



LEGENDScreen[™]

Lyophilized Antibody Array

Mouse PE Kit

Catalog Number:	700005
Size:	1 kit, ready to use 1 test per antibody/well
Reactivity:	Mouse
Antibody Format:	PE-conjugated, lyophilized
Configuration:	266 pre-titrated antibodies including 255 PE-conjugated antibodies and 11 isotype controls, arrayed on three 96-well plates, with one specificity per well

It is highly recommended that this manual be read in its entirety before using this product.

Do not use this kit beyond the expiration date.

For research use only

BioLegend, Inc
biolegend.com

LEGENDScreen™ Mouse PE Kit

For Research Purposes Only. Not for use in diagnostic or therapeutic procedures. Purchase does not include or carry the right to resell or transfer this product either as a stand-alone product or as a component of another product. Any use of this product other than the permitted use without the express written authorization of BioLegend is strictly prohibited.

Table of Contents

Description.....	1
Materials Provided.....	1
Materials to be Provided by the End-User.....	2
Storage and Handling.....	2
Preparation of Cells for Staining.....	2
Plate Preparation.....	3
Cell Staining Procedure.....	4
Tips for Successful Staining.....	5
Data analysis.....	6
Frequently Asked Questions.....	7
Product Performance.....	8
Representative data.....	9
Plate Maps.....	10
Antibody Information Tables.....	13

LEGENDScreen™ Mouse PE Kit

Description

The LEGENDScreen™ Mouse PE Kit contains 255 PE-conjugated monoclonal antibodies to cell surface markers as well as 11 mouse, rat, or hamster Ig isotype controls in a convenient 96 well, ready to use format. The kit can be used for screening cell lines, primary cells such as splenocytes, bone marrow-derived cells and cells isolated from tissues. If other co-stains are needed, they can be added to the wells following reconstitution. Positive “hits” from the screening can be quickly identified based on the plate map and the catalog number and clone information can be obtained using the tables on pages 12 through 17. Analysis with individual fluorochrome-conjugated antibodies should be performed to confirm the screening results. The LEGENDScreen™ Mouse PE Kit provides a convenient, easy to use, and powerful tool for immunology, stem cell, and cancer research.

Materials Provided

Description	Quantity	Size	Part#
LEGENDScreen™ Mouse PE Kit Plate #1	1 Plate	96-well	79489
LEGENDScreen™ Mouse PE Kit Plate #2	1 Plate	96-well	79490
LEGENDScreen™ Mouse PE Kit Plate #3	1 Plate	96-well	79491
Cell Staining Buffer	1 Bottle	500 mL	420201
Fixation Buffer	1 Bottle	100 mL	420801
Plate Sealers	8 Sheets	79.4mm x 141mm	78101

Materials to be Provided by the End User

1. Adjustable multichannel pipettes for measuring volumes ranging from 25 μ L to 1,000 μ L
2. Centrifuge with a rotor and adaptors for 96-well plates
3. Cell culture medium (for cell culture)
4. Cell dissociation buffer (for adherent cells)
5. 1X PBS (Phosphate-Buffered Saline): 8.0 g NaCl, 1.16 g Na_2HPO_4 , 0.2 g KH_2PO_4 , 0.2 g KCl, add deionized water to 1 liter; pH to 7.4, 0.2 μ m filtered
6. Plastic reservoirs for pipetting deionized water or Cell Staining or Fixation Buffer with a multichannel pipette
7. A flow cytometer, preferably compatible with reading 96-well plates

Storage and Handling

1. Store unopened kit components at 4°C. Do not use this kit beyond its expiration date.
2. Once opened, reconstitute plates with distilled, deionized water. The reconstituted plates can be used immediately or sealed for storage at 4°C in the dark for up to one month.
3. Keep the buffer/s at 4°C and use within one month after opening.

Preparation of Cells for Staining

1. Obtain desired tissue (*e.g. spleen, lymph node, thymus, bone marrow*) and prepare a single cell suspension. Wash cells in 1X PBS or cell culture medium of choice, and resuspend in Cell Staining Buffer at a density between 4×10^6 and 1×10^7 cells/mL.
2. For cultured cells in suspension, spin and resuspend cells in Cell Staining Buffer at a density between 4×10^6 and 1×10^7 cells/mL.
3. For cultured adherent cells, dissociate cells using a mild enzyme or non-enzymatic dissociation buffer. Wash cells in 1X PBS or cell culture medium of choice, and resuspend in Cell Staining Buffer at a density between 4×10^6 and 1×10^7 cells/mL.
4. Filter the cells through a 40 μ m cell strainer to remove any clumps. Keep the cells on ice before use. Approximately 22.5 mL of cells at a density between 4×10^6 and 1×10^7 cells/mL are needed for all

LEGENDScreen™ Mouse PE Kit

three plates. Lower cell density (e.g. 1.5×10^6 cells/mL) might be used depending on the application.

5. Optional: Reagents that block Fc receptors (*TruStain fcX™ (anti-mouse CD16/32) Antibody, Cat. No. 101320, BioLegend*) may be useful for reducing nonspecific immunofluorescent staining.

Plate Preparation

One hour before the staining, perform the following steps to prepare the plates.

1. Remove the lyophilized plates from the aluminum pouches.
2. Centrifuge the plates at $600 \times g$ for 10 minutes.
3. Make sure the lyophilized cakes settle to the bottom of the plates. Keep the plates upright at all times from this point forward. Handle plates with care so that the cakes are not agitated at any time.
4. Gently settle the plates on the bench away from drafts and air conditioning vents. If possible, discharge any static electricity on yourself prior to removing the cover. **Hold the plate firmly and carefully remove the plate cover from one corner, one well at a time, for easy opening and to prevent cross-well contamination.** You may need to wiggle the cover a little while pulling back slowly.
5. **Discard the cover. Do not move the plate, reseal with the original cover or apply any sealer until the lyophilized antibodies have been reconstituted.**
6. Reconstitute the lyophilized antibodies **immediately**, with 25 μ L/well of deionized water using a multichannel pipette. To avoid cross-contamination, do not let the pipette tips touch the wells. Make sure that all the cakes are dissolved in water. If a cake is stuck on the side of the well, aspirate the 25 μ L water added to the well and rinse the cake down to the bottom.
7. Seal the plates with the plate sealers provided.
8. Wait at least 15 minutes before the staining procedure. Keep the plates in the dark.

LEGENDScreen™ Mouse PE Kit

Note:

- PE-conjugated antibodies are light-sensitive. Try to minimize the exposure of the plates to light as much as practically possible.
- Do not open the pouches until the day you are ready to run the experiment. Once the plates are removed from the pouches, the antibodies must be reconstituted immediately.
- If an experiment is not performed after reconstitution, plates can be sealed and stored in the dark at 4°C for up to one month.

Cell Staining Procedure

1. Using a multichannel pipette, add 75 μL of cells ($\sim 3 - 7.5 \times 10^5$ cells/well) to each well of the plates.
2. Set up extra tubes to stain cells for flow cytometer setup and compensation, if needed.
3. Using the multichannel pipette, gently mix the cell suspensions by pipetting up and down 2 - 3 times. **Be sure to change tips between each row or column.** Avoid creating bubbles while pipetting.
4. Incubate for 20 - 30 minutes at 4°C in the dark.
5. Spin the plates at $500 \times g$ for 10 minutes to pellet cells in each well. Immediately after centrifugation, dump the supernatant into the sink by quickly inverting and flicking the plate. Gently blot the plate on a clean paper towel, being careful not to disturb the cell pellet.
6. Using a multi-channel pipette add 200 μL of Cell Staining Buffer to each well. Gently mix up and down to resuspend cells. **Be sure to change tips between each row.**
7. Repeat step 5.
8. To fix the cells, using a multichannel pipette, aliquot 100 μL of Fixation Buffer into each well. Gently mix up and down to resuspend cells. **Be sure to change tips between each row.**
9. Incubate for 10 minutes at room temperature in the dark.
10. Repeat steps 5 - 6.
11. Repeat step 5 one more time.

LEGENDScreen™ Mouse PE Kit

12. Resuspend cells completely in 160 μ L of Cell Staining Buffer per well and analyze using a flow cytometer. We recommend acquiring 70 μ L of sample and collecting 5-10, 000 events. Users should determine the optimal number of events to be collected based on specific application they are testing. While the first plate is being acquired, store the other plates at 4°C in the dark.

Tips for a Successful Staining

1. Read the entire manual carefully before the experiment.
2. Plan the experiment in advance. Designate a full day for this experiment. Do not rush any step.
3. Make sure that the flow cytometer's autosampler is well maintained and working well before the experiment. If the flow cytometer does not have an autosampler, the samples from each well of the plates should be transferred to individual FACS tubes and tube adjust the sample volume to approximately 300 μ l or more to avoid running dry. Alternatively, a Cluster Tube System (such as Corning catalog # 4410) and acquired manually. The cluster tube can then be transferred to a regular FACS tube for acquisition.
4. Make sure that enough cells have been prepared for the staining. If there are not enough cells, you may choose to divide the staining into two separate experiments.
5. Depending on the application, the cell number needed for staining can be decreased. Successful staining has been done with 1×10^5 cells/well.
6. Make sure to prepare cells for machine setup and compensation. These cells should be treated the same way as the cells for staining in the plates.
7. Handle the plates with care. Keep the plates upright at all times and be careful not to knock the plates over.
8. Protect the plates from exposure to light as much as possible.
9. Use extra caution when opening the plate cover (**see Plate Preparation for instructions**). After the cover is removed, do not move or apply any sealers on the plates until the cakes are reconstituted. Applying plate sealers or any cover onto the plate before the cakes are reconstituted can result in cakes being dislodged and stuck to the cover or escaping from the wells.

LEGENDScreen™ Mouse PE Kit

10. Some cell surface markers are sensitive to enzymatic digestion. If adherent cells are being used for staining, a mild enzyme or non-enzymatic dissociation buffer should be used when possible.
11. Make sure cells are in a single cell suspension. DNase treatment is recommended to avoid clumps caused by dead cells.
12. Acquire only 70 μ L of the 160 μ L total volume so that a second run can be performed if necessary.

Data Analysis

We recommend labeling acquisition files as Plate 1_A1, Plate 1_A2 and so on. After sample acquisition is complete, transfer the files into FlowJo™, or equivalent analysis software. Using the “Create Group” function, group samples based on the isotype. This will help set gates for all samples of the same isotype. Antibodies are arranged by isotypes (isotype control followed by all antibodies of that isotype) – for example, Plate 1, A2 is Armenian hamster IgG isotype control and antibodies A3 through E5 are all of the Armenian hamster IgG isotype. Plate 1, E6 is mouse IgG2a isotype control and antibodies E7 through E12 are of the mouse IgG2a isotype control.

Create 11 groups based on the different isotype control antibodies as described below –

1. Armenian Hamster IgG isotype – Plate 1_A2-E5
2. Mouse IgG2a, κ isotype – Plate 1_E6-E12
3. Mouse IgG1, κ isotype – Plate 1_F1-G1
4. Mouse IgG2b, κ isotype – Plate 1_G2-G5
5. Mouse IgM, κ isotype – Plate 1_G6-G7
6. Rat IgG1, κ isotype – Plate 1_G8-H12 and Plate 2_A2-A11
7. Rat IgG2a, κ isotype –Plate 2_A12-H12 and Plate 3_A2-B8
8. Rat IgG2b, κ isotype –Plate 3_B8-F4
9. Rat IgG2c, κ isotype –Plate 3_F6-F8
10. Rat IgM, κ isotype –Plate 3_F9-F11
11. Syrian Hamster IgG κ isotype –Plate 3_F12-G6

Once the groups are created in FlowJo™, go to the first group and gate around the population(s) of interest and then make a histogram plot for PE. Using the isotype control sample, set a gate on the positive

LEGENDScreen™ Mouse PE Kit

population and then apply it to all files in that group. Repeat this process for the remaining 10 groups. Add statistics for percentage of positive, median fluorescence intensity (MFI), event count and any other statistic of interest. Import into an excel file for further analysis using the table editor function.

If comparing control and test samples, then the histograms can be overlaid for visual comparison. If using software other than FlowJo™ please follow a similar strategy.

Frequently Asked Questions

Q: What is the level of variability from one experiment to the other?

A: If the protocol is followed the variability should be minimal. The variability should be similar to single vial antibody staining.

Q: How should the kit be stored?

A: The kit should be stored at 2 - 8°C upon receipt. Once opened, the plates must be reconstituted immediately. Reconstituted plates can be used or stored at 2 - 8°C sealed in the dark and used within a month.

Q: How do I request a custom LEGENDScreen™ product with only my specificities of interest?

A: For more info, visit: biolegend.com/custom_solutions

Q: What are the guarantees regarding the lyophilized plate compared to the reconstituted plate?

A: Lyophilized product has a guaranteed shelf life of 6 months unopened. Reconstituted plates can be used or stored at 2 - 8°C sealed in the dark and used within a month. Be sure to properly seal the plates to prevent evaporation and shield the antibodies from light.

Q: I have added my own antibody solution to the lyophilized product, will the lyophilized antibody work?

A: Yes, as long as the fluorophores on these antibodies are compatible and proper compensation has been applied during acquisition and analysis.

LEGENDScreen™ Mouse PE Kit

Q: I am not going to use all the reconstituted antibody solution. Can I keep the leftover for later or re-dry the solution in the dark?

A: The antibody is in a one test per well format. There will not be any antibody left if the full test is used. Customers may decide to use less than the recommended volume per test, but this is not recommended and the performance is not guaranteed. Customers may also selectively transfer certain antibodies from the original plate to a new plate and use after reconstitution. If any antibody is not used after reconstitution, the plate can be sealed and store at 2 - 8°C for a month in the dark. Once reconstituted, re-drying is not recommended, as this may result in a loss of signal.

Q: If I don't have enough cells and use less than 4×10^6 /ml (3×10^5 cells/well), will it still work?

A: The kit may work with lower numbers of total cells, but we recommend trying to keep higher concentrations of cells for faster analysis. Of course, how many cells are needed depends on the specific application. Successful staining has been done with 1×10^5 cells/well.

Q: Are these plates made under sterile conditions?

A: The plates are not sterile. Handle them as you would handle a typical flow cytometry staining protocol or reagent.

Q: Can I use half or less of the plate and keep the rest for later?

A: Yes. Customers can use half of the plate or whatever specificities they are interested in. However, the whole plate should be reconstituted. The half plate of antibodies must be transferred to another empty plate for the staining. The remaining half must be sealed and stored at 2 - 8°C in the dark and used within a month.

Product Performance

The LEGENDScreen™ Mouse PE Kit was tested and compared with BioLegend's cataloged single vial liquid antibody reagents. For cell staining, cells were isolated from spleen and lymph nodes, and 3×10^5 cells were added to each well after the lyophilized antibodies were reconstituted. The cells were then stained for 20 minutes at 4°C, washed, and fixed with Fixation Buffer. The cells were then washed,

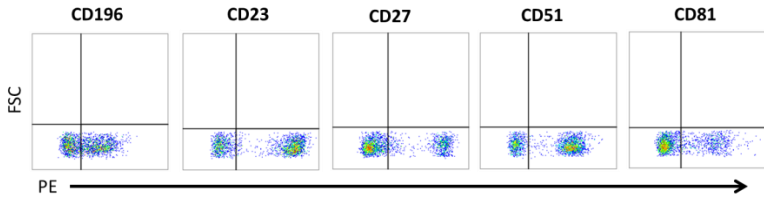
LEGENDScreen™ Mouse PE Kit

resuspended in 160 µL of Cell Staining Buffer, acquired using a BD FACSCanto™ II instrument, and analyzed using FlowJo™.

There was no significant difference in the staining patterns or median fluorescence intensity between the lyophilized product and liquid antibodies. Below are some representative data obtained when comparing the LEGENDScreen™ Mouse PE Kit vs. equivalent liquid antibodies.

Representative data

Liquid Reagents



LegendScreen Reagents

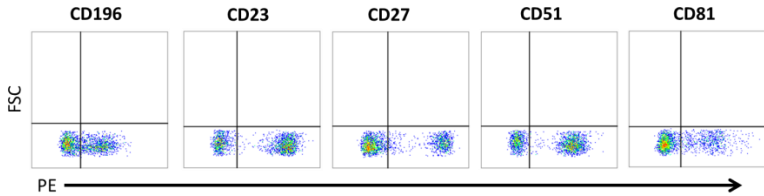


Plate Maps

Plate 1

	1	2	3	4	5	6	7	8	9	10	11	12
A	Blank	Armenian Hamster IgG isotype Ctrl	CD3ε	CD80	CD81	CD154	Notch 1	CD30	CD178	CD103	DLL4	CD195
B	Notch 4	CD229 (Ly-9)	CD69	Notch 3	JAML	Notch 2	CD194	CD152	CD120a	CD11c	DLL1	CD196
C	CD29	CD55	Jagged 2	CD79b	IFN-γ R β chain	CD61	CD121a	TCR β chain	FcεR1α	DcTRAIL -R2	CD36	DcTRAIL -R1
D	CD84	CD48	CD49b	CD120b	CD183	CD262	HVEM	TCR Vγ1.1 + Vγ1.2	B7-H4	CD339	CD49a	PD-1H
E	CD85k	Plexin B2	CD27	DR3	TCR γ/δ	Mouse IgG2a, k Isotype Ctrl	CD45.1	CD45.2	NK-1.1	Ly108	CD207	CX3CR1
F	Mouse IgG1, k Isotype Ctrl	CD66a	IFNAR-1	Tim-2	CD272	CD64	CD351	LAP	TIGIT	Trem-like 4	CD59a	Ly49H
G	CD90.1	Mouse IgG2b, k Isotype Ctrl	CD157	CD159a	XCR1	Mouse IgM, k Isotype Ctrl	SSEA-1	Rat IgG1, k Isotype Ctrl	Ig light chain k	Siglec H	CD255	CD202b
H	GITR Ligand	CD147	CD73	CD51	NKG2D	CD96	Integrin β7	CD210	CD83	Mac-3	CD223	CD134

LEGENDScreen™ Mouse PE Kit

Plate 2

	1	2	3	4	5	6	7	8	9	10	11	12
A	Blank	CD41	CD268	CD144	CD370	CD369	PIR-A/B	CD22	CD324	CD172a	CD319	Rat IgG2a, K Isotype Ctrl
B	MAIR-V	CD146	VISTA	CD8a	CD275	CD34	Ly-6A/E	CD40	CD45R/B220	CD197	CD47	CD98
C	CD14	CD107a	CD18	Ly-6G	CD21, CD35	Mac-2	CD199	Ly-51	IgD	Tim-4	CD71	H-2
D	CD45RB	CD326	IgM	CD155	CD200R	CD254	IL-21R	CD276	CD9	CD105	CD366	4-1BB Ligand
E	CD265	TLR4 (CD284)/MD2 Complex	CD19	LPAM-1	CD62L	CD23	CD5	CD273	CD31	F4/80	CD94	CD267
F	Ly-49A	CD180	CD11a	LTβR	CD122	CD106	CD365	CD115	CD140a	PDC-TREM	CD135	CD127
G	CD140b	ESAM	CD200	CD309	TLT-2	CD253	CD335	CD205	Galectin ₉	CD200R ₃	MAIR-IV	Ly49D
H	CD123	CD355	CD169	CD138	CD160	CD39	GARP	CD179a (VpreB)	CD371	CD63	CD49e	CD193

LEGENDScreen™ Mouse PE Kit

Plate 3

	1	2	3	4	5	6	7	8	9	10	11	12
A	Blank	CD300L G	CD301	IL-33R α	CD304	CD6	CD100	CD104	CD182	Mad CAM-1	MERTK (Mer)	CD226
B	Ly6K	CD16/3 2	CD150	CD25	CD38	CD133	CD301b	CD34	Rat IgG2b,k Isotype Ctrl	CD43	FR4	CD1d
C	CD70	CD4	I-A/I-E	CD153	CD54	33D1	CD90.2	TER-119	CD49d	CD24	Ly- 6G/Ly- 6C	CD86
D	CD11b	CD45	CD279	RAE-1 γ	CD8b	CD44	CD126	CD317	CD132	CD3	CD274	CD117
E	CD88	CD93	CD252	MD-1	CD357	CD185	CD37	CD300c/ d	CD186 (CXCR6)	CD130	CD198	CD20
F	CD124	IL-23R	CD184	CD2	Rat IgG2c,k Isotype Ctrl	Ly-6C	Ly-6D	Rat IgM, k Isotype Ctrl	CD49b	GL7	Syrian Hamster IgG Isotype Ctrl	CD28
G	Podopla nin	CD137	CD278	KLRG1	Ly-49 C/F//H	Blank	Blank	Blank	Blank	Blank	Blank	Blank
H	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank

Antibody information Tables
Plate 1

Well	Cat #	Clone	Description	Isotype
A1		Blank		
A2	400908	HTK888	Armenian Hamster IgG Isotype control	Armenian Hamster IgG
A3	100308	145-2C11	anti-mouse CD3 ϵ	Armenian Hamster IgG
A4	104708	16-10A1	anti-mouse CD80	Armenian Hamster IgG
A5	104906	Eat-2	anti-mouse/rat CD81	Armenian Hamster IgG
A6	106506	MR1	anti-mouse CD154	Armenian Hamster IgG
A7	130608	HMN1-12	anti-mouse Notch 1	Armenian Hamster IgG
A8	102306	mCD30.1	anti-mouse CD30	Armenian Hamster IgG
A9	106606	MFL3	anti-mouse CD178 (FasL)	Armenian Hamster IgG
A10	121405	2E7	anti-mouse CD103	Armenian Hamster IgG
A11	130808	HMD4-1	anti-mouse DLL4	Armenian Hamster IgG
A12	107006	HM-CCR5	anti-mouse CD195 (CCR5)	Armenian Hamster IgG
B1	128407	HMN4-14	anti-mouse Notch 4	Armenian Hamster IgG
B2	122905	Ly9ab3	anti-mouse CD229 (Ly-9)	Armenian Hamster IgG
B3	104508	H1.2F3	anti-mouse CD69	Armenian Hamster IgG
B4	130507	HMN3-133	anti-mouse Notch 3	Armenian Hamster IgG
B5	128503	4E10	anti-mouse JAML	Armenian Hamster IgG

Well	Cat #	Clone	Description	Isotype
E1	144904	H1.1	anti-mouse CD85k (gp49 Receptor)	Armenian Hamster IgG
E2	145904	3E7	anti-mouse Plexin B2	Armenian Hamster IgG
E3	124210	LG.3A10	anti-mouse/rat/human CD27	Armenian Hamster IgG1
E4	144406	4C12	anti-mouse DR3 (TNFRSF25)	Armenian Hamster IgG1
E5	118108	GL3	anti-mouse TCR γ/δ	Hamster IgG
E6	400212	MOPC-173	Mouse IgG2a, k isotype control	Mouse IgG2a, k
E7	110708	A20	anti-mouse CD45.1	Mouse (A.5W) IgG2a, k
E8	109808	104	anti-mouse CD45.2	Mouse (SJL) IgG2a, k
E9	108708	Pk136	anti-mouse NK-1.1	Mouse IgG2a, k
E10	134606	330-AJ	anti-mouse Ly108	Mouse IgG2a, k
E11	144204	4C7	anti-mouse/human CD207	Mouse IgG2a, k
E12	149006	SA011F11	anti-mouse CX3CR1	Mouse IgG2a, k
F1	400112	MOPC-21	Mouse IgG1 Isotype control	Mouse IgG1, k
F2	134506	MAB-CC1	anti-mouse CD66a (CEACAM1a)	Mouse IgG1, k
F3	127312	MAR1-5A3	anti-mouse IFNAR-1	Mouse IgG1, k
F4	129006	F37-2C4	anti-mouse Tim-2	Mouse IgG1, k
F5	134804	8F4	anti-mouse CD272 (BTLA)	Mouse IgG1, k

LEGENDScreen™ Mouse PE Kit

Well	Cat #	Clone	Description	Isotype
B6	130707	HMN2-35	anti-mouse Notch 2	Armenian Hamster IgG
B7	131204	2G12	anti-mouse CD194 (CCR4)	Armenian Hamster IgG
B8	106306	UC10-489	anti-mouse CD152	Armenian Hamster IgG
B9	113004	55R-286	anti-mouse CD120a (TNF R Type I/p55)	Armenian Hamster IgG
B10	117308	N418	anti-mouse CD11c	Armenian Hamster IgG
B11	128307	HMD1-3	anti-mouse DLL1	Armenian Hamster IgG
B12	129804	29-2L17	anti-mouse CD196 (CCR6)	Armenian Hamster IgG
C1	102208	HM81-1	anti-mouse/rat CD29	Armenian Hamster IgG
C2	131804	RiKO-3	anti-mouse CD55 (DAF)	Armenian Hamster IgG
C3	131007	HMI2-1	anti-mouse Jagged 2	Armenian Hamster IgG
C4	132804	HM79-12	anti-mouse CD79b (Igf3)	Armenian Hamster IgG
C5	113604	MOB-47	anti-mouse IFN- γ R β chain	Armenian Hamster IgG
C6	104308	2C9.G2 (HM83-1)	anti-mouse/rat CD61	Armenian Hamster IgG
C7	113505	JAMA-147	anti-mouse CD121a (IL-1 R, Type I/p80)	Armenian Hamster IgG
C8	109208	H57-597	anti-mouse TCR β chain	Armenian Hamster IgG
C9	134308	MAR-1	anti-mouse Fc ϵ R1 α	Armenian Hamster IgG
C10	133703	9E9	anti-mouse CD16.2	Armenian Hamster IgG
C11	102606	HM36	anti-mouse CD36	Armenian Hamster IgG
C12	133804	mDcR1-3	anti-mouse DcTRAIL-R1	Armenian Hamster IgG

Well	Cat #	Clone	Description	Isotype
F6	139304	X54-5/7.1	anti-mouse CD64 (Fc γ R1)	Mouse IgG1, k
F7	137306	TX61	anti-mouse/human CD351 (Fc α/μ α/μ receptor)	Mouse IgG1, k
F8	141306	TW7-20B9	anti-mouse LAP (TGF- β 1)	Mouse IgG1, k
F9	142104	1G9	anti-mouse TIGIT (Vstm3)	Mouse IgG1, k
F10	143304	16E5	anti-mouse Trem-like 4	Mouse IgG1, k
F11	143104	mCD59.3	anti-mouse CD59a	Mouse IgG1, k
F12	144706	3D10	anti-mouse Ly49H	Mouse IgG1, k
G1	202523	OX-7	anti-rat CD90/mouse CD90.1	Mouse IgG1, k
G2	400312	MPC-11	Mouse IgG2b, k isotype control	Mouse IgG2b, k
G3	140204	BP-3	anti-mouse CD157 (BST-1)	Mouse IgG2b, k
G4	142804	16A11	anti-mouse CD159a (NKG2AB6)	Mouse IgG2b, k
G5	148204	ZET	anti-mouse/rat XCR1	Mouse IgG2b, k
G6	401611	MM-30	Mouse IgM isotype control	Mouse IgM, k
G7	125606	MC-480	anti-mouse/human CD15 (SSEA-1)	Mouse IgM, k
G8	400408	RTK2071	Rat IgG1, k isotype control	Rat IgG1, k
G9	409506	RMK-45	anti-mouse IgG light chain k	Rat IgG
G10	129606	551	anti-mouse Siglec H	Rat IgG1, k
G11	120005	MTW-1	anti-mouse CD255 (TWEAK)	Rat IgG1, k
G12	124008	TEK4	anti-mouse CD202b (Tie-2)	Rat IgG1, k

LEGENDScreen™ Mouse PE Kit

Well	Cat #	Clone	Description	Isotype
D1	122806	mCD84.7	anti-mouse CD84	Armenian Hamster IgG
D2	103406	HM48-1	anti-mouse CD48	Armenian Hamster IgG
D3	103506	HM α 2	anti-mouse CD49b	Armenian Hamster IgG
D4	113406	TR75-89	anti-mouse CD120b (TNF R Type II/p75)	Armenian Hamster IgG
D5	126506	CXCR3-173	anti-mouse CD183 (CXCR3)	Armenian Hamster IgG
D6	119906	MD5-1	anti-mouse CD262 (DR5, TRAIL-R2)	Armenian Hamster IgG
D7	136304	HMHV-1818	anti-mouse CD270 (HVEM)	Armenian Hamster IgG
D8	142704	4B2.9	anti-mouse TCR V γ 1.1/V γ 1.2	Armenian Hamster IgG
D9	139406	HMH4-5G1	anti-mouse B7-H4 (B751, B7X)	Armenian Hamster IgG
D10	130908	HMJ1-29	anti-mouse CD339 (Jagged 1)	Armenian Hamster IgG
D11	142604	HM α 1	anti-mouse CD49a	Armenian Hamster IgG
D12	143708	MH5A	anti-mouse PD-1H	Armenian Hamster IgG

Well	Cat #	Clone	Description	Isotype
H1	120306	YGL 386	anti-mouse GITR Ligand	Rat IgG1, k
H2	123707	OX-114	anti-mouse CD147	Rat IgG1, k
H3	127206	TY/11.8	anti-mouse CD73	Rat IgG1, k
H4	104106	RMV-7	anti-mouse CD51	Rat IgG1, k
H5	130208	CX5	anti-mouse CD314 (NKG2D)	Rat IgG1, k
H6	131705	3.3	anti-mouse CD96 (TACTILE)	Rat IgG1, k
H7	121006	FIB27	anti-human/mouse integrin β 7	Rat IgG1, k
H8	112706	1B1.3a	anti-mouse CD210 (IL-10 R)	Rat IgG1, k
H9	121508	Michel-19	anti-mouse CD83	Rat IgG1, k
H10	108506	M3/84	anti-mouse CD107b (Mac-3)	Rat IgG1, k
H11	125208	C9B7W	anti-mouse CD223 (LAG-3)	Rat IgG1, k
H12	119410	OX-86	anti-mouse CD134 (OX-40)	Rat IgG1, k

LEGENDScreen™ Mouse PE Kit

Plate 2

Well	Cat #	Clone	Description	Isotype
A1			Blank	
A2	133906	MWReg30	anti-mouse CD41	Rat IgG1, k
A3	134104	7H22-E16	anti-mouse CD268 (BAFF-R)	Rat IgG1, k
A4	138010	BV13	anti-mouse CD144 (VE-cadherin)	Rat IgG1, k
A5	143504	7H11	anti-mouse CLECSA	Rat IgG1, k
A6	144304	RH1	anti-mouse CD369 (Dectin-1, CLECTA)	Rat IgG1, k
A7	144104	6C1	anti-mouse PIR-A/B	Rat IgG1, k
A8	126112	OX-97	anti-mouse CD22	Rat IgG1, k
A9	147304	DECMA-1	anti-mouse/human CD324 (E-Cadherin)	Rat IgG1, k
A10	144012	P84	anti-mouse CD172a (SIRP α)	Rat IgG1, k
A11	152006	4G2	anti-mouse CD319	Rat IgG1, k
A12	400508	RTK2758	Rat IgG2a, k isotype control	Rat IgG2a, k
B1	132704	TX70	anti-mouse MAIR-V	Rat IgG2a, k
B2	134704	ME-9F1	anti-mouse CD146	Rat IgG2a, k
B3	150204	MIH63	anti-mouse VISTA	Rat IgG2a, k
B4	100708	53-6.7	anti-mouse CD8a	Rat IgG2a, k
B5	107405	HK5.3	anti-mouse CD275 (B7-H2, B7-RP1, ICOS Ligand)	Rat IgG2a, k
B6	119308	MEC14.7	anti-mouse CD34	Rat IgG2a, k
B7	108108	D7	anti-mouse Ly-6A/E (Sca-1)	Rat IgG2a, k
B8	124610	3/23	anti-mouse CD40	Rat IgG2a, k

Well	Cat #	Clone	Description	Isotype
E1	119806	R12-31	anti-mouse CD265 (RANK)	Rat IgG2a, k
E2	117606	MT5510	anti-mouse TLR4 (CD284)/MD2 Complex	Rat IgG2a, k
E3	115508	6D5	anti-mouse CD19	Rat IgG2a, k
E4	120606	DATK32	anti-mouse LPAM-1 (Integrin α 4 β 7)	Rat IgG2a, k
E5	104408	MEL-14	anti-mouse CD62L	Rat IgG2a, k
E6	101608	B3B4	anti-mouse CD23	Rat IgG2a, k
E7	100608	53-7.3	anti-mouse CD5	Rat IgG2a, k
E8	107206	TY25	anti-mouse CD273 (B7-DC, PD-L2)	Rat IgG2a, k
E9	102408	390	anti-mouse CD31	Rat IgG2a, k
E10	123110	BMB8	anti-mouse F4/80	Rat IgG2a, k
E11	105508	18d3	anti-mouse CD94	Rat IgG2a, k
E12	133404	8F10	anti-mouse CD267 (TAC1)	Rat IgG2a, k
F1	116808	YE1/48.1 0.6	anti-mouse Ly-49A	Rat IgG2a, k
F2	117706	RP/14	anti-mouse CD180 (RP105)	Rat IgG2a, k
F3	101107	M17/4	anti-mouse CD11a	Rat IgG2a, k
F4	134403	5G11	anti-mouse lymphotoxin beta receptor (LTBR)	Rat IgG2a, k
F5	105906	5H4	anti-mouse CD122 (IL-2R β)	Rat IgG2a, k
F6	105714	429 (MVCAM.A)	anti-mouse CD106	Rat IgG2a, k
F7	119506	RMT1-4	anti-mouse CD365 (Tim-1)	Rat IgG2a, k
F8	135506	AFS98	anti-mouse CD115 (CSF-1R)	Rat IgG2a, k

LEGENDScreen™ Mouse PE Kit

Well	Cat #	Clone	Description	Isotype
B9	103208	RA3-6B2	anti-mouse/human CD45R/B220	Rat IgG2a, k
B10	120106	4B12	anti-mouse CD197 (CCR7)	Rat IgG2a, k
B11	127508	miap301	anti-mouse CD47	Rat IgG2a, k
B12	128208	RL388	anti-mouse CD98 (4F2)	Rat IgG2a, k
C1	123310	Sa14-2	anti-mouse CD14	Rat IgG2a, k
C2	121612	1D4B	anti-mouse CD107a (LAMP-1)	Rat IgG2a, k
C3	101408	M18/2	anti-mouse CD18	Rat IgG2a, k
C4	127608	1A8	anti-mouse Ly-6G	Rat IgG2a, k
C5	123410	7E9	anti-mouse CD21/CD35 (CR2/CR1)	Rat IgG2a, k
C6	125405	M3/38	anti-mouse/human Mac-2 (Galectin-3)	Rat IgG2a, k
C7	129708	9B1	anti-mouse CD199 (CCR9)	Rat IgG2a, k
C8	108308	6C3	anti-mouse Ly-51	Rat IgG2a, k
C9	405706	11-26c.2a	anti-mouse IgD	Rat IgG2a, k
C10	130006	RMT4-54	anti-mouse Tim-4	Rat IgG2a, k
C11	113808	RI7217	anti-mouse CD71	Rat IgG2a, k
C12	125506	M1/42	anti-mouse H-2	Rat IgG2a, k
D1	103308	C363-16A	anti-mouse CD45RB	Rat IgG2a, k
D2	118206	G8.8	anti-mouse CD326 (EPCAM)	Rat IgG2a, k
D3	406508	RMM-1	anti-mouse IgM	Rat IgG2a, k
D4	131508	TX56	anti-mouse CD155 (PVR)	Rat IgG2a, k
D5	123908	OX-110	anti-mouse CD200 R (OX2R)	Rat IgG2a, k
D6	510006	Ik22/5	anti-mouse CD254 (TRANSC, RANKL)	Rat IgG2a, k
D7	131906	4A9	anti-mouse IL-21 Receptor	Rat IgG2a, k
D8	124508	RTAA15	anti-mouse CD276 (B7-H3)	Rat IgG2a, k

Well	Cat #	Clone	Description	Isotype
F9	135906	APA5	anti-mouse CD140a	Rat IgG2a, k
F10	139204	4A6	anti-mouse PDC-TREM	Rat IgG2a, k
F11	135306	A2F10	anti-mouse CD135	Rat IgG2a, k
F12	135010	A7R34	anti-mouse CD127 (IL-7R α)	Rat IgG2a, k
G1	136006	APB5	anti-mouse CD140b	Rat IgG2a, k
G2	136204	1G8/ESAM	anti-mouse ESAM	Rat IgG2a, k
G3	123808	OX-90	anti-mouse CD200 (OX2)	Rat IgG2a, k
G4	136404	Avas12	anti-mouse CD309 (VEGFR2, Flk-1)	Rat IgG2a, k
G5	136603	MIH47	anti-mouse TLT-2	Rat IgG2a, k
G6	109306	N2B2	anti-mouse CD253 (TRAIL)	Rat IgG2a, k
G7	137604	29A1.4	anti-mouse CD335 (Nkp46)	Rat IgG2a, k
G8	138214	NLDC-145	anti-mouse CD205 (DEC-205)	Rat IgG2a, k
G9	137904	108A2	anti-mouse Galectin-9	Rat IgG2a, k
G10	142206	Ba13	anti-mouse CD200R3	Rat IgG2a, k
G11	139605	TX69	anti-mouse MAIR-IV (CLM-5)	Rat IgG2a, k
G12	138308	4E5	anti-mouse Ly49D	Rat IgG2a, k
H1	106005	5B11	anti-mouse CD123	Rat IgG2a, k
H2	142006	11-5/CRTAM	anti-mouse CD355 (CRTAM)	Rat IgG2a, k
H3	142404	3D6.112	anti-mouse CD169/Siglec-1	Rat IgG2a, k
H4	142504	281-2	anti-mouse CD138	Rat IgG2a, k
H5	143004	7H1	anti-mouse CD160	Rat IgG2a, k
H6	143804	Duha59	anti-mouse CD39	Rat IgG2a, k
H7	142904	F011-5	anti-mouse GARP	Rat IgG2a, k
H8	143604	R3	anti-mouse CD179a/VpreB	Rat IgG2a, k

LEGENDScreen™ Mouse PE Kit

Well	Cat #	Clone	Description	Isotype
D9	124806	MZ3	anti-mouse CD9	Rat IgG2a, k
D10	120408	MJ7/18	anti-mouse CD105	Rat IgG2a, k
D11	119704	RMT3-23	anti-mouse Tim-3	Rat IgG2a, k
D12	107105	Tks-1	anti-mouse 4-1BB Ligand (CD137L)	Rat IgG2a, k

Well	Cat #	Clone	Description	Isotype
H9	143404	5D3/CLEC12A	anti-mouse CLEC12A	Rat IgG2a, k
H10	143904	NVG-2	anti-mouse CD63	Rat IgG2a, k
H11	103805	5H10-27(MFR5)	anti-mouse CD49e	Rat IgG2a, k
H12	144506	J073E5	anti-mouse CCR3 (CD193)	Rat IgG2a, k

LEGENDScreen™ Mouse PE Kit

Plate 3

Well	Cat #	Clone	Description	Isotype
A1			Blank	
A2	147104	ZAQ5	anti-mouse CD300LG (Nepmucin)	Rat IgG2a, k
A3	145606	LOM-8.7	anti-mouse CD301a	Rat IgG2a, k
A4	145304	DIH9	anti-mouse IL-33R α (ST2/T1)	Rat IgG2a, k
A5	145204	3E12	anti-mouse CD304	Rat IgG2a, k
A6	146404	OX-129	anti-mouse CD6	Rat IgG2a, k
A7	147604	BMA-12	anti-mouse CD100/Sema4 D	Rat IgG2a, k
A8	123610	346-11A	anti-mouse CD104	Rat IgG2a, k
A9	149304	SA044G4	anti-mouse CD182 (CXCR2)	Rat IgG2a, k
A10	120710	MECA-367	anti-mouse MadCAM-1	Rat IgG2a, k
A11	151506	2B10C42	anti-mouse MERTK (Mer)	Rat IgG2a, k
A12	133603	TX42.1	anti-mouse CD226 (DNAM-1)	Rat IgG2a, k
B1	151304	Mk34	anti-mouse LY6K	Rat IgG2a, k
B2	101308	93	anti-mouse CD16/32	Rat IgG2a, λ
B3	115904	TC15-12F12.2	anti-mouse CD150 (SLAM)	Rat IgG2a, λ
B4	102008	PC61	anti-mouse CD25	Rat IgG2a, λ
B5	102708	90	anti-mouse CD38	Rat IgG2a, λ
B6	141204	315-2C11	anti-mouse CD133	Rat IgG2a, λ
B7	146804	URA-1	anti-mouse CD301b	Rat IgG2a, λ
B8	152204	SA376A4	anti-mouse CD34	Rat IgG2a, λ
B9	400608	RTK4530	Rat IgG2b, k isotype control	Rat IgG2b, k
B10	143206	S11	anti-mouse CD43	Rat IgG2b, k
B11	125007	12A5	anti-mouse FR4 (Folate Receptor 4)	Rat IgG2b, k

Well	Cat #	Clone	Description	Isotype
E1	135806	20/70	anti-mouse CD88 (CSaR)	Rat IgG2b, k
E2	136504	AA4.1	anti-mouse CD93 (AA4.1, early B lineage)	Rat IgG2b, k
E3	108806	RM134L	anti-mouse CD252 (OX40 Ligand)	Rat IgG2b, k
E4	117805	MD-113	anti-mouse MD-1	Rat IgG2b, k
E5	120208	YGITR 765	anti-mouse CD357 (GITR)	Rat IgG2b, k
E6	145504	L138D7	anti-mouse CXCR5	Rat IgG2b, k
E7	146204	Duno85	anti-mouse CD37	Rat IgG2b, k
E8	148004	TX52	anti-mouse CD300c/CD300d/MAIR-II	Rat IgG2b, k
E9	151104	SA051D1	anti-mouse CXCR6	Rat IgG2b, k
E10	149404	4H1B35	anti-mouse CD130 (gp130)	Rat IgG2b, k
E11	150312	SA214G2	anti-mouse CD198 (CCR8)	Rat IgG2b, k
E12	150410	SA275A11	anti-mouse CD20	Rat IgG2b, k
F1	144804	I015F8	anti-mouse CD124 (IL4Ra)	Rat IgG2b, k
F2	150904	12B2B64	anti-mouse IL-23R	Rat IgG2b, k
F3	146506	L276F12	anti-mouse CXCR4	Rat IgG2b, k
F4	100108	RM2-5	anti-mouse CD2	Rat IgG2b, λ
F5	400707	RTK4174	Rat IgG2c, k isotype control	Rat IgG2c, k
F6	128008	HK1.4	anti-mouse Ly-6C	Rat IgG2c, k
F7	138604	49-H4	anti-mouse Ly-6D	Rat IgG2c, k
F8	400808	RTK2118	Rat IgM, k isotype control	Rat IgM, k
F9	108908	DX5	anti-mouse CD49b (pan-Nk cells)	Rat IgM, k
F10	144608	GL7	anti-mouse GL7 antigen (Ly77)	Rat IgM, k
F11	402008	SHG-1	Syrian hamster IgG isotype control	Syrian Hamster IgG

LEGENDScreen™ Mouse PE Kit

Well	Cat #	Clone	Description	Isotype
B12	123510	1B1	anti-mouse CD1d (CD1.1, Ly-38)	Rat IgG2b, k
C1	104606	FR70	anti-mouse CD70	Rat IgG2b, k
C2	100408	GK1.5	anti-mouse CD4	Rat IgG2b, k
C3	107608	M5/114.15.2	anti-mouse I-A/I-E	Rat IgG2b, k
C4	106405	RM153	anti-mouse CD153	Rat IgG2b, k
C5	116108	YN1/1.7.4	anti-mouse CD54	Rat IgG2b, k
C6	124905	33D1	anti-mouse DC Marker (33D1)	Rat IgG2b, k
C7	105308	30-H12	anti-mouse CD90.2	Rat IgG2b, k
C8	116208	TER-119	anti-mouse TER-119/Erythroid Cells	Rat IgG2b, k
C9	103608	R1-2	anti-mouse CD49d	Rat IgG2b, k
C10	101808	M1/69	anti-mouse CD24	Rat IgG2b, k
C11	108408	RB6-8C5	anti-mouse Ly-6G/Ly-6C (Gr-1)	Rat IgG2b, k
C12	105106	PO3	anti-mouse CD86	Rat IgG2b, k
D1	101208	M1/70	anti-mouse/human CD11b	Rat IgG2b, k
D2	103106	30-F11	anti-mouse CD45	Rat IgG2b, k
D3	109104	RMP1-30	anti-mouse CD279 (PD-1)	Rat IgG2b, k
D4	130107	CX1	anti-mouse RAE-1 γ	Rat IgG2b, k
D5	126608	YTS156.7.7	anti-mouse CD8b (Ly-3)	Rat IgG2b, k
D6	103008	IM7	anti-mouse/human CD44	Rat IgG2b, k
D7	115806	D7715A7	anti-mouse/rat CD126 (IL-6 R α chain)	Rat IgG2b, k
D8	127010	927	anti-mouse CD317 (BST2, PDCA-1)	Rat IgG2b, k
D9	132306	TUGm2	anti-mouse CD132 (common γ chain)	Rat IgG2b, k

Well	Cat #	Clone	Description	Isotype
F12	102106	37.51	anti-mouse CD28	Syrian Hamster IgG
G1	127408	8.1.1	anti-mouse Podoplanin	Syrian Hamster IgG
G2	106106	17B5	anti-mouse CD137	Syrian Hamster IgG
G3	107706	15F9	anti-mouse CD278 (ICOS)	Syrian Hamster IgG
G4	138408	2F1/KLRG1	anti-mouse/human KLRG1 (MAFA)	Syrian Hamster IgG
G5	108208	14B11	anti-mouse Ly-49C/F/I/H	Syrian Hamster IgG
G6			Blank	
G7			Blank	
G8			Blank	
G9			Blank	
G10			Blank	
G11			Blank	
G12			Blank	
H1			Blank	
H2			Blank	
H3			Blank	
H4			Blank	
H5			Blank	
H6			Blank	
H7			Blank	
H8			Blank	
H9			Blank	

LEGENDScreen™ Mouse PE Kit

Well	Cat #	Clone	Description	Isotype
D10	100206	17A2	anti-mouse CD3	Rat IgG2b, k
D11	124308	10F.9G2	anti-mouse CD274 (B7-H1, PD-L1)	Rat IgG2b, k
D12	105808	2B8	anti-mouse CD117 (c-kit)	Rat IgG2b, k

Well	Cat #	Clone	Description	Isotype
H10			Blank	
H11			Blank	
H12			Blank	



LEGENDScreen™ Kits are manufactured by **BioLegend Inc.**
9727 Pacific Heights Blvd.
San Diego, CA 92121
Tel: 1.858.768.5800
Tel: US & Canada Toll-Free: 1.877.Bio-Legend
(1.877.246.5343)
Fax: 1.877.455.9587
Email: info@biolegend.com
biolegend.com

For a complete list of world-wide BioLegend offices and distributors, please visit our website at: biolegend.com

700005 V1