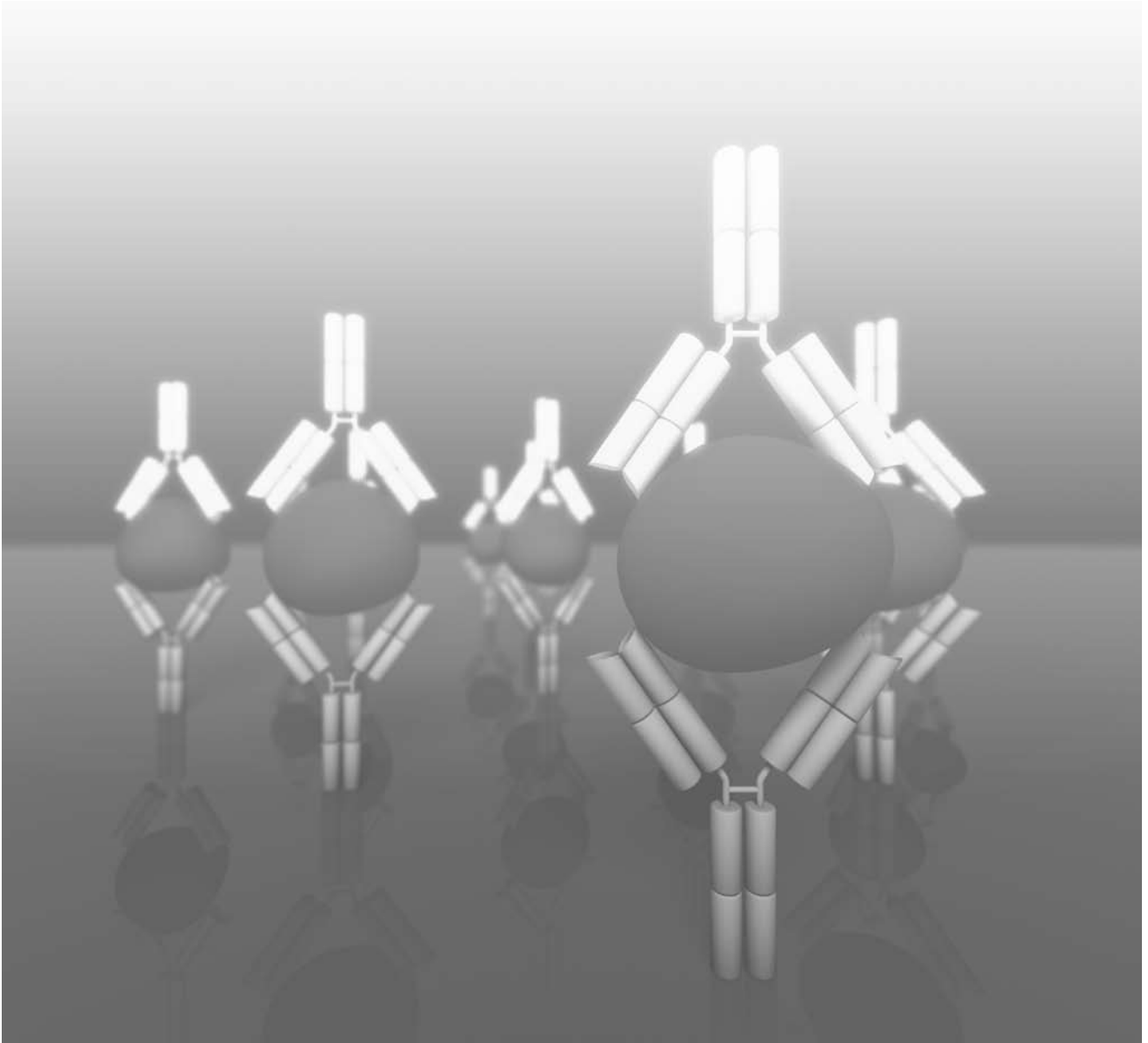

BD Biosciences
ELISA and ELISPOT Reagents

Product List



ELISA Reagents

ELISA Kits

DESCRIPTION	REG	SIZE	CAT. NO.
Canine C-reactive protein (CRP) ELISA Kit	RUO	1 plate	557826
Human C3a ELISA Kit	RUO	1 plate	550499
Human C4a ELISA Kit	RUO	1 plate	550947
Human C5a ELISA Kit II	RUO	1 plate	557965
Human IFN- γ ELISA Kit II	RUO	2 plates	550612
Human IL-1 β ELISA Kit II	RUO	2 plates	557966
Human IL-2 ELISA Kit II	RUO	2 plates	550611
Human IL-4 ELISA Kit II	RUO	2 plates	550614
Human IL-6 ELISA Kit II	RUO	2 plates	550799
Human IL-8 ELISA Kit II	RUO	2 plates	550999
Human IL-10 ELISA Kit II	RUO	2 plates	550613
Human IL-12 (p40) ELISA Kit II	RUO	2 plates	551116
Human IL-12 (p70) ELISA Kit	RUO	2 plates	559258
Human MCP-1 ELISA Kit	RUO	2 plates	559017
Human TNF ELISA Kit II	RUO	2 plates	550610
Mouse IFN- γ ELISA Kit II	RUO	2 plates	558258
Mouse IL-6 ELISA Kit	RUO	2 plates	550950
Mouse Immunoglobulin Isotyping ELISA Kit	RUO	80 tests	550487
Mouse TNF ELISA Kit	RUO	2 plates	560478
Rat C-reactive protein (CRP) ELISA Kit	RUO	1 plate	557825
Rat Immunoglobulin Isotyping ELISA Kit	RUO	80 tests	557081
Rat TNF ELISA Kit	RUO	2 plates	560479

ELISA Reagents

ELISA Sets

DESCRIPTION	REG	SIZE	CAT. NO.
Caspase-3, Active Form, Human ELISA Set (CPP32)	RUO	Reagents for 5 plates	550930
Human C5b-9 ELISA Set	RUO	Reagents for 5 plates	558315
Human CD54 (sICAM-1) ELISA Set	RUO	Reagents for 20 plates	551424
Human CD253 ELISA Set (TRAIL)	RUO	Reagents for 20 plates	550948
Human Cleaved PARP ELISA Set	RUO	Reagents for 5 plates	552592
Human Eotaxin ELISA Set	RUO	Reagents for 20 plates	555175
Human GM-CSF ELISA Set	RUO	Reagents for 20 plates	555126
Human IFN- γ ELISA Set	RUO	Reagents for 20 plates	555142
Human IL-1 β ELISA Set II	RUO	Reagents for 20 plates	557953
Human IL-2 ELISA Set	RUO	Reagents for 20 plates	555190
Human IL-2 sR α ELISA Set	RUO	Reagents for 20 plates	559104
Human IL-4 ELISA Set	RUO	Reagents for 20 plates	555194
Human IL-5 ELISA Set	RUO	Reagents for 20 plates	555202
Human IL-6 ELISA Set	RUO	Reagents for 20 plates	555220
Human IL-8 ELISA Set	RUO	Reagents for 20 plates	555244
Human IL-10 ELISA Set	RUO	Reagents for 20 plates	555157
Human IL-12 (p40) ELISA Set	RUO	Reagents for 20 plates	555171
Human IL-12 (p70) ELISA Set	RUO	Reagents for 20 plates	555183
Human IL-15 ELISA Set	RUO	Reagents for 20 plates	559268
Human IP-10 ELISA Set	RUO	Reagents for 20 plates	550926
Human LT- α ELISA Set	RUO	Reagents for 20 plates	550995
Human MCP-1 ELISA Set	RUO	Reagents for 20 plates	555179
Human MIG ELISA Set	RUO	Reagents for 20 plates	550998
Human TGF- β 1 ELISA Set	RUO	Reagents for 20 plates	559119
Human TNF ELISA Set	RUO	Reagents for 20 plates	555212
Monkey IFN- γ ELISA Set	RUO	Reagents for 20 plates	551492
Monkey IL-2 ELISA Set	RUO	Reagents for 20 plates	551494

ELISA Reagents

ELISA Sets

DESCRIPTION	REG	SIZE	CAT. NO.
Mouse GM-CSF ELISA Set	RUO	Reagents for 20 plates	555167
Mouse IFN- γ ELISA Set	RUO	Reagents for 20 plates	555138
Mouse IFN- γ ELISA Set (AN-18)	RUO	Reagents for 20 plates	551866
Mouse IgE ELISA Set	RUO	Reagents for 20 plates	555248
Mouse IgG _{2a} ELISA Set	RUO	Reagents for 20 plates	552576
Mouse IL-1 α ELISA Set	RUO	Reagents for 20 plates	550347
Mouse IL-1 β ELISA Set	RUO	Reagents for 20 plates	559603
Mouse IL-2 ELISA Set	RUO	Reagents for 20 plates	555148
Mouse IL-3 ELISA Set	RUO	Reagents for 20 plates	555228
Mouse IL-4 ELISA Set	RUO	Reagents for 20 plates	555232
Mouse IL-5 ELISA Set	RUO	Reagents for 20 plates	555236
Mouse IL-6 ELISA Set	RUO	Reagents for 20 plates	555240
Mouse IL-10 ELISA Set	RUO	Reagents for 20 plates	555252
Mouse IL-12 (p40) ELISA Set	RUO	Reagents for 20 plates	555165
Mouse IL-12 (p70) ELISA Set	RUO	Reagents for 20 plates	555256
Mouse MCP-1 ELISA Set	RUO	Reagents for 20 plates	555260
Mouse TNF (Mono/Mono) ELISA Set	RUO	Reagents for 20 plates	555268
Mouse TNF (Mono/Mono) ELISA Set II	RUO	Reagents for 20 plates	558534
Rat IFN- γ ELISA Set	RUO	Reagents for 20 plates	558861
Rat IL-4 ELISA Set	RUO	Reagents for 20 plates	555198
Rat IL-6 ELISA Set	RUO	Reagents for 20 plates	550319
Rat IL-10 ELISA Set	RUO	Reagents for 20 plates	555134
Rat MCP-1 ELISA Set	RUO	Reagents for 20 plates	555130
Rat TNF ELISA Set	RUO	Reagents for 20 plates	558535

ELISA Reagents

In Vivo Capture Assays

DESCRIPTION	REG	SIZE	CAT. NO.
IFN- γ , Mouse <i>In Vivo</i> Capture Assay	RUO	10 plates	558491
IL-2, Mouse <i>In Vivo</i> Capture Assay	RUO	10 plates	557937
IL-4, Mouse <i>In Vivo</i> Capture Assay	RUO	10 plates	557781
IL-6, Mouse <i>In Vivo</i> Capture Assay	RUO	10 plates	558075
IL-10, Mouse <i>In Vivo</i> Capture Assay	RUO	10 plates	558072
TNF, Mouse <i>In Vivo</i> Capture Assay	RUO	10 plates	558568

ELISA Reagents

ELISA Pairs - Human

DESCRIPTION	REACT	CLONE	ISOTYPE	REG	FORMAT	SIZE	CAT. NO.
Blys (BAFF)	Hu	H32-406	Mouse IgG ₁ , κ	RUO	Purified	0.1 mg	558289
		H32-544	Mouse IgG ₁ , κ	RUO	Biotin	0.1 mg	558290
Caspase-3, Active Form, ELISA Pair	Hu	19, C92-605		RUO	Purified, Purified	Reagents for 10 plates	550578
CD120a (TNF Receptor Type I)	Hu	MABTNFR1-A1	Mouse IgG ₁	RUO	Purified	0.5 mg	552535
		MABTNFR1-B1	Mouse IgG _{2a}	RUO	Biotin	0.5 mg	552536
CD126 (IL-6 Receptor α chain)	Hu	M182	Mouse IgG ₁ , κ	RUO	Biotin	0.5 mg	552503
		M5	Mouse IgG ₁ , κ	RUO	Purified	0.5 mg	551462
CD137 (4-1BB)	Hu	4B4-1	Mouse IgG ₁ , κ	RUO	Biotin	0.5 mg	552533
		H41BB-M127	Mouse IgG ₁ , κ	RUO	Purified	0.5 mg	552532
CD154 (CD40L)	Hu	hCD40L-M90.1	Mouse IgG ₁ , κ	RUO	Purified	0.5 mg	552559
		hCD40L-M91.2	Mouse IgG ₁ , κ	RUO	Biotin	0.5 mg	552560
Eotaxin	Hu	10C11	Mouse IgG _{2b}	RUO	Biotin	0.5 mg	555060
		3C7	Mouse IgG ₁	RUO	Purified	0.5 mg	555035
Eotaxin, Recombinant human				RUO		5 µg	555102
G-CSF	Hu	BVD11-37G10	Rat IgG _{2a}	RUO	Biotin	0.5 mg	554670
		BVD13-3A5	Rat IgG ₁	RUO	Purified	1.0 mg	551342
GM-CSF	Hu	BVD2-21C11	Rat IgG _{2a}	RUO	Biotin	0.5 mg	554505
		BVD2-23B6	Rat IgG _{2a}	RUO	Purified	0.5 mg	554502
GM-CSF, Recombinant human				RUO		10 µg	550068
IFN-γ	Hu	4S.B3	Mouse IgG ₁ , κ	RUO	Biotin	0.5 mg	554550
		NIB42	Mouse IgG ₁	RUO	Purified	1.0 mg	551221
IFN-γ, Recombinant human				RUO		25 µg	554616
				RUO		50 µg	554617
IL-2	Hu	5344.111	Mouse IgG ₁ , κ	RUO	Purified	0.5 mg	555051
		B33-2	Mouse IgG ₁	RUO	Biotin	0.5 mg	555040
IL-2, Recombinant human				RUO		10 µg	554603
IL-3	Hu	BVD3-1F9	Rat IgG ₁	RUO	Biotin	0.5 mg	554674
		BVD8-3G11	Rat IgG ₁	RUO	Purified	0.5 mg	554672
IL-3, Recombinant human				RUO		10 µg	554604
IL-4	Hu	8D4-8	Mouse IgG ₁ , κ	RUO	Purified	0.5 mg	554515
		MP4-25D2	Rat IgG ₁	RUO	Biotin	0.5 mg	554483
IL-4, Recombinant human				RUO		5 µg	554605
IL-5	Hu	JES1-39D10	Rat IgG _{2a}	RUO	Purified	0.5 mg	554488
		JES1-5A10	Rat IgG _{2a}	RUO	Biotin	0.5 mg	554491
IL-5, Recombinant human				RUO		5 µg	554606
IL-6	Hu	MQ2-13A5	Rat IgG ₁	RUO	Purified	0.5 mg	554543
		MQ2-39C3	Rat IgG _{2a}	RUO	Biotin	0.5 mg	554546
IL-6, Recombinant human				RUO		10 µg	550071
IL-7	Hu	BVD10-11C10	Rat IgG _{2a}	RUO	Biotin	0.5 mg	554494
		BVD10-40F6	Rat IgG ₁ , κ	RUO	Purified	0.5 mg	554493
IL-7, Recombinant human				RUO		5 µg	554608
IL-8	Hu	G265-5	Mouse IgG _{2b} , κ	RUO	Purified	0.5 mg	554716
		G265-8	Mouse IgG _{2b}	RUO	Biotin	0.5 mg	554718
IL-8, Recombinant human				RUO		20 µg	554609
IL-10	Hu, Vir	JES3-12G8	Rat IgG _{2a}	RUO	Biotin	0.5 mg	554499
		JES3-19F1	Rat IgG _{2a}	RUO	Purified	0.5 mg	554705
IL-10, Recombinant human				RUO		5 µg	554611
IL-12 (p40)	Hu	C8.3	Mouse IgG ₁	RUO	Purified	1.0 mg	551227
IL-12 (p40/p70)	Hu	C8.6	Mouse IgG ₁	RUO	Biotin	0.5 mg	554660
IL-12 (p40), Recombinant human				RUO		5 µg	554633
IL-12 (p70)	Hu	20C2	Rat IgG ₁	RUO	Purified	0.5 mg	555065
IL-12 (p40/p70)	Hu	C8.6	Mouse IgG ₁	RUO	Biotin	0.5 mg	554660
IL-12 (p70), Recombinant human				RUO		5 µg	554613

ELISA Reagents

ELISA Pairs - Human

DESCRIPTION	REACT	CLONE	ISOTYPE	REG	FORMAT	SIZE	CAT. NO.
IL-13	Hu	JES10-5A2	Rat IgG ₁	RUO	Purified	0.5 mg	554570
		B69-2	Mouse IgG ₁	RUO	Biotin	0.5 mg	555054
IL-15	Hu	G243-935	Mouse IgG ₁	RUO	Purified	0.5 mg	554712
		G243-886	Mouse IgG ₁	RUO	Biotin	0.5 mg	554713
IL-15, Recombinant human						5 µg	560254
IP-10	Hu	4D5	Mouse IgG _{2a}	RUO	Purified	0.5 mg	555046
		6D4/D6/G2	Mouse IgG _{2a}	RUO	Biotin	0.5 mg	555048
IP-10, Recombinant human						10 µg	551130
LT-α	Hu	359-238-8	Mouse IgG ₁	RUO	Purified	1.0 mg	551222
		359-81-11	Mouse IgG ₁	RUO	Biotin	0.5 mg	554555
LT-α, Recombinant human						10 µg	554619
MCP-1	Hu	10F7	Mouse IgG ₁	RUO	Purified	0.5 mg	555055
		5D3-F7	Mouse IgG ₁ , κ	RUO	Biotin	0.5 mg	554664
MCP-1, Recombinant human						10 µg	554620
MCP-3	Hu	9H11	Mouse IgG ₁	RUO	Purified	1.0 mg	551228
		6H5	Mouse IgG ₁	RUO	Biotin	0.5 mg	555034
MCP-3, Recombinant human						10 µg	554629
MIG	Hu	B8-11	Mouse IgG ₁	RUO	Purified	0.5 mg	555038
		B8-6	Mouse IgG ₁	RUO	Biotin	0.5 mg	555037
MIG, Recombinant human						5 µg	554636
sCD14 (Soluble CD14)	Hu	55-3	Rat IgG _{2a} , κ	RUO	Purified	0.5 mg	551403
		3-C39	Mouse IgG _{2a} , κ	RUO	Biotin	0.5 mg	551405
TGF-β1	Hu, Ms, Pig	A75-2	Rat IgG _{2a}	RUO	Purified	0.5 mg	555052
		A75-3	Rat IgG _{2a}	RUO	Biotin	0.5 mg	555053
TNF	Hu	MAB1	Mouse IgG ₁	RUO	Purified	1.0 mg	551220
		MAB11	Mouse IgG ₁ , κ	RUO	Biotin	0.5 mg	554511
TNF, Recombinant human						10 µg	554618

ELISA Reagents

ELISA Pairs - Mouse

DESCRIPTION	REACT	CLONE	ISOTYPE	REG	FORMAT	SIZE	CAT. NO.
C3a	Ms	I87-1162	Rat IgG ₁ , κ	RUO	Purified	0.1 mg	558250
		I87-419	Rat IgG ₁ , κ	RUO	Biotin	0.1 mg	558251
C3a Protein (native), mouse						10 tests	558618
C5a	Ms	I52-1486	Rat IgG ₁ , κ	RUO	Purified	0.1 mg	558027
		I52-278	Rat IgG _{2a} , κ	RUO	Biotin	0.1 mg	558028
C5a, Recombinant mouse						10 tests	622597
GM-CSF	Ms	MP1-22E9	Rat IgG _{2a}	RUO	Purified	0.5 mg	554404
		MP1-31G6	Rat IgG ₁	RUO	Biotin	0.5 mg	554407
GM-CSF, Recombinant mouse						10 µg	554586
IFN-γ	Ms	R4-6A2	Rat IgG ₁	RUO	Purified	1.0 mg	551216
		XMG1.2	Rat IgG ₁ , κ	RUO	Biotin	0.5 mg	554410
IFN-γ, Recombinant mouse						10 µg	554587
IgA	Ms	C10-3	Rat (LOU) IgG ₁ , κ	RUO	Purified	0.5 mg	556969
		C10-1	Rat (LOU) IgG ₁ , κ	RUO	Biotin	0.5 mg	556978
Mouse IgA, κ						0.5 mg	553476
IgE	Ms	R35-118	Rat (LOU) IgG ₁ , κ	RUO	Biotin	0.5 mg	553419
		R35-72	Rat (LOU) IgG ₁ , κ	RUO	Purified	0.5 mg	553413
Mouse IgE, κ						0.5 mg	557079
IgG ₁	Ms	A85-3	Rat (LOU) IgG _{2a} , κ	RUO	Purified	0.5 mg	553445
		A85-1	Rat (LOU) IgG ₁ , κ	RUO	Biotin	0.5 mg	553441
Mouse IgG ₁ , κ						0.5 mg	557273

ELISA Reagents

ELISA Pairs - Mouse

DESCRIPTION	REACT	CLONE	ISOTYPE	REG	FORMAT	SIZE	CAT. NO.
IgG _{2a}	Ms	R11-89	Rat (LOU) IgG ₁ , κ	RUO	Purified	0.5 mg	553446
		R19-15	Rat (LOU) IgG ₁ , κ	RUO	Biotin	0.5 mg	553388
Mouse IgG _{2a} , κ		G155-178	Mouse (BALB/c) IgG _{2a} , κ	RUO	Purified	0.5 mg	553454
IgG _{2b}	Ms	R9-91	Rat (LOU) IgG ₁	RUO	Purified	0.5 mg	553396
		R12-3	Rat (LOU) IgG _{2a} , κ	RUO	Biotin	0.5 mg	553393
Mouse IgG _{2b} , κ		MPC-11	Mouse IgG _{2b} , κ	RUO	Purified	0.5 mg	557351
IgG ₃	Ms	R2-38	Rat IgG ₁ , κ	RUO	Purified	0.5 mg	553404
		R40-82	Rat (LOU) IgG _{2a} , κ	RUO	Biotin	0.5 mg	553401
Mouse IgG ₃ , κ		A112-3	Mouse (BALB/c) IgG ₃ , κ	RUO	Purified	0.5 mg	553486
IgM	Ms	II/41	Rat IgG _{2a} , κ	RUO	Purified	0.5 mg	553435
		R6-60.2	Rat (LOU) IgG _{2a} , κ	RUO	Biotin	0.5 mg	553406
Mouse IgM, κ		G155-228	Mouse (BALB/c) IgM, κ	RUO	Purified	0.5 mg	553472
IL-1α	Ms	ALF-161	Armenian Hamster IgG1, λ 1	RUO	Purified	0.5 mg	550604
		C1150-27	Rabbit Ig	RUO	Biotin	0.5 mg	550606
IL-2	Ms	JES6-1A12	Rat IgG _{2a}	RUO	Purified	0.5 mg	554424
		JES6-5H4	Rat IgG _{2b}	RUO	Biotin	0.5 mg	554426
IL-2, Recombinant mouse				RUO		20 µg	550069
IL-3	Ms	MP2-43D1	Rat IgG _{2a}	RUO	Biotin	0.5 mg	554384
		MP2-8F8	Rat IgG ₁	RUO	Purified	0.5 mg	554381
IL-3, Recombinant mouse				RUO		10 µg	554579
IL-4	Ms	11B11	Rat IgG ₁	RUO	Purified	0.5 mg	554434
		BVD4-1D11	Rat IgG _{2b}	RUO	Purified	0.5 mg	554387
		BVD6-24G2	Rat IgG ₁	RUO	Biotin	0.5 mg	554390
IL-4, Recombinant mouse				RUO		10 µg	550067
IL-5	Hu, Ms	TRFK5	Rat IgG ₁	RUO	Purified	0.1 mg	554392
			RUO	Purified	0.5 mg	554393	
	Ms	TRFK4	Rat IgG _{2a}	RUO	Biotin	0.5 mg	554397
IL-5, Recombinant mouse				RUO		5 µg	554581
IL-6	Ms	MP5-20F3	Rat IgG ₁	RUO	Purified	0.5 mg	554400
		MP5-32C11	Rat IgG _{2a}	RUO	Biotin	0.5 mg	554402
IL-6, Recombinant mouse				RUO		5 µg	554582
IL-9	Ms	D8402E8	Rat IgG _{2a}	RUO	Purified	1.0 mg	551218
		D9302C12	Armenian Hamster IgG2, κ	RUO	Biotin	0.5 mg	554473
IL-10	Ms	JES5-2A5	Rat IgG ₁ , κ	RUO	Purified	1.0 mg	551215
		JES5-16E3	Rat IgG _{2b}	RUO	Biotin	0.5 mg	554465
		SXC-1	Rat IgM	RUO	Biotin	0.5 mg	554423
IL-10, Recombinant mouse				RUO		10 µg	550070
IL-12 (p40/p70)	Ms	C15.6	Rat IgG ₁	RUO	Purified	1.0 mg	551219
		C17.8	Rat IgG _{2a}	RUO	Biotin	0.5 mg	554476
IL-12 (p40), Recombinant mouse				RUO		2 µg	554594
IL-12 (p70)	Ms	9A5	Rat IgG _{2b}	RUO	Purified	0.5 mg	554658
		C17.8	Rat IgG _{2a}	RUO	Biotin	0.5 mg	554476
IL-12 (p70), Recombinant mouse				RUO		5 µg	554592
IL-17A	Ms	TC11-18H10	Rat IgG ₁ , κ	RUO	Purified	0.1 mg	559501
			RUO	Purified	0.5 mg	555068	
			Rat IgG ₁	RUO	Biotin	0.5 mg	555067
MCP-1	Ms	2H5	Armenian Hamster IgG1, κ	RUO	Purified	1.0 mg	551217
		4E2/MCP	Armenian Hamster IgG1, λ 1	RUO	Biotin	0.5 mg	554444
MCP-1, Recombinant mouse				RUO		5 µg	554590
TNF	Ms, Rat	TN3-19.12	Armenian Hamster IgG1, λ 1	RUO	Purified	0.5 mg	557516
		C1150-14	Rabbit Ig	RUO	Biotin	0.5 mg	557432
		516D1A1	Human Ig	RUO	Biotin	0.5 mg	558415
		G281-2626	Rat IgG ₁	RUO	Purified	1.0 mg	551225
	Ms	MP6-XT3	Rat IgG ₁	RUO	Biotin	0.5 mg	554415
TNF, Recombinant mouse				RUO		10 µg	554589

ELISA Reagents

ELISA Pairs - Rat

DESCRIPTION	REACT	CLONE	ISOTYPE	REG	FORMAT	SIZE	CAT. NO.
IgA	Rat	A93-3	Mouse (BALB/c) IgG ₁ , κ	RUO	Purified	0.5 mg	553913
				RUO	Biotin	0.5 mg	553912
Rat IgA, κ		R3-30	Rat IgA, κ	RUO	Purified	0.5 mg	553945
IgE	Rat	B41-1	Mouse (BALB/c) IgG ₁ , κ	RUO	Purified	0.5 mg	553914
		B41-3	Mouse (BALB/c) IgG ₁ , κ	RUO	Biotin	0.5 mg	553916
IgG_{2a}	Rat	B46-7	Mouse (BALB/c) IgG ₁ , κ	RUO	Purified	0.5 mg	553918
		RG7/1.30	Mouse (SJL) IgG _{2b} , κ	RUO	Biotin	0.5 mg	553894
Rat IgG_{2a}, κ (anti-KLH)		A110-2	Rat (LOU) IgG _{2a} , κ	RUO	Purified	0.5 mg	553992
IgG_{2b}	Rat	G15-337	Mouse IgG _{2b} , κ	RUO	Purified	0.5 mg	553882
		RG7/11.1	Mouse (SJL) IgG _{2b} , κ	RUO	Biotin	0.5 mg	553898
Rat IgG_{2b}, κ		A95-1	Rat (LOU) IgG _{2b} , κ	RUO	Purified	0.5 mg	553986
IgG₂	Rat	A92-3	Mouse (BALB/c) IgG ₁ , κ	RUO	Purified	0.5 mg	553910
		A92-1	Mouse (BALB/c) IgG ₁ , κ	RUO	Biotin	0.5 mg	553909
Rat IgG₂, κ		A23-1	Rat (LOU) IgG ₂ , κ	RUO	Purified	0.5 mg	553982
IgM	Rat	G53-238	Mouse IgG ₁ , κ	RUO	Purified	0.5 mg	553885
				RUO	Biotin	0.5 mg	553886
Rat IgM, κ		R4-22	Rat IgM, κ	RUO	Purified	0.5 mg	553940
IL-4	Rat	OX-81	Mouse IgG ₁ , κ	RUO	Purified	0.5 mg	555080
		B11-3	Mouse IgG ₁	RUO	Biotin	0.5 mg	555090
IL-4, Recombinant rat				RUO		5 µg	555107
IL-6	Rat	C3-4	Mouse IgG ₁	RUO	Purified	0.5 mg	550644
		G307-2	Mouse IgG ₁	RUO	Biotin	0.5 mg	550642
IL-6, Recombinant rat				RUO		5 µg	557008
IL-10	Rat	A5-6	Mouse IgG ₁	RUO	Biotin	0.5 mg	555084
		A5-7	Mouse IgG ₁	RUO	Purified	0.5 mg	555083
IL-10, Recombinant rat				RUO		5 µg	555113
MCP-1	Rat	C4	Mouse IgG ₁	RUO	Purified	0.5 mg	555072
		B4	Mouse IgG ₁	RUO	Biotin	0.5 mg	555074
MCP-1, Recombinant rat				RUO		5 µg	555110
TNF	Ms, Rat	TN3-19.12	Armenian Hamster IgG1, λ 1	RUO	Purified	0.5 mg	557516
		C1150-14	Rabbit Ig	RUO	Biotin	0.5 mg	557432
		516D1A1	Human Ig	RUO	Biotin	0.5 mg	558415
TNF, Recombinant rat				RUO		5 µg	555109

ELISA Reagents

ELISA Pairs - Other Species

DESCRIPTION	REACT	CLONE	ISOTYPE	REG	FORMAT	SIZE	CAT. NO.
IgG1	Ar Ham, Syr Ham	HIG-632	Mouse (BALB/c) IgG _{2b} , κ	RUO	Purified	0.5 mg	550637
		G94-56	Mouse (BALB/c) IgG _{2b} , κ	RUO	Biotin	0.5 mg	554007
Hamster IgG1, k		A19-3	Armenian Hamster IgG1, κ	RUO	Purified	0.5 mg	553969
IgG2	Ar Ham	HIG-65	Mouse (BALB/c) IgG ₁ , κ	RUO	Purified	0.5 mg	550638
IgG	Ar Ham, Syr Ham	G192-1	Mouse (BALB/c) IgG ₁ , κ	RUO	Biotin	0.5 mg	554025
		B81-3	Armenian Hamster IgG2, κ	RUO	Purified	0.5 mg	559277
IgM	Ar Ham, Rat	G188-2	Mouse (BALB/c) IgG ₁ , κ	RUO	Purified	0.5 mg	554031
		G188-9	Mouse (BALB/c) IgG ₁ , κ	RUO	Biotin	0.5 mg	554035
Hamster IgM		G235-1	Armenian Hamster IgM, λ1	RUO	Purified	0.5 mg	553958
IFN-γ	Pig	P2C11	Mouse IgG _{2a}	RUO	Biotin	0.5 mg	559958
		P2G10	Mouse IgG ₁ , κ	RUO	Purified	0.5 mg	559961
IgM	Pig	F008-1629	Mouse (BALB/c) IgG ₁ , κ	RUO	Purified	0.5 mg	552551
		F008-2005	Mouse (BALB/c) IgG ₁ , κ	RUO	Biotin	0.2 mg	552550
IgG	Rab	C101-359	Mouse (BALB/c) IgG ₁ , κ	RUO	Purified	0.5 mg	550326
		C101-167	Mouse (BALB/c) IgG ₁ , κ	RUO	Biotin	0.5 mg	550346
Rabbit IgG		Polyclonal	Rabbit Ig	RUO	Purified	0.1 mg	550875
TNF	Rab	Polyclonal	Goat IgG	RUO	Purified	1.0 mg	551214
		23H1.1	Mouse IgG ₁ , κ	RUO	Biotin	0.5 mg	552470
TNF, Recombinant Rabbit				RUO		10 tests	558135
TNP	TNP	107.3	Mouse (BALB/c) IgG ₁ , κ	RUO	Purified	0.5 mg	554054
		49.2	Mouse (BALB/c) IgG _{2b} , κ	RUO	Biotin	0.5 mg	554055

ELISA Reagents

Ancillary Products

DESCRIPTION	REACT	CLONE	ISOTYPE	REG	FORMAT	SIZE	CAT. NO.
96-well ELISA Microplates, Flat-bottom				Class I Medical		100 plates	353279
Assay Diluent				RUO		500 mL	555213
BD OptEIA™ Reagent Set A (pH 6.5 buffer)				RUO	Coating Buffer (pH 6.5), Assay Diluent, Wash Concentrate, Substrate, and Stop Solution	Reagents for 20 plates	550536
BD OptEIA Reagent Set B (pH 9.5 buffer)				RUO	Coating Buffer (pH 9.5), Assay Diluent, Wash Concentrate, Substrate, and Stop Solution	Reagents for 20 plates	550534
Ig (human, mouse, and rat adsorbed)	Rab	Polyclonal	Ig	RUO	HRP	1.0 mL	554021
Ig (multiple adsorption)	Rat	Polyclonal	Goat Ig	RUO	Biotin	0.5 mg	554014
				RUO	HRP	1.0 mL	554017
Ig (multiple adsorption)	Ms	Polyclonal	Goat Ig	RUO	HRP	1.0 mL	554002
TMB Substrate Reagent Set				RUO	buffered solution/ organic solvent for BD OptEIA™	1 Set (300 mL each of 2 reagents)	555214

ELISPOT Reagents

BD ELISPOT Kits

DESCRIPTION	REG	SIZE	CAT. NO.
Human Granzyme B ELISPOT Kit	RUO	2 plates	552573
Human IFN- γ ELISPOT Kit	RUO	2 plates	552138
Mouse IFN- γ ELISPOT Kit	RUO	2 plates	552569

ELISPOT Reagents

BD ELISPOT Sets

DESCRIPTION	REG	SIZE	CAT. NO.
Human Granzyme B ELISPOT Set	RUO	10 plates	552572
Human IFN- γ ELISPOT Set	RUO	10 plates	551849
Human IL-2 ELISPOT Set	RUO	10 plates	551282
Human IL-4 ELISPOT Set	RUO	10 plates	551084
Human IL-5 ELISPOT Set	RUO	10 plates	551085
Human IL-10 ELISPOT Set	RUO	10 plates	551018
Human IL-12p70 ELISPOT Set	RUO	10 plates	552574
Human TNF ELISPOT Set	RUO	10 plates	551446
Mouse IFN- γ ELISPOT Set	RUO	10 plates	551083
Mouse IL-2 ELISPOT Set	RUO	10 plates	551076
Mouse IL-4 ELISPOT Set	RUO	10 plates	551017
Mouse IL-5 ELISPOT Set	RUO	10 plates	551075
Mouse IL-6 ELISPOT Set	RUO	10 plates	552567
Mouse IL-12p70 ELISPOT Set	RUO	10 plates	558671
Mouse TNF ELISPOT Set	RUO	10 plates	551491
Rat IFN- γ ELISPOT Set	RUO	10 plates	557819

ELISPOT Reagents

BD ELISPOT Pairs

DESCRIPTION	REG	SIZE	CAT. NO.
Human IFN- γ ELISPOT Pair	RUO	5 plates	551873
Human IL-2 ELISPOT Pair	RUO	5 plates	551884
Human IL-4 ELISPOT Pair	RUO	5 plates	551885
Human IL-5 ELISPOT Pair	RUO	5 plates	551886
Human IL-10 ELISPOT Pair	RUO	5 plates	551883
Human TNF ELISPOT Pair	RUO	5 plates	551882
Mouse IFN- γ ELISPOT Pair	RUO	5 plates	551881
Mouse IL-2 ELISPOT Pair	RUO	5 plates	551876
Mouse IL-4 ELISPOT Pair	RUO	5 plates	551878
Mouse IL-5 ELISPOT Pair	RUO	5 plates	551880
Mouse IL-10 ELISPOT Pair	RUO	5 plates	557982
Mouse TNF ELISPOT Pair	RUO	5 plates	551875

ELISPOT Reagents

Ancillary reagents

DESCRIPTION	REG	SIZE	CAT. NO.
AEC Substrate Set	RUO	1 set	551951
Streptavidin-HRP ELISPOT	RUO	1.0 mL	557630

