

# BD® RUO/GMP Reagents

Manufactured in accordance with Good Manufacturing Practices (GMP) for devices under 21 CFR Part 820

Marker	Violet laser			Blue laser					Red laser				Biotin	Pure
	Emission max	448 nm	491 nm	602 nm	519 nm	578 nm	378 nm	695 nm	785 nm	660 nm	704 nm	785 nm		
	BD Horizon™ Reagent V450	AmCyan	BD Horizon™ Reagent BV605	FITC	PE	PerCP	PerCP-Cy5.5	PE-Cy7	APC	APC-R700	APC-Cy7	APC-H7		
CD1a			663805		341651									
CD2	644485			347593	347597			335786	341022					347590
CD3	652355	339186			347347	347344	340949	341091	340440	659119	341090	641397		347340
CD4 v4				340422	340419									
CD4	651849	339187			347327	347324	341654	348789	340443		341095	641398	347321	346320
CD5	644487			347303	347307		341089	348790	340583				347301	346300
CD7	642916			688220	340581				653311	659124				
CD8		339188		347313		347314	341051	335787	340584		348793	641400		346310
CD9	658167			341646	341647		341649		341648			655409		
CD10				688428	999287			341092	340923	659120		655404		347500
CD11a				347983										
CD11b					347557				340937					
CD11c			663807		347637		658330	652358	340544					
CD13					347837			338425		657698				347830
CD14				347493	347497	340585			340436		333945	641394		347490
CD15	642917			347423										347420
CD16	644489			1003669	347617			338426	335788					347530
CD18				347953										
CD19	644491	339190		687952	340364	687965	340951	341093	340437	659121	348794	641395		347540
CD20	642274			347673	346595	347674	340955	335793	340941		335794	641396		347670
CD21	658169													
CD22				347573	347577		658329		340933					
CD23				656148	341007				340935	659125				347790
CD24												658331		
CD25				347643	341009			335789	340939	659126				347640
CD26				340426	340423									
CD27				340424	340425		656643		337169					
CD28					348047	337181								689015
CD30				341644	341645									
CD31					340297									
CD33				340533	347787		341650	333946	340474	659122				347780
CD34				348053	348057	340430	347203	348791	340441	659123				695710
CD36				656151										
CD38				340927	347687			335790	340439	659127		656646		347680
CD41a				340929			340931							
CD42a				348083		340537								
CD43												655407		
CD44				347943										
CD45RA								337167						
CD45RO					347967			337168	340438					
CD45	642275	339192		695186		347464	340953		340943		348795	641399		347460
CD49d					340296							658332		340976
CD54					347977									
CD55					341030									
CD56				340410	688110			335791	341025	657886				347740
CD57				347393									347391	347390
CD58					340295									
CD61				348093		340506								
CD62L				347443	341012									
CD62P					348107									
CD64					644385					657700				
CD69				347823	341652	340548		335792	340560					347820
CD71				347513					341028			655408		347510
CD79a					340579		341643		340578					
CD79b				335798	335799		656644		335800	657702				
CD80					340294									
CD81												656647		
CD95				340479	340480				340481					
CD103				340945	340998									
CD107a									641581					
CD117					340529		333944	339195	341096					
CD122					340254									
CD123					340545									
CD134					340420				658171					
CD138				347191	347192		341087		347193					
CD154					340477				648887					
CD158a				340531										
CD161				340536	340530									
CD200									655406	657868				
Avidin				349011										
β2-Microglobulin							656645							347580
Bcl-2				340575	340576									
BrdU				347583										
Cytokeratin				347653	347204									
EpCAM				347197	347198		347199		347200					
FMCF7	644493			340919										
Glycophorin A					340947									
Goat Anti-Ms Ig				349031										
HER-2/neu				340553	340552				340554					
HLA-DQ				347453										
HLA-DR	642276			347363	347367	347364	339194	335795	340549		335796	641393	347361	347360
IFN-γ				340449	340452				341117					
IgG														345140
IgM														345160
IL-1				340513										
IL-1α					340514									
IL-1β				340515	340516									
IL-1ra					340525									
IL-2				340448	340450				341116					
IL-4					340451									
IL-6				340526	340527									
IL-8				340509	340510									
IL-13					340508									
IREM-2									656158					
Kappa	658286			643773	346601				341088					345120
Kappa F(ab)2				348063										
KIR				340483	340484									
Lambda				346600	642919							656648		345130
Lambda F(ab)2				348073										
Myeloperoxidase (MPO)				340580	341642									
Multiclonal CD4				347413										
PAC-1				340507										340535
P-glycoprotein					340555									
Streptavidin					349023	340130			349024					
TCR-α/β				347773										347770
TCRγ/δ				347903	347907			655410		657706				347900
Terminal-Deoxynucleotidyl Transferase (TdT)				347194	347195				347196					
TNF-α				340511	340512				340534					
ZAP-70				344931	344635									
Mouse IgG1 Isotype Control	642268			349041	349043 340013	349044	347202	348788	340442		348792	641401		349040
Mouse IgG2a Isotype Control				349051	349053	349054			340473					349050
Ms IgG1 Isotype Control		339185	657720							657708				
Ms IgG2a Isotype Control					340459									
Rat Anti-Ms Ig Kpa Isotype Control					349073									
Rat Anti-Ms IgG1 Isotype Control					340270	340272								
Rat Anti-Ms IgG2a/b Isotype Control					340269	340271								

Visit [bdbiosciences.com/RUOGMP](https://bdbiosciences.com/RUOGMP) for more information.

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

BD Life Sciences, San Jose, CA, 95131, U.S.

[bdbiosciences.com](https://bdbiosciences.com)

BD, the BD Logo and Horizon are trademarks of Becton, Dickinson and Company. All other trademarks are the property of their respective owners. © 2023 BD.

All rights reserved. BD-83159 (v1.0) 0223

Cy is a trademark of Global Life Sciences Solutions Germany GmbH or an affiliate doing business as Cytiva.

