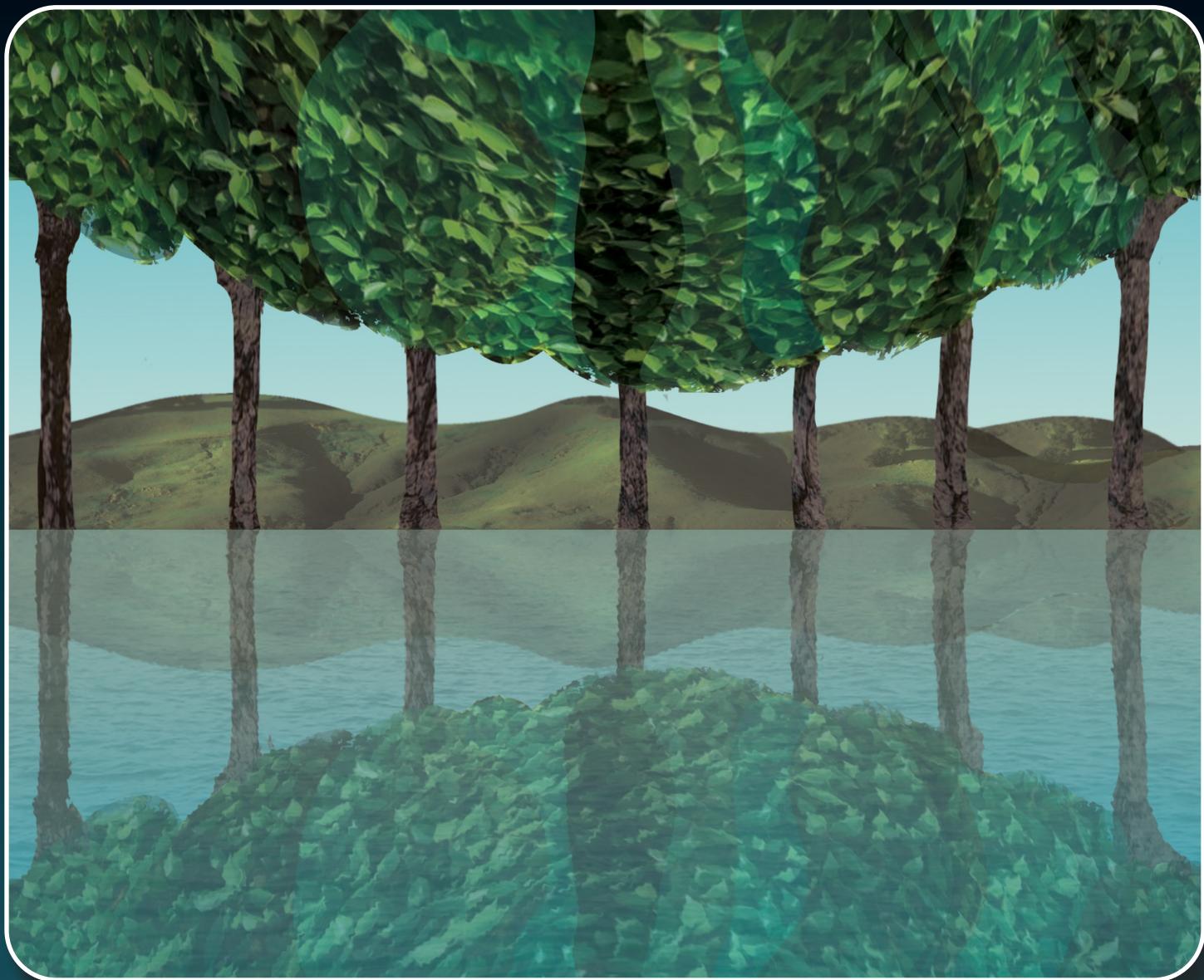




IMMUNOTHERAPY

TOOLS FOR DRUG DISCOVERY

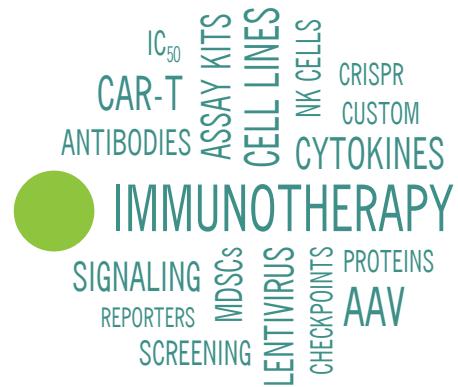
Bioactive Proteins & Antibodies | Biochemical Assay Kits | Cell Lines
Virus-Based Tools | Screening & Profiling



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 IC_{50} 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)

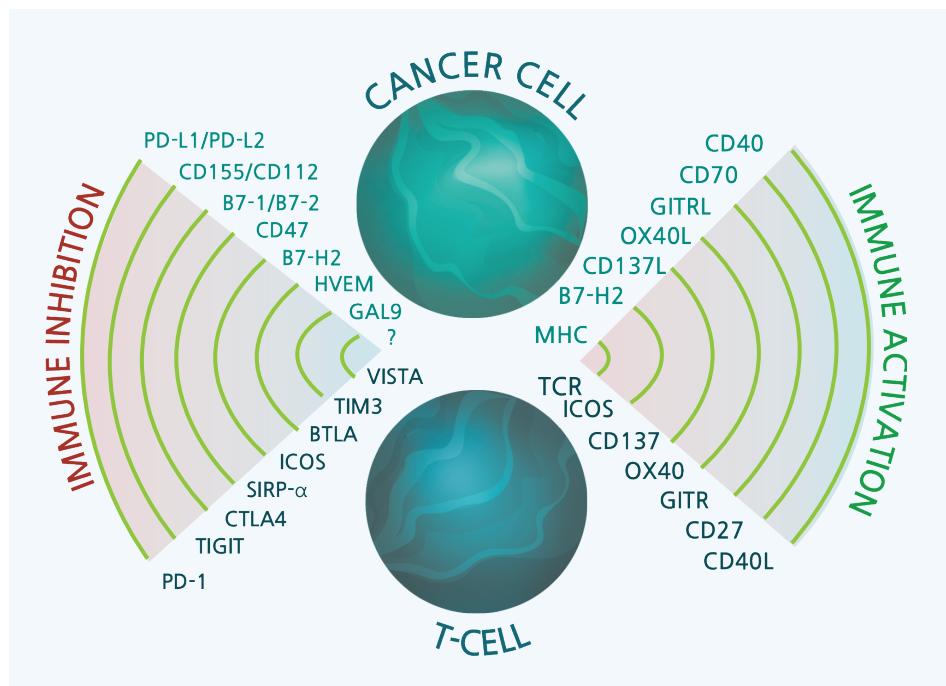
- >90% purity
- Verified via HPLC to guarantee <10% aggregation
- Ideal for studying enzyme activity, ligand binding, protein folding, and more

Bulk Production and Customization

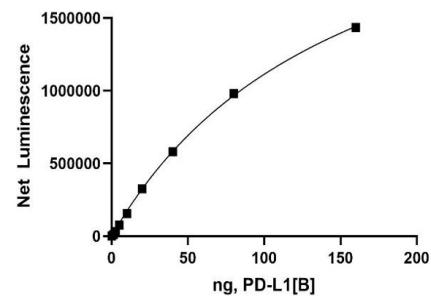
Product sizing, formulations, tags, mutations, and activity testing can be customized to meet your research needs.

Activity Testing

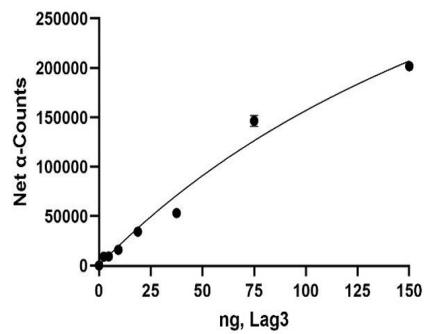
- Our proteins and antibodies are validated on each lot to confirm activity, when relevant.
- In-house developed assay kits use our quality manufactured proteins for consistent performance.



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.

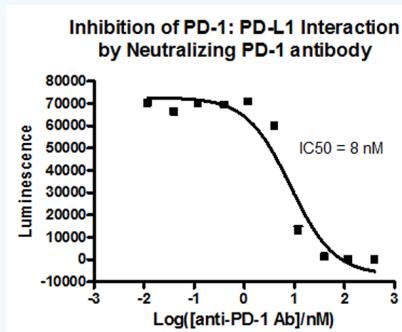
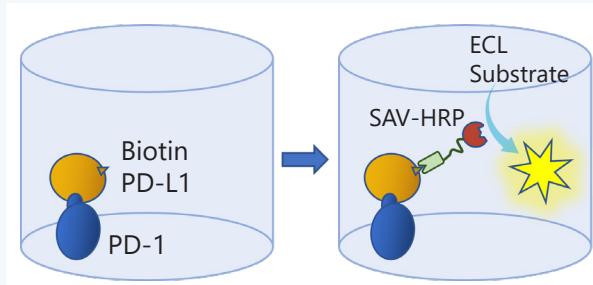
Assay kit formats include:

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

Featured Products

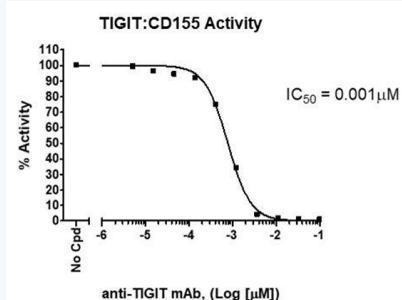
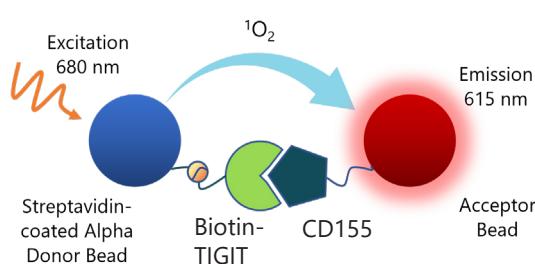
PD-1:PD-L1[Biotinylated] Inhibitor Screening Assay Kit (#72003)

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



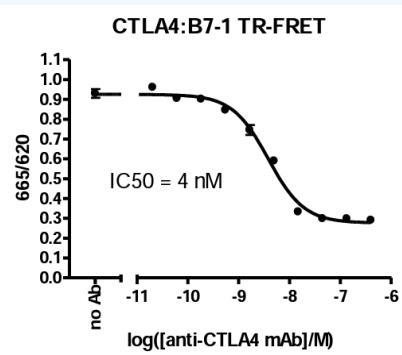
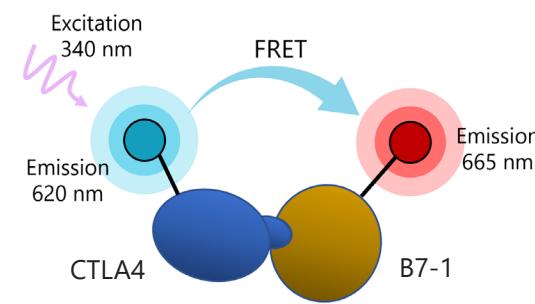
TIGIT:CD155 Homogeneous Assay Kit (#72029)

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)

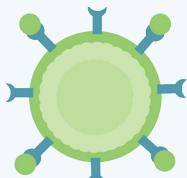
Designed to screen for inhibitors of CTLA4 binding to B7-1 in a homogeneous 384 reaction format.



Cell Lines

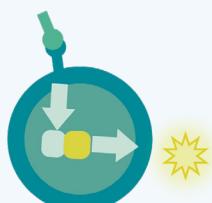
Recombinant cell lines can serve as excellent model systems in a variety of applications for immunotherapy research. Our broad selection of many novel cell line tools offers an excellent set of options for your research.

● Overexpression/Knockout Cell Lines



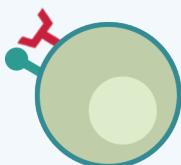
- Antibody screening
- Study ion channels, GPCR screens, and metabolite-based assays
- Flow cytometry analysis
- Binding studies

● 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



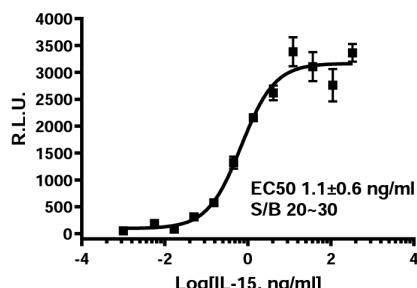
- Inhibitor screening
- Antibody affinity
- Co-culture studies

● CAR-T Cells

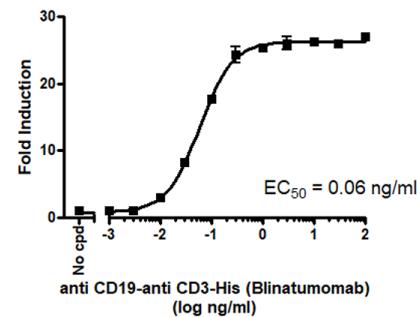


- For design of co-culture studies
- Positive controls for developing new CARs

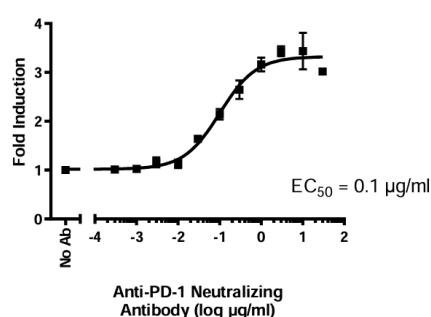
IL-15 agonist stimulation of luciferase activity in IL-15 responsive reporter cell line (#78402)



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

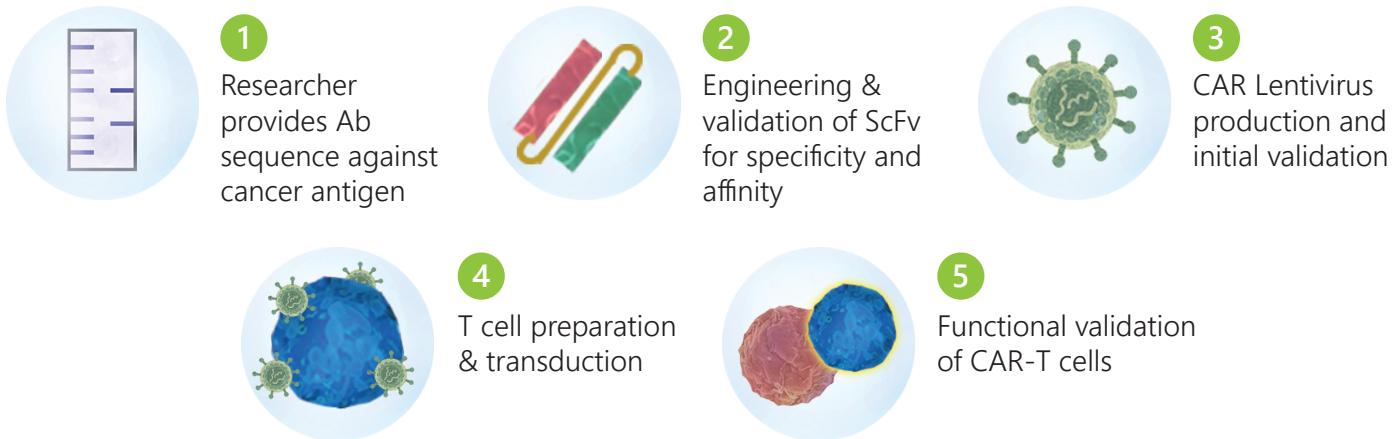


PD-L1/TCR activator cells to screen for PD-1 blocking antibodies (#60536)



CAR-T Cell Therapy Solutions

Custom CAR-T Cell Development



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



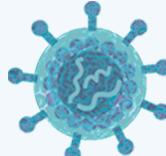
- Bispecific T cell Engager antibodies, including Anti-CD19-Anti-CD3
- CAR-T targets, such as Anti-BCMA, anti-CD19, and more
- T cell agonist antibodies, such as Anti-CD3 and Anti-CD28
- Human Ig isotype controls

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

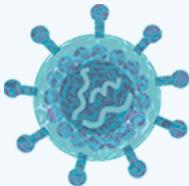


- Capable of transducing primary cells
- For long-term stable expression of a transgene
- CAR lentiviruses for Anti-CD19, Anti-CD20, Anti-BCMA and more
- CRISPR/Cas9 lentiviruses for target protein knockout

Virus-Based Tools

Virus-based tools such 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

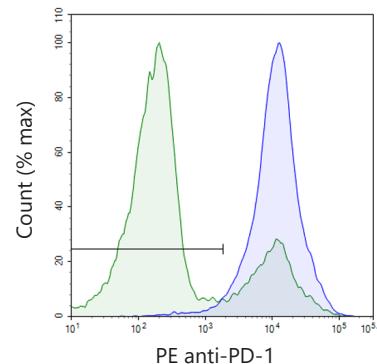
- Stable cell line generation
- Protein expression
- CRISPR/Cas9 knockout
- Generating cellular reporter assays (GFP, luciferase)
- Screen for neutralizing antibodies

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.

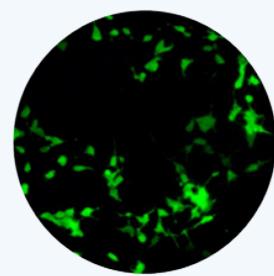


AAV Reporter Vectors



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.



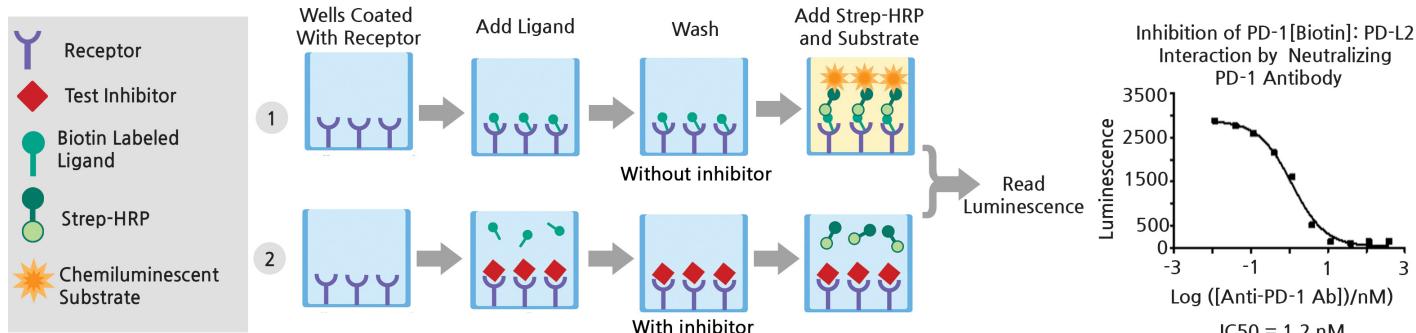
Biochemical & Cell-Based Screening & Profiling

BPS Bioscience offers rapid and dependable Immunotherapy Screening Services to determine IC₅₀ 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 IC₅₀ 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

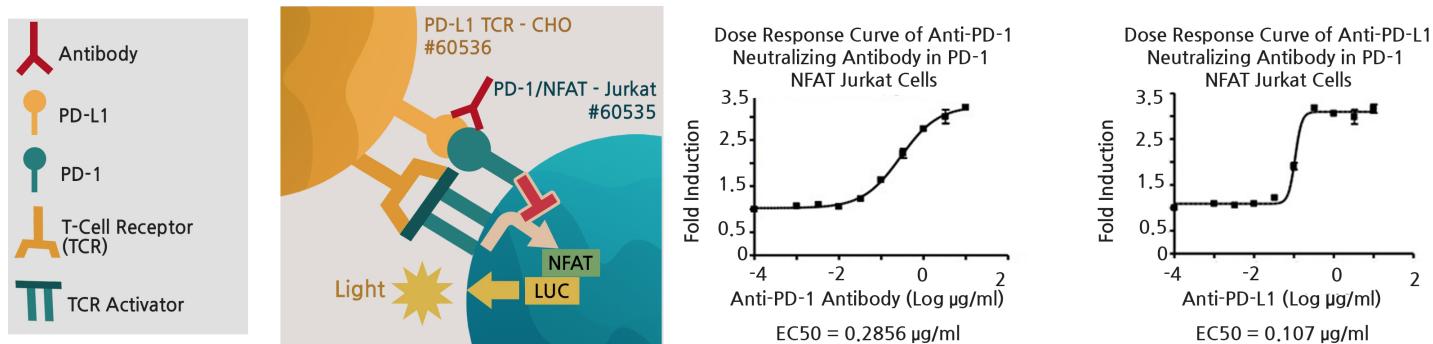
Biochemical Assays

- 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, GITR:GITR-Ligand and many more.



Antibodies	Catalog #	Antibodies	Catalog #
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-CD123 IgG Antibody, Biotin-Labeled	101140	Anti-CD8 Antibody, PE-Labeled	102011
Anti-CD137 Agonist Antibody	79097	Anti-Claudin-18 Isoform 2 Antibody, FITC-Labeled	101866
Anti-CD14 Antibody, FITC-Labeled	101884	Anti-Claudin-18 Isoform 2 Antibody, PE-Labeled	101676
Anti-CD14 Antibody, PE-Labeled	101883	Anti-Claudin-18 Isoform 2 IgG Antibody	101564
Anti-CD16a Antibody, FITC-Labeled	101979	Anti-Claudin-18 Isoform 2 IgG Antibody, Biotin-Labeled	101565
Anti-CD16a Antibody, PE-Labeled	101978	Anti-Claudin-18 Isoform2-Anti-CD3 IgG Bispecific Antibody	101541
Anti-CD19 Antibody, FITC-Labeled	101863	Anti-CTLA4 Neutralizing Antibody	71212
Anti-CD19 Antibody, PE-Labeled	101625	Anti-FcGR3A (CD16A) IgG Antibody, Biotin-Labeled	101225
Anti-CD19 IgG Antibody	100981	Anti-FcGR3A-Anti-EGFR Bispecific Antibody	101535
Anti-CD19 IgG Antibody, Biotin-Labeled	101093	Anti-GITR Agonist Antibody	79053
Anti-CD19-Anti-CD3 Bispecific Molecule	100441	Anti-GITR Antibody, PE-labeled	71295-1
Anti-CD19-Anti-CD3 IgG format Bispecific Antibody	101076	Anti-GITR Antibody, PE-labeled	71295-2
Anti-CD2 IgG Antibody	101213	Anti-HLA-DR Biotin-Labeled, Antibody	101769
Anti-CD20 Antibody, FITC-Labeled	101864	Anti-Human IgG, Unconjugated Antibody	100736
Anti-CD20 Antibody, PE-Labeled	101672	Anti-Human IgM, Unconjugated Antibody	100737
Anti-CD20 Functional Antibody	71209	Anti-IL-17A Neutralizing Antibody	91015
Anti-CD20 IgG Antibody, Biotin-labeled	101207	Anti-IL-2RA (CD25) Neutralizing Antibody	101593
Anti-CD20-Anti-CD3 Bispecific Antibody	100836	Anti-IL23 α Neutralizing Antibody	101453
Anti-CD20-Anti-CD3 IgM format Bispecific Antibody	100860	Anti-LAG3 Antibody, PE-Labeled	71226-1
Anti-CD235a Antibody, PE-Labeled	101675	Anti-LAG3 Antibody, PE-Labeled	71226-2
Anti-CD235a IgG1 Antibody, Biotin-Labeled	101098	Anti-LAG3, Neutralizing Antibody	71219
Anti-CD27 Agonist Antibody	100111	Anti-Mesothelin-Anti-CD3 Bispecific Antibody	101621
Anti-CD28 Agonist Antibody	100182	Anti-NCAM1 (CD56) IgG Antibody, Biotin-labeled	101112
Anti-CD28 Agonist Antibody (Humanized)	100186	Anti-NCAM1 Antibody, FITC-Labeled	101865
Anti-CD3 Agonist Antibody	71274-1	Anti-NCAM1 Antibody, PE-Labeled	101673
Anti-CD3 Agonist Antibody	71274-2	Anti-OX40 Antagonist Antibody	72063-1
Anti-CD3 Antibody, FITC-Labeled	102008	Anti-OX40 Antagonist Antibody	72063-2
Anti-CD3 Antibody, PE-Labeled	102009	Anti-OX40 Antibody, PE-Labeled	72064-2
Anti-CD38 Antibody	79120	Anti-OX40 Antibody, PE-Labeled	72064-1

Antibodies	Catalog #	Biochemical Assay Kits	Catalog #
Anti-OX40 Competitive Antibody	72558-1	Cathepsin D Inhibitor Screening Assay Kit	82141
Anti-OX40 Competitive Antibody	72558-2	CD112R:CD112[Biotinylated] Inhibitor Screening Assay Kit	79732
Anti-PD-1 (S228P) Neutralizing Antibody	71326	CD137:CD137L[Biotinylated] Inhibitor Screening Assay Kit	78590
Anti-PD-1 Agonist Antibody	101178	CD137[Biotinylated]:CD137L Inhibitor Screening Assay Kit	72025
Anti-PD-1 Neutralizing Antibody	71120	CD226:CD112 Homogeneous Assay Kit	72051
Anti-PD-1 Neutralizing Antibody, Biotin-labeled	71215	CD226:CD155 Homogeneous Assay Kit	72052
Anti-PD-1 Neutralizing Antibody, PE-labeled	71290	CD27:CD70[Biotinylated] Inhibitor Screening Assay Kit	79695
Anti-PD-L1 (CD274) Neutralizing Antibody	71213	CD28:B7-1[Biotinylated] Inhibitor Screening Assay Kit	72007
Anti-PD-L1 Antibody, PE-labeled	71128	CD28:B7-2[Biotinylated] Inhibitor Screening Assay Kit	72062
Anti-PD-L1 Neutralizing Antibody, Biotin-Labeled	71214	CD38 (Dog) Inhibitor Screening Assay Kit (Hydrolase Activity)	78108
Anti-PD-L1-Fab Avi-Tag, Biotin-Labeled	11032	CD38 (Mouse) Fluorogenic Assay Kit (Cyclase Activity)	78285
Anti-PSMA Antibody	101695	CD38 (Mouse) Inhibitor Screening Assay Kit (Hydrolase Activity)	79682
Anti-PSMA Antibody, FITC-Labeled	101977	CD38 (Pig) Fluorogenic Assay Kit (Hydrolase Activity)	78178
Anti-PSMA Antibody, PE-Labeled	101976	CD38 (Rat) Inhibitor Screening Assay Kit (Hydrolase Activity)	79690
Anti-PSMA, Biotin Label Antibody	101757	CD38 Inhibitor Screening Assay Kit (Cyclase Activity)	71275
Anti-PSMA-Anti-CD3 IgG format Bispecific Antibody	101242	CD38 Inhibitor Screening Assay Kit (Hydrolase Activity)	79287
Anti-PVRIG Neutralizing Antibody	101712	CD38 Inhibitor Screening Assay Kit (Hydrolase Activity)	79672
Anti-RANKL Neutralizing Antibody	100874	CD39 Inhibitor Screening Assay Kit	79278
Anti-TFR Antibody, Biotin-Labeled	101596	CD40:CD40L TR-FRET Assay	79258
Anti-TIGIT Antibody, PE-Labeled	71228-1	CD40:CD40L[Biotinylated] Inhibitor Screening Assay Kit	79257
Anti-TIGIT Antibody, PE-Labeled	71228-2	CD47:SIRP- α [Biotinylated] Inhibitor Screening Assay Kit	72044
Anti-TIGIT Inhibitor Antibody	71218	CD47:SIRP- γ [Biotinylated] Inhibitor Screening Assay Kit	72059
Anti-TIGIT Neutralizing Antibody	71340	CD73 Inhibitor Screening Assay Kit	72055
Anti-TIM-3 Antibody, PE-Labeled	71121	CD73 Inhibitor Screening Assay Kit	72058
Anti-Tim3 Antibody	79098	Chemi-Verse™ BLK Kinase Assay Kit	82229
Anti-TL1A Neutralizing Antibody	101729	Chemi-Verse™ CDK1/CyclinA2 Kinase Assay Kit	82242
Anti-Transferrin Receptor Antibody, FITC-Labeled	101876	Chemi-Verse™ CDK2/CyclinA2 Kinase Assay Kit	82247
Anti-Transferrin Receptor Antibody, PE-Labeled	101875	Chemi-Verse™ CDK2/CyclinE1 Kinase Assay Kit	82228
Anti-VEGF Neutralizing Antibody	79478-1	Chemi-Verse™ CDK9/CyclinT Kinase Assay Kit	82233
Anti-VEGF Neutralizing Antibody	79478-2	Chemi-Verse™ LYN A Kinase Assay Kit	82224
Anti-VISTA Neutralizing Antibody	101293	Chemi-Verse™ MAPKAPK5 Kinase Assay Kit	82236
		Chemi-Verse™ NUAK1 Kinase Assay Kit	82231
		Chemi-Verse™ NUAK2 Kinase Assay Kit	82232
ADAR1:RNA TR-FRET Assay Kit	82252	Chemi-Verse™ PAK6 Kinase Assay Kit	82182
ARG1 Inhibitor Screening Assay Kit	72048	Chemi-Verse™ PAK7 Kinase Assay Kit	82183
ARG2 Inhibitor Screening Assay Kit	72043	Chemi-Verse™ VRK1 Kinase Assay Kit	82230
BAFF:BCMA[Biotinylated] Inhibitor Screening Assay Kit	79667	CTLA4 : B7-2[Biotinylated] Inhibitor Screening Assay Kit	79658
BCMA:APRIL[Biotinylated] Inhibitor Screening Assay Kit	79722	CTLA4 [Biotinylated]:B7-1 Inhibitor Screening Assay Kit	79973
BTLA:HVEM[Biotinylated] Inhibitor Screening Assay Kit	72008	CTLA4:B7-1 TR-FRET Assay Kit	72120

Biochemical Assay Kits	Catalog #	Biochemical Assay Kits	Catalog #
CTLA4:B7-1[Biotinylated] Inhibitor Screening Assay Kit	72009	Mouse ITK Kinase Assay Kit	78520
CTLA4[Biotinylated]:B7-2 Inhibitor Screening Assay Kit	72024	Mouse PD-1[Biotinylated]:PD-L1 Inhibitor Screening Assay Kit	72027
DcR3:TL1A Inhibitor Screening Assay Kit	82160	NAMPT Inhibitor Screening Assay Kit	71276
DR3:TL1A Inhibitor Screening Assay Kit	82241	NEU2 Fluorogenic Assay Kit	82139
ERK1 Kinase Assay Kit	82282	OX40[Biotinylated]:OX40L Inhibitor Screening Assay Kit	72045
Fc (IgG1): FcRn Inhibitor Screening Colorimetric Assay Kit	78501	PADI4 Fluorogenic Assay Kit	82111
FGL1:LAG3 Inhibitor Screening Assay Kit	78824	PD-1:PD-L1 Cell-Based Inhibitor Screening Assay Kit	79377
FGL1:LAG3 TR-FRET Assay Kit	79739	PD-1:PD-L1 Homogeneous Assay Kit	72014
GDF15:GFRAL[Biotinylated] Inhibitor Screening Assay Kit	78807	PD-1:PD-L1 TR-FRET Assay	72032
GITR:GITRL TR-FRET Assay Kit	79054	PD-1:PD-L1 TR-FRET Assay	72038
GITRL:GITR[Biotinylated] Inhibitor Screening Assay Kit	72061	PD-1:PD-L1/PD-L2 Cell-Based Inhibitor Screening Assay Kit	60800
GPX4 Fluorogenic Assay Kit	82165	PD-1:PD-L1[Biotinylated] Inhibitor Screening Assay Kit	72003
HSA:FcRn Neutralizing Antibody Screening Chemiluminescence Assay Kit	82151	PD-1:PD-L1[Biotinylated] Inhibitor Screening Colorimetric Assay Kit	72016
ICOSL:ICOS[Biotinylated] Inhibitor Screening Assay Kit	79673	PD-1:PD-L2 Cell-Based Inhibitor Screening Assay Kit	79378
IDO1 Fluorogenic Inhibitor Screening Assay Kit	72037	PD-1:PD-L2 Homogeneous Assay Kit	72015
IDO1 Fluorogenic Inhibitor Screening Assay Kit -384	72047	PD-1:PD-L2 TR-FRET Assay Kit	72012
IDO1 Inhibitor Mechanism of Action Assay Kit	72050	PD-1:PD-L2[Biotinylated] Inhibitor Screening Assay Kit	72004
IDO1 Inhibitor Screening Assay Kit	72021	PD-1:PD-L2[Biotinylated] Inhibitor Screening Colorimetric Assay Kit	72017
IDO1 Inhibitor Screening Assay Kit -384	72034	PD-1[Biotinylated]: PD-L1 Inhibitor Screening Assay Kit	79523
IDO2 Inhibitor Screening Assay Kit	72022	PD-1[Biotinylated]:PD-L1 Homogeneous Assay Kit	72028
IL-17RA:IL-17A[Biotin] Inhibitor Screening Assay Kit	79891	PD-1[Biotinylated]:PD-L1 Inhibitor Screening Assay Kit	72005
IL-17RA[Biotin]:IL-17A Inhibitor Screening Assay Kit	72060	PD-1[Biotinylated]:PD-L1 Inhibitor Screening Colorimetric Assay Kit	72018
IL-6:IL-6R Inhibitor Screening Assay Kit	78027	PD-1[Biotinylated]:PD-L2 Inhibitor Screening Assay Kit	72006
KMO Inhibitor Screening Assay Kit	79513	PD-1[Biotinylated]:PD-L2 Inhibitor Screening Colorimetric Assay Kit	72019
Kynureninase Inhibitor Screening Assay Kit	79514	PD-1[Biotinylated]:PD-L2 Inhibitor Screening Assay Kit	72019
LAG3:FGL1[Biotinylated] Inhibitor Screening Assay Kit	79789	PD-L1:B7-1 [Biotinylated] Inhibitor Screening Assay Kit (Mouse)	79821
LIGHT:HVEM [Biotinylated] Inhibitor Screening Assay Kit	79684	PD-L1:B7-1[Biotinylated] Inhibitor Screening Assay Kit	72026
LT β R:LIGHT[Biotinylated] Inhibitor Screening Assay Kit	79808	RANK:RANKL TR-FRET Assay	79101
LysA™ Universal PARylation Assay Kit	82123	TDO Fluorogenic Inhibitor Screening Assay Kit -384 (Human)	72049
MAT2A Inhibitor Screening Assay Kit	71402	TDO Fluorogenic Inhibitor Screening Assay Kit (Human)	72039
MDM2-Driven p53 Ubiquitination Assay Kit	82179	TDO Inhibitor Screening Assay Kit	72023
MMP7 Fluorogenic Assay Kit	78864	TDO Inhibitor Screening Assay Kit -384	72036
Molecular Glue/PROTAC® Optimization Kit for RBM39-DCAF15	82251	Tie2 Kinase Assay Kit	78573
Monkey PD-1:PD-L1[Biotinylated] Inhibitor Screening Assay Kit	79863	TIGIT (Mouse): CD155 (Mouse) Homogeneous Assay Kit	78141
Monkey PD-1[Biotinylated]:PD-L1 Inhibitor Screening Assay Kit	79865	TIGIT:CD112 Homogeneous Assay Kit	72030
Mouse CTLA4:B7-1 TR-FRET Assay	79508	TIGIT:CD155 Homogeneous Assay Kit	72029
Mouse CTLA4[Biotin]:B7-1 Inhibitor Screening Assay Kit	79515	TNFR2:TNF-alpha[Biotinylated] Inhibitor Screening Assay Kit	79756
Mouse IDO1 Inhibitor Screening Assay Kit	72041	TPH1 Inhibitor Screening Assay Kit	72053
Mouse IDO2 Inhibitor Screening Assay Kit	72042	TPH1 Inhibitor Screening Assay Kit	72056
		TPH2 Inhibitor Screening Assay Kit	72054

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

Biochemical Assay Kits	Catalog #	Cell Lines	Catalog #
TPH2 Inhibitor Screening Assay Kit	72057	CD137 (4-1BB) HEK293 Recombinant Cell Line	60691
Universal IDO1/IDO2/TDO Inhibitor Screening Assay Kit	72035	CD137L (4-1BBL) CHO-K1 Recombinant Cell Line	60523
VSIG-3:VISTA [Biotinylated] Inhibitor Screening Assay Kit	79782	CD14 CHO Recombinant Cell Line	78335
WRN Helicase Activity Assay Kit	78852	CD155 (PVR) - HEK293 Recombinant Cell Line	60537
		CD155 / TCR Activator - CHO Recombinant Cell line	60548
Cell Isolation Kits/Components	Catalog #		
10x Cell Isolation Buffer	78563	CD160/NFAT - Luciferase Reporter - Jurkat Recombinant Cell Line	79594
CD14 Positive Cell Isolation Kit	78897	CD19 / BCMA / Firefly Luciferase - CHO Recombinant Cell Line	78030
CD19 Positive Cell Isolation Kit	78564	CD19 / CD20 / Firefly Luciferase CHO Cell Line	78186
CD20 Positive Cell Isolation Kit	82132	CD19 / Firefly Luciferase - CHO Recombinant Cell Line	79714
CD235a Positive Cell Isolation Kit	78896	CD19 CHO Recombinant Cell Line (Low, Medium and High Expression)	79561
Cell Isolation Magnetic Tube Rack	78571	CD19 Knockout Raji Cell Line	82166
NCAM1/CD56 Positive Cell Isolation Kit	78808	CD2 CHO Cell Line	78330
		CD20 CHO Recombinant Cell Line (High or Medium Expression)	79624
Cell Lines	Catalog #		
ADAR1 Responsive Luciferase Reporter HEK293 Cell Line	82238	CD20/Firefly Luciferase CHO Cell Line	78620
ADCC Bioassay Effector Cell (Mouse) Jurkat Cell Line	79733	CD22 / Luciferase - CHO Recombinant Cell Line	79715
ADCC Bioassay Effector Cell F Variant (Low Affinity)/ NFAT Luciferase Reporter Jurkat Cell Line	60540	CD22 CHO Recombinant Cell Line (Medium and High Expression)	79557
ADCC Bioassay Effector Cell V Variant (High Affinity)/ NFAT Luciferase Reporter Jurkat Cell Line	60541	CD27 CHO-K1 Stable Recombinant Cell Line	60624
ADCP Bioassay Effector Cell FcγRIIa (H Variant)/ NFAT Luciferase Reporter Jurkat Cell Line	71273	CD27/NF-κB Reporter-Jurkat Recombinant Cell Line	79509
Adenosine A2A Receptor Functional HEK293 Cell Line	79381	CD33- CHO K1 Recombinant Cell Line (High, Medium and Low Expression)	79935
Anti-BCMA CAR /NFAT (Luciferase) Reporter Jurkat Cell Line	79694	CD34 CHO Recombinant Cell Line	78336
Anti-CD19 CAR / NFAT (Luciferase) Reporter Jurkat Cell Line (CD19 SCFV-CD28-4-1BB-CD3ζ)	79853	CD36 HEK293 Cell Line (High or Medium Expression)	78526
Anti-CD19 CAR Negative Control/NFAT (Luciferase) Reporter Jurkat Cell Line (CD19 SCFV-CD28 Transmembrane Motif)	79854	CD36 Knockout THP-1 Cell Line	82190
B2M Knockout Jurkat Cell Line	78342	CD37 CHO Recombinant Cell Line (High or Low Expression)	79607
B2M Knockout NFAT Luciferase Reporter Jurkat Cell Line	78363	CD38 / BCMA / Firefly Luciferase CHO Recombinant Cell Line	78148
B2M Knockout THP-1 Cell Line	78389	CD38 / CD19 / Firefly Luciferase CHO Recombinant Cell Line	78149
B2M/CIITA Double Knockout THP-1 Cell Line	78391	CD38 CHO Recombinant Cell Line (High, Medium or Low Expression)	79615
B7-H3 CHO Cell Line (High, Medium, or Low Expression)	78541	CD40 - HEK293 Cell Line	71257
B7-H4 CHO Cell Line (Low, Medium or High Expression)	78586	CD40/NF-κB Luciferase Reporter HEK293 Cell Line	60626
B7-H7 (HHLA2)/TCR Activator CHO Cell Line	78321	CD40L K562 Cell Line (Medium and High Expression)	78942
BAFF-R - CHO K1 Recombinant Cell Line (High, Medium or Low Expression)	79921	CD43 CHO Cell Line	78334
BCMA / CD20 / Firefly Luciferase CHO Cell Line	78185	CD47 - HEK293 Cell Line	71249
BCMA / GLuc - CHO Recombinant Cell Line	79830	CD5- CHO K1 Recombinant Cell Line (High, Medium, or Low Expression)	78290
BCMA / Luciferase - CHO Recombinant Cell Line	79724	CD7 CHO Cell Line (Medium or High Expression)	78324
BCMA / NF-κB - Reporter HEK293 Recombinant Cell Line	79755	CD70-CHO Recombinant Cell line	79510
BCMA CHO Recombinant Cell Line (High or Low Expression)	79500	CD8+ TCR Knockout NFAT-Luciferase Reporter Jurkat Cell Line	78757
CD123 (IL3Ra) - CHO Recombinant Cell Line (High, Medium, or Low Expression)	79640	CD8a/CD8b Jurkat Cell Line	78917
CD137 (4-1BB) /NF-κB Reporter - HEK293 Recombinant Cell Line	79289	CD95 CHO Cell Line (Medium or High Expression)	78499
		CEACAM5 CHO Cell Line (High, Medium, or Low Expression)	78704
		CEACAM6 CHO Cell Line	78550

Cell Lines	Catalog #	Cell Lines	Catalog #
Claudin-1 CHO Cell Line	82097	FOLR1 CHO Cell Line	78360
Claudin-18 Isoform 1 CHO Cell Line	78361	Full Length Notch1/CSL Luciferase Reporter HEK293 Cell Line	82201
Claudin-18 Isoform 2 CHO Cell Line (High, Medium, or Low Expression)	78533	FUT4-CD15 CHO Cell Line	78417
Claudin-3 CHO Cell Line	78566	GITR / NF-κB Luciferase Reporter - Jurkat Cell Line	60546
Claudin-4 CHO Cell Line	78567	GITR - HEK293 Recombinant Cell Line	79092
Claudin-6 CHO Cell Line (High, Medium, or Low) Expressions	78527	GITRL CHO-K1 Recombinant Cell Line	60547
Claudin-9 CHO Cell Line	78687	Glycophorin A (CD235a) CHO Cell Line	78353
CTLA4 - HEK293 Recombinant Cell Line	60681	GPC3 - CHO-K1 Recombinant Cell Line	78100
CTLA4 / IL-2 Reporter - Jurkat Recombinant Cell Line	79525	GPRC5D (Cynomolgus) CHO Cell Line	78338
DLL3 CHO Cell Line	78882	GPRC5D (Cynomolgus) HEK293 Cell Line	78346
DLL4 CHO Cell Line	82217	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	HLA-C*08:02 K562 Cell Line	78974
FAP- CHO K1 Recombinant Cell Line (High, Medium or Low Expression)	79947	Human EGFRvIII CHO K1 Cell Line (High or Low Expression)	78145
FcGR1a (CD64) Knockout THP-1 Cell Line	82191	Human Mesothelin - CHO-K1 Recombinant Cell Line	78132
FCGR2A (CD32A) Knockout Jurkat Cell Line	78549	HVEM - CHO Recombinant Cell Line	79297
FcGR2B CHO K1 Recombinant Cell Line	79511	HVEM - HEK293 Recombinant Cell Line	79313
FcGR3A (CD16A) CHO Cell Line	78332	HVEM/NF-κB Reporter Jurkat Recombinant Cell Line	79310
FcGR3B (CD16B) CHO Cell Line	78333	ICOS/NFAT Reporter-Jurkat Recombinant Cell Line	79668
FcRL5 CHO Cell Line	78375	ICOSL-CHO Recombinant Cell Line	79635
FcRL5 HEK293 Cell Line	78374	IDO1-HEK293 Recombinant Cell Line (Human)	60532
FcRL5/Firefly Luciferase RPMI 8226 Cell Line	78378	IDO2 - HEK293 Recombinant Cell Line (Human)	60533
FcRn: IgG Recycling HMEC-1 Cell Pool	82163	IL-15 Responsive Luciferase Reporter Cell Line	78402
Firefly Luciferase CCRF-CEM Cell Line	78495	IL-2 Luciferase Reporter Jurkat Cell Line	60481
Firefly Luciferase CD19 Knockout NALM6 Cell Line	82168	IL-4/IL-13 Responsive STAT6 Luciferase Reporter HEK293 Cell Line	78941
Firefly Luciferase HL-60 Cell Line	78410	IL-6 Responsive Luciferase Reporter HEK293 Cell Line	82235
Firefly Luciferase Jurkat Cell Line	78373	KIR3DL3/IL-2 Luciferase Reporter Jurkat Cell Line	78322
Firefly Luciferase K562 Cell Line	78621	LAG3 / IL-2 Reporter - Jurkat Recombinant Cell Line	79813
Firefly Luciferase KG1 Cell Line	78493	LAG3 / NFAT Reporter - Jurkat Recombinant Cell Line	71278
Firefly Luciferase MC38 Cell Line	78383	LAIR1 CHO Cell Line	78913
Firefly Luciferase Molm13 Cell Line	78372	LIGHT-CHO Recombinant Cell Line	79262
Firefly Luciferase NALM6 Cell Line	78494	LILRB4 NFAT-Luciferase Reporter Jurkat Recombinant Cell Line	79750
Firefly Luciferase Raji Cell Line	78622	MUC16 (CA125), variant 4 (region 13785-14507) CHO Cell Line	78848
Firefly Luciferase SKOV-3 Cell Line	78425	NCAM1 (CD56) CHO Cell Line (High, Medium, or Low Expression)	78352
Firefly Luciferase THP-1 Cell Line	78409	Nectin4 - CHO K1 Recombinant Cell Line (High, Medium, or Low Expression)	78097
Firefly Luciferase - RPMI 8226 Recombinant Cell Line	79834	NF-κB Luciferase Reporter Jurkat Cell Line	60651

Cell Lines	Catalog #	Cell Lines	Catalog #
NFAT Luciferase Reporter Jurkat Cell Line	60621	TIGIT / NFAT Reporter - Jurkat Cell Line	60538
NFAT Luciferase-eGFP Reporter Jurkat Cell Line	78662	Transferrin Receptor Protein-1 (TFRC) HEK293 Cell Line	78308
NFAT Reporter (eGFP) Jurkat Recombinant Cell Line	78384	TROP2 - CHO-K1 Recombinant Cell Line	78099
NKp46 CHO Cell Line (High, Medium, or Low Expression)	78915	VSIG4 CHO Cell Line	78914
NY-ESO-1 TCR (1G4) CD8+ NFAT-Luciferase Reporter Jurkat Cell Line	78769		
		Cell-Based Assays and Expression Kits	Catalog #
NY-ESO-1 TCR (c259) CD8+ NFAT-Luciferase Reporter Jurkat Cell Line	78771	CSF1R / SRE Reporter Kit (MAPK/ERK Signaling Pathway)	79379
OX40 - HEK293 Recombinant Cell Line	60682	IDO1 Cell-Based Assay Kit	72031
OX40/NF-κB Luciferase Reporter HEK293 Cell Line	60482	Mouse PD-L1 / TCR Activator Mammalian Expression Kit	79778
p53 Luciferase Reporter HCT116 Cell Line	78681	PBMC Cytotoxicity Bioassay Kit (CFSE, 7-ADD)	82173
PAI-1 Reporter (Luc) - Mv1 Lu Cell Line	60544	PBMC Cytotoxicity Luciferase Assay Kit	82214
PARG Knockout HeLa Cell Line	82171	PBMC Cytotoxicity Luciferase Assay Kit (NALM6)	82174
PARP1 Knockout HeLa Cell Line	82169	PD-L1 (Woodchuck) / TCR Activator Mammalian Expression Kit	79455
PD-1 (Mouse) / NFAT - Reporter - Jurkat Recombinant Cell Line	79762	PD-L1 / TCR Activator Mammalian Expression Kit	60610
PD-1 (Woodchuck) / NFAT - Reporter - Jurkat Recombinant Cell Line	79456	PD-L2/TCR Activator Mammalian Expression Kit	60620
PD-1 - HEK293 Recombinant Cell Line	60680	TDO Cell-Based Assay Kit	72033
PD-1 / NFAT Reporter - Jurkat Recombinant Cell Line	60535		
		Inhibitors/Activators	Catalog #
PD-1 CHO Cell Line	78530	AMP (500 μM)	79496
PD-1/IL-2 Luciferase Reporter Jurkat Cell Line	82117	Epacadostat (INCBO24360)	27339-1
PD-L1 (Mouse)/TCR Activator CHO Cell Line	79763	Epacadostat (INCBO24360)	27339-2
PD-L1 (Woodchuck) / TCR activator - CHO Recombinant Cell line	79457	INCBO24360 Analog	27338-1
PD-L1 - CHO Recombinant Cell Line	60543	INCBO24360 Analog	27338-2
PD-L1 / TCR Activator - CHO Recombinant Cell Line	60536	NLG919	27337-1
PD-L2 / TCR activator - CHO Recombinant Cell Line	79632	NLG919	27337-2
PSMA (FOLH1) - CHO Recombinant Cell Line (High, Medium, or Low Expression)	79641	Telotristat	27340
RANK/NF-κB Luciferase Reporter HEK293 Cell Line	78102	Telotristat	27341
ROR1 CHO Recombinant Cell Line (High, Medium, or Low Expression)	79609		
		Lentiviruses	Catalog #
SIRP-α / HEK293 Recombinant Cell Line	60689	Anti-BCMA CAR Lentivirus (Clone C11D5.3 ScFv-CD8-4-1BB-CD3ζ)	78655
SLAMF7 (CS1) CHO Recombinant Cell Line (High, Medium, or Low Expression)	79608	Anti-BCMA CAR Lentivirus (VHH1/VHH2 ScFv-CD8-4-1BB-CD3ζ)	78783
STAT3 Luciferase Reporter THP-1 Cell Line	78498	Anti-CD19 CAR Lentivirus (CD19 ScFv-CD8-4-1BB-CD3ζ)	78600
STAT3 Reporter Jurkat Cell Line	78497	Anti-CD19 CAR Lentivirus (CD19 ScFv-CD8-4-1BB-CD3ζ, eGFP)	78775
TCR Activator CHO Recombinant Cell line	60539	Anti-CD19 CAR Lentivirus (CD19 ScFv-CD8-4-1BB-CD3ζ, PuroR)	78602
TCR Activator Raji Cell Line	60556	Anti-CD19 CAR Lentivirus (CD19 ScFv-CD8-4-1BB-CD3ζ; SIN Vector)	78601
TCR Activator/FcGR2B CHO Cell Line	78436	Anti-CD19/CD22 Bispecific CAR Lentivirus (Clones FMC63/m971 ScFv-CD8-4-1BB-CD3ζ)	78609
TCR Knockout Jurkat Cell Line	78539	Anti-CD20 CAR Lentivirus (Clone Leu-16 ScFv-CD8-4-1BB-CD3ζ)	78606
TCR Knockout NFAT-Luciferase Reporter Jurkat Cell Line	78556	Anti-CD22 CAR Lentivirus (Clone m971 ScFv-CD8-4-1BB-CD3ζ)	78608
TCR/B2M Knockout Jurkat Cell Line	78552	Anti-Mesothelin CAR Lentivirus (P4 ScFv-CD8-4-1BB-CD3ζ)	78703
TCR/B2M Knockout NFAT Luciferase Reporter Jurkat Cell Line	78557		
TDO-HEK293 Recombinant Cell line	60534		
TIGIT - HEK293 Recombinant Cell Line	79332		

Lentiviruses	Catalog #	Lentiviruses	Catalog #
ATF6 Luciferase Reporter Lentivirus (ATF6 Pathway)	78667	Nectin-4 Lentivirus	78712
B7-H4 Lentivirus	78727	NKp46 Lentivirus (Macaca fascicularis/Cynomolgus)	78779
BCMA CRISPR/Cas9 Lentivirus (Integrating)	78893	NLRP3 CRISPR/Cas9 Lentivirus (Integrating)	78545
BCMA CRISPR/Cas9 Lentivirus (Non-Integrating)	78894	NLRP3 CRISPR/Cas9 Lentivirus (Non-Integrating)	78546
BCMA Lentivirus	78714	PD-1 (Human) sgRNA-MS2 Lentivirus (Integrating)	78190
CD40 Ligand (CD40L) Lentivirus	78931	PD-1 CRISPR/Cas9 Lentivirus (Integrating)	78052
CD47 CRISPR/Cas9 Lentivirus (Integrating)	78056	PD-1 CRISPR/Cas9 Lentivirus (Non-Integrating)	78059
CD47 CRISPR/Cas9 Lentivirus (Non-Integrating)	78063	PD-L1 CRISPR/Cas9 Lentivirus (Integrating)	78057
CD5 (Human) CRISPR/Cas9 Lentivirus (Integrating)	78119	PD-L1 CRISPR/Cas9 Lentivirus (Non-Integrating)	78064
CD5 (Human) CRISPR/Cas9 Lentivirus (Non-Integrating)	78198	PD-L1 Lentivirus	78925
CD8a Lentivirus	78648	PSMA Lentivirus	78726
CD8a/CD8b Lentivirus	78650	STAT6 Luciferase Reporter Lentivirus (STAT6 Signaling Pathway)	78799
CEACAM5 Lentivirus	78719	TCR CRISPR/Cas9 Lentivirus (Non-Integrating)	78062
CEACAM6 Lentivirus	78720	TGFBR2 CRISPR/Cas9 Lentivirus (Integrating)	78535
Claudin-3 Lentivirus	78722	TGFBR2 CRISPR/Cas9 Lentivirus (Non-Integrating)	78536
Claudin-4 Lentivirus	78723	TIGIT CRISPR/Cas9 Lentivirus (Integrating)	78058
Claudin-9 Lentivirus	78721	TIGIT CRISPR/Cas9 Lentivirus (Non-Integrating)	78065
CSL (CBF1/RBP-Jk) Luciferase Reporter Lentivirus (Notch Signaling Pathway)	78746	TNFR2 Lentivirus	78765
CTLA4 CRISPR/Cas9 Lentivirus (Integrating)	78054	Trop2 Lentivirus (Macaca fascicularis/Cynomolgus)	78776
CTLA4 CRISPR/Cas9 Lentivirus (Non-Integrating)	78061	ULBP2 Lentivirus	78744
DLL3 Lentivirus	78909	ULBP2 Lentivirus (Macaca fascicularis/Cynomolgus)	78777
Dominant Negative TGF-β Receptor Type II (TGF-βRII) Lentivirus	78928	VSIG4 Lentivirus	78902
EGR1 Promoter Luciferase Reporter Lentivirus	78664	Primary Cells	Catalog #
FCGR2A CRISPR/Cas9 Lentivirus (Integrating)	78537	Anti-Mesothelin CAR-T Cells	78729
FCGR2A CRISPR/Cas9 Lentivirus (Non-Integrating)	78538	Anti-BCMA CAR-T Cells	78660
FcRL5 Lentivirus (Macaca fascicularis/Cynomolgus)	78781	Anti-CD19 CAR-T Cells	78171
FOLR1 Lentivirus (Macaca fascicularis/Cynomolgus)	78778	Anti-CD20 CAR-T Cells	78611
GPRC5D Lentivirus	78716	CD4+ T cells, Negatively Selected (Human)	79752
GPRC5D Lentivirus (Macaca fascicularis/Cynomolgus)	78780	CD8+ T cells, Negatively Selected (Human)	79753
HLA-C*08:02 Lentivirus	78930	Expanded Human Peripheral Blood NK Cells, Frozen	78798
HLA-E Lentivirus	78929	Normal Human Peripheral Blood Mononuclear Cells, Frozen	79059
HSE Luciferase Reporter Lentivirus (Heat Shock Response)	78669	Proteins	Catalog #
IL15/IL15Ra Lentivirus	78938	ADAR1, FLAG-Tag	100472
KRAS G12D-Specific TCR Lentivirus (Clone 10)	78937	ADAR2 (ADARB1), FLAG-Tag	101164
KRAS G12D-Specific TCR Lentivirus (Clone 9c)	78936	APRIL, His-Avi-Tag	100254
LAG3 CRISPR/Cas9 Lentivirus (Integrating)	78053	APRIL, His-Avi-Tag, Biotin-Labeled	100262
LAG3 CRISPR/Cas9 Lentivirus (Non-Integrating)	78060	ARG1, His-Tag	71658
LAIR1 Lentivirus	78903		
MAGE-A1-Specific TCR Lentivirus (Clone 1367)	78934		

Proteins	Catalog #	Proteins	Catalog #
ARG2, His-Tag	71659	CBL-B (Y363F), His-tag, Biotin-labeled (Human)	80413
B7-1 (CD80) (Mouse), Fc fusion (Human), Avi-Tag	79058	CBL-B, His-Avi-Tag, Biotin-labeled (Human)	80412
B7-1 (CD80), Fc fusion (Human) HiP™	71125	CD111, Fc-Fusion, Avi-Tag	77001
B7-1 (CD80), Fc fusion, Biotin-labeled (Human) HiP™	71114	CD111, Fc-Fusion, Avi-Tag, Biotin-Labeled HiP™	77002
B7-1 (CD80), FLAG-Avi-His-Tag (Mouse)	79002	CD112, Fc Fusion, Avi-Tag	11079
B7-1 (CD80), FLAG-Avi-His-Tag	71261	CD112, Fc-Fusion, Avi-Tag, Biotin-labeled (Human)	72231
B7-1 (Woodchuck), Fc-fusion (Human), Avi-Tag, Biotin-labeled	100040	CD112, His-tag (Human) HiP™	71197
B7-1(CD80) (Mouse), Fc fusion, Avi-Tag, Biotin-labeled, HiP™	100478	CD112, His-Tag, Biotin-Labeled (Human) HiP™	71234
B7-2 (CD86), Fc fusion (Human)	71150	CD112R (PVRIG), Fc-Fusion, Avi-Tag (Human) HiP™	79116
B7-2 (CD86), Fc fusion, Biotin labeled (Human) HiP™	71159	CD112R (PVRIG), Fc-Fusion, Avi-Tag, Biotin-Labeled (Human) HiP™	79270
B7-2 (CD86), FLAG-Avi-His-Tag, Biotin Labeled	71263	CD113, Fc-Fusion (IgG1), Avi-Tag	72510
B7-H2 (CD275, ICOSL), Fc fusion (Human) HiP™	71130	CD123, Avi-His-Tag	101035
B7-H2, Avi-His-Tag	79119	CD123, Fc-Fusion (IgG1) Avi-Tag	100058
B7-H2, Avi-His-Tag, Biotin-Labeled	79300	CD123, Fc-Fusion (IgG1) Avi-Tag, Biotin-Labeled	100068
B7-H3, Avi-His-Tag HiP™	79337	CD125, Fc Fusion (IgG1), Avi-Tag	101050
B7-H3, Avi-His-Tag, Biotin-Labeled HiP™	79367	CD125, Fc-Fusion (IgG1), Avi-Tag, Biotin-Labeled	101051
B7-H4, His-Tag (Human)	71144	CD137 (4-1BB), Fc fusion (Human) HiP™	71170
B7-H4, His-Tag, Biotin-Labeled (Human) HiP™	71129	CD137 (4-1BB), Fc Fusion, Avi-tag Biotin-Labeled HiP™	71171
B7-H5 (VISTA), Fc fusion (Human) HiP™	71148	CD137, Fc fusion (hlgG1) (Mouse)	71250
B7-H5 Fc Fusion, Avi-Tag, Biotin-Labeled (Human)	71327	CD137, Fc fusion (mlgG2a), Avi-tag (Mouse) HiP™	71254
B7-H5, Avi-His-Tag	100358	CD137, Fc fusion (mlgG2a), Biotin-labeled (Mouse) HiP™	71255
B7-H5, Avi-His-Tag, Biotin Labeled, HiP™	100359	CD137L (4-1BB Ligand), His-Avi-Tag	100238
B7-H6, Fc-Fusion-Avi-Tag (Human)	79383	CD137L, His-Tag (E. Coli-Derived)	100562
B7-H6, Fc-Fusion-Avi-Tag, Biotin-Labeled (Human)	79384	CD137L, His-Tag (Mouse)	72517
B7-H7, Fc-Fusion, Avi-Tag	79365	CD14, Avi-His-Tag HiP™	101187
B7-H7, Fc-Fusion, Avi-Tag, Biotin-Labeled	79366	CD14, Avi-His-Tag, Biotin-Label	101188
BAFF-R, Fc-fusion (IgG1), Avi-Tag	100286	CD150, Fc Fusion (IgG1), Biotin Labeled (Human)	72307
BAFF-R, Fc-fusion (IgG1), Avi-Tag, Biotin-labeled	100287	CD150, Fc Fusion, Avi-Tag (Human)	72304
BCMA, Fc-fusion (IgG1), Avi-Tag, Biotin-Labeled	79467	CD155 (PVR), His-tag (Human)	71181
BCMA, Fc-Fusion, Avi-Tag	79465	CD155 (PVR), His-tag (Mouse) HiP™	71167
BCMA, Fc-Fusion, Avi-Tag, PE-Labeled	100733	CD155, Fc-Avi-Tag	79063
BTLA(CD272), Fc fusion (Human)	71141	CD155, His-tag, Biotin-labeled (Mouse) HiP™	71168
BTLA, Fc Fusion, Avi-Tag HiP™	100033	CD160, His-tag (Human) HiP™	71173
BTLA, Fc Fusion, Avi-Tag, Biotin-Labeled HiP™	100047	CD160, His-tag, Biotin-labeled (Human) HiP™	71135
BTN1A1 Avi-His-Tag	100469	CD161, Fc Fusion (IgG1), Avi-Tag HiP™	101211
BTN2A1, Fc Fusion, Avi-Tag	102083	CD19, Avi-His-Tag	101015
BTNL8, Fc-fusion (IgG1), Avi-Tag	100467	CD19, Fc-Fusion (IgG1), Avi-Tag	79472
BTNL8, Fc-fusion (IgG1), Avi-Tag, Biotin-labeled	100468	CD19, Fc-Fusion (IgG1), Avi-Tag, Biotin-Labeled	79475
c-Mer, GST-tag (Human)	40254	CD19, Fc-Fusion (IgG1), Avi-Tag, PE-labeled	100732

Proteins	Catalog #	Proteins	Catalog #
CD2, Avi-His-Tag HiP™	101238	CD36, Fc Fusion, Avi-Tag	101813
CD2, Avi-His-Tag, Biotin-Labeled HiP™	101239	CD37, Avi-His-Tag	101330
CD2, Fc fusion, Avi-Tag, (Human) HiP™	71264	CD37, Avi-His-Tag, Biotin-Labeled	101331
CD2, Fc fusion, Biotin-labeled (Human) HiP™	71270	CD38, Avi-His-Tag	100346
CD20, FLAG-Tag	101572	CD38, Avi-His-Tag, Biotin-Labeled HiP™	100352
CD200, Avi-Tag, His-Tag	102030	CD38, FLAG-Tag (Pig), HiP™	101019
CD200, Avi-Tag, His-Tag, Biotin-Labeled	102031	CD38, His-Tag (Dog)	100955
CD200R1, Fc Fusion, Avi-Tag	102032	CD38, His-Tag (Human), HiP™	71277
CD200R1, Fc Fusion, Avi-Tag, Biotin-Labeled	102033	CD38, His-Tag (Mouse), HiP™	79070
CD22, Fc Fusion, Avi-Tag, PE-labeled	101028	CD38, His-Tag, PE-labeled	71882
CD22, Fc-fusion, Avi-Tag HiP™	79464	CD38-APC, His-Tag	71883
CD22, Fc-fusion, Avi-Tag, Biotin-labeled HiP™	79466	CD39, His-Tag (Human)	71284
CD226, Fc fusion (Human) HiP™	71252	CD4, Avi-Tag, His-Tag, HiP™	101862
CD226, Fc fusion, Biotin-labeled (Human) HiP™	71253	CD40, Fc fusion (Human)	71174
CD235a, Fc Fusion (IgG1), Avi-Tag HiP™	101196	CD40, Fc Fusion (IgG1) Avi-Tag, Biotin Labeled	79102
CD235a, Fc Fusion (IgG1), Avi-Tag, Biotin-Labeled HiP™	101197	CD40L (CD154), His-tag (Human)	71191
CD24, Fc Fusion, Avi-tag	100527	CD44, Fc fusion (Human)	71187
CD24, Fc Fusion, Avi-tag, Biotin-labeled, HiP™	100528	CD44, Fc fusion, Biotin-labeled (Human) HiP™	71133
CD244, Fc Fusion (IgG1), Avi-Tag (Human)	72305	CD45(PTPRC), His-tag	30044
CD244, Fc Fusion (IgG1), Biotin-Labeled (Human)	72306	CD46-BC, His-tag (Human)	71230
CD27(Mouse), Fc-Fusion (Human), Avi-Tag, Biotin	79107	CD46-BC, His-Tag (Monkey) HiP™	71232
CD27, Fc fusion (Human)	71176	CD46-C, His-tag (Human)	71231
CD27, Fc Fusion (IgG1) Avi-Tag, Biotin Labeled	100250	CD46-C, His-tag (Monkey)	71233
CD27, Fc-Fusion (IgG1), Avi-Tag (Mouse) HiP™	79067	CD47 (Monkey), Fc Fusion (Human), Avi-Tag HiP™	79118
CD277, Fc-Fusion (IgG1) Avi-Tag	100073	CD47 (Monkey), Fc Fusion (Human), Avi-Tag, Biotin HiP™	79302
CD277, Fc-Fusion (IgG1) Avi-Tag, Biotin-Labeled	100074	CD47, Fc Fusion (IgG1)	71177
CD28, Fc fusion (Human)	71113	CD47, Fc fusion, Avi-Tag (Human) HiP™	79051
CD28, Fc Fusion, Avi-Tag, Biotin Labeled	101580	CD47, Fc Fusion, Avi-Tag, Biotin-Labeled (Mouse)	72514
CD28H, Fc Fusion (IgG1), Avi-Tag	11083	CD47, Fc fusion, Biotin-labeled (Human) HiP™	71169
CD28H, Fc fusion, Biotin-labeled (Human) HiP™	71239	CD47, Fc-Fusion, Streptavidin-Labeled	71292
CD28H, His-Tag	79375	CD47, His-Tag (Human)	71127
CD30, Fc-Fusion (IgG1), Avi-Tag, Biotin-Labeled	72511	CD48, Fc fusion (IgG1), Avi-Tag™ HiP™ (Human)	72309
CD33, Fc Fusion, Avi-tag	100155	CD48, Fc-Avi, Biotin-Labeled	72308
CD33, Fc Fusion, Avi-tag, Biotin Labeled, HiP™	100482	CD5, Fc Fusion, Avi-Tag, Biotin-labeled, HiP™	101006
CD33, Fc Fusion, Avi-Tag, PE-labeled	101029	CD5, Fc-Fusion, Avi-Tag, HiP™	101005
CD34, Avi-His-Tag HiP™	101190	CD55, Avi-His-Tag, Biotin-labeled	100943
CD34, Avi-His-Tag, Biotin-Labeled	101191	CD55, Avi-His-Tag, HiP™	100942
CD34, Fc Fusion, Avi-Tag	101812	CD58, Fc fusion, Biotin-labeled HiP™	71269
CD36, Avi-His-Tag	101184	CD5L, Avi-His-Tag	101007

Proteins	Catalog #	Proteins	Catalog #
CD66b, Fc Fusion, Avi-Tag	102029	CRTAM, Fc-Fusion (IgG1), Avi-Tag, Biotin-Labeled	40687
CD70 His-Avi-Tag, Biotin-Labeled	100228	CSF1R, Fc-Fusion, Avi-Tag (Human) HiP™	72550
CD70, His-Avi-Tag	100240	CSF1R, Fc-Fusion, Avi-Tag, Biotin-Labeled	79343
CD70, His-tag (Human)	71178	CTLA4 (CD152), Fc fusion, Biotin-labeled (Human)	71152
CD70, His-Tag (Mouse)	79066	CTLA4 (CD152), Fc-fusion (IgG1)	71149
CD73, Avi-His-Tag (Mouse)	72523	CTLA4 (Mouse), Fc-Fusion (Human), Avi-Tag	79062
CD73, Avi-His-Tag, Biotin Labeled	79127	CTLA4 (Mouse), Fc-Fusion (Human), Avi-Tag, Biotin Labeled	79001
CD73, Avi-His-Tag, Biotin-Labeled (Mouse)	79416	CTLA4 (Woodchuck), Fc-Fusion (Human), Avi-tag HiP™	100037
CD73, His-Tag (Human)	71184	DDX58, FLAG-Tag	100070
CD83, Avi-His-Tag HiP™	100224	Diablo, His-Tag	101118
CD83, Avi-His-Tag, Biotin-Labeled HiP™	100225	DPP4, His-tag	80040
CD83, Fc-fusion (IgG1), Avi-Tag	100347	DRAK2	71225
CD83, Fc-fusion (IgG1), Avi-Tag, Biotin labeled	100348	DRAK2, GST-tag (Mouse)	100444
CD8a, Avi-Tag, His-Tag	101872	DRAK2, GST-Tag	71224
CD8a, Avi-Tag, His-Tag, Biotin-Labeled	101873	EGFR (del746-750, T790M, C797S), GST-Tag, His-Tag	100187
CD8a, Fc Fusion, Avi-Tag, Biotin-Labeled, HiP™	101871	ELANE (Elastase), Avi-His-Tag HiP™	101141
CD8a, Fc Fusion, Avi-Tag, HiP™	101870	ELANE (Elastase), Avi-His-Tag, Biotin-Labeled HiP™	101142
CD8B, Avi-His-Tag	101843	EpCAM, Avi-His-Tag	100461
CD8b, Avi-Tag, His-Tag, Biotin-Labeled	101844	EpCAM, Avi-His-Tag, Biotin Labeled	100462
CD8b, Fc Fusion, Avi-Tag	101841	EPOR, Fc fusion, Avi-Tag, Biotin-labeled HiP™	100612
CD8B, Fc Fusion, Avi-Tag, Biotin-Labeled	101842	ERAP1, Avi-His-Tag, Biotin-labeled	100577
CD93, Fc Fusion (IgG1), Avi-Tag	101543	FCGR2A, Avi-His-Tag	100094
CD93, Fc Fusion (IgG1), Avi-Tag, Biotin-Labeled	101544	FCGR2B, Avi-His-Tag	100089
CD96, Fc-Fusion, Avi-Tag (Human)	71265	FCGR2B, Avi-His-Tag, Biotin-Labeled	100474
CD96, Fc-Fusion, Avi-Tag, Biotin Labeled (Human)	71280	FCGR3A (F176V), Avi-His-Tag	100079
CEACAM, Avi-His-Tag, Biotin Labeled	70201	FCGR3A, Avi-His-Tag	79013
CEACAM1, Avi-His-Tag	70206	FCGR3B, Avi-His-Tag	79016
CEACAM5, Avi-His-Tag	100509	FcRn (FCGRT/B2M) Blocker	101468
CERK, FLAG-Tag	102043	FcRn (FCGRT/B2M), His-Avi-Tag, Biotin-Labeled (Mouse)	71286
Claudin-18 Isoform 2, FLAG-Tag	101570	FcRn (FCGRT/B2M), His-Tag (Mouse) HiP™	11349
Claudin-6, FLAG-Tag	101599	FcRn (FCGRT/B2M), His-Tag	71285
CLEC1A, His-Avi-Tag, Biotin-Labeled	101011	FcRn Complex (FCGRT/B2M), His-Avi-Tag, Biotin-Labeled, HiP™	71283
CLEC1A, His-Tag	101010	FGF19, Avi-Tag, His-Tag	101886
CLEC1B, His-Avi-Tag	101021	FGF19, Avi-Tag, His-Tag, Biotin-Labeled	101888
CLEC1B, His-Avi-Tag, Biotin-Labeled	101022	FGFR1a (IIIc), Fc Fusion, Avi-Tag	101836
CLEC2D, Fc Fusion (IgG1), Avi-Tag	101186	FGFR2C, Fc Fusion, Avi-Tag	101877
CLEC2D, Fc Fusion (IgG1), Avi-Tag, Biotin-Labeled HiP™	101189	FGFR3, Avi-Tag, His-Tag, Biotin-Labeled, HiP™	101818
CLEC2D, His-Avi-Tag, Biotin-Labeled	101470	FGFR3, Fc Fusion, Avi-Tag	101819
CRTAM, Fc-Fusion (IgG1), Avi-Tag	40686	FGFR3, Fc Fusion, Avi-Tag, Biotin-Labeled	101820

Proteins	Catalog #	Proteins	Catalog #
GITR (CD357), Fc fusion (Human) HiP™	71172	IL-2RG, Avi-His-Tag, Biotin-Labeled	101150
GITR (CD357), Fc Fusion, Biotin-labeled (Human)	71256	IL12RB1, Avi-His-Tag	101336
GITRL, His-tag (Human)	71190	IL18R1, Fc Fusion (IgG1), Avi-Tag HiP™	101283
GPC3, His-tag [Monkey (M. fascicularis)]	71260	IL18RAP, Fc Fusion (IgG1), Avi-Tag HiP™	101367
HMGB1, Avi-His-Tag	79082	IL23R, Fc Fusion (IgG1), Avi-Tag, Biotin-Labeled HiP™	101322
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

Proteins	Catalog #	Proteins	Catalog #
LILRB2, Avi-His-Tag, Biotin-labeled	100335	OX40L, His-Avi-Tag	100808
LILRB4, Avi-His-Tag, Biotin-Labeled HiP™	100336	p21, GST-Tag	101584
LILRB4, Avi-His-Tag, HiP™	100236	PD-1 (CD279), Fc fusion (Human) HiP™	71106
Lin28A, FLAG-Avi-Tag	79254	PD-1 (CD279), Fc fusion (Mouse) HiP™	71116
Lin28A, FLAG-Avi-Tag, Biotin-Labeled	79072	PD-1 (CD279), Fc fusion, Biotin-labeled (Human) HiP™	71109
Lin28B, FLAG-Avi-Tag	79255	PD-1 (CD279), Fc fusion, Biotin-labeled (Mouse) HiP™	71118
Lin28B, FLAG-Avi-Tag, Biotin-Labeled	79073	PD-1 (E136A), Fc fusion, Biotin-labeled (Human) HiP™	71247
MAT2A, His-Avi-Tag, Biotin-Labeled	79308	PD-1 (Human), Fc Fusion (Mouse IgG2a) HiP™	71997
MAT2A, His-Tag	71401	PD-1 (I126A), Fc fusion, Biotin-labeled (Human) HiP™	71238
Mesothelin, Avi-His-Tag, Biotin-Labeled, HiP™	100291	PD-1 (I134A), Fc fusion, Biotin-labeled (Human) HiP™	71245
Mesothelin, Avi-His-Tag, HiP™	100290	PD-1 (K78A), Fc fusion, Biotin-labeled (Human) HiP™	71244
MICA, Fc fusion, Avi-Tag	100295	PD-1 (L128A), Fc fusion, Biotin-labeled (Human) HiP™	71248
MICA, Fc fusion, Avi-Tag, Biotin-labeled	100300	PD-1 (Monkey), Fc fusion (Human IgG1) HiP™	71158
MICB, Avi-His-Tag HiP™	100505	PD-1 (Monkey), Fc fusion (Human IgG1), Biotin-labeled HiP™	71235
MMP8, His-tag	100552	PD-1 (Mouse), Fc fusion (Human), PE-labeled	71267
Mouse LTBR, Fc Fusion HiP™	71122	PD-1 (P130A), Fc fusion, Biotin-labeled (Human) HiP™	71246
mTOR/Raptor/MLST8 Complex, FLAG-Tag, His-Tag (Human)	40300	PD-1, Avi-His-Tag (Mouse)	79288
NCAM1, Avi-His-Tag, Biotin-Labeled	101044	PD-1, Avi-His-Tag, Biotin-Labeled (Mouse)	79189
NCAM1, Avi-His-Tag, HiP™	101043	PD-1, Fc Fusion (IgG1), Eu-labeled	79011
Nectin4, His-Avi-Tag HiP™	100674	PD-1, Fc Fusion (Woodchuck)	79314
Nectin4, His-Avi-Tag, Biotin-labeled	100675	PD-1, Fc fusion, Avi-Tag	100406
NKG2A, His-Avi-Tag	100638	PD-1, Fc Fusion, PE-Labeled (Human)	71243
NKG2D, Avi-Tag, Fc fusion	100252	PD-1, FLAG-Avi-His-tag (Human) HiP™	71198
NKG2D, Avi-Tag, Fc fusion, Biotin-labeled	100313	PD-1, FLAG-Avi-His-Tag, Biotin-Labeled (Human) HiP™	71325
Nkp30, Fc Fusion, Avi-tag	100546	PD-L1 (A121Q), Fc fusion, Biotin-labeled (Human) HiP™	71320
NKp46, Fc-fusion (IgG1), Avi-Tag	100465	PD-L1 (CD274)	100443
NKp46, Fc-fusion (IgG1), Avi-Tag, Biotin-Labeled	100466	PD-L1 (CD274), Fc fusion (Human) HiP™	71104
NLRP3 (delPYD) / NEK7 Complex	100477	PD-L1 (CD274), Fc fusion (Mouse) HiP™	71117
NLRP3 (delPYD), FLAG-Tag	100454	PD-L1 (CD274), Fc fusion, Biotin-labeled (Human) HiP™	71105
NLRP3 (K232A), His-FLAG-Tag (Sf9-derived)	40741	PD-L1 (CD274), Fc fusion, Biotin-labeled (Mouse) HiP™	71119
NLRP3 (NALP3) (GKT(231-233)AAA), His-Flag-Tags (HEK293-derived)	100190	PD-L1 (CD274), FLAG-tag (Human) HiP™	71183
NLRP3 (NALP3), His-Flag-Tags	100189	PD-L1 (D122A), Fc fusion, Biotin-labeled (Human) HiP™	71321
NPM1, His-Tag	102057	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 (I154A), Fc fusion, Biotin-labeled (Human) HiP™	71314
OX40L (CD252), His-tag (Human)	71185	PD-L1 (K124A), Fc fusion, Avi-Tag (Human) HiP™	71993
OX40L (CD252), His-Tag (Mouse)	72516	PD-L1 (K124A), Fc fusion, Biotin-labeled (Human) HiP™	71323

Proteins	Catalog #	Proteins	Catalog #
PD-L1 (M115A), Fc fusion, Biotin-labeled (Human) HiP™	71318	Siglec7, Fc Fusion, Avi-tag, Biotin-labeled	100636
PD-L1 (Monkey), Fc fusion (Human IgG1) HiP™	71153	Siglec8, Fc Fusion, Avi-Tag	100421
PD-L1 (Monkey), Fc fusion (Human IgG1), Biotin-labeled HiP™	71154	Siglec9, Fc Fusion, Avi-tag	100558
PD-L1 (Q66A), Fc fusion, Biotin-labeled (Human) HiP™	71316	Siglec9, Fc Fusion, Avi-tag, Biotin-labeled	100559
PD-L1 (R113A), Fc fusion, Biotin-labeled (Human) HiP™	71317	SIRP-α (CD172a), Avi-His-Tag (Monkey) HiP™	79117
PD-L1 (R125A), Fc fusion, Biotin-labeled (Human) HiP™	71324	SIRP-α (CD172a), Avi-His-Tag, Biotin (Monkey) HiP™	79303
PD-L1 (S117Q), Fc fusion, Biotin-labeled (Human) HiP™	71319	SIRP-α (CD172a), His-tag (Human)	71145
PD-L1 (T20Q), Fc fusion, Biotin-labeled (Human) HiP™	71312	SIRP-α (SIRP alpha), His-tag, Biotin-labeled (Human) HiP™	71138
PD-L1 (Y123A), Fc fusion, Avi-Tag (Human)	71992	SIRP-γ (CD172g), Fc fusion, Biotin-labeled (Human) HiP™	71236
PD-L1 (Y123A), Fc fusion, Biotin-labeled (Human) HiP™	71322	SLAMF7, Avi-His-Tag HiP™	100271
PD-L1 (Y56A), Fc fusion, Biotin-labeled (Human) HiP™	71315	SLAMF7, Avi-His-Tag, Biotin-Labeled	100272
PD-L1(CD274), FLAG-Avi-His-Tag, (Human) HiP™	71279	TACI, Fc-Fusion (IgG1), Avi-Tag	100283
PD-L1, Fc Fusion (Human IgG1) PE-Labeled	71881	TDO, His-Tag (Human)	71195
PD-L1, Fc Fusion, PE-Labeled (Mouse)	71268	TDO2, His-tag (Mouse)	71241
PD-L2 (CD273), Fc fusion (Human) HiP™	71107	TFRC (Transferrin Receptor), Avi-Tag, Fc Fusion, Biotinylated	101315
PD-L2 (CD273), Fc fusion, Biotin-labeled (Human)	71108	TFRC, Avi-Tag, Fc Fusion (IgG1)	101287
PDPN, Fc-Fusion, Avi-Tag HiP™	79341	TGFβR2, Fc Fusion, Avi-tag, Biotin-labeled HiP™	100629
PDPN, Fc-Fusion, Avi-Tag, Biotin-Labeled HiP™	79342	TIGIT, Fc Fusion (Human)	71186
PSGL1 (CD162), Fc Fusion, Avi-tag	100631	TIGIT, Fc Fusion (IgG1) Avi-Tag, Biotin (Mouse) HiP™	79269
PSGL1 (CD162), Fc Fusion, Avi-tag, Biotin-labeled	100632	TIGIT, Fc Fusion, Avi-Tag, Biotin-Labeled (Human) HiP™	71251
PSMA, His-Avi-Tag	100463	TIM-1, Avi-His-Tag	72512
PSMA, His-Avi-Tag, Biotin-Labeled	100464	TIM-3 (Mouse) Fc-Fusion (Human)-Avi-Tag HiP™	72549
RANK, Fc fusion (IgG1), Avi-tag (Human)	70823	TIM-3 (Mouse), Fc-Fusion (Human)-Avi-Tag, Biotin-Labeled HiP™	77000
RANK, Fc fusion (IgG1), Biotin Labeled (Human)	70822	TIM-3, Avi-His-Tag	71262
RANKL, His-Tag (Human)	71051	TIM-3, Avi-Tag, Biotin Labeled (Human)	79096
ROR1, Fc-Fusion (IgG1)-Avi-Tag, PE-labeled	100995	TIM-3, Biotin-labeled, His-tag (Human) HiP™	71237
Sestrin 1, FLAG-tag	50284	TIM-3, Fc fusion (Human)	71151
Sestrin 2, FLAG-Tag	50286	TIM-3, FLAG-tag (Human)	71180
Siglec10, Avi-His-tag	100529	TIM-4, Avi-His-Tag	72513
Siglec11, Fc Fusion, Avi-Tag	100662	TNFR2, Fc-Fusion (IgG1), His-Avi-Tag	79363
Siglec11, Fc-Fusion (IgG1), Avi-Tag, Biotin-Labeled	100663	TNFR2, Fc-Fusion (IgG1), His-Avi-Tag, Biotin-labeled	100205
Siglec15, Fc Fusion, Avi-tag	100423	TPBG, Avi-His-Tag	101514
Siglec15, Fc Fusion, Avi-tag, Biotin-labeled	100424	TPBG, Avi-His-Tag, Biotin Labeled,	101515
Siglec4, Fc Fusion, Avi-tag, Biotin-labeled	100532	TPBG, Fc Fusion (IgG), Avi-Tag	101516
Siglec5, Fc Fusion, Avi-tag	100533	TPBG, Fc Fusion (IgG1), Avi-Tag, Biotin Labeled	101517
Siglec5, Fc Fusion, Avi-tag, Biotin-labeled	100534	TPH1, His-tag (Human)	71192
Siglec6, Fc Fusion, Avi-tag	100535	TPH2, His-tag (Human)	71193
Siglec6, Fc Fusion, Avi-tag, Biotin-labeled	100536	TREM2, Fc-fusion (IgG1), Avi-Tag (Mouse)	79458
Siglec7, Fc Fusion, Avi-tag	100589	TREM2, Fc-fusion (IgG1), Avi-Tag (Mouse), Biotin	79459

Proteins	Catalog #
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
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

