

MaxPar® Ready

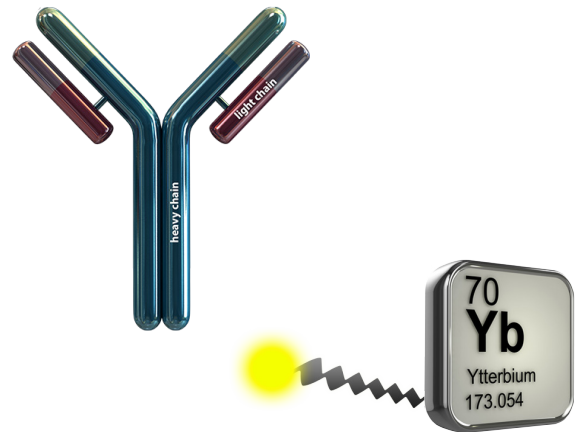
CyTOF® Validated Antibodies

So you have a new CyTOF® instrument? Then get CyTOF®-validated purified antibodies for metal labeling from BioLegend. MaxPar® Ready purified antibodies are optimized and ready-to-use with MaxPar® metal labeling kits from DVS Sciences.

Why choose MaxPar® Ready purified antibodies?

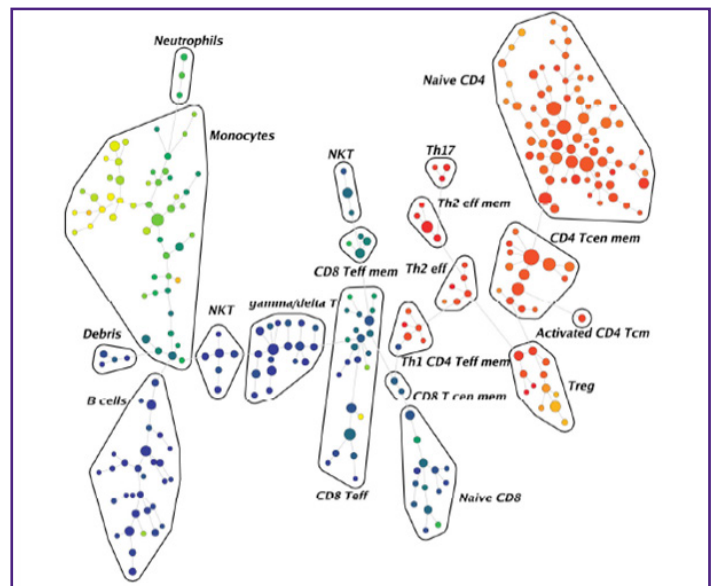
- High concentration (1.0 mg/ml), optimal for conjugation
- No additional proteins in the buffer, ready-to-use
- Ultra-low protein binding tubes guarantees accuracy of antibody amount
- High quality, low cost

Request a sample or bulk today: sales@biolegend.com



CyTOF® mass cytometry is an emerging and exciting application for research

CyTOF® mass cytometry utilizes rare metals tagged to antibodies, allowing researchers to examine over 30 different proteins on and within a single cell without concern of overlapping emission spectra. This single cell proteomics platform enables simultaneous deep phenotyping with functional profiling and has already been used in a diverse range of research areas, including immunology, oncology and hematopoietic stem cell research.



SPADE diagram of PBMCs analyzed by CyTOF® mass cytometry. Node size represents the number of cells and the color indicates the expression of CD4.



Toll-Free Tel: (US & Canada): 1.877.BIOLEGEND (246.5343)

Tel: 858.768.5800

biolegend.com

07-0073-00

MaxPar[®] Ready Purified Antibodies

For the latest product updates and product information, visit: [biolegend.com maxpar_ready](http://biolegend.com/maxpar_ready)

Human

Specificity	Clone	Cat. No.
CD3	UCHT1	300443
CD4	RPA-T4	300541
CD8a	RPA-T8	301053
CD10	HI10a	312223
CD11b	M1/70	101249
CD11b	ICRF44	301337
CD11c	3.9	301639
CD11c	Bu15	337221
CD14	M5E2	301843
CD15	W6D3	323035
CD16	3G8	302051
CD19	HIB19	302247
CD20	2H7	302343
CD27	O323	302839
CD33	WMS3	303419
CD34	581	343531
CD38	HIT2	303535
CD41	HIP8	303721
CD44	IM7	103051
CD44	BJ18	338811

Specificity	Clone	Cat. No.
CD45	HI30	304045
CD45R	RA3-6B2	103249
CD45RA	HI100	304143
CD56 (NCAM)	HCD56	318345
CD61	VI-PL2	336413
CD69	FN50	310939
CD90 (Thy1)	5E10	328129
CD117 (c-kit)	104D2	313223
CD123	6H6	306027
CD161	HP-3G10	339919
CD184 (CXCR4)	12G5	306523
CD235ab	HIR2	306615
HLA-DR	L243	307651
IFN- γ	B27	506521
IL-2	MQ1-17H12	500339
IL-4	MP4-25D2	500829
IL-5	TRFK5	504309
IL-6	MQ2-13A5	501115
Ki-67	Ki-67	350523
TNF- α	MAB11	502941

Mouse

Specificity	Clone	Cat. No.
CD4	RM4-5	100561
CD5	53-7.3	100619
CD8a	53-6.7	100755
CD11b	M1/70	101249
CD11c	N418	117341
CD19	6D5	115547
CD25	3C7	101913
CD44	IM7	103051
CD45	30-F11	103141
CD45R	RA3-6B2	103249

Specificity	Clone	Cat. No.
CD62L	MEL-14	104443
IFN- γ	XMG1.2	505843
IL-4	11B11	504129
IL-5	TRFK5	504309
IL-6	MP5-20F3	504509
IL-10	JES5-16E3	505029
IL-17A	TC11-18H10.1	506935
NK-1.1	PK136	108743
TCR- β chain	H57-597	109235
TNF- α	MP6-XT22	506335

MaxPar[®] and CyTOF[®] are registered trademarks of DVS Sciences



Listen to The LEGENDeers new song:

Heavy Metal Cytometry.

This is the first song ever written about mass cytometry.

Enjoy: biolegend.com/maxpar_ready

Download it for free at: cdbaby.com/cd/thelegendeers2



Toll-Free Tel: (US & Canada): 1.877.BIOLEGEND (246.5343)

Tel: 858.768.5800

biolegend.com

07-0073-00