

Catalog #	Protein Name	Total Lots	Total Remaining Amount (mg)
100317	GYG/GYS2	1	0.11
100638	NKG2A	1	0.12
100410	FXR	1	0.15
100663	Siglec11 (biotinylated)	1	0.15
102342	PTK7(biotinylated)	1	0.15
101332	HDAC6 (rat)	2	0.16
50007	HDAC7	1	0.21
100201	PDGFRa (D842Y)	2	0.22
101834	FGFR1	1	0.25
40403	BRK	1	0.26
102419	IL33	1	0.26
40092	MAPK10	2	0.27
101091	Spike(B.1.351 delta242-244)(SARS-CoV-2)	1	0.27
102167	KAT8	1	0.27
100466	NKp46 (biotinylated)	1	0.28
100662	Siglec11	1	0.29
100465	NKp46	1	0.36
50260	AHCY	1	0.39
40630	p85a	1	0.42
40267	Ret	1	0.43
100294	CUL2	1	0.43
100645	TGFbR1 (biotinylated)	1	0.5
100264	PDE3A (dog)	2	0.51
60099	PDE10A1	1	0.55
102360	RNF43	1	0.56
100078	PRDM9	3	0.59
40595	RB	2	0.62
101495	CUL4B/ Rbx1(M5I)/ DDB1/DCAF11	2	0.66
50015	SIRT4	4	0.73
71256	GITR	1	0.77
79472	CD19	2	0.77
102826	USP30	1	0.78
40371	RIPK1	2	0.82
102421	IL33(biotinylated)	1	0.83
60042	PDE4B2	1	0.95
100071	GPC3	2	0.96
100644	TGFbR1	1	1.06
71180	TIM-3	1	1.1
102249	NcoR2, GST-Tag Recombinant	1	1.12
50009	HDAC9	1	1.34
101755	Cul1 /Skp1/Skp2	1	1.38
101336	IL12RB1	1	1.48
71258	TUT4(biotinylated)	4	1.5
79366	B7-H7(biotinylated)	1	1.56
80401	XIAP	3	1.6
71099	IDH1(R132H)	1	1.63

101872	CD8a	1	1.68
101991	ACC1	3	1.74
100426	FNDC5 (biotinylated)	1	1.76
102823	NEU1(48-415e)	1	1.8
102163	TSLPR	2	1.84
102198	HER2	2	1.96
70020	Peroxiredoxin I	1	2.02
50014	SIRT3	5	2.17
80506	PARP6	2	2.18
101470	CLEC2D (biotinylated)	1	2.23
102231	HER3 (biotinylated)	1	2.37
101813	CD36	1	2.51
100284	TACI (biotinylated)	1	2.53
102043	CERK	1	2.68
71077	Fucokinase	1	2.72
101543	CD93	1	2.74
100421	Siglec8	1	2.75
101414	KIR3DL3	1	2.82
100657	FGFR3 (PTP1B coexpressed)(V555F,dephosphorylated)	2	2.89
102516	IL31Ra (20-519-Fc (IgG1)-AviBiotin)	1	2.92
100068	cd123 (biotinylated)	1	2.93
101886	FGF19	1	3.43
102238	HER4, Fc Fusion, Avi-Tag Recombinant	1	3.47
101725	PCSK9	2	3.56
101469	CLEC2D (LLT1), His-Tag, Avi-Tag Recombinant	2	3.62
50270	ALAS2	1	3.63
50012	SIRT1	1	3.64
51001	G9a	2	3.73
101377	DDR2 (biotinylated)	1	3.81
79016	FCGR3b	1	3.87
70823	RANK	1	3.94
100349	VSIG3	2	3.96
102535	ADAR1 (E1008Q)	2	4.03
80213	MMP2	1	4.09
102035	CD200R1(biotinylated)	1	4.14
100677	Tgfbr2 (mouse,biotinylated)	1	4.26
79384	B7-H6 (biotinylated)	1	4.47
100467	BTNL8	1	4.48
100695	Spike(RBD) (SARS-CoV-2,biotinylated)	1	4.57
101882	TL1A(TNFSF15)(biotinylated)	1	4.57
101888	FGF19(biotinylated)	1	4.65
100628	TGFBR2	2	4.71
101812	CD34	1	4.84
40195	EGFR (mouse)	3	5.28
102538	IL4	1	5.31
102539	IL4 (Biotin)	1	5.4
101944	FGF21 (His-Avi-29-209e)	1	5.43

101820	FGFR3 (biotinylated)	1	5.54
102468	IL4Ra (Biotin)	1	5.58
102570	ACVR2B(biotin)	1	5.58
71093	Carbonic Anhydrase 1	1	5.68
100546	NKp30	1	5.81
100468	BTNL8 (biotinylated)	1	5.85
102675	EGFR	1	5.96
100527	CD24	1	6.09
102495	MADCAM1 (biotin)	1	6.12
50201	ACC2	4	6.36
102507	IL5 (His-Avi-20-134eBiotin)	1	6.48
102467	IL4Ra	1	6.53
100545	ULBP2 (biotinylated)	3	6.54
100424	Siglec15 (biotinylated)	1	6.57
102202	TSLP(R127A, R130A,biotinylated)	1	6.58
101517	TPBG (biotinylated)	1	6.6
100676	Tgfr2 (mouse)	1	6.63
101516	TPBG	1	7.01
102292	Caspase3(rat)	4	7.14
101597	TRAF6	2	7.59
40510	EIF4EBP1	2	7.66
100155	CD33	1	7.69
100423	Siglec15	1	8.05
102515	IL31Ra (20-519-Fc (IgG1)-Avi)	1	8.13
40254	c-Mer	2	8.24
40640	PI3K (p110α(E545K)/p85α)(E545K)	2	8.52
102506	IL5 (His-Avi-20-134e)	1	8.61
71240	IDO2 (mouse)	2	8.64
102494	MADCAM1	1	8.76
101044	CD56(NCAM1)(biotinylated)	1	9.53
101602	MGMT	2	9.67
79481	ROR1	1	10.15
100698	Spike(RBD) (SARS-CoV-2)	1	11.53
50013	SIRT2	2	11.79
101667	USP1 (G670A, G671A)/UAF1 Complex Recombinant	2	12.12
100469	BTN1A1	1	12.3
102057	NPM1	1	12.43
100509	CEACAM5	2	12.76
102596	PARP1	2	12.8
100425	FNDC5	3	13.71
102451	IL1RL1	1	14.25
101945	DNA Polymerase Theta	5	14.47
101669	3C-like Proteasetagless(SARS-CoV-2,T21I, S144A)	1	15.54
102676	EGFR (biotin)	1	17.37
102293	CD157	1	18.25
101781	DLK1 (biotinylated)	1	18.38

80503	PARP3	3	18.68
79124	PRMT9	2	18.69
101690	PARP4	5	18.83
102101	VEGFR3	2	18.86
40449	Jak1	3	19.17
71098	NAMPT	5	20.42
101913	PI3K(p110α(E542K)/p85α)	1	20.69
101811	FGFR3	2	21.62
101726	PARG	3	21.83
40051	DAPK1	3	22.93
40122	MEKK2	1	23.81
101836	FGFR1	1	25.08
101515	TPBG (biotinylated)	2	25.1
40612	Sphingosine Kinase 2	1	33.16
101877	FGFR2	2	33.62
80500	Caspase 3	4	39.98
52013	Histone H3	1	43.38
40089	MKK6	2	82.16
100490	COMT	3	272.48
101880	TL1A(TNFSF15)	2	274.36
11005	ACHLY (B. pertussis)	4	292.49