783
Claudin-18 Isoform 1 CHO Cell Line	78361	Anti-CD19 CAR Lentivirus (CD19 ScFv-CD8-4-1BB-CD3ζ)	78600
Claudin-18 Isoform 2 CHO Cell Line (High, Medium, or Low Expression)	78533	Anti-CD19 CAR Lentivirus (CD19 ScFv-CD8-4-1BB-CD3ζ, eGFP)	78775
eGFP/Firefly Luciferase K562 Cell Line	78911	Anti-CD19 CAR Lentivirus (CD19 ScFv-CD8-4-1BB-CD3ζ, PuroR)	78602
eGFP/Firefly Luciferase MIA PaCa-2 Cell Line	78766	Anti-CD19 CAR Lentivirus (CD19 ScFv-CD8-4-1BB-CD3ζ; SIN Vector)	78601
eGFP/Firefly Luciferase OVCAR3 Cell Line	78953	Anti-CD19/CD22 Bispecific CAR Lentivirus (Clones FMC63/m971 ScFv-CD8-4-1BB-CD3ζ)	78609
eGFP/Firefly Luciferase Ramos (RA 1) Cell Line	82149	Anti-CD20 CAR Lentivirus (Clone Leu-16 ScFv-CD8-4-1BB-CD3ζ)	78606
eGFP/Firefly Luciferase RS4;11 Cell Line	78926	Anti-CD22 CAR Lentivirus (Clone m971 ScFv-CD8-4-1BB-CD3ζ)	78608
FAP- CHO K1 Recombinant Cell Line (High, Medium or Low Expression)	79947	Anti-Mesothelin CAR Lentivirus (P4 ScFv-CD8-4-1BB-CD3ζ)	78703
FcGR1a (CD64) Knockout THP-1 Cell Line	82191	B2M (Human) CRISPR/Cas9 Lentivirus (Integrating)	78340
Firefly Luciferase - CHO Recombinant Cell Line	79725	B2M (Human) CRISPR/Cas9 Lentivirus (Non-Integrating)	78341
Firefly Luciferase CD19 Knockout NALM6 Cell Line	82168	BCMA Lentivirus	78714
GPRC5D (Cynomolgus) CHO Cell Line	78338	CD47 CRISPR/Cas9 Lentivirus (Integrating)	78056
GPRC5D (Cynomolgus) HEK293 Cell Line	78346	CD47 CRISPR/Cas9 Lentivirus (Non-Integrating)	78063
GPRC5D CHO Cell Line	78337	CIITA (Human) CRISPR/Cas9 Lentivirus (Integrating)	78435
GPRC5D HEK293 Cell Line	78345	CIITA (Human) CRISPR/Cas9 Lentivirus (Non-integrating)	78434
HER2 (ERBB2) CHO Recombinant Cell Line (High, Medium, or Low Expression)	79612	Dominant Negative TGF-β Receptor Type II (TGF-βRII) Lentivirus	78928
Human Mesothelin - CHO-K1 Recombinant Cell Line	78132	GPRC5D Lentivirus	78716
IL-2 Luciferase Reporter Jurkat Cell Line	60481	GPRC5D Lentivirus (Macaca fascicularis/Cynomolgus)	78780
MART-1 TCR (DMF4) CD8+ NFAT-Luciferase Reporter Jurkat Cell Line	78772	HLA-C*08:02 Lentivirus	78930
MUC16 (CA125), variant 4 (region 13785-14507) CHO Cell Line	78848	HLA-E Lentivirus	78929
NY-ESO-1 TCR (1G4) CD8+ NFAT-Luciferase Reporter Jurkat Cell Line	78769	MART-1-Specific TCR Lentivirus (Clone DMF4)	78678
NY-ESO-1 TCR (c259) CD8+ NFAT-Luciferase Reporter Jurkat Cell Line	78771	MART-1-Specific TCR Lentivirus (Clone DMF5)	78679
PSMA (FOLH1) - CHO Recombinant Cell Line (High, Medium, or Low Expression)	79641	NY-ESO-1-Specific TCR Lentivirus (Clone 1G4)	78675
TCR Activator CHO Recombinant Cell line	60539	NY-ESO-1-Specific TCR Lentivirus (Clone c259)	78676
TCR Activator Raji Cell Line	60556	PSMA Lentivirus	78726
TCR Activator/FcGR2B CHO Cell Line	78436	TCR CRISPR/Cas9 Lentivirus (Integrating)	78055
TCR Knockout Jurkat Cell Line	78539	Trop2 Lentivirus (Macaca fascicularis/Cynomolgus)	78776
TCR Knockout NFAT-Luciferase Reporter Jurkat Cell Line	78556		
TCR/B2M Knockout Jurkat Cell Line	78552		
TCR/B2M Knockout NFAT Luciferase Reporter Jurkat Cell Line	78557		
TROP2 - CHO-K1 Recombinant Cell Line	78099		

Lentiviruses	Catalog#
Anti-BCMA CAR Lentivirus (Clone C11D5.3 ScFv-CD8-4-1BB-CD3ζ)	78655

Peptides	Catalog#
MART-1 Peptide (26-35)	78759
MART-1 Peptide (26-35, Leu27)	78760
MART-1 Peptide (27-35)	78761
NY-ESO-1 Peptide (157-165)	78758

Primary Cells	Catalog#	Primary Cells	Catalog#	Proteins	Catalog#
Anti-Mesothelin CAR-T Cells	78729	Anti-BCMA CAR-T Cells	78660	CD47, Fc fusion, Biotin-labeled (Human) HiP™	71169
		Anti-CD19 CAR-T Cells	78171	CD47, Fc-Fusion, Streptavidin-Labeled	71292
		Anti-CD20 CAR-T Cells	78611	CD47, His-Tag (Human)	71127
		Untransduced T Cells	78170	Claudin-18 Isoform 2, FLAG-Tag	101570
Proteins	Catalog#	Proteins	Catalog#	EGFR, His-tag, GST-tag	40187
BCMA, Fc-fusion (IgG1), Avi-Tag, Biotin-Labeled	79467	IL-12 (p40/p35) Fc Fusion (IgG1), Avi-Tag	101431	GPC3, Avi-His-Tag	100071
BCMA, Fc-Fusion, Avi-Tag	79465	IL-12 (p40/p35) Fc Fusion (IgG1), Avi-Tag, Biotin-Labeled	101432	HER2, GST-Tag	40230
BCMA, Fc-Fusion, Avi-Tag, PE-Labeled	100733	IL-2 (C145A)	100159	IL-12 (p40/p35) Fc Fusion (IgG1), Avi-Tag	101431
c-Met, GST-tag	40255	IL-2 (R58A, F62A, Y65A, E82A, C145A) (Woodchuck)	100156	IL-12 (p40/p35) Fc Fusion (IgG1), Avi-Tag, Biotin-Labeled	101432
Carbonic Anhydrase 9 (CA9), His-tag	71101	IL-2 (R58D, K63E, E81R, C146A) (Woodchuck)	100157	IL-2 (C145A)	100159
CD19, Avi-His-Tag	101015	IL-2, Fc Fusion (IgG1), Avi-Tag, Biotin-Labeled	101381	IL-2 (R58A, F62A, Y65A, E82A, C145A) (Woodchuck)	100156
CD19, Fc-Fusion (IgG1), Avi-Tag	79472	IL-2RB (CD122)	79655	IL-2 (R58D, K63E, E81R, C146A) (Woodchuck)	100157
CD19, Fc-Fusion (IgG1), Avi-Tag, Biotin-Labeled	79475	IL-2RB, Avi-His-Tag	100427	IL-2, Fc Fusion (IgG1), Avi-Tag, Biotin-Labeled	101381
CD19, Fc-Fusion (IgG1), Avi-Tag, PE-labeled	100732	IL-2RB, Avi-His-Tag	100427	IL-2RB (CD122)	79655
CD20, FLAG-Tag	101572	IL-2RG, Avi-His-Tag	101149	IL-2RB, Avi-FLAG-Tag, Biotin-Labeled HiP™	101314
CD22, Fc Fusion, Avi-Tag, PE-labeled	101028	IL-2RG, Avi-His-Tag, Biotin-Labeled	101150	IL2RB, Avi-FLAG-Tag, HiP™	101313
CD22, Fc-fusion, Avi-Tag HiP™	79464	IL2RB, Avi-FLAG-Tag, Biotin-Labeled HiP™	101314	IL2RB, Avi-FLAG-Tag, HiP™	101313
CD22, Fc-fusion, Avi-Tag, Biotin-labeled HiP™	79466	IL2RB, Avi-His-Tag, Biotin-Labeled	100428	IL2RB, Avi-His-Tag, Biotin-Labeled	100428
CD277, Fc-Fusion (IgG1) Avi-Tag	100073	Mesothelin, Avi-His-Tag, Biotin-Labeled, HiP™	100291	Mesothelin, Avi-His-Tag, Biotin-Labeled, HiP™	100291
CD38, Avi-His-Tag	100346	Mesothelin, Avi-His-Tag, HiP™	100290	Mesothelin, Avi-His-Tag, HiP™	100290
CD38, Avi-His-Tag, Biotin-Labeled HiP™	100352	PDPN, Fc-Fusion, Avi-Tag HiP™	79341	PDPN, Fc-Fusion, Avi-Tag HiP™	79341
CD38, FLAG-Tag (Pig), HiP™	101019	PDPN, Fc-Fusion, Avi-Tag, Biotin-Labeled HiP™	79342	PDPN, Fc-Fusion, Avi-Tag, Biotin-Labeled HiP™	79342
CD38, His-Tag (Dog)	100955	PSMA, His-Avi-Tag	100463	PSMA, His-Avi-Tag	100463
CD38, His-Tag (Human), HiP™	71277	ROR1, Fc-Fusion (IgG1), Avi-Tag	79481	ROR1, Fc-Fusion (IgG1), Avi-Tag	79481
CD38, His-Tag (Mouse), HiP™	79070	ROR1, Fc-Fusion (IgG1), Avi-Tag, Biotin-Labeled	79482	ROR1, Fc-Fusion (IgG1), Avi-Tag, Biotin-Labeled	79482
CD38, His-Tag, PE-labeled	71882	ROR1, GST-tag	40396	ROR1, GST-tag	40396
CD38-APC, His-Tag	71883	ROR2, Fc-Fusion (IgG1), Avi-Tag HiP™	100029	ROR2, Fc-Fusion (IgG1), Avi-Tag HiP™	100029
CD47 (Monkey), Fc Fusion (Human), Avi-Tag HiP™	79118	ROR2, Fc-Fusion (IgG1), Avi-Tag, Biotin-Labeled HiP™	100046	ROR2, Fc-Fusion (IgG1), Avi-Tag, Biotin-Labeled HiP™	100046
CD47 (Monkey), Fc Fusion (Human), Avi-Tag, Biotin HiP™	79302	ROR2, GST-tag	40296	ROR2, GST-tag	40296
CD47, Fc Fusion (IgG1)	71177	Trop2 (88-274), Fc Fusion (IgG1), Avi-Tag	101346	Trop2 (88-274), Fc Fusion (IgG1), Avi-Tag	101346
CD47, Fc fusion, Avi-Tag (Human) HiP™	79051	Trop2 (88-274), Fc Fusion (IgG1), Avi-Tag, Biotin-Labeled	101347	Trop2 (88-274), Fc Fusion (IgG1), Avi-Tag, Biotin-Labeled	101347
CD47, Fc Fusion, Avi-Tag, Biotin-Labeled (Mouse)	72514	Trop2, Fc Fusion (IgG1), Avi-Tag	101344	Trop2, Fc Fusion (IgG1), Avi-Tag	101344



FIND A TRUSTED PARTNER FOR YOUR NEXT PROJECT



Ordering Information
orders@bpsbioscience.com



Technical Support
support@bpsbioscience.com



US Sales Support
sales-team@bpsbioscience.com



International Sales Support
international@bpsbioscience.com

BPS Bioscience, Inc.
6405 Mira Mesa Blvd, Suite 100
San Diego, CA 92121
Tel: 858-202-1401

bpsbioscience.com

