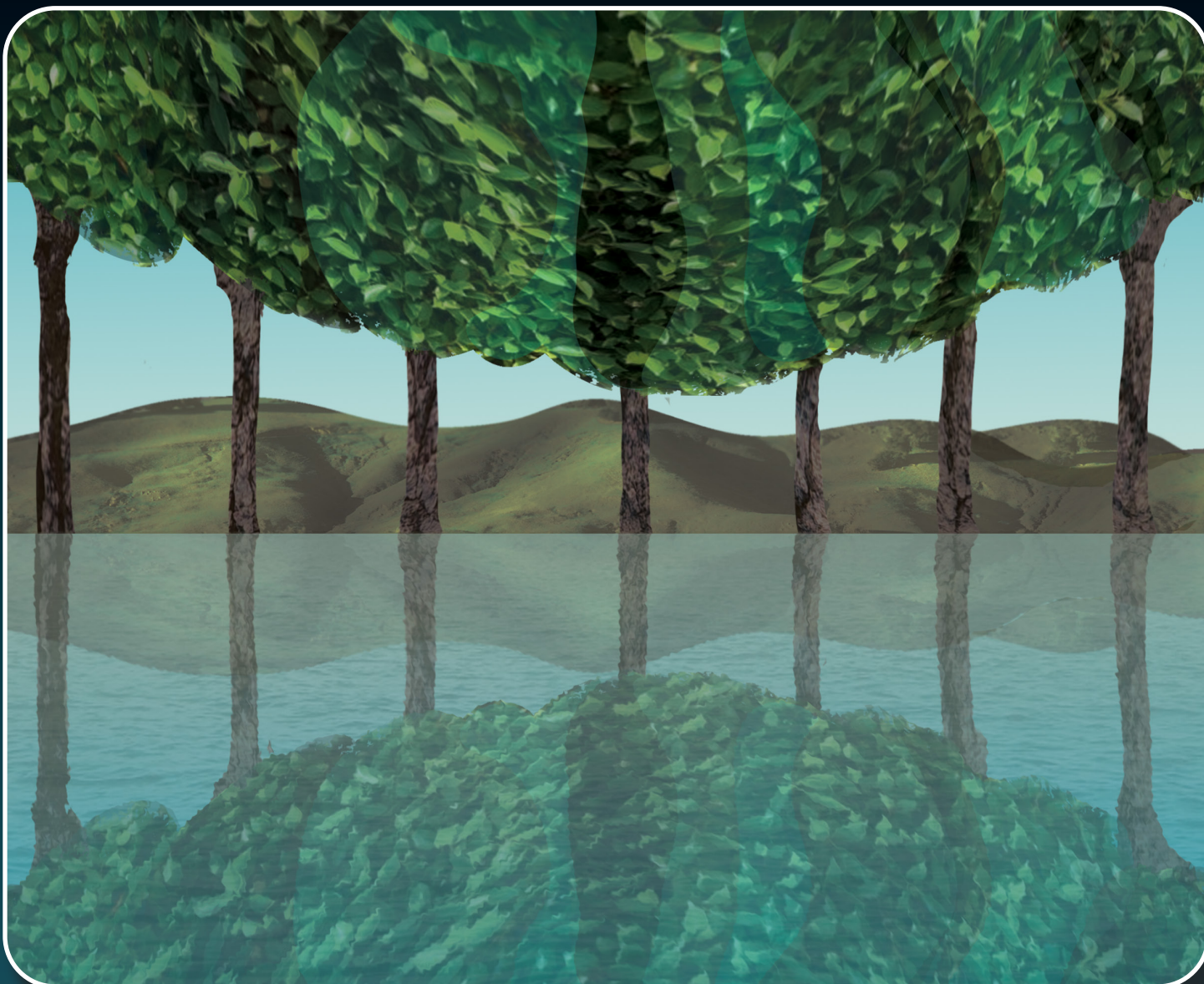


IMMUNOTHERAPY

TOOLS FOR DRUG DISCOVERY

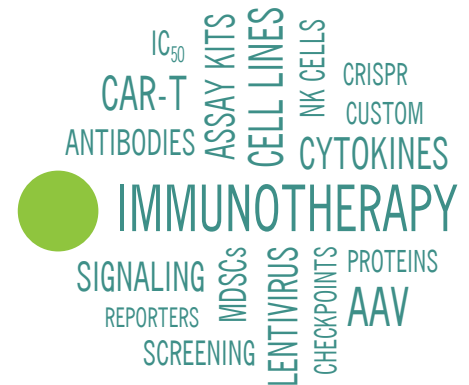
Bioactive Proteins & Antibodies | Biochemical Assay Kits | Cell Lines
Virus-Based Tools | Screening & Profiling



Expanding Immunology Solutions for Cancer

Immunotherapeutic approaches to treat cancer or other diseases holds exceptional promise for eradicating or controlling diseases. Some approaches, including immune checkpoint blockade, CAR-T cells, CAR-NK cells, cytokines, and antibodies have already proven effective in the clinic. Yet more work needs to be done to improve therapeutic effectiveness across a broader range of cancers, reduce harmful side effects, and expand immunotherapy to other diseases.

BPS Bioscience offers a comprehensive selection of immunotherapy products, featuring novel tools, such as our bispecific and trispecific antibodies, bioactive proteins, biochemical and cellular bioassay systems, recombinant cell lines, and virus-based tools, as well as screening and profiling services to test your compounds in our assay systems.



Our Advantages



Produced In-House

- Made in the USA at our San Diego, CA laboratory
- 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 400+ proteins and antibodies, 100+ cell lines, 90+ assay kits, and other quality tools
- Immunology areas: immune checkpoint proteins, CAR-T cells, cytokines, adoptive T cell therapy, MDSCs, NK cells, immune signaling, tumor microenvironment



Immunotherapy Screening and Profiling

- Unique panel of 50+ assays for compound or antibody screening
- Select from IC₅₀ determination and single point concentrations
- Perform follow-up studies with the same proteins manufactured in-house

Bioactive Recombinant Proteins and Antibodies

BPS Bioscience specializes in manufacturing bioactive enzymes and protein complexes, offering over 2000 protein products critical for drug discovery. Our immunotherapy portfolio consists of immune checkpoints, cytokines, activating cell surface receptors, CD molecules, adhesion molecules, and intracellular signaling proteins.

Our antibody portfolio consists of recombinantly produced, fit-for-purpose antibodies for highly specific target identification, protein/receptor neutralization, or activity induction.

Protein options include:



HiP™ Proteins (High Purity, Low Aggregation)

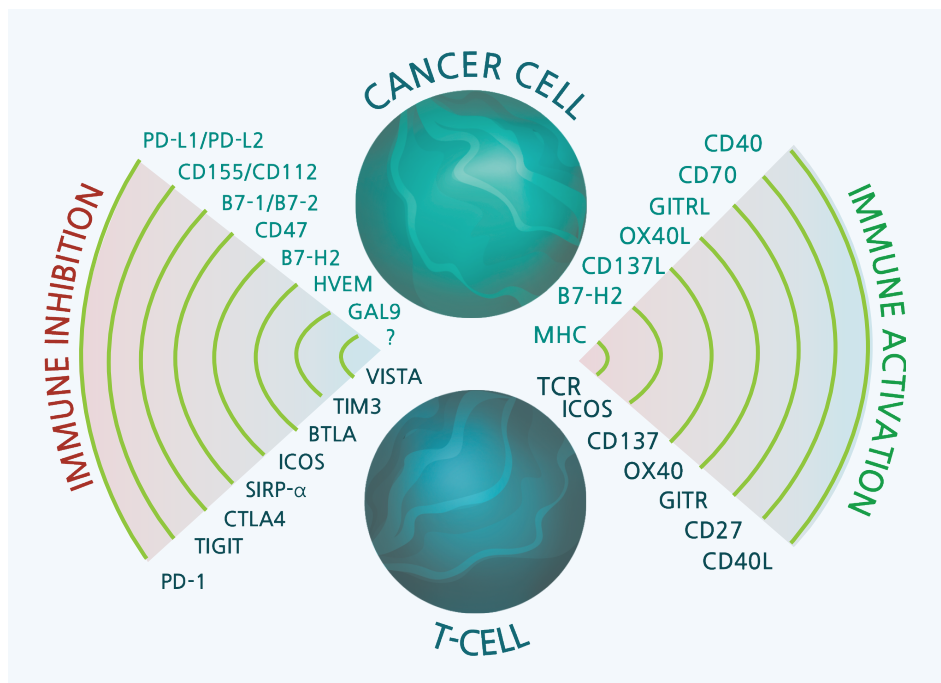
- >90% purity
- Verified via HPLC to guarantee <10% aggregation
- Ideal for studying enzyme activity, ligand binding, protein folding, and more

Bulk Production and Customization

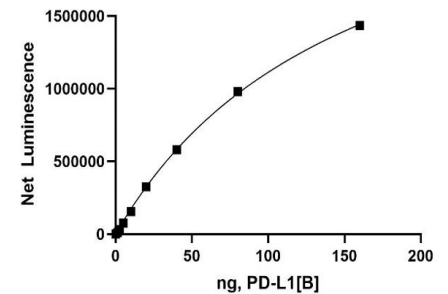
Product sizing, formulations, tags, mutations, and activity testing can be customized to meet your research needs.

Activity Testing

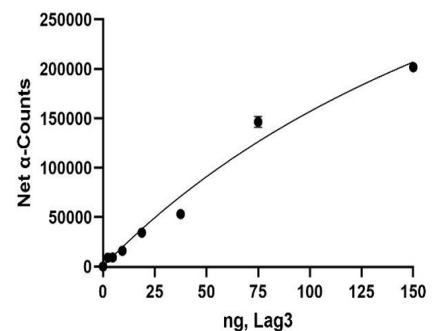
- Our proteins and antibodies are validated on each lot to confirm activity, when relevant.
- In-house developed assay kits use our quality manufactured proteins for consistent performance.



PD-1:PD-L1 Binding Activity



LAG3:FGL1 Binding Activity



Biochemical and Cell-Based Assay Kits

Our 90+ biochemical and cellular assay kits for immunotherapy enable your compound and biotherapeutic drug discovery by providing specific, highly sensitive analysis in a variety of formats. Many of our kits are provided as homogenous assays, with no wash steps, for efficient high throughput screening.

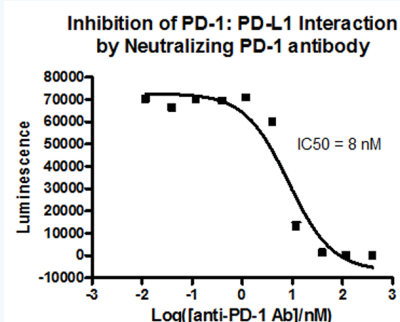
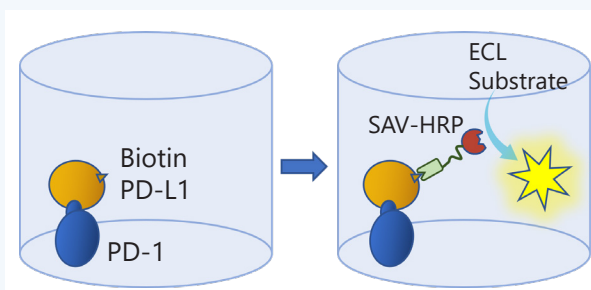
Assay kit formats include:

- AlphaLISA®
- Chemiluminescent
- Colorimetric
- Fluorogenic
- Fluorescence Polarization
- Luminescent
- TR-FRET
- UV Spectrophotometric

Featured Products

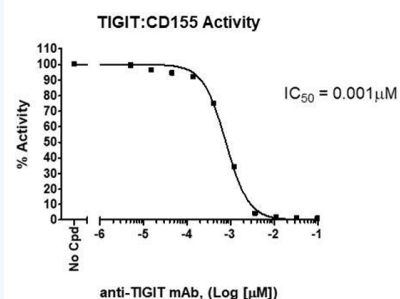
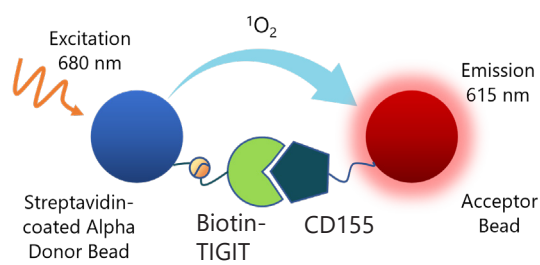
PD-1:PD-L1[Biotinylated] Inhibitor Screening Assay Kit (#72003)

Designed to screen for inhibitors of PD-1:PD-L1 interaction using a chemiluminescent readout



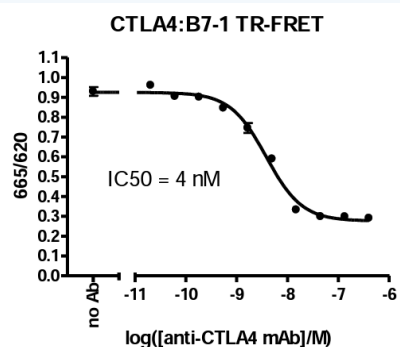
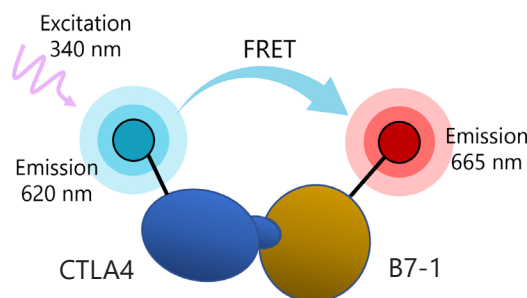
TIGIT:CD155 Homogeneous Assay Kit (#72029)

Designed to screen for inhibitors of TIGIT:CD155 interaction using a homogeneous, no-wash AlphaLISA® 384-well assay system.



CTLA4:B7-1 TR-FRET Assay Kit (#72120)

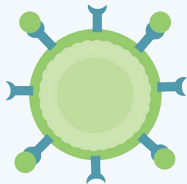
Designed to screen for inhibitors of CTLA4 binding to B7-1 in a homogeneous 384 reaction format.



Cell Lines

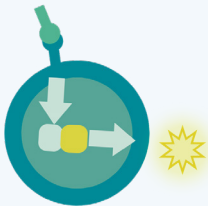
Recombinant cell lines can serve as excellent model systems in a variety of applications for immunotherapy research. Our broad selection of many novel cell line tools offers an excellent set of options for your research.

● Overexpression/Knockout Cell Lines



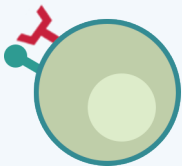
- Antibody screening
- Study ion channels, GPCR screens, and metabolite-based assays
- Flow cytometry analysis
- Binding studies

● Reporter Cell Lines



- Measure signaling activity
- Measure promoter activity
- Measure transcription factor activity
- Use as a control cell line
- Use with our ONE-Step™ Luciferase Assay System

● Co-stimulatory Cell Lines



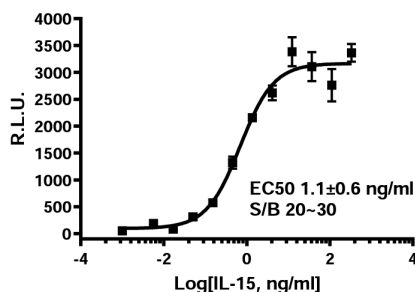
- Inhibitor screening
- Antibody affinity
- Co-culture studies

● CAR-T Cells

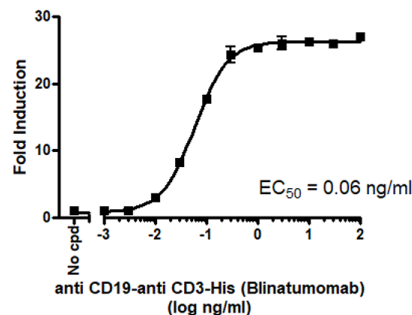


- For design of co-culture studies
- Positive controls for developing new CARs

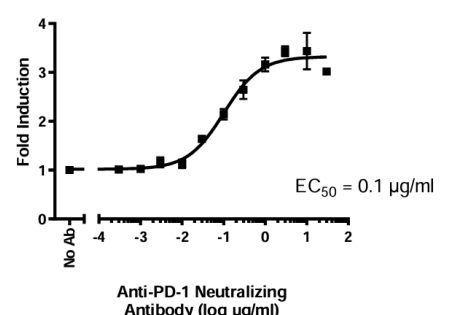
IL-15 agonist stimulation of luciferase activity in IL-15 responsive reporter cell line (#78402)



Bispecific antibody stimulation of NFAT luciferase reporter Jurkat cells (#60621) in the presence of CD19+ Raji cells



PD-L1/TCR activator cells to screen for PD-1 blocking antibodies (#60536)

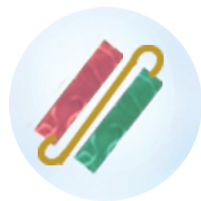


CAR-T Cell Therapy Solutions

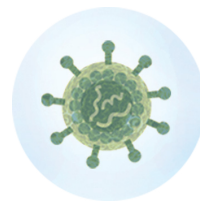
Custom CAR-T Cell Development



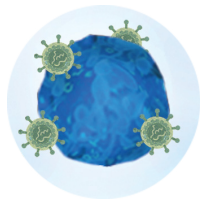
1
Researcher provides Ab sequence against cancer antigen



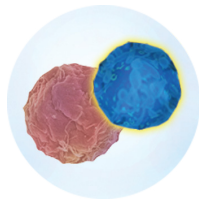
2
Engineering & validation of ScFv for specificity and affinity



3
CAR Lentivirus production and initial validation



4
T cell preparation & transduction



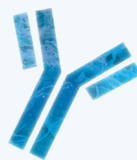
5
Functional validation of CAR-T cells

● Proteins



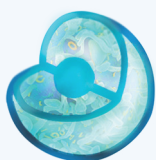
- Functional proteins for binding and blocking assays
- CAR-T target proteins
- Immune checkpoint inhibitor proteins
- Cell activating and co-stimulatory molecules
- Epitope tagged or biotin-labeled

● Antibodies



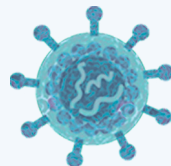
- Bispecific T cell Engager antibodies, including Anti-CD19-Anti-CD3
- CAR-T targets, such as Anti-BCMA, anti-CD19, and more
- T cell agonist antibodies, such as Anti-CD3 and Anti-CD28
- Human Ig isotype controls

● Cell Lines & Primary Cells



- Overexpressing CAR-target cell lines, including BCMA, CD19, and more
- Luciferase reporter cell lines
- TCR, B2M, and CIITA knockout cell lines to model universal CAR-T cells
- Primary CD4+ and CD8+ T cells and PBMCs

● Lentiviruses

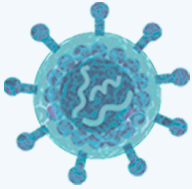


- Capable of transducing primary cells
- For long-term stable expression of a transgene
- CAR lentiviruses for Anti-CD19, Anti-CD20, Anti-BCMA and more
- CRISPR/Cas9 lentiviruses for target protein knockout

Virus-Based Tools

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 viral tools for research in immunotherapy.

Lentiviruses



Lentiviruses are enveloped retroviruses that fuse with the target cell membrane, delivering genetic material into the cytoplasm of the cell. 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.

Applications

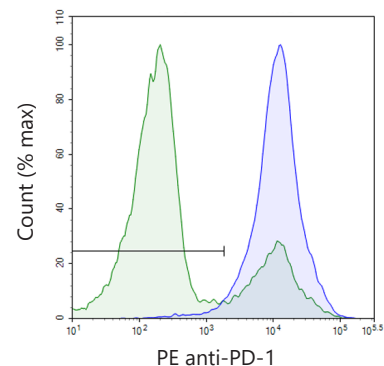
- Stable cell line generation
- Protein expression
- CRISPR/Cas9 knockout
- Generating cellular reporter assays (GFP, luciferase)
- Screen for neutralizing antibodies

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

Knockdown of PD-1 in PD-1 Overexpressing Jurkat cells using PD-1 CRISPR/Cas9 Lentivirus (Non-Integrating) (#78059)

PD-1 NFAT-Reporter Jurkat cells (BPS Bioscience, Cat #60535) were transduced via spinoculation with 5,000,000 TU/well of PD-1 CRISPR/Cas9 lentivirus. 72 hours after transduction, cells were stained with PE anti-human PD-1 antibody and analyzed by flow cytometry. Parental PD-1 NFAT-Reporter Jurkat cells are shown in blue, and the transduced cells are shown in green.

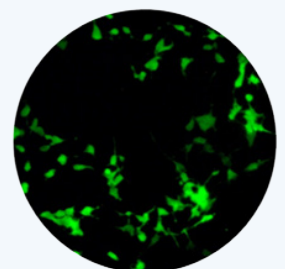


AAV Reporter Vectors



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.



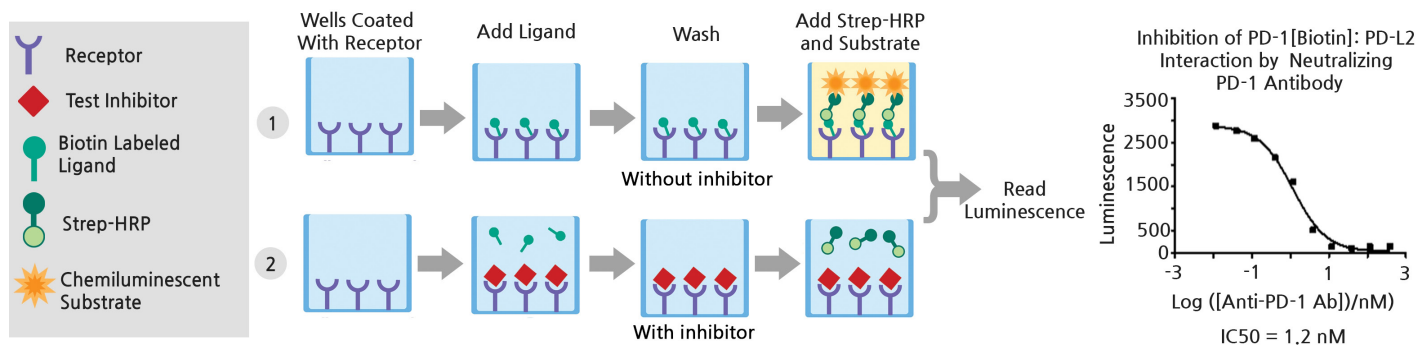
Biochemical & Cell-Based Screening & Profiling

BPS Bioscience offers rapid and dependable Immunotherapy Screening Services to determine IC_{50} values of small molecules, peptides, and antibodies against requested targets. Our team of experts along with our broad services portfolio make it easy to:

- Screen for inhibitors/targets
- Select from IC_{50} determination and single point concentrations
- Receive data within weeks of compound submission
- Perform follow-up studies using the same soluble receptors manufactured in-house
- Get questions answered or project guidance in a time-efficient manner

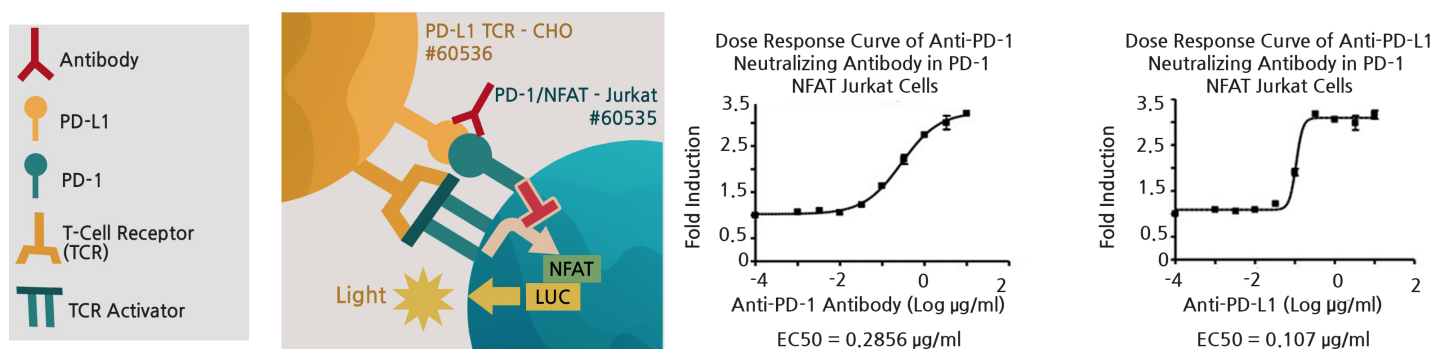
Biochemical Assays

- Over 50 assays to choose from. Featuring CBL-B, PD-1:PD-L2, CD40:CD40L, and many more.



Cell-Based Assays

- Cell research is facilitated by the engineering of target-expressing cells, the design of cell-based assays using reporter target or effector cells, or the design of co-culture assays. These research tools are especially useful in the field of immunotherapeutics.
- Over 15 assays to choose from. Featuring A2AR, BCMA, IDO1, G1TR:G1TR-Ligand and many more.



Antibodies	Catalog#	Antibodies	Catalog#
Anti-RARA polyclonal antibody	25310	Anti-CD38 Antibody	79120
Anti-BCMA Antibody	101219	Anti-CD4 Antibody, Biotin-Labeled	101761
Anti-BCMA Antibody (Single-Chain Variable Fragment), His-Tag	100173	Anti-CD4 Antibody, PE-Labeled	102010
Anti-BCMA-Anti-CD19-Anti-CD3 Trispecific Molecule	100761	Anti-CD47 Antagonist Antibody	79065-1
Anti-BCMA-Anti-CD3 Bispecific Molecule	100689	Anti-CD47 Antagonist Antibody	79065-2
Anti-BCMA-Anti-CD3 IgG Bispecific Antibody	101968	Anti-CD47 Antibody, PE-Labeled	79105-2
Anti-BTLA Neutralizing Antibody	100244	Anti-CD47 Antibody, PE-Labeled	79105-1
Anti-CD123 Antibody, Biotin-Labeled	101733	Anti-CD66b Antibody, Biotin-Labeled	102078
Anti-CD123 Antibody, FITC-Labeled	101955	Anti-CD73 Antagonist Ab (clone #7F3)	100231
Anti-CD123 Antibody, PE-Labeled	101954	Anti-CD73 Antagonist Antibody	79421
Anti-CD123 IgG Antibody	101036	Anti-CD8 Antibody, Biotin-Labeled	101779
Anti-CD137 Agonist Antibody	79097	Anti-CD8 Antibody, FITC-Labeled	102224
Anti-CD14 Antibody, FITC-Labeled	101884	Anti-CD8 Antibody, PE-Labeled	102011
Anti-CD14 Antibody, PE-Labeled	101883	Anti-Claudin-18 Isoform 2 Antibody, FITC-Labeled	101866
Anti-CD16a Antibody, FITC-Labeled	101979	Anti-Claudin-18 Isoform 2 Antibody, PE-Labeled	101676
Anti-CD16a Antibody, PE-Labeled	101978	Anti-Claudin-18 Isoform 2 IgG Antibody	101564
Anti-CD19 Antibody, FITC-Labeled	101863	Anti-Claudin-18 Isoform 2 IgG Antibody, Biotin-Labeled	101565
Anti-CD19 Antibody, PE-Labeled	101625	Anti-Claudin-18 Isoform2-Anti-CD3 IgG Bispecific Antibody	101541
Anti-CD19 IgG Antibody	100981	Anti-CTLA4 Neutralizing Antibody	71212
Anti-CD19 IgG Antibody, Biotin-Labeled	101093	Anti-FcGR3A (CD16A) IgG Antibody, Biotin-Labeled	101225
Anti-CD19-Anti-CD3 Bispecific Molecule	100441	Anti-FcGR3A-Anti-EGFR Bispecific Antibody	101535
Anti-CD19-Anti-CD3 IgG format Bispecific Antibody	101076	Anti-GITR Agonist Antibody	79053
Anti-CD2 IgG Antibody	101213	Anti-GITR Antibody, PE-labeled	71295-1
Anti-CD20 Antibody, FITC-Labeled	101864	Anti-GITR Antibody, PE-labeled	71295-2
Anti-CD20 Antibody, PE-Labeled	101672	Anti-HLA-DR Biotin-Labeled, Antibody	101769
Anti-CD20 Functional Antibody	71209	Anti-Human IgG, Unconjugated Antibody	100736
Anti-CD20 IgG Antibody, Biotin-labeled	101207	Anti-Human IgM, Unconjugated Antibody	100737
Anti-CD20-Anti-CD3 Bispecific Antibody	100836	Anti-IL-17A Neutralizing Antibody	91015
Anti-CD20-Anti-CD3 IgM format Bispecific Antibody	100860	Anti-IL-2RA (CD25) Neutralizing Antibody	101593
Anti-CD200 Neutralizing Antibody	102062	Anti-IL23 α Neutralizing Antibody	101453
Anti-CD235a Antibody, PE-Labeled	101675	Anti-LAG3 Antibody, PE-Labeled	71226-1
Anti-CD235a IgG1 Antibody, Biotin-Labeled	101098	Anti-LAG3 Antibody, PE-Labeled	71226-2
Anti-CD27 Agonist Antibody	100111	Anti-LAG3, Neutralizing Antibody	71219
Anti-CD28 Agonist Antibody	100182	Anti-Mesothelin-Anti-CD3 Bispecific Antibody	101621
Anti-CD28 Agonist Antibody (Humanized)	100186	Anti-NCAM1 (CD56) IgG Antibody, Biotin-labeled	101112
Anti-CD3 Agonist Antibody	71274-1	Anti-NCAM1 Antibody, FITC-Labeled	101865
Anti-CD3 Agonist Antibody	71274-2	Anti-NCAM1 Antibody, PE-Labeled	101673
Anti-CD3 Antibody, FITC-Labeled	102008	Anti-Notch Neutralizing Antibody	102065
Anti-CD3 Antibody, PE-Labeled	102009	Anti-OX40 Antagonist Antibody	72063-1

Antibodies	Catalog#	Biochemical Assay Kits	Catalog#
Anti-OX40 Antagonist Antibody	72063-2	ARG1 Inhibitor Screening Assay Kit	72048
Anti-OX40 Antibody, PE-Labeled	72064-2	ARG2 Inhibitor Screening Assay Kit	72043
Anti-OX40 Antibody, PE-Labeled	72064-1	BAFF:BCMA[Biotinylated] Inhibitor Screening Assay Kit	79667
Anti-OX40 Competitive Antibody	72558-1	BCMA:APRIL[Biotinylated] Inhibitor Screening Assay Kit	79722
Anti-OX40 Competitive Antibody	72558-2	BLM Helicase Activity Assay Kit	82512
Anti-PD-1 (S228P) Neutralizing Antibody	71326	BTLA:HVEM[Biotinylated] Inhibitor Screening Assay Kit	72008
Anti-PD-1 Agonist Antibody	101178	Cathepsin D Inhibitor Screening Assay Kit	82141
Anti-PD-1 Neutralizing Antibody	71120	CD112R:CD112[Biotinylated] Inhibitor Screening Assay Kit	79732
Anti-PD-1 Neutralizing Antibody, Biotin-labeled	71215	CD137:CD137L[Biotinylated] Inhibitor Screening Assay Kit	78590
Anti-PD-1 Neutralizing Antibody, PE-labeled	71290	CD137[Biotinylated]:CD137L Inhibitor Screening Assay Kit	72025
Anti-PD-L1 (CD274) Neutralizing Antibody	71213	CD155:TIGIT [Biotinylated] Inhibitor Screening Chemilumescence Assay Kit	82526
Anti-PD-L1 Antibody, PE-labeled	71128	CD200:CD200R1 [Biotinylated] Inhibitor Screening Chemiluminescence Assay Kit	82285
Anti-PD-L1 Neutralizing Antibody, Biotin-Labeled	71214	CD226:CD112 Homogeneous Assay Kit	72051
Anti-PD-L1-Fab Avi-Tag, Biotin-Labeled	11032	CD226:CD155 Homogeneous Assay Kit	72052
Anti-PSMA Antibody	101695	CD27:CD70[Biotinylated] Inhibitor Screening Assay Kit	79695
Anti-PSMA Antibody, FITC-Labeled	101977	CD28:B7-1[Biotinylated] Inhibitor Screening Assay Kit	72007
Anti-PSMA Antibody, PE-Labeled	101976	CD28:B7-2[Biotinylated] Inhibitor Screening Assay Kit	72062
Anti-PSMA, Biotin Label Antibody	101757	CD38 (Dog) Inhibitor Screening Assay Kit (Hydrolase Activity)	78108
Anti-PSMA-Anti-CD3 IgG format Bispecific Antibody	101242	CD38 (Mouse) Fluorogenic Assay Kit (Cyclase Activity)	78285
Anti-PVRIG Neutralizing Antibody	101712	CD38 (Mouse) Inhibitor Screening Assay Kit (Hydrolase Activity)	79682
Anti-RANKL Neutralizing Antibody	100874	CD38 (Pig) Fluorogenic Assay Kit (Hydrolase Activity)	78178
Anti-TFR Antibody, Biotin-Labeled	101596	CD38 (Rat) Inhibitor Screening Assay Kit (Hydrolase Activity)	79690
Anti-TIGIT Antibody, PE-Labeled	71228-1	CD38 Inhibitor Screening Assay Kit (Cyclase Activity)	71275
Anti-TIGIT Antibody, PE-Labeled	71228-2	CD38 Inhibitor Screening Assay Kit (Hydrolase Activity)	79287
Anti-TIGIT Inhibitor Antibody	71218	CD38 Inhibitor Screening Assay Kit (Hydrolase Activity)	79672
Anti-TIGIT Neutralizing Antibody	71340	CD39 Inhibitor Screening Assay Kit	79278
Anti-TIM-3 Antibody, PE-Labeled	71121	CD40:CD40L TR-FRET Assay	79258
Anti-Tim3 Antibody	79098	CD40:CD40L[Biotinylated] Inhibitor Screening Assay Kit	79257
Anti-TL1A Neutralizing Antibody	101729	CD47:SIRP- α [Biotinylated] Inhibitor Screening Assay Kit	72044
Anti-Transferrin Receptor Antibody, FITC-Labeled	101876	CD47:SIRP- γ [Biotinylated] Inhibitor Screening Assay Kit	72059
Anti-Transferrin Receptor Antibody, PE-Labeled	101875	CD73 Inhibitor Screening Assay Kit	72055
Anti-TSLP Neutralizing Antibody	102138	CD73 Inhibitor Screening Assay Kit	72058
Anti-VEGF Neutralizing Antibody	79478-1	Chemi-Verse™ BLK Kinase Assay Kit	82229
Anti-VEGF Neutralizing Antibody	79478-2	Chemi-Verse™ CDK1/CyclinA2 Kinase Assay Kit	82242
Anti-VISTA Neutralizing Antibody	101293	Chemi-Verse™ CDK12/CyclinK Kinase Assay Kit	82295
		Chemi-Verse™ CDK2/CyclinA2 Kinase Assay Kit	82247
		Chemi-Verse™ CDK2/CyclinE1 Kinase Assay Kit	82228
		Chemi-Verse™ CDK9/CyclinT Kinase Assay Kit	82233
		Chemi-Verse™ LYN A Kinase Assay Kit	82224
Biochemical Assay Kits	Catalog#		
Activin A: Activin RII[Biotinylated] Inhibitor Screening Assay Kit	82531		
ADAR1:RNA TR-FRET Assay Kit	82252		

Biochemical Assay Kits	Catalog#	Biochemical Assay Kits	Catalog#
Chemi-Verse™ MAPKAPK5 Kinase Assay Kit	82236	LysA™ Universal PARYlation Assay Kit	82123
Chemi-Verse™ NUA1 Kinase Assay Kit	82231	MAT2A Inhibitor Screening Assay Kit	71402
Chemi-Verse™ NUA2 Kinase Assay Kit	82232	MDM2-Driven p53 Ubiquitination Assay Kit	82179
Chemi-Verse™ PAK6 Kinase Assay Kit	82182	MLL1 (KMT2A): WDR5 Binding Chemiluminescent Assay Kit	82502
Chemi-Verse™ PAK7 Kinase Assay Kit	82183	MLL4 (KMT2B): WDR5 Binding Chemiluminescent Assay Kit	82503
Chemi-Verse™ PIM1 Kinase Assay Kit	82525	MMP7 Fluorogenic Assay Kit	78864
Chemi-Verse™ VRK1 Kinase Assay Kit	82230	Molecular Glue/PROTAC® Optimization Kit for RBM39-DCAF15	82251
CTLA4 : B7-2[Biotinylated] Inhibitor Screening Assay Kit	79658	Monkey PD-1:PD-L1[Biotinylated] Inhibitor Screening Assay Kit	79863
CTLA4 [Biotinylated]:B7-1 Inhibitor Screening Assay Kit	79973	Monkey PD-1[Biotinylated]:PD-L1 Inhibitor Screening Assay Kit	79865
CTLA4:B7-1 TR-FRET Assay Kit	72120	Mouse CTLA4:B7-1 TR-FRET Assay	79508
CTLA4:B7-1[Biotinylated] Inhibitor Screening Assay Kit	72009	Mouse CTLA4[Biotin]:B7-1 Inhibitor Screening Assay Kit	79515
CTLA4[Biotinylated]:B7-2 Inhibitor Screening Assay Kit	72024	Mouse IDO1 Inhibitor Screening Assay Kit	72041
Cytotoxicity Dye Kit (CFSE, 7-ADD)	82296	Mouse IDO2 Inhibitor Screening Assay Kit	72042
DcR3:TL1A Inhibitor Screening Assay Kit	82160	Mouse ITK Kinase Assay Kit	78520
DR3:TL1A Inhibitor Screening Assay Kit	82241	Mouse PD-1[Biotinylated]:PD-L1 Inhibitor Screening Assay Kit	72027
ERK1 Kinase Assay Kit	82282	NAMPT Inhibitor Screening Assay Kit	71276
Fc (IgG1): FcRn Inhibitor Screening Colorimetric Assay Kit	78501	NEU2 Fluorogenic Assay Kit	82139
FGL1:LAG3 Inhibitor Screening Assay Kit	78824	Notch1:DLL4 [Biotinylated] Inhibitor Screening Chemiluminescence Assay Kit	82284
FGL1:LAG3 TR-FRET Assay Kit	79739	OX40[Biotinylated]:OX40L Inhibitor Screening Assay Kit	72045
GDF15:GFRAL[Biotinylated] Inhibitor Screening Assay Kit	78807	PADI4 Fluorogenic Assay Kit	82111
GITR:GITRL TR-FRET Assay Kit	79054	PARP1 Olaparib Competitive Inhibitor Assay Kit	82293
GITRL:GITR[Biotinylated] Inhibitor Screening Assay Kit	72061	PARP2 Olaparib Competitive Inhibitor Assay Kit	82294
GPX4 Fluorogenic Assay Kit	82165	PD-1:PD-L1 Cell-Based Inhibitor Screening Assay Kit	79377
HSA:FcRn Neutralizing Antibody Screening Chemiluminescence Assay Kit	82151	PD-1:PD-L1 Homogeneous Assay Kit	72014
ICOSL:ICOS[Biotinylated] Inhibitor Screening Assay Kit	79673	PD-1:PD-L1 TR-FRET Assay	72032
IDO1 Fluorogenic Inhibitor Screening Assay Kit	72037	PD-1:PD-L1 TR-FRET Assay	72038
IDO1 Fluorogenic Inhibitor Screening Assay Kit -384	72047	PD-1:PD-L1/PD-L2 Cell-Based Inhibitor Screening Assay Kit	60800
IDO1 Inhibitor Mechanism of Action Assay Kit	72050	PD-1:PD-L1[Biotinylated] Inhibitor Screening Assay Kit	72003
IDO1 Inhibitor Screening Assay Kit	72021	PD-1:PD-L1[Biotinylated] Inhibitor Screening Colorimetric Assay Kit	72016
IDO1 Inhibitor Screening Assay Kit -384	72034	PD-1:PD-L2 Cell-Based Inhibitor Screening Assay Kit	79378
IDO2 Inhibitor Screening Assay Kit	72022	PD-1:PD-L2 Homogeneous Assay Kit	72015
IL-17RA:IL-17A[Biotin] Inhibitor Screening Assay Kit	79891	PD-1:PD-L2 TR-FRET Assay Kit	72012
IL-17RA[Biotin]:IL-17A Inhibitor Screening Assay Kit	72060	PD-1:PD-L2[Biotinylated] Inhibitor Screening Assay Kit	72004
IL-6:IL-6R Inhibitor Screening Assay Kit	78027	PD-1:PD-L2[Biotinylated] Inhibitor Screening Colorimetric Assay Kit	72017
KMO Inhibitor Screening Assay Kit	79513	PD-1[Biotinylated]: PD-L1 Inhibitor Screening Assay Kit	79523
Kynureninase Inhibitor Screening Assay Kit	79514	PD-1[Biotinylated]:PD-L1 Homogeneous Assay Kit	72028
LAG3:FGL1[Biotinylated] Inhibitor Screening Assay Kit	79789	PD-1[Biotinylated]:PD-L1 Inhibitor Screening Assay Kit	72005
LIGHT:HVEM [Biotinylated] Inhibitor Screening Assay Kit	79684	PD-1[Biotinylated]:PD-L1 Inhibitor Screening Colorimetric Assay Kit	72018
LTβR:LIGHT[Biotinylated] Inhibitor Screening Assay Kit	79808	PD-1[Biotinylated]:PD-L2 Inhibitor Screening Assay Kit	72006

Biochemical Assay Kits, Cell Isolation Kits/Components, | Product Listing

Cell Lines

Biochemical Assay Kits	Catalog#	Cell Lines	Catalog#
PD-1[Biotinylated]:PD-L2 Inhibitor Screening Colorimetric Assay Kit	72019	ADCC Bioassay Effector Cell F Variant (Low Affinity)/ NFAT Luciferase Reporter Jurkat Cell Line	60540
PD-L1:B7-1 [Biotinylated] Inhibitor Screening Assay Kit (Mouse)	79821	ADCC Bioassay Effector Cell V Variant (High Affinity)/ NFAT Luciferase Reporter Jurkat Cell Line	60541
PD-L1:B7-1[Biotinylated] Inhibitor Screening Assay Kit	72026	ADCP Bioassay Effector Cell FcγRIIa (H Variant)/ NFAT Luciferase Reporter Jurkat Cell Line	71273
RANK:RANKL TR-FRET Assay	79101	Adenosine A2A Receptor Functional HEK293 Cell Line	79381
Sialidase Fluorogenic Assay Kit	82154	Androgen Luciferase Reporter 22RV1 Cell Line	78972
TDO Fluorogenic Inhibitor Screening Assay Kit -384 (Human)	72049	Anti-BCMA CAR /NFAT (Luciferase) Reporter Jurkat Cell Line	79694
TDO Fluorogenic Inhibitor Screening Assay Kit (Human)	72039	Anti-CD19 CAR / NFAT (Luciferase) Reporter Jurkat Cell Line (CD19 SCFV-CD28-4-1BB-CD3ζ)	79853
TDO Inhibitor Screening Assay Kit	72023	Anti-CD19 CAR Negative Control/NFAT (Luciferase) Reporter Jurkat Cell Line (CD19 SCFV-CD28 Transmembrane Motif)	79854
TDO Inhibitor Screening Assay Kit -384	72036	B2M Knockout Jurkat Cell Line	78342
Tie2 Kinase Assay Kit	78573	B2M Knockout NFAT Luciferase Reporter Jurkat Cell Line	78363
TIGIT (Mouse): CD155 (Mouse) Homogeneous Assay Kit	78141	B2M Knockout THP-1 Cell Line	78389
TIGIT:CD112 Homogeneous Assay Kit	72030	B2M/CIITA Double Knockout THP-1 Cell Line	78391
TIGIT:CD155 Homogeneous Assay Kit	72029	B7-H3 CHO Cell Line (High, Medium, or Low Expression)	78541
TNFR2:TNF-alpha[Biotinylated] Inhibitor Screening Assay Kit	79756	B7-H4 CHO Cell Line (Low, Medium or High Expression)	78586
TPH1 Inhibitor Screening Assay Kit	72053	B7-H7 (HLA2)/TCR Activator CHO Cell Line	78321
TPH1 Inhibitor Screening Assay Kit	72056	BAFF-R – CHO K1 Recombinant Cell Line (High, Medium or Low Expression)	79921
TPH2 Inhibitor Screening Assay Kit	72054	BCMA / CD20 / Firefly Luciferase CHO Cell Line	78185
TPH2 Inhibitor Screening Assay Kit	72057	BCMA / GLuc - CHO Recombinant Cell Line	79830
TSLPR:TSLP [Biotinylated] Inhibitor Screening Chemiluminescence Assay Kit	82507	BCMA / Luciferase - CHO Recombinant Cell Line	79724
Universal IDO1/IDO2/TDO Inhibitor Screening Assay Kit	72035	BCMA / NF-κB – Reporter HEK293 Recombinant Cell Line	79755
VSIG-3:VISTA [Biotinylated] Inhibitor Screening Assay Kit	79782	BCMA CHO Recombinant Cell Line (High or Low Expression)	79500
WRN Helicase Activity Assay Kit	78852	CD123 (IL3Ra) - CHO Recombinant Cell Line (High, Medium, or Low Expression)	79640
		CD137 (4-1BB) /NF-κB Reporter - HEK293 Recombinant Cell Line	79289
		CD137 (4-1BB) HEK293 Recombinant Cell Line	60691
		CD137L (4-1BBL) CHO-K1 Recombinant Cell Line	60523
		CD14 CHO Recombinant Cell Line	78335
		CD155 (PVR) - HEK293 Recombinant Cell Line	60537
		CD155 / TCR Activator - CHO Recombinant Cell line	60548
		CD160/NFAT - Luciferase Reporter - Jurkat Recombinant Cell Line	79594
		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
		CD2 CHO Cell Line	78330
		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
		CD27 CHO-K1 Stable Recombinant Cell Line	60624

Cell Lines	Catalog#	Cell Lines	Catalog#
CD27/NF-κB Reporter-Jurkat Recombinant Cell Line	79509	FAP- CHO K1 Recombinant Cell Line (High, Medium or Low Expression)	79947
CD33- CHO K1 Recombinant Cell Line (High, Medium and Low Expression)	79935	FcGR1a (CD64) Knockout THP-1 Cell Line	82191
CD34 CHO Recombinant Cell Line	78336	FCGR2A (CD32A) Knockout Jurkat Cell Line	78549
CD36 HEK293 Cell Line (High or Medium Expression)	78526	FcGR2B CHO K1 Recombinant Cell Line	79511
CD36 Knockout THP-1 Cell Line	82190	FcGR3A (CD16A) CHO Cell Line	78332
CD37 CHO Recombinant Cell Line (High or Low Expression)	79607	FcGR3B (CD16B) CHO Cell Line	78333
CD38 / BCMA / Firefly Luciferase CHO Recombinant Cell Line	78148	FcRL5 CHO Cell Line	78375
CD38 / CD19 / Firefly Luciferase CHO Recombinant Cell Line	78149	FcRL5 HEK293 Cell Line	78374
CD38 CHO Recombinant Cell Line (High, Medium or Low Expression)	79615	FcRL5/Firefly Luciferase RPMI 8226 Cell Line	78378
CD40 - HEK293 Cell Line	71257	FcRn: IgG Recycling HMEC-1 Cell Pool	82163
CD40/NF-κB Luciferase Reporter HEK293 Cell Line	60626	Firefly Luciferase CCRF-CEM Cell Line	78495
CD40L K562 Cell Line (Medium and High Expression)	78942	Firefly Luciferase CD19 Knockout NALM6 Cell Line	82168
CD43 CHO Cell Line	78334	Firefly Luciferase HL-60 Cell Line	78410
CD47 - HEK293 Cell Line	71249	Firefly Luciferase Jurkat Cell Line	78373
CD5- CHO K1 Recombinant Cell Line (High, Medium, or Low Expression)	78290	Firefly Luciferase K562 Cell Line	78621
CD7 CHO Cell Line (Medium or High Expression)	78324	Firefly Luciferase KG1 Cell Line	78493
CD70-CHO Recombinant Cell line	79510	Firefly Luciferase MC38 Cell Line	78383
CD8+ TCR Knockout NFAT-Luciferase Reporter Jurkat Cell Line	78757	Firefly Luciferase Molm13 Cell Line	78372
CD8a/CD8b Jurkat Cell Line	78917	Firefly Luciferase NALM6 Cell Line	78494
CD95 CHO Cell Line (Medium or High Expression)	78499	Firefly Luciferase Raji Cell Line	78622
CEACAM5 CHO Cell Line (High, Medium, or Low Expression)	78704	Firefly Luciferase SKOV-3 Cell Line	78425
CEACAM6 CHO Cell Line	78550	Firefly Luciferase THP-1 Cell Line	78409
Claudin-1 CHO Cell Line	82097	Firefly Luciferase - RPMI 8226 Recombinant Cell Line	79834
Claudin-18 Isoform 1 CHO Cell Line	78361	FOLR1 CHO Cell Line	78360
Claudin-18 Isoform 2 CHO Cell Line (High, Medium, or Low Expression)	78533	Full Length Notch1/CSL Luciferase Reporter HEK293 Cell Line	82201
Claudin-3 CHO Cell Line	78566	FUT4-CD15 CHO Cell Line	78417
Claudin-4 CHO Cell Line	78567	GITR / NF-κB Luciferase Reporter - Jurkat Cell Line	60546
Claudin-6 CHO Cell Line (High, Medium, or Low) Expressions	78527	GITR - HEK293 Recombinant Cell Line	79092
Claudin-9 CHO Cell Line	78687	GITRL CHO-K1 Recombinant Cell Line	60547
CTLA4 - HEK293 Recombinant Cell Line	60681	Glycophorin A (CD235a) CHO Cell Line	78353
CTLA4 / IL-2 Reporter - Jurkat Recombinant Cell Line	79525	GPC3 - CHO-K1 Recombinant Cell Line	78100
DLL3 CHO Cell Line	78882	GPRC5D (Cynomolgus) CHO Cell Line	78338
DLL4 CHO Cell Line	82217	GPRC5D (Cynomolgus) HEK293 Cell Line	78346
eGFP Daudi Cell Line	82297	GPRC5D CHO Cell Line	78337
eGFP/Firefly Luciferase MM.1S Cell Line	78376	GPRC5D HEK293 Cell Line	78345
eGFP/Firefly Luciferase OVCAR3 Cell Line	78953	Growth Arrested PD-1 / NFAT Reporter Jurkat Cell Line	79687
eGFP/Firefly Luciferase Raji Cell Line	78916	Growth-Arrested NK Feeder Cells	78912
Empty vector control - CHO-K1 Recombinant Cell line	60545	HER2 (ERBB2) CHO Recombinant Cell Line (High, Medium, or Low Expression)	79612
EpCAM CHO Cell Line	78683	HER3 CHO Cell Line	82299

Cell Lines	Catalog#	Cell Lines	Catalog#
HLA-C*08:02 K562 Cell Line	78974	p53 Luciferase Reporter HCT116 Cell Line	78681
Human EGFRvIII CHO K1 Cell Line (High or Low Expression)	78145	PAI-1 Reporter (Luc) - Mv1 Lu Cell Line	60544
Human Mesothelin - CHO-K1 Recombinant Cell Line	78132	PARG Knockout HeLa Cell Line	82171
HVEM - CHO Recombinant Cell Line	79297	PARP1 Knockout HeLa Cell Line	82169
HVEM - HEK293 Recombinant Cell Line	79313	PD-1 (Mouse) / NFAT - Reporter - Jurkat Recombinant Cell Line	79762
ICOS/NFAT Reporter-Jurkat Recombinant Cell Line	79668	PD-1 (Woodchuck) / NFAT - Reporter - Jurkat Recombinant Cell Line	79456
ICOSL-CHO Recombinant Cell Line	79635	PD-1 - HEK293 Recombinant Cell Line	60680
IDO1-HEK293 Recombinant Cell Line (Human)	60532	PD-1 / NFAT Reporter - Jurkat Recombinant Cell Line	60535
IDO2 - HEK293 Recombinant Cell Line (Human)	60533	PD-1 CHO Cell Line	78530
IL-15 Responsive Luciferase Reporter Cell Line	78402	PD-1/IL-2 Luciferase Reporter Jurkat Cell Line	82117
IL-2 Luciferase Reporter Jurkat Cell Line	60481	PD-L1 (Mouse)/TCR Activator CHO Cell Line	79763
IL-4/IL-13 Responsive STAT6 Luciferase Reporter HEK293 Cell Line	78941	PD-L1 (Woodchuck) / TCR activator - CHO Recombinant Cell line	79457
IL-6 Responsive Luciferase Reporter HEK293 Cell Line	82235	PD-L1 - CHO Recombinant Cell Line	60543
KIR3DL3/IL-2 Luciferase Reporter Jurkat Cell Line	78322	PD-L1 / TCR Activator - CHO Recombinant Cell Line	60536
KRAS G12D TCR (Clone 10) CD8+ NFAT-Luciferase Reporter Jurkat Cell Line	82303	PD-L2 / TCR activator - CHO Recombinant Cell Line	79632
KRAS G12D TCR (Clone 9c) CD8+ NFAT-Luciferase Reporter Jurkat Cell Line	82304	PSMA (FOLH1) - CHO Recombinant Cell Line (High, Medium, or Low Expression)	79641
LAG3 / IL-2 Reporter - Jurkat Recombinant Cell Line	79813	RANK/NF-κB Luciferase Reporter HEK293 Cell Line	78102
LAG3 / NFAT Reporter - Jurkat Recombinant Cell Line	71278	ROR1 CHO Recombinant Cell Line (High, Medium, or Low Expression)	79609
LAIR1 CHO Cell Line	78913	SIRP-α / HEK293 Recombinant Cell Line	60689
LIGHT-CHO Recombinant Cell Line	79262	SLAMF7 (CS1) CHO Recombinant Cell Line (High, Medium, or Low Expression)	79608
LIGHT-Responsive HVEM/NF-κB Reporter Jurkat Cell Line	79310	STAT3 Luciferase Reporter THP-1 Cell Line	78498
LILRB4 NFAT-Luciferase Reporter Jurkat Recombinant Cell Line	79750	STAT3 Reporter Jurkat Cell Line	78497
MAGE-A1 TCR CD8+ NFAT-Luciferase Reporter Jurkat Cell Line	78993	TCR Activator CHO Recombinant Cell line	60539
MAGE-A4 TCR CD8+ NFAT-Luciferase Reporter Jurkat Cell Line	78984	TCR Activator Raji Cell Line	60556
MART-1 TCR (DMF5) CD8+ NFAT-Luciferase Reporter Jurkat Cell Line	78773	TCR Activator/FcGR2B CHO Cell Line	78436
Membrane-Bound TNFα CHO Cell Line	78971	TCR Knockout Jurkat Cell Line	78539
MUC16 (CA125), variant 4 (region 13785-14507) CHO Cell Line	78848	TCR Knockout NFAT-Luciferase Reporter Jurkat Cell Line	78556
NCAM1 (CD56) CHO Cell Line (High, Medium, or Low Expression)	78352	TCR/B2M Knockout Jurkat Cell Line	78552
Nectin4 - CHO K1 Recombinant Cell Line (High, Medium, or Low Expression)	78097	TCR/B2M Knockout NFAT Luciferase Reporter Jurkat Cell Line	78557
NF-κB Luciferase Reporter Jurkat Cell Line	60651	TDO-HEK293 Recombinant Cell line	60534
NFAT Luciferase Reporter Jurkat Cell Line	60621	TIGIT - HEK293 Recombinant Cell Line	79332
NFAT Luciferase-eGFP Reporter Jurkat Cell Line	78662	TIGIT / NFAT Reporter - Jurkat Cell Line	60538
NFAT Reporter (eGFP) Jurkat Recombinant Cell Line	78384	TL1A-Responsive Luciferase Reporter Jurkat Cell Line	78811
NKp46 CHO Cell Line (High, Medium, or Low Expression)	78915	TNFR2 HEK293 Cell Line	78828
NLRP3 Knockdown THP-1 Cell Pool	82121	Transferrin Receptor Protein-1 (TFRC) HEK293 Cell Line	78308
NY-ESO-1 TCR (1G4) CD8+ NFAT-Luciferase Reporter Jurkat Cell Line	78769	TROP2 - CHO-K1 Recombinant Cell Line	78099
NY-ESO-1 TCR (c259) CD8+ NFAT-Luciferase Reporter Jurkat Cell Line	78771	VSIG4 CHO Cell Line	78914
OX40 - HEK293 Recombinant Cell Line	60682		
OX40/NF-κB Luciferase Reporter HEK293 Cell Line	60482		

Cell-Based Assays and Expression Kits	Catalog#	Lentiviruses	Catalog#
CSF1R / SRE Reporter Kit (MAPK/ERK Signaling Pathway)	79379	BCMA CRISPR/Cas9 Lentivirus (Integrating)	78893
IDO1 Cell-Based Assay Kit	72031	BCMA CRISPR/Cas9 Lentivirus (Non-Integrating)	78894
Mouse PD-L1 / TCR Activator Mammalian Expression Kit	79778	BCMA Lentivirus	78714
PBMC Cytotoxicity Bioassay Kit (CFSE, 7-ADD)	82173	CD19 Lentivirus	78657
PBMC Cytotoxicity Luciferase Assay Kit	82214	CD20 Lentivirus	78658
PBMC Cytotoxicity Luciferase Assay Kit (NALM6)	82174	CD22 Lentivirus	78659
PD-L1 (Woodchuck) / TCR Activator Mammalian Expression Kit	79455	CD40 Ligand (CD40L) Lentivirus	78931
PD-L1 / TCR Activator Mammalian Expression Kit	60610	CD47 CRISPR/Cas9 Lentivirus (Integrating)	78056
PD-L2/TCR Activator Mammalian Expression Kit	60620	CD47 CRISPR/Cas9 Lentivirus (Non-Integrating)	78063
TDO Cell-Based Assay Kit	72033	CD5 (Human) CRISPR/Cas9 Lentivirus (Integrating)	78119
		CD5 (Human) CRISPR/Cas9 Lentivirus (Non-Integrating)	78198
		CD8a Lentivirus	78648
		CD8a/CD8b Lentivirus	78650
		CEACAM5 Lentivirus	78719
		CEACAM6 Lentivirus	78720
		Claudin-3 Lentivirus	78722
		Claudin-4 Lentivirus	78723
		Claudin-9 Lentivirus	78721
		CSL (CBF1/RBP-Jk) Luciferase Reporter Lentivirus (Notch Signaling Pathway)	78746
		CTLA4 CRISPR/Cas9 Lentivirus (Integrating)	78054
		CTLA4 CRISPR/Cas9 Lentivirus (Non-Integrating)	78061
		Cyno EpCAM Lentivirus	78978
		DLL3 Lentivirus	78909
		Dominant Negative TGF- β Receptor Type II (TGF- β RII) Lentivirus	78928
		EGR1 Promoter Luciferase Reporter Lentivirus	78664
		FCGR2A CRISPR/Cas9 Lentivirus (Integrating)	78537
		FCGR2A CRISPR/Cas9 Lentivirus (Non-Integrating)	78538
		FcRL5 Lentivirus (Macaca fascicularis/Cynomolgus)	78781
		FOLR1 Lentivirus (Macaca fascicularis/Cynomolgus)	78778
		GPRC5D Lentivirus	78716
		GPRC5D Lentivirus (Macaca fascicularis/Cynomolgus)	78780
		HLA-C*08:02 Lentivirus	78930
		HLA-E Lentivirus	78929
		HSE Luciferase Reporter Lentivirus (Heat Shock Response)	78669
		IL15/IL15Ra Lentivirus	78938
		KRAS G12D-Specific TCR Lentivirus (Clone 10)	78937
		KRAS G12D-Specific TCR Lentivirus (Clone 9c)	78936
		LAG3 CRISPR/Cas9 Lentivirus (Integrating)	78053
		LAG3 CRISPR/Cas9 Lentivirus (Non-Integrating)	78060
Inhibitors/Activators	Catalog#		
AMP (500 μ M)	79496		
Cyclopamine, Free Base	27013		
Epacadostat (INCB024360)	27339-1		
Epacadostat (INCB024360)	27339-2		
Fludarabine-Cl	82534		
INCB024360 Analog	27338-1		
INCB024360 Analog	27338-2		
NLG919	27337-1		
NLG919	27337-2		
Telotristat	27340		
Telotristat	27341		
Vismodegib (GDC-04491)	27015		
Vismodegib (GDC-04491)	27010		
Lentiviruses	Catalog#		
Anti-BCMA CAR Lentivirus (Clone C11D5.3 ScFv-CD8-4-1BB-CD3 ζ)	78655		
Anti-BCMA CAR Lentivirus (VHH1/VHH2 ScFv-CD8-4-1BB-CD3 ζ)	78783		
Anti-CD19 CAR Lentivirus (CD19 ScFv-CD8-4-1BB-CD3 ζ)	78600		
Anti-CD19 CAR Lentivirus (CD19 ScFv-CD8-4-1BB-CD3 ζ , eGFP)	78775		
Anti-CD19 CAR Lentivirus (CD19 ScFv-CD8-4-1BB-CD3 ζ , PuroR)	78602		
Anti-CD19 CAR Lentivirus (CD19 ScFv-CD8-4-1BB-CD3 ζ ; SIN Vector)	78601		
Anti-CD19/CD22 Bispecific CAR Lentivirus (Clones FMC63/m971 ScFv-CD8-4-1BB-CD3 ζ)	78609		
Anti-CD20 CAR Lentivirus (Clone Leu-16 ScFv-CD8-4-1BB-CD3 ζ)	78606		
Anti-CD22 CAR Lentivirus (Clone m971 ScFv-CD8-4-1BB-CD3 ζ)	78608		
Anti-Mesothelin CAR Lentivirus (P4 ScFv-CD8-4-1BB-CD3 ζ)	78703		
ATF6 Luciferase Reporter Lentivirus (ATF6 Pathway)	78667		
B7-H4 Lentivirus	78727		

Lentiviruses	Catalog#	Primary Cells	Catalog#
LAIR1 Lentivirus	78903	Anti-CD20 CAR-T Cells	78611
MAGE-A1-Specific TCR Lentivirus (Clone 1367)	78934	CD4+ T cells, Negatively Selected (Human)	79752
MAGE-A4 Specific TCR Lentivirus	78935	CD8+ T cells, Negatively Selected (Human)	79753
Membrane-Bound TNF α (mTNF α) Lentivirus	78955	Expanded Human Peripheral Blood NK Cells, Frozen	78798
Nectin-4 Lentivirus	78712	Normal Human Peripheral Blood Mononuclear Cells, Frozen	79059
NKp46 Lentivirus (Macaca fascicularis/Cynomolgus)	78779		
NLRP3 CRISPR/Cas9 Lentivirus (Integrating)	78545		
NLRP3 CRISPR/Cas9 Lentivirus (Non-Integrating)	78546		
NLRP3 Human shRNA Lentivirus	82122		
Notch1dE Lentivirus	78747		
PD-1 (Human) sgRNA-MS2 Lentivirus (Integrating)	78190		
PD-1 CRISPR/Cas9 Lentivirus (Integrating)	78052		
PD-1 CRISPR/Cas9 Lentivirus (Non-Integrating)	78059		
PD-L1 CRISPR/Cas9 Lentivirus (Integrating)	78057		
PD-L1 CRISPR/Cas9 Lentivirus (Non-Integrating)	78064		
PD-L1 Lentivirus	78925		
PSMA Lentivirus	78726		
STAT6 Luciferase Reporter Lentivirus (STAT6 Signaling Pathway)	78799		
TCR CRISPR/Cas9 Lentivirus (Non-Integrating)	78062		
TGFB2 CRISPR/Cas9 Lentivirus (Integrating)	78535		
TGFB2 CRISPR/Cas9 Lentivirus (Non-Integrating)	78536		
TIGIT CRISPR/Cas9 Lentivirus (Integrating)	78058		
TIGIT CRISPR/Cas9 Lentivirus (Non-Integrating)	78065		
TNFR2 Lentivirus	78765		
Trop2 Lentivirus (Macaca fascicularis/Cynomolgus)	78776		
ULBP2 Lentivirus	78744		
ULBP2 Lentivirus (Macaca fascicularis/Cynomolgus)	78777		
VSIG4 Lentivirus	78902		
Vy4V61 TCR Lentivirus	78986		
Vy9V62 TCR Lentivirus	78985		
Primary Cells	Catalog#	Proteins	Catalog#
Anti-Mesothelin CAR-T Cells	78729	AC-HLY, His-FLAG-Tags	11005
Dual Epitope Anti-BCMA CAR-T Cells	78790	Activin Blocker	102121
NY-ESO-1 (c259) TCR-T Cells	78990	ADAR1, FLAG-Tag	100472
Untransduced T Cells (NY-ESO-1 TCR-T Negative Control)	78989	ADAR2 (ADARB1), FLAG-Tag	101164
Anti-BCMA CAR-T Cells	78660	APRIL, His-Avi-Tag	100254
Anti-CD19 CAR-T Cells	78171	APRIL, His-Avi-Tag, Biotin-Labeled	100262
Anti-CD19 CAR-T Cells (eGFP)	78789	ARG1, His-Tag	71658
		ARG2, His-Tag	71659
		B7-1 (CD80) (Mouse), Fc fusion (Human), Avi-Tag	79058
		B7-1 (CD80), Fc fusion (Human) HiP™	71125
		B7-1 (CD80), Fc fusion, Biotin-labeled (Human) HiP™	71114
		B7-1 (CD80), FLAG-Avi-His-Tag (Mouse)	79002
		B7-1 (CD80), FLAG-Avi-His-Tag	71261
		B7-1 (Woodchuck), Fc-fusion (Human), Avi-Tag, Biotin-labeled	100040
		B7-1(CD80) (Mouse), Fc fusion, Avi-Tag, Biotin-labeled, HiP™	100478
		B7-2 (CD86), Fc fusion (Human)	71150
		B7-2 (CD86), Fc fusion, Biotin labeled (Human) HiP™	71159
		B7-2 (CD86), FLAG-Avi-His-Tag, Biotin Labeled	71263
		B7-H2 (CD275, ICOSL), Fc fusion (Human) HiP™	71130
		B7-H2, Avi-His-Tag	79119
		B7-H2, Avi-His-Tag, Biotin-Labeled	79300
		B7-H3, Avi-His-Tag HiP™	79337
		B7-H3, Avi-His-Tag, Biotin-Labeled HiP™	79367
		B7-H4, His-Tag (Human)	71144
		B7-H4, His-Tag, Biotin-Labeled (Human) HiP™	71129
		B7-H5 (VISTA), Fc fusion (Human) HiP™	71148
		B7-H5 Fc Fusion, Avi-Tag, Biotin-Labeled (Human)	71327
		B7-H5, Avi-His-Tag	100358
		B7-H5, Avi-His-Tag, Biotin Labeled, HiP™	100359
		B7-H6, Fc-Fusion-Avi-Tag (Human)	79383
		B7-H6, Fc-Fusion-Avi-Tag, Biotin-Labeled (Human)	79384
		B7-H7, Fc-Fusion, Avi-Tag	79365

Proteins	Catalog#	Proteins	Catalog#
B7-H7, Fc-Fusion, Avi-Tag, Biotin-Labeled	79366	CD137L, His-Tag (Mouse)	72517
BAFF-R, Fc-fusion (IgG1), Avi-Tag	100286	CD14, Avi-His-Tag HiP™	101187
BAFF-R, Fc-fusion (IgG1), Avi-Tag, Biotin-labeled	100287	CD14, Avi-His-Tag, Biotin-Label	101188
BCMA, Fc-fusion (IgG1), Avi-Tag, Biotin-Labeled	79467	CD150, Fc Fusion (IgG1), Biotin Labeled (Human)	72307
BCMA, Fc-Fusion, Avi-Tag	79465	CD150, Fc Fusion, Avi-Tag (Human)	72304
BCMA, Fc-Fusion, Avi-Tag, PE-Labeled	100733	CD155 (PVR), His-tag (Human)	71181
BTLA(CD272), Fc fusion (Human)	71141	CD155 (PVR), His-tag (Mouse) HiP™	71167
BTLA, Fc Fusion, Avi-Tag HiP™	100033	CD155, Fc Fusion, Avi-Tag, Biotin-Labeled	79064
BTLA, Fc Fusion, Avi-Tag, Biotin-Labeled HiP™	100047	CD155, Fc-Avi-Tag	79063
BTN1A1 Avi-His-Tag	100469	CD155, His-tag, Biotin-labeled (Mouse) HiP™	71168
BTN2A1, Avi-Tag, His-Tag Biotin Labeled	102086	CD160, His-tag (Human) HiP™	71173
BTN2A1, Avi-Tag, His-Tag	102085	CD160, His-tag, Biotin-labeled (Human) HiP™	71135
BTN2A1, Fc Fusion, Avi-Tag	102083	CD161, Fc Fusion (IgG1), Avi-Tag HiP™	101211
BTNL8, Fc-fusion (IgG1), Avi-Tag	100467	CD19, Avi-His-Tag	101015
BTNL8, Fc-fusion (IgG1), Avi-Tag, Biotin-labeled	100468	CD19, Fc-Fusion (IgG1), Avi-Tag	79472
c-Mer, GST-tag (Human)	40254	CD19, Fc-Fusion (IgG1), Avi-Tag, Biotin-Labeled	79475
CBL-B (Y363F), His-tag, Biotin-labeled (Human)	80413	CD19, Fc-Fusion (IgG1), Avi-Tag, PE-labeled	100732
CBL-B, His-Avi-Tag, Biotin-labeled (Human)	80412	CD2, Avi-His-Tag HiP™	101238
CD111, Fc-Fusion, Avi-Tag	77001	CD2, Avi-His-Tag, Biotin-Labeled HiP™	101239
CD111, Fc-Fusion, Avi-Tag, Biotin-Labeled HiP™	77002	CD2, Fc fusion, Avi-Tag, (Human) HiP™	71264
CD112, Fc Fusion, Avi-Tag	11079	CD2, Fc fusion, Biotin-labeled (Human) HiP™	71270
CD112, Fc-Fusion, Avi-Tag, Biotin-labeled (Human)	72231	CD20, FLAG-Tag	101572
CD112, His-tag (Human) HiP™	71197	CD200, Avi-Tag, His-Tag	102030
CD112, His-Tag, Biotin-Labeled (Human) HiP™	71234	CD200, Avi-Tag, His-Tag, Biotin-Labeled	102031
CD112R (PVRIG), Fc-Fusion, Avi-Tag (Human) HiP™	79116	CD200, Fc Fusion, Avi-Tag	102013
CD112R (PVRIG), Fc-Fusion, Avi-Tag, Biotin-Labeled (Human) HiP™	79270	CD200, Fc Fusion, Avi-Tag, Biotin-Labeled	102014
CD113, Fc-Fusion (IgG1), Avi-Tag	72510	CD200R1, Fc Fusion, Avi-Tag	102032
CD123, Avi-His-Tag	101035	CD200R1, Fc Fusion, Avi-Tag, Biotin-Labeled	102033
CD123, Fc-Fusion (IgG1) Avi-Tag	100058	CD22, Fc Fusion, Avi-Tag, PE-labeled	101028
CD123, Fc-Fusion (IgG1) Avi-Tag, Biotin-Labeled	100068	CD22, Fc-fusion, Avi-Tag HiP™	79464
CD125, Fc Fusion (IgG1), Avi-Tag	101050	CD22, Fc-fusion, Avi-Tag, Biotin-labeled HiP™	79466
CD125, Fc-Fusion (IgG1), Avi-Tag, Biotin-Labeled	101051	CD226, Fc fusion (Human) HiP™	71252
CD137 (4-1BB), Fc fusion (Human) HiP™	71170	CD226, Fc fusion, Biotin-labeled (Human) HiP™	71253
CD137 (4-1BB), Fc Fusion, Avi-tag Biotin-Labeled HiP™	71171	CD235a, Fc Fusion (IgG1), Avi-Tag HiP™	101196
CD137, Fc fusion (hlgG1) (Mouse)	71250	CD235a, Fc Fusion (IgG1), Avi-Tag, Biotin-Labeled HiP™	101197
CD137, Fc fusion (mlgG2a), Avi-tag (Mouse) HiP™	71254	CD24, Fc Fusion, Avi-tag	100527
CD137, Fc fusion (mlgG2a), Biotin-labeled (Mouse) HiP™	71255	CD24, Fc Fusion, Avi-tag, Biotin-labeled, HiP™	100528
CD137L (4-1BB Ligand), His-Avi-Tag	100238	CD244, Fc Fusion (IgG1), Avi-Tag (Human)	72305
CD137L, His-Tag (E. Coli-Derived)	100562	CD244, Fc Fusion (IgG1), Biotin-Labeled (Human)	72306

Proteins	Catalog#	Proteins	Catalog#
CD27(Mouse), Fc-Fusion (Human), Avi-Tag, Biotin	79107	CD45(PTPRC), His-tag	30044
CD27, Fc fusion (Human)	71176	CD46-BC, His-tag (Human)	71230
CD27, Fc Fusion (IgG1) Avi-Tag, Biotin Labeled	100250	CD46-BC, His-Tag (Monkey) HiP™	71232
CD27, Fc-Fusion (IgG1), Avi-Tag (Mouse) HiP™	79067	CD46-C, His-tag (Human)	71231
CD277, Avi-Tag, His-Tag	102113	CD46-C, His-tag (Monkey)	71233
CD277, Fc-Fusion (IgG1) Avi-Tag	100073	CD47 (Monkey), Fc Fusion (Human), Avi-Tag HiP™	79118
CD277, Fc-Fusion (IgG1) Avi-Tag, Biotin-Labeled	100074	CD47 (Monkey), Fc Fusion (Human), Avi-Tag, Biotin HiP™	79302
CD28, Fc fusion (Human)	71113	CD47, Fc Fusion (IgG1)	71177
CD28, Fc Fusion, Avi-Tag, Biotin Labeled	101580	CD47, Fc fusion, Avi-Tag (Human) HiP™	79051
CD28H, Fc Fusion (IgG1), Avi-Tag	11083	CD47, Fc Fusion, Avi-Tag, Biotin-Labeled (Mouse)	72514
CD28H, Fc fusion, Biotin-labeled (Human) HiP™	71239	CD47, Fc fusion, Biotin-labeled (Human) HiP™	71169
CD28H, His-Tag	79375	CD47, Fc-Fusion, Streptavidin-Labeled	71292
CD30, Fc-Fusion (IgG1), Avi-Tag, Biotin-Labeled	72511	CD47, His-Tag (Human)	71127
CD33, Fc Fusion, Avi-tag	100155	CD48, Fc fusion (IgG1), Avi-Tag™ HiP™ (Human)	72309
CD33, Fc Fusion, Avi-tag, Biotin Labeled, HiP™	100482	CD48, Fc-Avi, Biotin-Labeled	72308
CD33, Fc Fusion, Avi-Tag, PE-labeled	101029	CD5, Fc Fusion, Avi-Tag, Biotin-labeled, HiP™	101006
CD34, Avi-His-Tag HiP™	101190	CD5, Fc-Fusion, Avi-Tag, HiP™	101005
CD34, Avi-His-Tag, Biotin-Labeled	101191	CD55, Avi-His-Tag, Biotin-labeled	100943
CD34, Fc Fusion, Avi-Tag	101812	CD55, Avi-His-Tag, HiP™	100942
CD36, Avi-His-Tag	101184	CD58, Fc fusion, Biotin-labeled HiP™	71269
CD36, Fc Fusion, Avi-Tag	101813	CD5L, Avi-His-Tag	101007
CD37, Avi-His-Tag	101330	CD66b, Fc Fusion, Avi-Tag	102029
CD37, Avi-His-Tag, Biotin-Labeled	101331	CD70 His-Avi-Tag, Biotin-Labeled	100228
CD38, Avi-His-Tag	100346	CD70, His-Avi-Tag	100240
CD38, Avi-His-Tag, Biotin-Labeled HiP™	100352	CD70, His-tag (Human)	71178
CD38, FLAG-Tag (Pig), HiP™	101019	CD70, His-Tag (Mouse)	79066
CD38, His-Tag (Dog)	100955	CD73, Avi-His-Tag (Mouse)	72523
CD38, His-Tag (Human), HiP™	71277	CD73, Avi-His-Tag, Biotin Labeled	79127
CD38, His-Tag (Mouse), HiP™	79070	CD73, Avi-His-Tag, Biotin-Labeled (Mouse)	79416
CD38, His-Tag, PE-labeled	71882	CD73, His-Tag (Human)	71184
CD38-APC, His-Tag	71883	CD83, Avi-His-Tag HiP™	100224
CD39, His-Tag (Human)	71284	CD83, Avi-His-Tag, Biotin-Labeled HiP™	100225
CD4, Avi-Tag, His-Tag, Biotin-Labeled	101869	CD83, Fc-fusion (IgG1), Avi-Tag	100347
CD4, Avi-Tag, His-Tag, HiP™	101862	CD83, Fc-fusion (IgG1), Avi-Tag, Biotin labeled	100348
CD40, Fc fusion (Human)	71174	CD8a, Avi-Tag, His-Tag	101872
CD40, Fc Fusion (IgG1) Avi-Tag, Biotin Labeled	79102	CD8a, Avi-Tag, His-Tag, Biotin-Labeled	101873
CD40L (CD154), His-tag (Human)	71191	CD8a, Fc Fusion, Avi-Tag, Biotin-Labeled, HiP™	101871
CD44, Fc fusion (Human)	71187	CD8a, Fc Fusion, Avi-Tag, HiP™	101870
CD44, Fc fusion, Biotin-labeled (Human) HiP™	71133	CD8B, Avi-His-Tag	101843

Proteins	Catalog#	Proteins	Catalog#
CD8b, Avi-Tag, His-Tag, Biotin-Labeled	101844	DRAK2	71225
CD8b, Fc Fusion, Avi-Tag	101841	DRAK2, GST-tag (Mouse)	100444
CD8B, Fc Fusion, Avi-Tag, Biotin-Labeled	101842	DRAK2, GST-Tag	71224
CD93, Fc Fusion (IgG1), Avi-Tag	101543	EGFR (del746-750, T790M, C797S), GST-Tag, His-Tag	100187
CD93, Fc Fusion (IgG1), Avi-Tag, Biotin-Labeled	101544	eIF4E, GST-tag	40530
CD96, Fc-Fusion, Avi-Tag (Human)	71265	EIF4EBP1, His-tag	40510
CD96, Fc-Fusion, Avi-Tag, Biotin Labeled (Human)	71280	ELANE (Elastase), Avi-His-Tag HiP™	101141
CEACAM, Avi-His-Tag, Biotin Labeled	70201	ELANE (Elastase), Avi-His-Tag, Biotin-Labeled HiP™	101142
CEACAM1, Avi-His-Tag	70206	EpCAM, Avi-His-Tag	100461
CEACAM5, Avi-His-Tag	100509	EpCAM, Avi-His-Tag, Biotin Labeled	100462
CERK (K298G, K303G), FLAG-Tag	102044	EPOR, Fc fusion, Avi-Tag	100611
CERK, FLAG-Tag	102043	EPOR, Fc fusion, Avi-Tag, Biotin-labeled HiP™	100612
Claudin-18 Isoform 2, FLAG-Tag	101570	ERAP1, Avi-His-Tag, Biotin-labeled	100577
Claudin-6, FLAG-Tag	101599	FCGR2A, Avi-His-Tag	100094
CLEC1A, His-Avi-Tag, Biotin-Labeled	101011	FCGR2B, Avi-His-Tag	100089
CLEC1A, His-Tag	101010	FCGR2B, Avi-His-Tag, Biotin-Labeled	100474
CLEC1B, His-Avi-Tag	101021	FCGR3A (F176V), Avi-His-Tag	100079
CLEC1B, His-Avi-Tag, Biotin-Labeled	101022	FCGR3A, Avi-His-Tag	79013
CLEC2D, Fc Fusion (IgG1), Avi-Tag	101186	FCGR3B, Avi-His-Tag	79016
CLEC2D, Fc Fusion (IgG1), Avi-Tag, Biotin-Labeled HiP™	101189	FcRn (FCGRT/B2M) Blocker	101468
CLEC2D, His-Avi-Tag, Biotin-Labeled	101470	FcRn (FCGRT/B2M), His-Avi-Tag, Biotin-Labeled (Mouse)	71286
CRTAM, Fc-Fusion (IgG1), Avi-Tag	40686	FcRn (FCGRT/B2M), His-Tag (Mouse) HiP™	11349
CRTAM, Fc-Fusion (IgG1), Avi-Tag, Biotin-Labeled	40687	FcRn (FCGRT/B2M), His-Tag	71285
CSF1R, Fc-Fusion, Avi-Tag (Human) HiP™	72550	FcRn Complex (FCGRT/B2M), His-Avi-Tag, Biotin-Labeled, HiP™	71283
CSF1R, Fc-Fusion, Avi-Tag, Biotin-Labeled	79343	FGF19, Avi-Tag, His-Tag	101886
CTLA4 (CD152), Fc fusion, Biotin-labeled (Human)	71152	FGF19, Avi-Tag, His-Tag, Biotin-Labeled	101888
CTLA4 (CD152), Fc-fusion (IgG1)	71149	FGFR1a (IIIc), Fc Fusion, Avi-Tag	101836
CTLA4 (Mouse), Fc-Fusion (Human), Avi-Tag	79062	FGFR2C, Avi-His-Tag	101830
CTLA4 (Mouse), Fc-Fusion (Human), Avi-Tag, Biotin Labeled	79001	FGFR2C, Fc Fusion, Avi-Tag	101877
CTLA4 (Woodchuck), Fc-Fusion (Human), Avi-tag HiP™	100037	FGFR3, Avi-Tag, His-Tag, Biotin-Labeled, HiP™	101818
DDX58, FLAG-Tag	100070	FGFR3, Fc Fusion, Avi-Tag	101819
Diablo, His-Tag	101118	FGFR3, Fc Fusion, Avi-Tag, Biotin-Labeled	101820
DLK1, Avi-His-Tag HiP™	101745	GITR (CD357), Fc fusion (Human) HiP™	71172
DLK1, Avi-Tag, His-Tag, Biotin-Label	101781	GITR (CD357), Fc Fusion, Biotin-labeled (Human)	71256
DLL1, Avi-Tag, His-Tag, Biotin-Labeled	101810	GITRL, His-tag (Human)	71190
DLL1, Avi-Tag, His-Tag,	101805	GPC3, His-tag [Monkey (M. fascicularis)]	71260
DLL4, Fc Fusion, Avi-Tag	101903	HER3, Avi-Tag, His-Tag	102230
DLL4, Fc Fusion, Avi-Tag, Biotin-Labeled	101904	HER3, Fc Fusion, Avi-Tag	102228
DPP4, His-tag	80040	HMGB1, Avi-His-Tag	79082

Proteins	Catalog#	Proteins	Catalog#
HMGB1, Avi-His-Tag, Biotin-labeled	79007	IL23R, Fc-Fusion (IgG1), Avi-Tag HiP™	101321
HPK1, GST-Tag, His-Avi-Tag (Human)	40398	IL2RA, Avi-His-Tag	101203
Human IL-8	90258-A	IL2RA, Avi-His-Tag, Biotin-Labeled	101204
Human IL-8	90258-B	IL2RB, Avi-FLAG-Tag, Biotin-Labeled HiP™	101314
Human Interleukin-15	90180	IL2RB, Avi-FLAG-Tag, HiP™	101313
Human Interleukin-18	90183-A	IL2RB, Avi-His-Tag, Biotin-Labeled	100428
Human Interleukin-18	90183-B	IL6R, Fc-Fusion (IgG1)	100916
Human Interleukin-21	90186-A	KIR2DL1 Fc-fusion, Biotin-labeled HiP™	71296
Human Interleukin-21	90186-B	KIR2DL1, Fc-fusion, Avi-Tag HiP™	79649
Human Interleukin-8 (72AA)	90199-A	KIR2DL2 Fc-fusion, Avi-Tag HiP™	79650
Human Interleukin-8 (72AA)	90199-B	KIR2DL2 Fc-fusion, Biotin-labeled HiP™	71297
Human LTBR, Fc Fusion HiP™	71112	KIR2DL3, Fc fusion, Avi-Tag HiP™	71299
HVEM, Fc fusion (Human)	71142	KIR2DL3, Fc fusion, Biotin-labeled (Human)	71298
HVEM, Fc fusion, Biotin-labeled (Human)	71143	KIR3DL3, Avi-His-Tag	101408
ICOS (CD278), Fc fusion (Human)	71179	KIR3DL3, Fc Fusion (IgG1), Avi-Tag	101414
IDO1, His-Avi-Tag, Biotin Labeled	79081	KIR3DS1, Fc Fusion (IgG1), Avi-Tag, Biotin-Labeled	101248
IDO1, His-tag (Mouse)	71196	KIR3DS1, Fc-Fusion (IgG1), Avi-Tag	101247
IDO1, His-tag	71182	KMO, His-FLAG-Tags	11307
IDO2, His-tag (Mouse)	71240	KYNU, His-Tag	79485
IDO2, His-Tag	71194	LAG3 (CD223) (Mouse), Fc fusion (Human) Avi-Tag	79050
IgG1, Fc (Human) HiP™	71456	LAG3 (CD223) (Mouse), Fc(Human) Avi-Tag, Biotin HiP™	79003
IgG4, Fc (Human) HiP™	71457	LAG3 (CD223), Fc fusion (Human)	71146
IL-12 (p40/p35) Fc Fusion (IgG1), Avi-Tag	101431	LAG3 (CD223), Fc-Fusion, Avi-Tag, Biotin-Labeled (Human) HiP™	71147
IL-12 (p40/p35) Fc Fusion (IgG1), Avi-Tag, Biotin-Labeled	101432	LAG3, His-Tag, HiP™	71229
IL-17A, Avi-Tag	91014	LAIR1, Avi-His-Tag HiP™	79483
IL-17RA, Fc fusion, Avi-Tag	11264	LAT, His-tag (Human)	71188
IL-17RA, Fc fusion, Biotin-labeled (Human)	91013	LIGHT His-Avi-Tag, Biotin-labeled	100032
IL-2 (C145A)	100159	LIGHT, His-Tag (Human)	71266
IL-2 (R58A, F62A, Y65A, E82A, C145A) (Woodchuck)	100156	LIGHT, His-Tag (Mouse)	79068
IL-2 (R58D, K63E, E81R, C146A) (Woodchuck)	100157	LILRA1, Avi-His Tag	100245
IL-2, Fc Fusion (IgG1), Avi-Tag, Biotin-Labeled	101381	LILRA1, Avi-His Tag, Biotin-labeled	100246
IL-2RB (CD122)	79655	LILRB1, Fc-Fusion (IgG1), Avi-Tag HiP™	79473
IL-2RB, Avi-His-Tag	100427	LILRB1, Fc-Fusion (IgG1), Avi-Tag, Biotin-Labeled HiP™	79474
IL-2RG, Avi-His-Tag	101149	LILRB2, Avi-His-Tag HiP™	100234
IL-2RG, Avi-His-Tag, Biotin-Labeled	101150	LILRB2, Avi-His-Tag, Biotin-labeled	100335
IL12RB1, Avi-His-Tag	101336	LILRB4, Avi-His-Tag, Biotin-Labeled HiP™	100336
IL18R1, Fc Fusion (IgG1), Avi-Tag HiP™	101283	LILRB4, Avi-His-Tag, HiP™	100236
IL18RAP, Fc Fusion (IgG1), Avi-Tag HiP™	101367	Lin28A, FLAG-Avi-Tag	79254
IL23R, Fc Fusion (IgG1), Avi-Tag, Biotin-Labeled HiP™	101322	Lin28A, FLAG-Avi-Tag, Biotin-Labeled	79072

Proteins	Catalog#	Proteins	Catalog#
Lin28B, FLAG-Avi-Tag	79255	OX40L (CD252), His-Tag (Mouse)	72516
Lin28B, FLAG-Avi-Tag, Biotin-Labeled	79073	OX40L, His-Avi-Tag	100808
MAT2A, His-Avi-Tag, Biotin-Labeled	79308	p21, GST-Tag	101584
MAT2A, His-Tag	71401	PD-1 (CD279), Fc fusion (Human) HiP™	71106
Mesothelin, Avi-His-Tag, Biotin-Labeled, HiP™	100291	PD-1 (CD279), Fc fusion (Mouse) HiP™	71116
Mesothelin, Avi-His-Tag, HiP™	100290	PD-1 (CD279), Fc fusion, Biotin-labeled (Human) HiP™	71109
MICA, Fc fusion, Avi-Tag	100295	PD-1 (CD279), Fc fusion, Biotin-labeled (Mouse) HiP™	71118
MICA, Fc fusion, Avi-Tag, Biotin-labeled	100300	PD-1 (E136A), Fc fusion, Biotin-labeled (Human) HiP™	71247
MICB, Avi-His-Tag HiP™	100505	PD-1 (Human), Fc Fusion (Mouse IgG2a) HiP™	71997
MMP8, His-tag	100552	PD-1 (I126A), Fc fusion, Biotin-labeled (Human) HiP™	71238
Mouse LTBR, Fc Fusion HiP™	71122	PD-1 (I134A), Fc fusion, Biotin-labeled (Human) HiP™	71245
mTOR/MLST8/RPTOR Complex	102073	PD-1 (K78A), Fc fusion, Biotin-labeled (Human) HiP™	71244
mTOR/Raptor/MLST8 Complex, FLAG-Tag, His-Tag (Human)	40300	PD-1 (L128A), Fc fusion, Biotin-labeled (Human) HiP™	71248
NCAM1, Avi-His-Tag, Biotin-Labeled	101044	PD-1 (Monkey), Fc fusion (Human IgG1) HiP™	71158
NCAM1, Avi-His-Tag, HiP™	101043	PD-1 (Monkey), Fc fusion (Human IgG1), Biotin-labeled HiP™	71235
Nectin4, His-Avi-Tag HiP™	100674	PD-1 (Mouse), Fc fusion (Human), PE-labeled	71267
Nectin4, His-Avi-Tag, Biotin-labeled	100675	PD-1 (P130A), Fc fusion, Biotin-labeled (Human) HiP™	71246
NEU2, His-Tag	101722	PD-1, Avi-His-Tag (Mouse)	79288
NKG2A, His-Avi-Tag	100638	PD-1, Avi-His-Tag, Biotin-Labeled (Mouse)	79189
NKG2D, Avi-Tag, Fc fusion	100252	PD-1, Fc Fusion (IgG1), Eu-labeled	79011
NKG2D, Avi-Tag, Fc fusion, Biotin-labeled	100313	PD-1, Fc Fusion (Woodchuck)	79314
Nkp30, Fc Fusion, Avi-tag	100546	PD-1, Fc fusion, Avi-Tag	100406
NKp46, Fc-fusion (IgG1), Avi-Tag	100465	PD-1, Fc Fusion, PE-Labeled (Human)	71243
NKp46, Fc-fusion (IgG1), Avi-Tag, Biotin-Labeled	100466	PD-1, FLAG-Avi-His-tag (Human) HiP™	71198
NLRP3 (delPYD) / NEK7 Complex	100477	PD-1, FLAG-Avi-His-Tag, Biotin-Labeled (Human) HiP™	71325
NLRP3 (delPYD), FLAG-Tag	100454	PD-L1 (A121Q), Fc fusion, Biotin-labeled (Human) HiP™	71320
NLRP3 (K232A), His-FLAG-Tag (Sf9-derived)	40741	PD-L1 (CD274)	100443
NLRP3 (NALP3) (GKT(231-233)AAA), His-Flag-Tags (HEK293-derived)	100190	PD-L1 (CD274), Fc fusion (Human) HiP™	71104
NLRP3 (NALP3), His-Flag-Tags	100189	PD-L1 (CD274), Fc fusion (Mouse) HiP™	71117
Notch1, Avi-His-Tag	101897	PD-L1 (CD274), Fc fusion, Biotin-labeled (Human) HiP™	71105
Notch2, Fc Fusion, Avi-Tag	101986	PD-L1 (CD274), Fc fusion, Biotin-labeled (Mouse) HiP™	71119
Notch2, Fc Fusion, Avi-Tag, Biotin-Labeled	101987	PD-L1 (CD274), FLAG-tag (Human) HiP™	71183
NPM1, His-Tag	102057	PD-L1 (D122A), Fc fusion, Biotin-labeled (Human) HiP™	71321
OX40 (CD134), Avi-Tag, His-Tag	11372	PD-L1 (D26A), Fc fusion, Biotin-labeled (Human) HiP™	71313
OX40 (CD134), Biotin-labeled, His-tag (Human)	71310	PD-L1 (F19A), Fc fusion, Biotin-labeled (Human) HiP™	71311
OX40 (CD134), Fc fusion (Human)	71175	PD-L1 (Human), Fc fusion (Mouse IgG2a) HiP™	71999
OX40, Fc-Fusion, Avi-Tag (Mouse) HiP™	72546	PD-L1 (Human), Fc Fusion (Mouse IgG2a) PE-Labeled	71998
OX40, Fc-Fusion, Avi-Tag, Biotin-Labeled (Mouse) HiP™	79357	PD-L1 (I54A), Fc fusion, Biotin-labeled (Human) HiP™	71314
OX40L (CD252), His-tag (Human)	71185	PD-L1 (K124A), Fc fusion, Avi-Tag (Human) HiP™	71993

Proteins	Catalog#	Proteins	Catalog#
PD-L1 (K124A), Fc fusion, Biotin-labeled (Human) HiP™	71323	Siglec5, Fc Fusion, Avi-tag, Biotin-labeled	100534
PD-L1 (M115A), Fc fusion, Biotin-labeled (Human) HiP™	71318	Siglec6, Fc Fusion, Avi-tag	100535
PD-L1 (Monkey), Fc fusion (Human IgG1) HiP™	71153	Siglec6, Fc Fusion, Avi-tag, Biotin-labeled	100536
PD-L1 (Monkey), Fc fusion (Human IgG1), Biotin-labeled HiP™	71154	Siglec7, Fc Fusion, Avi-tag	100589
PD-L1 (Q66A), Fc fusion, Biotin-labeled (Human) HiP™	71316	Siglec7, Fc Fusion, Avi-tag, Biotin-labeled	100636
PD-L1 (R113A), Fc fusion, Biotin-labeled (Human) HiP™	71317	Siglec8, Fc Fusion, Avi-Tag	100421
PD-L1 (R125A), Fc fusion, Biotin-labeled (Human) HiP™	71324	Siglec9, Fc Fusion, Avi-tag	100558
PD-L1 (S117Q), Fc fusion, Biotin-labeled (Human) HiP™	71319	Siglec9, Fc Fusion, Avi-tag, Biotin-labeled	100559
PD-L1 (T20Q), Fc fusion, Biotin-labeled (Human) HiP™	71312	SIRP-α (CD172a), Avi-His-Tag (Monkey) HiP™	79117
PD-L1 (Y123A), Fc fusion, Avi-Tag (Human)	71992	SIRP-α (CD172a), Avi-His-Tag, Biotin (Monkey) HiP™	79303
PD-L1 (Y123A), Fc fusion, Biotin-labeled (Human) HiP™	71322	SIRP-α (CD172a), His-tag (Human)	71145
PD-L1 (Y56A), Fc fusion, Biotin-labeled (Human) HiP™	71315	SIRP-α (SIRP alpha), His-tag, Biotin-labeled (Human) HiP™	71138
PD-L1(CD274), FLAG-Avi-His-Tag, (Human) HiP™	71279	SIRP-γ (CD172g), Fc fusion, Biotin-labeled (Human) HiP™	71236
PD-L1, Fc Fusion (Human IgG1) PE-Labeled	71881	SLAMF7, Avi-His-Tag HiP™	100271
PD-L1, Fc Fusion, PE-Labeled (Mouse)	71268	SLAMF7, Avi-His-Tag, Biotin-Labeled	100272
PD-L2 (CD273), Fc fusion (Human) HiP™	71107	STAT3, GST-tag	75003
PD-L2 (CD273), Fc fusion, Biotin-labeled (Human)	71108	STING, GST-Tag, His-Tag	100043
PDPN, Fc-Fusion, Avi-Tag HiP™	79341	TACI, Fc-Fusion (IgG1), Avi-Tag	100283
PDPN, Fc-Fusion, Avi-Tag, Biotin-Labeled HiP™	79342	TDO, His-Tag (Human)	71195
PSGL1 (CD162), Fc Fusion, Avi-tag	100631	TDO2, His-tag (Mouse)	71241
PSGL1 (CD162), Fc Fusion, Avi-tag, Biotin-labeled	100632	TFRC (Transferrin Receptor), Avi-Tag, Fc Fusion, Biotinylated	101315
PSMA, His-Avi-Tag	100463	TFRC, Avi-Tag, Fc Fusion (IgG1)	101287
PSMA, His-Avi-Tag, Biotin-Labeled	100464	TGFβR2, Fc Fusion, Avi-tag, Biotin-labeled HiP™	100629
RANK, Fc fusion (IgG1), Avi-tag (Human)	70823	TIGIT, Fc Fusion (Human)	71186
RANK, Fc fusion (IgG1), Biotin Labeled (Human)	70822	TIGIT, Fc Fusion (IgG1) Avi-Tag, Biotin (Mouse) HiP™	79269
RANKL, His-Tag (Human)	71051	TIGIT, Fc Fusion, Avi-Tag, Biotin-Labeled (Human) HiP™	71251
RB, GST-tag	40595	TIM-1, Avi-His-Tag	72512
ROR1, Fc-Fusion (IgG1)-Avi-Tag, PE-labeled	100995	TIM-3 (Mouse) Fc-Fusion (Human)-Avi-Tag HiP™	72549
RPTOR, Strep-Tag	102070	TIM-3 (Mouse), Fc-Fusion (Human)-Avi-Tag, Biotin-Labeled HiP™	77000
Sestrin 1, FLAG-tag	50284	TIM-3, Avi-His-Tag	71262
Sestrin 2, FLAG-Tag	50286	TIM-3, Avi-Tag, Biotin Labeled (Human)	79096
Sialidase, His-Tag	101734	TIM-3, Biotin-labeled, His-tag (Human) HiP™	71237
Siglec10, Avi-His-tag	100529	TIM-3, Fc fusion (Human)	71151
Siglec11, Fc Fusion, Avi-Tag	100662	TIM-3, FLAG-tag (Human)	71180
Siglec11, Fc-Fusion (IgG1), Avi-Tag, Biotin-Labeled	100663	TIM-4, Avi-His-Tag	72513
Siglec15, Fc Fusion, Avi-tag	100423	TL1A, His-Tag, Avi-Tag	101880
Siglec15, Fc Fusion, Avi-tag, Biotin-labeled	100424	TL1A, His-Tag, Avi-Tag, Biotin-Labeled	101882
Siglec4, Fc Fusion, Avi-tag, Biotin-labeled	100532	TNFR2, Fc-Fusion (IgG1), His-Avi-Tag	79363
Siglec5, Fc Fusion, Avi-tag	100533	TNFR2, Fc-Fusion (IgG1), His-Avi-Tag, Biotin-labeled	100205

Proteins	Catalog#
TPBG, Avi-His-Tag	101514
TPBG, Avi-His-Tag, Biotin Labeled,	101515
TPBG, Fc Fusion (IgG), Avi-Tag	101516
TPBG, Fc Fusion (IgG1), Avi-Tag, Biotin Labeled	101517
TPH1, His-tag (Human)	71192
TPH2, His-tag (Human)	71193
TREM2, Fc-fusion (IgG1), Avi-Tag (Mouse)	79458
TREM2, Fc-fusion (IgG1), Avi-Tag (Mouse), Biotin	79459
TREM2, Fc-fusion (IgG1), Avi-Tag, Biotin-Labeled	79418
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
TSLP, Avi-Tag, His-Tag	102168
TSLP, Avi-Tag, His-Tag, Biotin-Labeled	102169
TSLPR, Fc Fusion	102163
TUT4, FLAG-tag (Human)	71259
TUT4, FLAG-tag, Biotin-Labeled (Human)	71258
ULBP1/NKG2DL1, Fc-Fusion (IgG1) Avi-Tag	100253
ULBP2, Avi-His-Tag	100544
ULBP2, Avi-His-Tag, Biotin-labeled	100545
ULBP3, Avi-His-Tag	100506
VEGFR3, Avi-Tag, His-Tag, Biotin-Labeled	102123
VEGFR3, Fc Fusion, Avi-Tag,	102101
VSIG3, Avi-His-Tag	100349
VSIG3, Fc-fusion (IgG1), Avi-Tag	79491
VSIG3, Fc-fusion (IgG1), Avi-Tag, Biotin-Labeled	100044
VSIG8, Fc Fusion, Avi-Tag	40743
WRN, GST-Tag	101264

Vectors	Catalog#
PD-1 sgRNA-MS2 for CRISPRa (Plasmid)	78091



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

