

ADOPTIVE CELL THERAPY

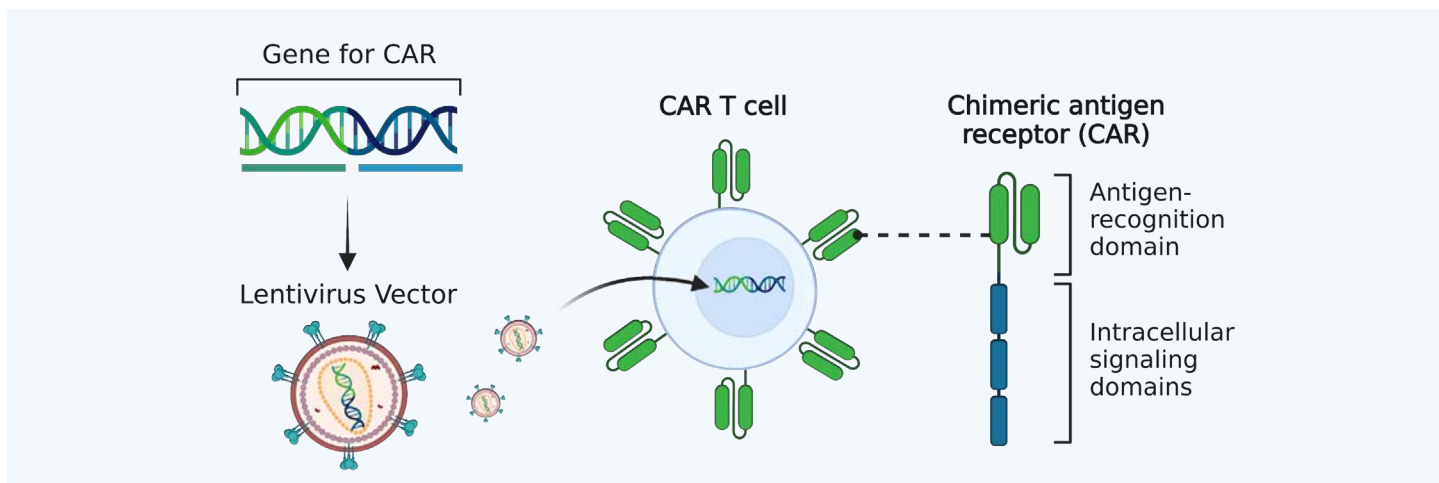
TOOLS FOR ENGINEERING THE IMMUNE SYSTEM

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



Engineering T & NK Cells for Adoptive Cell Therapy

The development of Chimeric Antigen Receptor (CAR)-expressing T cells has translated into significant progress in the treatment of some types of cancer, with potential for applications in transplantation and autoimmunity as well. Engineered CAR-NK cells also hold promise for improved therapies. In addition, cancer-targeting TCRs may provide expanded adoptive cell therapy options. The future is bright for CAR-T and TCR-T cell research, and BPS Bioscience continues to develop unique cell lines and other tools to help researchers create, evaluate, and enhance immune 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 cell lines, proteins/peptides, primary cells, antibodies, BiTE® (Bispecific T-cell Engager) molecules, lentivirus, and AAV
- Consistently launching new and novel products to meet research needs

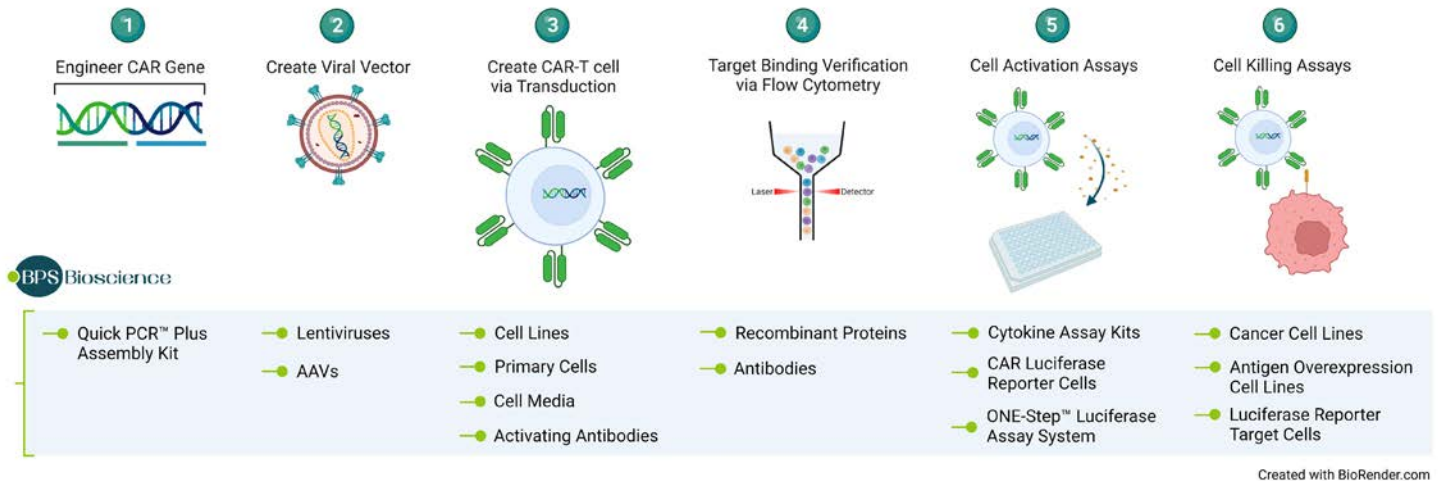


CAR-T Cell Service Platform

- Lentivirus vector design and construction
- Preparation and validation of functional CAR-T and TCR-T cells
- Cytotoxic (cell killing) assays
- Custom cell line development

Products to Enable the Cell Therapy Workflow

Our solutions support critical steps in CAR-T and CAR-NK cell therapy research and development spanning from creation of CAR-T cells through to functional testing and validation, according to the workflow below. Our off-the-shelf products are ready to ship upon order, and we also have custom capabilities to build unique reagents to your specifications and quantities. We can also help you design your CAR and generate CAR cells from start to finish.



Unique Solutions for TCR-T Research

Ready-to-use antigen-specific TCR-expressing reporter cells are available to support the design and optimization of co-culture bioassays, or to use as positive controls in bioassays. Robust signal-to-noise enables sensitive detection of TCR activation via cancer antigen peptides.

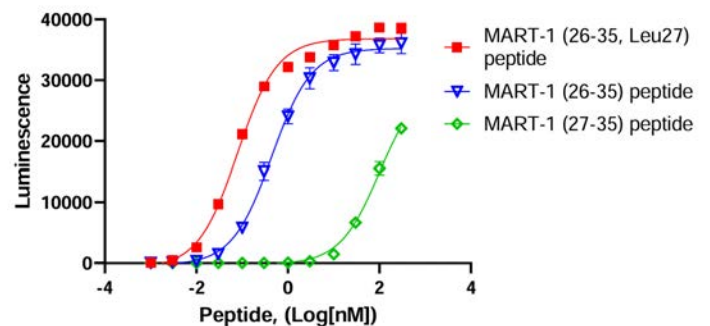
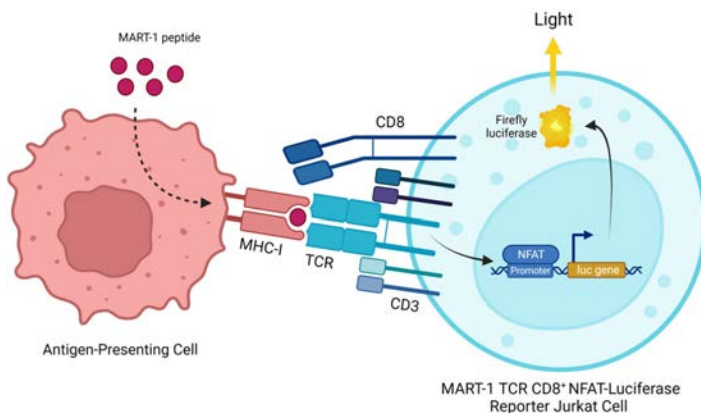
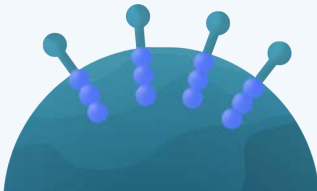


Illustration of the functional co-culture assay used to validate the MART-1 TCR (DMF4) CD8⁺ NFAT-Luciferase Reporter Jurkat cell line (#78772).

MART-1 TCR (DMF4) CD8⁺ NFAT Luciferase Reporter Jurkat cells were co-cultured for 6 hours with T2 cells loaded with various concentration of MART-1 peptides (#78759, #78760 and #78761). MART-1 peptide variants display differing affinities for MART-1. ONE-Step™ Luciferase Assay was performed and the results are shown as raw luminescence readings.

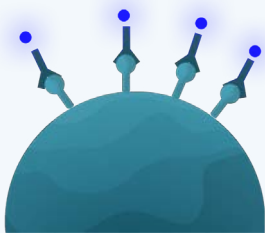
Cell Lines and Primary Cells

● CAR-T and TCR-T Cells



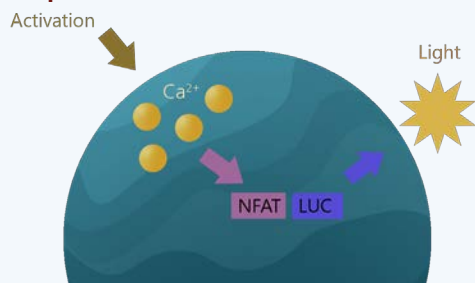
- For design of co-culture studies
- Positive controls for developing new CARs and TCRs
- Cell lines and primary cells

● 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 or re-engineering TCRs

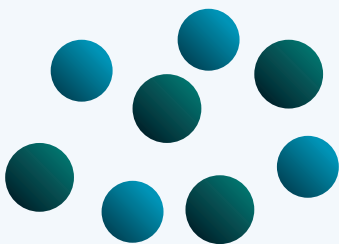
● Reporter Cell Lines



Measure:

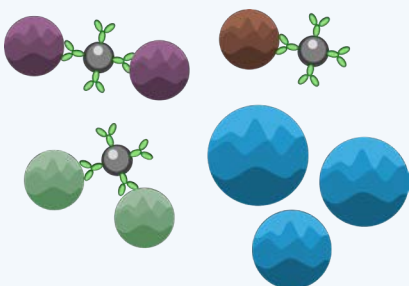
- TCR and CAR signaling activity
- Promoter activity
- Transcription factor activity

● Primary Cells



- Human CD4⁺ and CD8⁺ T cells, negatively selected
- CAR and TCR-expressing CD4⁺/CD8⁺ T cells
- PBMCs
- Isolated from peripheral blood of healthy donors

● Human T Cell Isolation Kit

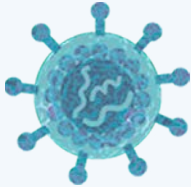


- Immuno-magnetic negative selection yields a highly purified T cell population
- Fast & efficient

Lentivirus and AAV Vectors

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

Lentiviruses

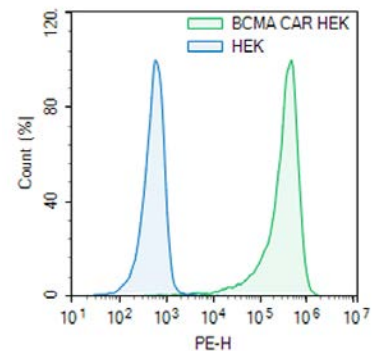
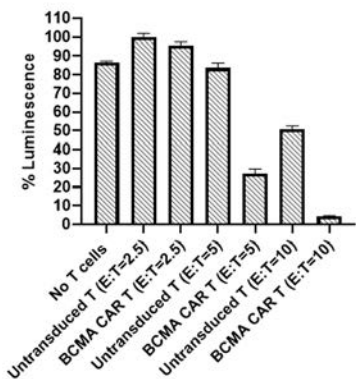


Lentiviruses are a popular tool for transducing CAR and TCR genes into primary T cells. Our replication-incompetent lentiviruses have been VSV-G pseudotyped, making these virus particles safe, stable and especially useful to target a wide range of cell types, particularly T cells in culture. Lentiviruses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

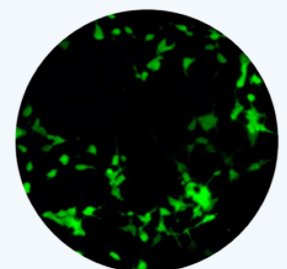
Anti-BCMA CAR Lentivirus-transduced HEK293 cells express CAR molecules that bind to biotinylated recombinant BCMA and PE-streptavidin.

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.



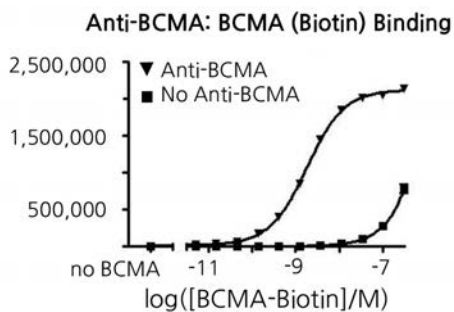
Proteins and Antibodies

● Recombinant Proteins

We specialize in the development and manufacture of bioactive enzymes and proteins, including a large selection related to CAR-T and TCR-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.

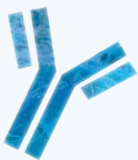


- Functional proteins for binding and blocking assays
- CAR-T target proteins
- Immune checkpoint inhibitor proteins
- Cell activating and co-stimulatory molecules, including TCR-specific peptides
- Epitope tagged, biotin-labeled, or fluorophore-labeled



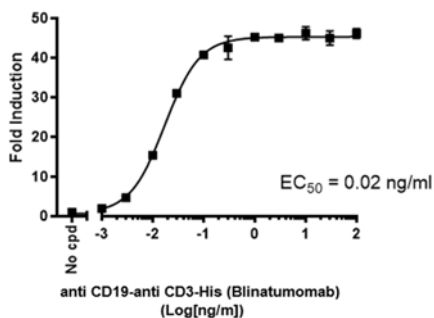
Anti-BCMA antibody (single-chain variable fragment) (#100173) was used to coat wells at 100 ng/well. Biotin-labeled BCMA recombinant protein (#79467) was titrated from 0-300 nM with a 3-fold dilution series and added to wells. Binding was detected with Streptavidin-HRP (#79742) and developed with a chemiluminescence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

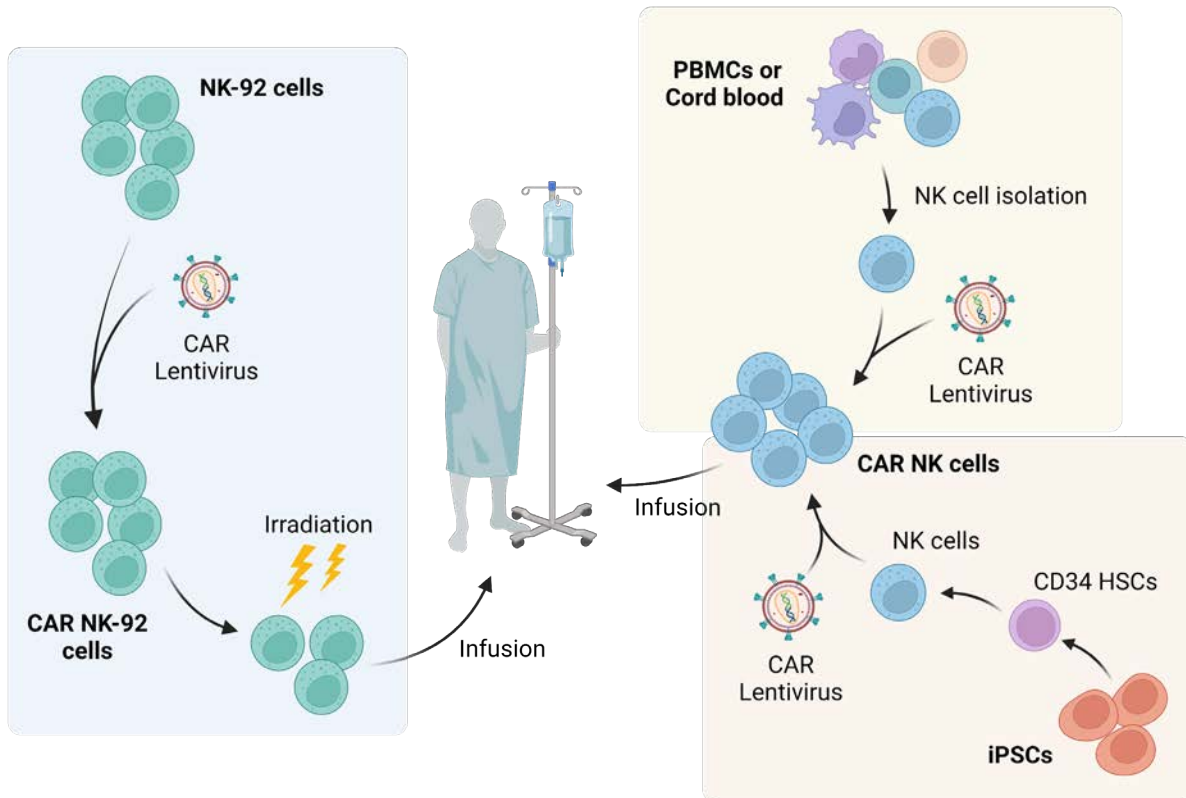
Activation of Jurkat Reporter Cells by Anti-CD19-Anti-CD3 BiTE® in Presence of CD19-positive Raji Cells



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).

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 Cells

- Magnetic NCAM1/CD56 Positive Cell Isolation Kit (#78808)
- NK Cell Expansion Kit (#78927)
- Expanded Human NK Cells (#78798)

● 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

- Growth-Arrested NK Feeder Cells (#78912)
- KIR3DL3/IL-2 Luciferase Reporter Jurkat Cell Line (#78322)
- FcGR3A (CD16A) (#78332) and FcGR3B (CD16B) (#78333) CHO Cell Lines
- NCAM1 (CD56) CHO Cell Line (High, Medium, & Low)

● Lentiviruses

- Ideal for introducing transgenes into primary cells and both dividing and non-dividing cells
- Stable integration for long term expression
- High transduction efficiency/low toxicity
- Custom production available

Custom CAR or TCR-T Cells

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



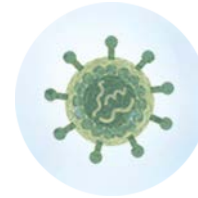
1

Researcher provides Ab sequence against antigen



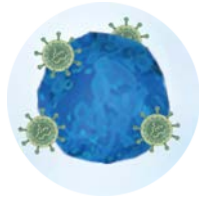
2

Engineering & validation of CAR or TCR for specificity and affinity



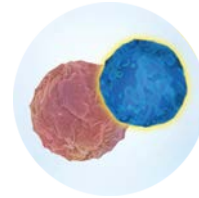
3

Lentivirus production and initial validation



4

T cell preparation & transduction

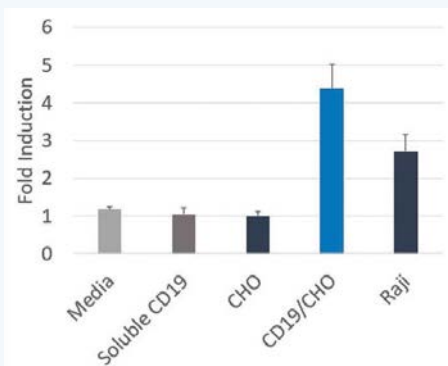


5

Functional validation of CAR or TCR-T cells

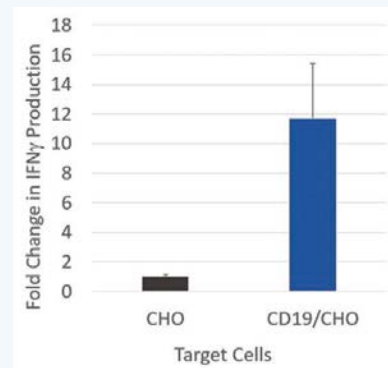
● Functional Validation

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



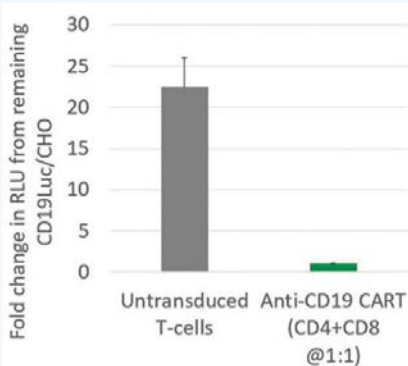
Luciferase activity in a stable cell pool of anti-CD19 CAR-expressing NFAT luciferase Jurkat cells co-cultured with the indicated targets and controls.

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).

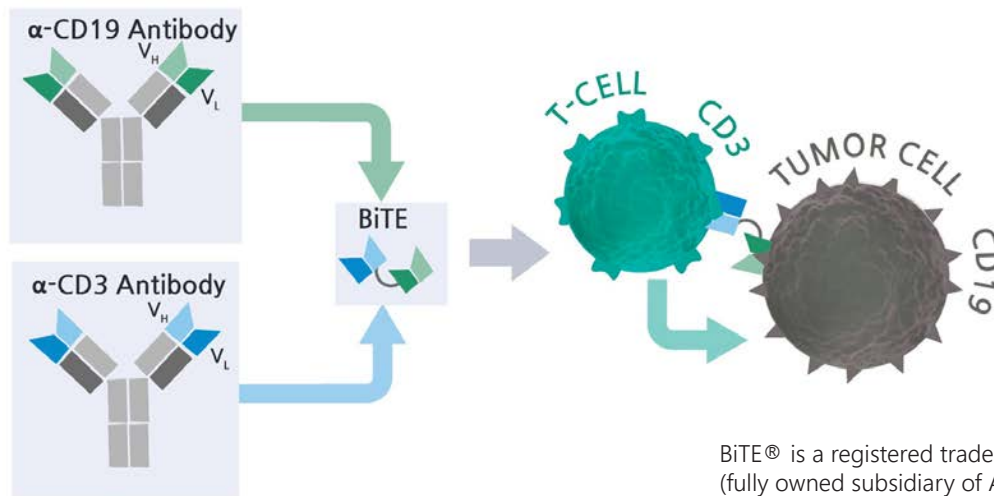
Additional Assays

- Flow cytometry to confirm CAR or TCR expression
- Mycoplasma testing

bpsbioscience.com/custom-car-t-cell-development

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
Human T Cell Isolation Kit	82288

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

Cell Lines	Catalog#
CD8+ TCR Knockout NFAT-Luciferase Reporter Jurkat Cell Line	78757
CIITA Knockout THP-1 Cell Line	78390
Claudin-18 Isoform 1 CHO Cell Line	78361
Claudin-18 Isoform 2 CHO Cell Line (High, Medium, or Low Expression)	78533
eGFP/Firefly Luciferase K562 Cell Line	78911
eGFP/Firefly Luciferase MIA PaCa-2 Cell Line	78766
eGFP/Firefly Luciferase OVCAR3 Cell Line	78953
eGFP/Firefly Luciferase Ramos (RA 1) Cell Line	82149
eGFP/Firefly Luciferase RS4;11 Cell Line	78926
FAP- CHO K1 Recombinant Cell Line (High, Medium or Low Expression)	79947
FcGR1a (CD64) Knockout THP-1 Cell Line	82191
Firefly Luciferase - CHO Recombinant Cell Line	79725
Firefly Luciferase CD19 Knockout NALM6 Cell Line	82168
GPRC5D (Cynomolgus) CHO Cell Line	78338
GPRC5D (Cynomolgus) HEK293 Cell Line	78346
GPRC5D CHO Cell Line	78337
GPRC5D HEK293 Cell Line	78345
HER2 (ERBB2) CHO Recombinant Cell Line (High, Medium, or Low Expression)	79612
Human Mesothelin - CHO-K1 Recombinant Cell Line	78132
IL-2 Luciferase Reporter Jurkat Cell Line	60481
KRAS G12D TCR (Clone 10) CD8+ NFAT-Luciferase Reporter Jurkat Cell Line	82303
KRAS G12D TCR (Clone 9c) CD8+ NFAT-Luciferase Reporter Jurkat Cell Line	82304
MAGE-A1 TCR CD8+ NFAT-Luciferase Reporter Jurkat Cell Line	78993
MAGE-A4 TCR CD8+ NFAT-Luciferase Reporter Jurkat Cell Line	78984
MART-1 TCR (DMF4) CD8+ NFAT-Luciferase Reporter Jurkat Cell Line	78772
MART-1 TCR (DMF5) CD8+ NFAT-Luciferase Reporter Jurkat Cell Line	78773
MUC16 (CA125), variant 4 (region 13785-14507) CHO Cell Line	78848
NY-ESO-1 TCR (1G4) CD8+ NFAT-Luciferase Reporter Jurkat Cell Line	78769
NY-ESO-1 TCR (c259) CD8+ NFAT-Luciferase Reporter Jurkat Cell Line	78771
PSMA (FOLH1) - CHO Recombinant Cell Line (High, Medium, or Low Expression)	79641
TCR Activator CHO Recombinant Cell line	60539
TCR Activator Raji Cell Line	60556
TCR Activator/FcGR2B CHO Cell Line	78436
TCR Knockout Jurkat Cell Line	78539
TCR Knockout NFAT-Luciferase Reporter Jurkat Cell Line	78556

Cell Lines	Catalog#
TCR/B2M Knockout Jurkat Cell Line	78552
TCR/B2M Knockout NFAT Luciferase Reporter Jurkat Cell Line	78557
TROP2 - CHO-K1 Recombinant Cell Line	78099

Cell-Based Assays and Expression Kits	Catalog#
PBMC Cytotoxicity Bioassay Kit (CFSE, 7-ADD)	82173
PBMC Cytotoxicity Luciferase Assay Kit	82214
PBMC Cytotoxicity Luciferase Assay Kit (NALM6)	82174

Lentiviruses	Catalog#
Anti-BCMA CAR Lentivirus (Clone C11D5.3 ScFv-CD8-4-1BB-CD3ζ)	78655
Anti-BCMA CAR Lentivirus (VHH1/VHH2 ScFv-CD8-4-1BB-CD3ζ)	78783
Anti-CD19 CAR Lentivirus (CD19 ScFv-CD8-4-1BB-CD3ζ)	78600
Anti-CD19 CAR Lentivirus (CD19 ScFv-CD8-4-1BB-CD3ζ, eGFP)	78775
Anti-CD19 CAR Lentivirus (CD19 ScFv-CD8-4-1BB-CD3ζ, PuroR)	78602
Anti-CD19 CAR Lentivirus (CD19 ScFv-CD8-4-1BB-CD3ζ; SIN Vector)	78601
Anti-CD19/CD22 Bispecific CAR Lentivirus (Clones FMC63/m971 ScFv-CD8-4-1BB-CD3ζ)	78609
Anti-CD20 CAR Lentivirus (Clone Leu-16 ScFv-CD8-4-1BB-CD3ζ)	78606
Anti-CD22 CAR Lentivirus (Clone m971 ScFv-CD8-4-1BB-CD3ζ)	78608
Anti-Mesothelin CAR Lentivirus (P4 ScFv-CD8-4-1BB-CD3ζ)	78703
B2M (Human) CRISPR/Cas9 Lentivirus (Integrating)	78340
B2M (Human) CRISPR/Cas9 Lentivirus (Non-Integrating)	78341
BCMA Lentivirus	78714
CD47 CRISPR/Cas9 Lentivirus (Integrating)	78056
CD47 CRISPR/Cas9 Lentivirus (Non-Integrating)	78063
CIITA (Human) CRISPR/Cas9 Lentivirus (Integrating)	78435
CIITA (Human) CRISPR/Cas9 Lentivirus (Non-integrating)	78434
Cyno EpCAM Lentivirus	78978
Dominant Negative TGF-β Receptor Type II (TGF-βRII) Lentivirus	78928
GPRC5D Lentivirus	78716
GPRC5D Lentivirus (Macaca fascicularis/Cynomolgus)	78780
HLA-C*08:02 Lentivirus	78930
HLA-E Lentivirus	78929
KRAS G12D-Specific TCR Lentivirus (Clone 10)	78937
KRAS G12D-Specific TCR Lentivirus (Clone 9c)	78936

Lentiviruses	Catalog#	Primary Cells	Catalog#
MAGE-A1-Specific TCR Lentivirus (Clone 1367)	78934	Anti-CD19 CAR-T Cells (eGFP)	78789
MAGE-A4 Specific TCR Lentivirus	78935	Anti-CD20 CAR-T Cells	78611
MART-1-Specific TCR Lentivirus (Clone DMF4)	78678	Human T Cell Activation Reagent	82283
MART-1-Specific TCR Lentivirus (Clone DMF5)	78679	Untransduced T Cells	78170
NY-ESO-1-Specific TCR Lentivirus (Clone 1G4)	78675		
NY-ESO-1-Specific TCR Lentivirus (Clone c259)	78676		
PSMA Lentivirus	78726		
TCR CRISPR/Cas9 Lentivirus (Integrating)	78055		
Trop2 Lentivirus (Macaca fascicularis/Cynomolgus)	78776		
Vy4V61 TCR Lentivirus	78986		
Vy9V62 TCR Lentivirus	78985		
Peptides	Catalog#	Proteins	Catalog#
KRAS G12D Peptide (10-18)	78967	BCMA, Fc-fusion (IgG1), Avi-Tag, Biotin-Labeled	79467
KRAS G12D Peptide (10-19)	78969	BCMA, Fc-Fusion, Avi-Tag	79465
KRAS Wild Type Peptide (10-18)	78968	BCMA, Fc-Fusion, Avi-Tag, PE-Labeled	100733
KRAS Wild Type Peptide (10-19)	78970	c-Met, GST-tag	40255
MAGE-A1 Peptide (278-286)	78965	Carbonic Anhydrase 9 (CA9), His-tag	71101
MAGE-A4 Peptide (230-239)	78966	CD19, Avi-His-Tag	101015
MAGE-A4 Peptide (286-294)	82305	CD19, Fc-Fusion (IgG1), Avi-Tag	79472
MART-1 Peptide (26-35)	78759	CD19, Fc-Fusion (IgG1), Avi-Tag, Biotin-Labeled	79475
MART-1 Peptide (26-35, Leu27)	78760	CD19, Fc-Fusion (IgG1), Avi-Tag, PE-labeled	100732
MART-1 Peptide (27-35)	78761	CD20, FLAG-Tag	101572
NY-ESO-1 Peptide (157-165)	78758	CD22, Fc Fusion, Avi-Tag, PE-labeled	101028
PRAME Peptide (394-402)	82307	CD22, Fc-fusion, Avi-Tag HiP™	79464
PRAME Peptide (425-433)	78991	CD22, Fc-fusion, Avi-Tag, Biotin-labeled HiP™	79466
PRAME Peptide (432-440)	82306	CD277, Fc-Fusion (IgG1) Avi-Tag	100073
		CD38, Avi-His-Tag	100346
		CD38, Avi-His-Tag, Biotin-Labeled HiP™	100352
		CD38, FLAG-Tag (Pig), HiP™	101019
		CD38, His-Tag (Dog)	100955
		CD38, His-Tag (Human), HiP™	71277
		CD38, His-Tag (Mouse), HiP™	79070
		CD38, His-Tag, PE-labeled	71882
		CD38-APC, His-Tag	71883
		CD47 (Monkey), Fc Fusion (Human), Avi-Tag HiP™	79118
		CD47 (Monkey), Fc Fusion (Human), Avi-Tag, Biotin HiP™	79302
		CD47, Fc Fusion (IgG1)	71177
		CD47, Fc fusion, Avi-Tag (Human) HiP™	79051
		CD47, Fc Fusion, Avi-Tag, Biotin-Labeled (Mouse)	72514
		CD47, Fc fusion, Biotin-labeled (Human) HiP™	71169
		CD47, Fc-Fusion, Streptavidin-Labeled	71292
Primary Cells	Catalog#		
Anti-Mesothelin CAR-T Cells	78729		
Dual Epitope Anti-BCMA CAR-T Cells	78790		
NY-ESO-1 (c259) TCR-T Cells	78990		
Untransduced T Cells (NY-ESO-1 TCR-T Negative Control)	78989		
Anti-BCMA CAR-T Cells	78660		
Anti-CD19 CAR-T Cells	78171		

Proteins	Catalog#
CD47, His-Tag (Human)	71127
Claudin-18 Isoform 2, FLAG-Tag	101570
EGFR, His-tag, GST-tag	40187
GPC3, Avi-His-Tag	100071
GPC3, Avi-His-Tagged, Biotin-Labeled	100072
HER2, GST-Tag	40230
IL-12 (p40/p35) Fc Fusion (IgG1), Avi-Tag	101431
IL-12 (p40/p35) Fc Fusion (IgG1), Avi-Tag, Biotin-Labeled	101432
IL-2 (C145A)	100159
IL-2 (R58A, F62A, Y65A, E82A, C145A) (Woodchuck)	100156
IL-2 (R58D, K63E, E81R, C146A) (Woodchuck)	100157
IL-2, Fc Fusion (IgG1), Avi-Tag, Biotin-Labeled	101381
IL-2RB (CD122)	79655
IL-2RB, Avi-His-Tag	100427
IL-2RG, Avi-His-Tag	101149
IL-2RG, Avi-His-Tag, Biotin-Labeled	101150
IL2RB, Avi-FLAG-Tag, Biotin-Labeled HiP™	101314
IL2RB, Avi-FLAG-Tag, HiP™	101313
IL2RB, Avi-His-Tag, Biotin-Labeled	100428
Mesothelin, Avi-His-Tag, Biotin-Labeled, HiP™	100291
Mesothelin, Avi-His-Tag, HiP™	100290
PDPN, Fc-Fusion, Avi-Tag HiP™	79341
PDPN, Fc-Fusion, Avi-Tag, Biotin-Labeled HiP™	79342
PSMA, His-Avi-Tag	100463
ROR1, Fc-Fusion (IgG1), Avi-Tag	79481
ROR1, Fc-Fusion (IgG1), Avi-Tag, Biotin-Labeled	79482
ROR1, GST-tag	40396
ROR2, Fc-Fusion (IgG1), Avi-Tag HiP™	100029
ROR2, Fc-Fusion (IgG1), Avi-Tag, Biotin-Labeled HiP™	100046
ROR2, GST-tag	40296
Trop2 (88-274), Fc Fusion (IgG1), Avi-Tag	101346
Trop2 (88-274), Fc Fusion (IgG1), Avi-Tag, Biotin-Labeled	101347
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

