



# Matched Antibody Pairs for ELISA

October 2017

Analyte	Reagent	Catalog #	Size
Human 2B4/CD244/SLAMF4	Capture	AF1039	100 µg
	Detection (Biotin)	BAF1039	50 µg
	Protein	1039-2B-050	50 µg
Human 4-1BB/TNFRSF9/CD137	Capture	AF838	100 µg
	Detection (Biotin)	BAF838	50 µg
	Protein	838-4B-100	100 µg
Mouse 4-1BB/TNFRSF9/CD137	Capture	MAB937	500 µg
	Detection (Biotin)	BAF937	50 µg
	Protein	937-4B-050	50 µg
Mouse ACE/CD143	Capture	MAB15132	500 µg
	Detection (Biotin)	BAF1513	50 µg
	Protein	1513-ZN-010	10 µg
Human/Mouse/Rat Activin A	Capture	MAB3381-500	500 µg
	Detection (Biotin)	BAM3381	100 µg
Human ADAM9	Capture	AF939	100 µg
	Detection (Biotin)	BAF939	50 µg
	Protein	939-AD-020	20 µg
Human Adiponectin/Acrp30	Capture	MAB10651-50	500 µg
	Capture	MAB10651R-1	100 µg
	Detection (Biotin)	BAM1065	250 µg
	Protein	1065-AP-050	50 µg
Rat Agrin	Capture	AF550	100 µg
	Detection (Biotin)	BAF550	50 µg
	Protein	550-AG-100	100 µg
Human AgRP/ART	Capture	MAB8312	500 µg
	Detection (Biotin)	BAF704	50 µg
	Protein	704-AG-002	2 µg
Human ALCAM/CD166	Capture	MAB6561-500	500 µg
	Detection (Biotin)	BAF656	50 µg
	Protein	656-AL-100	100 µg
Mouse ALK-1	Capture	AF770	100 µg
	Detection (Biotin)	BAF770	50 µg
	Protein	770-MA-100	100 µg
Human AMICA	Capture	AF3449	100 µg
	Detection (Biotin)	BAF3449	50 µg
	Protein	3449-AM-050	50 µg

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Human Amphiregulin	Capture	MAB262-500	500 µg
	Detection (Biotin)	BAF262	50 µg
	Protein	262-AR-100	100 µg
Mouse Amphiregulin	Capture	AF989	100 µg
	Detection (Biotin)	BAF989	50 µg
	Protein	989-AR-100	100 µg
Human Angiogenin	Capture	MAB265-500	500 µg
	Detection (Biotin)	BAF265	50 µg
	Protein	265-AN-050	50 µg
Human Angiopoietin-1	Capture	MAB9231-500	500 µg
	Detection (Biotin)	BAF923	50 µg
	Protein	923-AN-025	25 µg
Human Angiopoietin-2	Capture	MAB098-500	500 µg
	Detection (Biotin)	BAM0981	250 µg
	Protein	623-AN-025	25 µg
Human Angiopoietin-4	Capture	MAB9641	500 µg
	Detection (Biotin)	BAF964	50 µg
	Protein	964-AN-025	25 µg
Human Angiopoietin-like Protein 3/ANGPTL3	Capture	MAB38291	500 µg
	Detection (Biotin)	BAF3829	50 µg
	Protein	3829-AN-050	50 µg
Mouse Angiopoietin-like Protein 3/ANGPTL3	Capture	MAB13621	500 µg
	Detection (Biotin)	BAF136	50 µg
	Protein	136-AN-050	50 µg
Human/Primate Angiopoietin-like Protein 4/ANGPTL4	Capture	AF3485	100 µg
	Detection (Biotin)	BAF3485	50 µg
	Protein	4487-AN-050	50 µg
Human APRIL/TNFSF13	Capture	MAB8842	500 µg
	Detection (Biotin)	BAM8841	250 µg
	Protein	884-AP-010	10 µg
Human Axl	Capture	MAB154-500	500 µg
	Capture	MAB154R-100	100 µg
	Detection (Biotin)	BAF154	50 µg
	Protein	154-AL-100	100 µg
Mouse Axl	Capture	MAB854	500 µg
	Detection (Biotin)	BAF854	50 µg
	Protein	854-AX-100	100 µg
Human B7-1/CD80	Capture	MAB140-500	500 µg
	Detection (Biotin)	BAM1402	100 µg
	Protein	140-B1-100	100 µg
Mouse B7-1/CD80	Capture	AF740	100 µg
	Detection (Biotin)	BAF740	50 µg
	Protein	740-B1-100	100 µg
Rat B7-2/CD86	Capture	AF1340	100 µg
	Detection (Biotin)	BAF1340	50 µg
	Protein	1340-B2-050	50 µg
Mouse BAFF R/TNFRSF13C	Capture	AF1357	100 µg
	Detection (Biotin)	BAF1357	50 µg
	Protein	1357-BR-050	50 µg

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Human BCAM	Capture	MAB1481	500 µg
	Detection (Biotin)	BAF148	50 µg
	Protein	148-BC-100	100 µg
Human/Primate BCMA/TNFRSF17	Capture	AF193	100 µg
	Detection (Biotin)	BAF193	50 µg
	Protein	193-BC-050	50 µg
Human BDNF	Capture	MAB848-500	500 µg
	Detection (Biotin)	BAM648	50 µg
	Protein	248-BD-005	5 µg
Human Betacellulin/BTC	Capture	MAB261-500	500 µg
	Detection (Biotin)	BAF261	50 µg
	Protein	261-CE-010	10 µg
Mouse Betacellulin/BTC	Capture	MAB10251	500 µg
	Detection (Biotin)	BAF1025	50 µg
	Protein	1025-CE-025	25 µg
Human βIG-H3	Capture	MAB2935	500 µg
	Detection (Biotin)	BAF2935	50 µg
	Protein	3409-BG-050	50 µg
Mouse βIG-H3	Capture	AF2559	100 µg
	Detection (Biotin)	BAF2559	50 µg
	Protein	2559-BG-050	50 µg
Human BLAME/SLAMF8	Capture	MAB19073	500 µg
	Detection (Biotin)	BAM19074	250 µg
	Protein	1907-BL-050	50 µg
Human BMP-2	Capture	MAB3551-500	500 µg
	Detection (Biotin)	BAM3552	250 µg
	Protein	355-BM-010	10 µg
Human/Primate BMP-4	Capture	MAB7571-500	500 µg
	Detection (Biotin)	BAM7572	250 µg
	Protein	314-BP-010	10 µg
Human BMP-6	Capture	MAB507-500	500 µg
	Detection (Biotin)	BAF507	50 µg
	Protein	507-BP-020	20 µg
Human BMP-7	Capture	MAB3542-500	500 µg
	Detection (Biotin)	BAM354	250 µg
	Protein	354-BP-010	10 µg
Human/Primate BMP-9	Capture	MAB3209-500	500 µg
	Detection (Biotin)	BAF3209	50 µg
	Protein	3209-BP-010	10 µg
Mouse E-Cadherin	Capture	MAB7482	500 µg
	Detection (Biotin)	BAF748	50 µg
	Protein	748-EC-050	50 µg
Human P-Cadherin	Capture	MAB861-500	500 µg
	Detection (Biotin)	BAF861	50 µg
	Protein	861-PC-100	100 µg
Mouse P-Cadherin	Capture	AF761	100 µg
	Detection (Biotin)	BAF761	50 µg
	Protein	761-MP-050	50 µg
Human Carbonic Anhydrase IX	Capture	MAB2188-500	500 µg
	Detection (Biotin)	BAF2188	50 µg
	Protein	2188-CA-010	10 µg

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Mouse Cardiotrophin-1/CT-1	Capture	MAB438	500 µg
	Detection (Biotin)	BAF438	50 µg
	Protein	438-CT-010	10 µg
Human Cathepsin B	Capture	MAB2176	500 µg
	Detection (Biotin)	BAF953	50 µg
	Protein	953-CY-010	10 µg
Human Pro-Cathepsin B	Capture	MAB2176	500 µg
	Detection (Biotin)	BAM21761	250 µg
	Protein	953-CY-010	10 µg
Human Cathepsin L	Capture	AF952	100 µg
	Detection (Biotin)	BAF952	50 µg
	Protein	952-CY-010	10 µg
Human Cathepsin S	Capture	AF1183	100 µg
	Detection (Biotin)	BAF1183	50 µg
	Protein	1183-CY-010	10 µg
Human CCL1/I-309	Capture	MAB272-500	500 µg
	Detection (Biotin)	BAF272	50 µg
	Protein	272-I-010	10 µg
Mouse CCL1/TCA-3	Capture	MAB8451	500 µg
	Detection (Biotin)	BAF845	50 µg
	Protein	845-TC-025	25 µg
Canine CCL2/JE/MCP-1	Capture	MAB28171	500 µg
	Detection (Biotin)	BAF1774	50 µg
	Protein	1774-MC-025	25 µg
Human CCL2/JE/MCP-1	Capture	MAB679-500	500 µg
	Detection (Biotin)	BAF279	50 µg
	Protein	279-MC-010	10 µg
Mouse CCL2/JE/MCP-1	Capture	AF-479-NA	100 µg
	Detection (Biotin)	BAF479	50 µg
	Protein	479-JE-010	10 µg
Human CCL3/MIP-1α	Capture	AF-270-NA	100 µg
	Detection (Biotin)	BAF270	50 µg
	Protein	270-LD-010	10 µg
Mouse CCL3/MIP-1α	Capture	AF-450-NA	100 µg
	Detection (Biotin)	BAF450	50 µg
	Protein	450-MA-010	10 µg
Human CCL4/MIP-1β	Capture	MAB271-100	100 µg
	Detection (Biotin)	BAF271	50 µg
	Protein	271-BME-010	10 µg
Feline CCL5/RANTES	Capture	AF3819	100 µg
	Detection (Biotin)	BAF3819	50 µg
	Protein	3819-FR-025	25 µg
Human/Primate CCL5/RANTES	Capture	MAB678-500	500 µg
	Detection (Biotin)	BAF278	50 µg
	Protein	278-RN-010	10 µg
Mouse CCL5/RANTES	Capture	MAB4781-500	500 µg
	Detection (Biotin)	BAF478	50 µg
	Protein	478-MR-025	25 µg
Mouse CCL6/C10	Capture	MAB4871	500 µg
	Detection (Biotin)	BAM4872	250 µg
	Protein	487-C-010	10 µg

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Human CCL7/MCP-3	Capture	MAB282-500	500 µg
	Detection (Biotin)	BAF282	50 µg
	Protein	282-P3-010	10 µg
Human CCL8/MCP-2	Capture	MAB281-500	500 µg
	Detection (Biotin)	BAF281	50 µg
	Protein	281-CP-010	10 µg
Mouse CCL9/10/MIP-1γ	Capture	MAB463-500	500 µg
	Detection (Biotin)	BAF463	50 µg
	Protein	463-MG-025	25 µg
Human CCL11/Eotaxin	Capture	MAB320-500	500 µg
	Detection (Biotin)	BAF320	50 µg
	Protein	320-EO-020	20 µg
Mouse CCL11/Eotaxin	Capture	AF-420-NA	100 µg
	Detection (Biotin)	BAF420	50 µg
	Protein	420-ME-020	20 µg
Mouse CCL12/MCP-5	Capture	MAB428-500	500 µg
	Detection (Biotin)	BAF428	50 µg
	Protein	428-P5-025	25 µg
Human CCL13/MCP-4	Capture	MAB327	500 µg
	Detection (Biotin)	BAF327	50 µg
	Protein	327-P4-025	25 µg
Human CCL14a/HCC-1	Capture	MAB3241	500 µg
	Capture	MAB3241R-10	100 µg
	Detection (Biotin)	BAF324	50 µg
	Protein	324-HC-010	10 µg
Human CCL15/MIP-1δ	Capture	MAB363	500 µg
	Detection (Biotin)	BAF363	50 µg
	Protein	363-MG-025	25 µg
Human CCL16/HCC-4	Capture	MAB328	500 µg
	Detection (Biotin)	BAF802	50 µg
	Protein	802-HC-025	25 µg
Human CCL17/TARC	Capture	MAB364-500	500 µg
	Detection (Biotin)	BAF364	50 µg
	Protein	364-DN-025	25 µg
Mouse CCL17/TARC	Capture	MAB529-500	500 µg
	Detection (Biotin)	BAF529	50 µg
	Protein	529-TR-025	25 µg
Human CCL18/PARC	Capture	MAB394-500	500 µg
	Detection (Biotin)	BAF394	50 µg
	Protein	394-PA-010	10 µg
Human CCL19/MIP-3β	Capture	AF361	100 µg
	Detection (Biotin)	BAF361	50 µg
	Protein	361-MI-025	25 µg
Mouse CCL19/MIP-3β	Capture	MAB880-500	500 µg
	Detection (Biotin)	BAF880	50 µg
	Protein	440-M3-025	25 µg
Human CCL20/MIP-3α	Capture	MAB360-500	500 µg
	Detection (Biotin)	BAF360	50 µg
	Protein	360-MP-025	25 µg
Mouse CCL20/MIP-3α	Capture	MAB7601-500	500 µg
	Detection (Biotin)	BAF760	50 µg
	Protein	760-M3-025	25 µg

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Rat CCL20/MIP-3 $\alpha$	Capture	MAB540	500 $\mu$ g
	Detection (Biotin)	BAF540	50 $\mu$ g
	Protein	540-RM-025	25 $\mu$ g
Human CCL21/6Ckine	Capture	MAB3661	500 $\mu$ g
	Detection (Biotin)	BAF366	50 $\mu$ g
	Protein	366-6C-025	25 $\mu$ g
Mouse CCL21/6Ckine	Capture	MAB457-500	500 $\mu$ g
	Detection (Biotin)	BAF457	50 $\mu$ g
	Protein	457-6C-025	25 $\mu$ g
Human CCL22/MDC	Capture	MAB336-500	500 $\mu$ g
	Detection (Biotin)	BAF336	50 $\mu$ g
	Protein	336-MD-025	25 $\mu$ g
Mouse CCL22/MDC	Capture	MAB4391-500	500 $\mu$ g
	Detection (Biotin)	BAF439	50 $\mu$ g
	Protein	439-MD-025	25 $\mu$ g
Human CCL23/Ck $\beta$ 8-1	Capture	MAB508	500 $\mu$ g
	Detection (Biotin)	BAF508	50 $\mu$ g
	Protein	508-CK-025	25 $\mu$ g
Human CCL23/MPIF-1	Capture	MAB3711	500 $\mu$ g
	Detection (Biotin)	BAF371	50 $\mu$ g
	Protein	371-MP-025	25 $\mu$ g
Human CCL24/Eotaxin-2/MPIF-2	Capture	MAB343-500	500 $\mu$ g
	Detection (Biotin)	BAF343	50 $\mu$ g
	Protein	343-E2-025	25 $\mu$ g
Mouse CCL24/Eotaxin-2/MPIF-2	Capture	MAB528-500	500 $\mu$ g
	Detection (Biotin)	BAF528	50 $\mu$ g
	Protein	528-MP-025	25 $\mu$ g
Human CCL25/TECK	Capture	MAB3341	500 $\mu$ g
	Detection (Biotin)	BAF334	50 $\mu$ g
	Protein	9046-TK-025	25 $\mu$ g
Mouse CCL25/TECK	Capture	MAB4811	500 $\mu$ g
	Detection (Biotin)	BAF481	50 $\mu$ g
	Protein	481-TK-025	25 $\mu$ g
Human CCL26/Eotaxin-3	Capture	MAB653-500	500 $\mu$ g
	Detection (Biotin)	BAF346	50 $\mu$ g
	Protein	653-E3-025	25 $\mu$ g
Human CCL27/CTACK	Capture	MAB3761	500 $\mu$ g
	Detection (Biotin)	BAF376	50 $\mu$ g
	Protein	376-CT-025	25 $\mu$ g
Mouse CCL27/CTACK	Capture	MAB725	500 $\mu$ g
	Detection (Biotin)	BAF725	50 $\mu$ g
	Protein	725-CK-025	25 $\mu$ g
Human CCL28	Capture	MAB7171	500 $\mu$ g
	Detection (Biotin)	BAF717	50 $\mu$ g
	Protein	717-VC-025	25 $\mu$ g
Mouse CCL28	Capture	MAB5331	500 $\mu$ g
	Detection (Biotin)	BAF533	50 $\mu$ g
	Protein	533-VI-025	25 $\mu$ g
Mouse CD6	Capture	MAB7271	500 $\mu$ g
	Detection (Biotin)	BAM727	250 $\mu$ g
	Protein	727-CD-100	100 $\mu$ g

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Human CD14	Capture	MAB3833-500	500 µg
	Detection (Biotin)	BAF383	50 µg
	Protein	383-CD-050	50 µg
Human CD23/Fcε RII	Capture	MAB1231	500 µg
	Detection (Biotin)	BAF123	50 µg
	Protein	123-FE-050	50 µg
Mouse CD27/TNFRSF7	Capture	MAB5741	500 µg
	Detection (Biotin)	BAM5741	250 µg
	Protein	574-CD-050	50 µg
Mouse CD27 Ligand/TNFSF7	Capture	MAB783	500 µg
	Detection (Biotin)	BAF783	50 µg
	Protein	783-CL-050	50 µg
Mouse CD30/TNFRSF8	Capture	MAB8521	500 µg
	Detection (Biotin)	BAF852	50 µg
	Protein	852-CD-100	100 µg
Mouse CD30 Ligand/TNFSF8	Capture	MAB732	500 µg
	Detection (Biotin)	BAF732	50 µg
	Protein	732-CL-050	50 µg
Mouse CD36/SR-B3	Capture	AF2519	100 µg
	Detection (Biotin)	BAF2519	50 µg
	Protein	2519-CD-050	50 µg
Mouse CD40/TNFRSF5	Capture	MAB4401	500 µg
	Detection (Biotin)	BAF440	50 µg
	Protein	1215-CD-050	50 µg
Mouse CD48/SLAMF2	Capture	AF3327	100 µg
	Detection (Biotin)	BAF3327	50 µg
	Protein	3327-CD-050	50 µg
Human CD97	Capture	MAB25291	500 µg
	Detection (Biotin)	BAF2529	50 µg
	Protein	2529-CD-050	50 µg
Human CD117/c-kit	Capture	MAB332-500	500 µg
	Detection (Biotin)	BAF332	50 µg
	Protein	332-SR-050	50 µg
Human CD163	Capture	MAB16071-50	500 µg
	Detection (Biotin)	BAM16072	250 µg
	Protein	1607-CD-050	50 µg
Human CD200	Capture	MAB27243-50	500 µg
	Detection (Biotin)	BAM27242	250 µg
	Protein	2724-CD-050	50 µg
Human CD25/IL-2 Rα	Capture	MAB623-500	500 µg
	Detection (Biotin)	BAF223	50 µg
	Protein	223-2A-005	5 µg
Mouse CD25/IL-2 Rα	Capture	AF2438	100 µg
	Detection (Biotin)	BAF2438	50 µg
	Protein	2438-RM-050	50 µg
Human CEACAM-1/CD66a	Capture	MAB22441-50	500 µg
	Detection (Biotin)	BAF2244	50 µg
	Protein	2244-CM-050	50 µg
Human Chemerin	Capture	MAB23241-50	500 µg
	Detection (Biotin)	BAF2324	50 µg
	Protein	2324-CM-025	25 µg

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Mouse Chemerin	Capture	MAB23251	500 µg
	Detection (Biotin)	BAM2325	250 µg
	Protein	2325-CM-025	25 µg
Human/Primate Chitinase 3-like 1	Capture	MAB25991-50	500 µg
	Detection (Biotin)	BAF2599	50 µg
	Protein	2599-CH-050	50 µg
Mouse Chordin	Capture	MAB758	500 µg
	Detection (Biotin)	BAF758	50 µg
	Protein	758-CN-050	50 µg
Human Clusterin	Capture	MAB29372-50	500 µg
	Detection (Biotin)	BAM29373	250 µg
	Protein	2937-HS-050	50 µg
Mouse Clusterin	Capture	AF2747	100 µg
	Detection (Biotin)	BAF2747	50 µg
	Protein	2747-HS-050	50 µg
Human CNTF	Capture	MAB657	500 µg
	Detection (Biotin)	BAF257	50 µg
	Protein	257-NT-010	10 µg
Rat CNTF	Capture	MAB557	500 µg
	Detection (Biotin)	BAF557	50 µg
	Protein	557-NT-010	10 µg
Human Coagulation Factor III/Tissue Factor	Capture	MAB2339-500	500 µg
	Detection (Biotin)	BAF2339	50 µg
	Protein	2339-PA-010	10 µg
Human COCO	Capture	MAB30471	500 µg
	Detection (Biotin)	BAM3047	250 µg
	Protein	3047-CC-050	50 µg
Human Complement Component C5/C5a	Capture	MAB2037-500	500 µg
	Detection (Biotin)	BAM20371	250 µg
	Protein	2037-C5-025	25 µg
Mouse Complement Component C5a	Capture	MAB21501-50	500 µg
	Detection (Biotin)	BAF2150	50 µg
	Protein	2150-C5-025	25 µg
Human Complement Factor D/Adipsin	Capture	MAB18241-50	500 µg
	Detection (Biotin)	BAF1824	50 µg
	Protein	1824-SE-010	10 µg
Human C-Reactive Protein/CRP	Capture	MAB17071-50	500 µg
	Detection (Biotin)	BAM17072	50 µg
	Protein	1707-CR-200	200 µg
Mouse C-Reactive Protein/CRP	Capture	MAB1829	500 µg
	Detection (Biotin)	BAF1829	50 µg
	Protein	1829-CR-200	200 µg
Rat C-Reactive Protein/CRP	Capture	MAB17441	500 µg
	Detection (Biotin)	BAM17442	250 µg
	Protein	1744-CR-200	200 µg
Human/Primate Cripto-1	Capture	MAB2772-500	500 µg
	Detection (Biotin)	BAM2773	250 µg
	Protein	145-CR-010	10 µg
Human CRTAM	Capture	AF1695	100 µg
	Detection (Biotin)	BAF1695	50 µg
	Protein	1695-CR-050	50 µg

For research use only. Not for use in diagnostic procedures.



Analyte	Reagent	Catalog #	Size
Mouse CTLA-4	Capture	AF476	100 µg
	Detection (Biotin)	BAF476	50 µg
	Protein	434-CT-200	200 µg
Human CX3CL1/Fractalkine	Capture	MAB3652-500	500 µg
	Detection (Biotin)	BAM365	100 µg
	Protein	365-FR-025	25 µg
Mouse CX3CL1/Fractalkine	Capture	MAB571-500	500 µg
	Detection (Biotin)	BAF472	50 µg
	Protein	472-FF-025	25 µg
Rat CX3CL1/Fractalkine	Capture	AF537	100 µg
	Detection (Biotin)	BAF537	50 µg
	Protein	536-FR-025	25 µg
Human/Primate CXCL1/GRO $\alpha$	Capture	MAB275-500	500 µg
	Capture	MAB275R-100	100 µg
	Detection (Biotin)	BAF275	50 µg
	Protein	275-GR-010	10 µg
Mouse CXCL1/KC	Capture	MAB453-500	500 µg
	Detection (Biotin)	BAF453	50 µg
	Protein	453-KC-010	10 µg
Mouse CXCL2/MIP-2	Capture	MAB452-500	500 µg
	Detection (Biotin)	BAF452	50 µg
	Protein	452-M2-010	10 µg
Rat CXCL2/CINC-3	Capture	MAB525	500 µg
	Detection (Biotin)	BAF525	50 µg
	Protein	525-C3-025	25 µg
Rat CXCL3/CINC-2 $\alpha$ / $\beta$	Capture	MAB516	500 µg
	Detection (Biotin)	BAF516	50 µg
	Protein	516-CA-025	25 µg
	Protein	517-CB-025	25 µg
Human CXCL4/PF4	Capture	MAB7951-500	500 µg
	Detection (Biotin)	BAF795	50 µg
	Protein	795-P4-025	25 µg
Mouse CXCL4/PF4	Capture	MAB595-500	500 µg
	Detection (Biotin)	BAF595	50 µg
	Protein	595-P4-025	25 µg
Human CXCL5/ENA-78	Capture	MAB654-500	500 µg
	Detection (Biotin)	BAF254	50 µg
	Protein	254-XB-025	25 µg
Human CXCL6/GCP-2	Capture	MAB333	500 µg
	Detection (Biotin)	BAF333	50 µg
	Protein	333-GC-025	25 µg
Human CXCL7/NAP-2	Capture	MAB393	500 µg
	Detection (Biotin)	BAF393	50 µg
	Protein	393-NP-010	10 µg
Mouse CXCL7/Thymus Chemokine-1	Capture	MAB10911	500 µg
	Detection (Biotin)	BAF793	50 µg
	Protein	1091-CK-025	25 µg
Human CXCL9/MIG	Capture	MAB392-500	500 µg
	Detection (Biotin)	BAF392	50 µg
	Protein	392-MG-010	10 µg
Mouse CXCL9/MIG	Capture	AF-492-NA	100 µg
	Detection (Biotin)	BAF492	50 µg
	Protein	492-MM-010	10 µg

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Human CXCL10/IP-10	Capture	MAB266-500	500 µg
	Detection (Biotin)	BAF266	50 µg
	Protein	266-IP-010	10 µg
Mouse CXCL10/IP-10/CRG-2	Capture	MAB466-500	500 µg
	Detection (Biotin)	BAF466	50 µg
	Protein	466-CR-010	10 µg
Human CXCL11/I-TAC	Capture	MAB672-500	500 µg
	Detection (Biotin)	BAF672	50 µg
	Protein	672-IT-025	25 µg
Mouse CXCL11/I-TAC	Capture	AF572	100 µg
	Detection (Biotin)	BAF572	50 µg
	Protein	572-MC-025	25 µg
Human CXCL12/SDF-1 $\beta$	Capture	MAB350-500	500 µg
	Detection (Biotin)	BAF351	50 µg
	Protein	351-FS-010	10 µg
Human/Mouse CXCL12/SDF-1	Capture	MAB350-500	500 µg
	Detection (Biotin)	BAF310	50 µg
	Protein	350-NS-010	10 µg
	Protein	460-SD-010	10 µg
Human CXCL13/BLC/BCA-1	Capture	MAB801-500	500 µg
	Detection (Biotin)	BAF801	50 µg
	Protein	801-CX-025	25 µg
Mouse CXCL13/BLC/BCA-1	Capture	MAB470-500	500 µg
	Capture	MAB470R-100	100 µg
	Detection (Biotin)	BAF470	50 µg
	Protein	470-BC-025	25 µg
Human CXCL14/BRAK	Capture	MAB866	500 µg
	Detection (Biotin)	BAM8661	50 µg
	Protein	866-CX-025	25 µg
Mouse CXCL15/Lungkine	Capture	MAB442	500 µg
	Detection (Biotin)	BAF442	50 µg
	Protein	442-LK-002	2 µg
Human/Primate CXCL16	Capture	AF976	100 µg
	Detection (Biotin)	BAF976	50 µg
	Protein	1164-CX-025	25 µg
Mouse CXCL16	Capture	MAB503-500	500 µg
	Detection (Biotin)	BAF503	50 µg
	Protein	503-CX-025	25 µg
Human Cystatin C	Capture	MAB11962-50	500 µg
	Detection (Biotin)	BAM11961	100 µg
	Protein	1196-PI-010	10 µg
Mouse Cystatin C	Capture	AF1238	100 µg
	Detection (Biotin)	BAF1238	50 µg
	Protein	1238-PI-010	10 µg
Human DAN	Capture	MAB955	500 µg
	Detection (Biotin)	BAF955	50 µg
	Protein	955-DA-050	50 µg
Mouse DAN	Capture	MAB7552	500 µg
	Detection (Biotin)	BAF755	50 µg
	Protein	755-DA-050	50 µg
Human DcR3/TNFRSF6B	Capture	MAB21351	500 µg
	Detection (Biotin)	BAM2135	250 µg
	Protein	142-DC-100	100 µg

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Human Decorin	Capture	MAB1432	500 µg
	Detection (Biotin)	BAM1431	100 µg
	Protein	143-DE-100	100 µg
Mouse Decorin	Capture	AF1060	100 µg
	Detection (Biotin)	BAF1060	50 µg
Human Dkk-1	Capture	MAB10962-50	500 µg
	Detection (Biotin)	BAF1096	50 µg
	Protein	5439-DK-010	10 µg
Mouse Dkk-1	Capture	AF1765	100 µg
	Detection (Biotin)	BAF1765	50 µg
	Protein	1765-DK-010	10 µg
Human Dkk-3	Capture	MAB11181	500 µg
	Detection (Biotin)	BAF1118	50 µg
	Protein	1118-DK-050	50 µg
Human Dkk-4	Capture	MAB1269-500	500 µg
	Detection (Biotin)	BAF1269	50 µg
	Protein	1269-DK-010	10 µg
Human DLL1	Capture	MAB1818-500	500 µg
	Detection (Biotin)	BAF1818	50 µg
	Protein	1818-DL-050	50 µg
Mouse DLL4	Capture	MAB13891	500 µg
	Detection (Biotin)	BAM13892	250 µg
	Protein	1389-D4-050	50 µg
Human DMP-1	Capture	MAB4129	500 µg
	Detection (Biotin)	BAM41291	250 µg
	Protein	4129-DM-050	50 µg
Human DPPIV/CD26	Capture	MAB1180-500	500 µg
	Detection (Biotin)	BAF1180	50 µg
	Protein	1180-SE-010	10 µg
Mouse DPPIV/CD26	Capture	MAB954	500 µg
	Detection (Biotin)	BAF954	50 µg
	Protein	954-SE-010	10 µg
Human DR6/TNFRSF21	Capture	AF144	100 µg
	Detection (Biotin)	BAF144	50 µg
	Protein	144-DR-100	100 µg
Human Dtk	Capture	MAB859-500	500 µg
	Detection (Biotin)	BAF859	50 µg
	Protein	859-DK-100	100 µg
Mouse Dtk	Capture	MAB759	500 µg
	Detection (Biotin)	BAF759	50 µg
	Protein	759-DT-100	100 µg
Human EDA2R/TNFRSF27/XEDAR	Capture	MAB10931	500 µg
	Detection (Biotin)	BAF1093	50 µg
	Protein	1093-XD-050	50 µg
Human EDA-A2/Ectodysplasin A2	Capture	MAB9222	500 µg
	Detection (Biotin)	BAM9221	100 µg
	Protein	922-ED-010	10 µg
Mouse EDAR	Capture	MAB7451	500 µg
	Detection (Biotin)	BAM7452	250 µg
	Protein	745-ED-050	50 µg
Human/Primate EGF	Capture	MAB636-500	500 µg
	Detection (Biotin)	BAF236	50 µg
	Protein	236-EG-200	200 µg

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Mouse EGF	Capture	MAB2028-500	500 µg
	Detection (Biotin)	BAF2028	50 µg
	Protein	2028-EG-200	200 µg
Rat EGF	Capture	AF3214	100 µg
	Detection (Biotin)	BAF3214	50 µg
	Protein	3214-EG-100	100 µg
Human EGFR	Capture	AF231	100 µg
	Detection (Biotin)	BAF231	50 µg
	Protein	1095-ER-002	2 µg
Human EG-VEGF/PK1	Capture	MAB12091	500 µg
	Detection (Biotin)	BAF1209	50 µg
	Protein	1209-EV-025	25 µg
Mouse Endocan/ESM-1	Capture	MAB1999	500 µg
	Detection (Biotin)	BAF1999	50 µg
	Protein	1999-EC-050	50 µg
Human Endoglin/CD105	Capture	MAB10972-50	500 µg
	Detection (Biotin)	BAF1097	50 µg
	Protein	1097-EN-025	25 µg
Mouse Endoglin/CD105	Capture	MAB13201-50	500 µg
	Detection (Biotin)	BAF1320	50 µg
	Protein	1320-EN-025	25 µg
Human EpCAM/TROP-1	Capture	MAB9601-500	500 µg
	Detection (Biotin)	BAF960	50 µg
	Protein	960-EP-050	50 µg
Mouse Epigen	Capture	MAB11271	500 µg
	Detection (Biotin)	BAF1127	50 µg
	Protein	1127-EP-025	25 µg
Mouse Epiregulin	Capture	MAB1068	500 µg
	Detection (Biotin)	BAF1068	50 µg
	Protein	1068-EP-050	50 µg
Human ErbB2/Her2	Capture	MAB1129-500	500 µg
	Detection (Biotin)	BAF1129	50 µg
	Protein	1129-ER-050	50 µg
Human ErbB3/Her3	Capture	MAB3481-500	500 µg
	Detection (Biotin)	BAM348	250 µg
	Protein	348-RB-050	50 µg
Mouse Erythropoietin	Capture	MAB9591-500	500 µg
	Detection (Biotin)	BAM9592	250 µg
	Protein	959-ME-010	10 µg
Human Erythropoietin R	Capture	MAB3071-500	500 µg
	Detection (Biotin)	BAM3072	250 µg
	Protein	307-ER-050	50 µg
Human FABP2	Capture	MAB30781-50	500 µg
	Detection (Biotin)	BAF3078	50 µg
	Protein	2694-CL-100/C	100 µg
Human Fas/TNFRSF6/CD95	Capture	MAB144	500 µg
	Detection (Biotin)	BAF326	50 µg
	Protein	326-FS-050	50 µg
Mouse Fas/TNFRSF6/CD95	Capture	AF435	100 µg
	Detection (Biotin)	BAF435	50 µg
	Protein	435-FA-050	50 µg

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Human Fas Ligand/TNFSF6	Capture	MAB126-500	500 µg
	Detection (Biotin)	BAF126	50 µg
	Protein	126-FL-010	10 µg
Mouse Fas Ligand/TNFSF6	Capture	MAB5262	500 µg
	Detection (Biotin)	BAF526	50 µg
	Protein	526-SA-050	50 µg
Mouse Fcγ RIIA/CD32a	Capture	MAB14601-50	500 µg
	Detection (Biotin)	BAF1460	50 µg
	Protein	1460-CD-050	50 µg
Human Fcγ RIIb/CD32b	Capture	MAB18751-50	500 µg
	Detection (Biotin)	BAM1875	250 µg
	Protein	1875-CD-050	50 µg
Human Fetuin A/AHSG	Capture	MAB11841	500 µg
	Detection (Biotin)	BAF1184	50 µg
	Protein	1184-PI-050	50 µg
Mouse Fetuin A/AHSG	Capture	MAB15632	500 µg
	Detection (Biotin)	BAM15631	250 µg
	Protein	1563-PI-050	50 µg
Human Fetuin B	Capture	MAB17251	500 µg
	Detection (Biotin)	BAM17251	250 µg
	Protein	1725-PI-050	50 µg
Human FGF basic	Capture	MAB233-500	500 µg
	Detection (Biotin)	BAM233	250 µg
	Protein	233-FB-025	25 µg
Human FGF-4	Capture	MAB635	500 µg
	Detection (Biotin)	BAF235	50 µg
	Protein	235-F4-025	25 µg
Human FGF-6	Capture	MAB238	500 µg
	Detection (Biotin)	BAF238	50 µg
	Protein	238-F6-025	25 µg
Human KGF/FGF-7	Capture	MAB251-500	500 µg
	Detection (Biotin)	BAF251	50 µg
	Protein	251-KG-010	10 µg
Human FGF-9	Capture	MAB273-500	500 µg
	Detection (Biotin)	BAF273	50 µg
	Protein	273-F9-025	25 µg
Human FGF-19	Capture	MAB9691-500	500 µg
	Detection (Biotin)	BAF969	50 µg
	Protein	969-FG-025	25 µg
Human Fibroblast Activation Protein α/FAP	Capture	MAB3715-500	500 µg
	Detection (Biotin)	BAF3715	50 µg
	Protein	3715-SE-010	10 µg
Human Flt-3 Ligand	Capture	MAB608-500	500 µg
	Detection (Biotin)	BAF308	50 µg
	Protein	308-FKN-005	5 µg
Mouse Flt-3 Ligand	Capture	AF427	100 µg
	Detection (Biotin)	BAF427	50 µg
	Protein	427-FL-005	5 µg
Human Follistatin-related Gene Protein/FLRG	Capture	MAB1288	500 µg
	Detection (Biotin)	BAF1288	50 µg
	Protein	1288-F3-025	25 µg

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Human FOLR1	Capture	MAB5646	100 µg
	Detection (Biotin)	BAF5646	50 µg
	Protein	5646-FR-050	50 µg
Human Furin	Capture	MAB15032	500 µg
	Detection (Biotin)	BAF1503	50 µg
	Protein	1503-SE-010	10 µg
Mouse Galectin-1	Capture	AF1245	100 µg
	Detection (Biotin)	BAF1245	50 µg
	Protein	1245-GA-050	50 µg
Human Galectin-3	Capture	MAB11541-50	500 µg
	Detection (Biotin)	BAF1154	50 µg
	Protein	1154-GA-050	50 µg
Mouse Galectin-3	Capture	MAB1197	500 µg
	Detection (Biotin)	BAF1197	50 µg
	Protein	1197-GA-050	50 µg
Human Galectin-7	Capture	AF1339	100 µg
	Detection (Biotin)	BAF1339	50 µg
	Protein	1339-GA-050	50 µg
Mouse Galectin-7	Capture	MAB1304	500 µg
	Detection (Biotin)	BAF1304	50 µg
	Protein	1304-GA-050	50 µg
Human Gas1	Capture	MAB2636	500 µg
	Detection (Biotin)	BAF2636	50 µg
	Protein	2636-GS-050	50 µg
Mouse Gas1	Capture	AF2644	100 µg
	Detection (Biotin)	BAF2644	50 µg
	Protein	2644-GS-050	50 µg
Mouse Gas6	Capture	AF986	100 µg
	Detection (Biotin)	BAF986	50 µg
	Protein	986-GS-025	25 µg
Human GASP-1/WFIKKNRP	Capture	MAB20702	500 µg
	Detection (Biotin)	BAF2070	50 µg
	Protein	2070-GS-025	25 µg
Human GASP-2/WFIKKN	Capture	MAB21361	500 µg
	Detection (Biotin)	BAM21362	250 µg
	Protein	2136-GS-025	25 µg
Human G-CSF	Capture	MAB214-500	500 µg
	Capture	MAB214R-100	100 µg
	Detection (Biotin)	BAF214	50 µg
	Protein	214-CS-005	5 µg
Mouse G-CSF	Capture	MAB414-500	500 µg
	Detection (Biotin)	BAF414	50 µg
	Protein	414-CS-005	5 µg
Human G-CSF R/CD114	Capture	MAB3811	500 µg
	Detection (Biotin)	BAF381	50 µg
	Protein	381-GR-050	50 µg
Human/Primate GDF-15	Capture	MAB957-500	500 µg
	Detection (Biotin)	BAF940	50 µg
	Protein	957-GD-025	25 µg
Human GDNF	Capture	MAB212-500	500 µg
	Detection (Biotin)	BAF212	50 µg
	Protein	212-GD-010	10 µg

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Human GITR/TNFRSF18	Capture	MAB689-500	500 µg
	Detection (Biotin)	BAF689	50 µg
	Protein	689-GR-100	100 µg
Mouse GITR/TNFRSF18	Capture	MAB524	500 µg
	Detection (Biotin)	BAF524	50 µg
	Protein	524-GR-050	50 µg
Human GITR Ligand/TNFSF18	Capture	MAB6942-500	500 µg
	Detection (Biotin)	BAM6943	250 µg
	Protein	694-GL-025	25 µg
Mouse GITR Ligand/TNFSF18	Capture	MAB2177	500 µg
	Detection (Biotin)	BAF2177	50 µg
	Protein	2177-GL-025	25 µg
Canine GM-CSF	Capture	MAB1546	500 µg
	Detection (Biotin)	BAF1546	50 µg
	Protein	1546-GM-025	25 µg
Feline GM-CSF	Capture	MAB9872	500 µg
	Detection (Biotin)	BAM9873	250 µg
	Protein	987-FL-010	10 µg
Human GM-CSF	Capture	MAB615-500	500 µg
	Detection (Biotin)	BAM215	250 µg
	Protein	215-GM-010	10 µg
Mouse GM-CSF	Capture	MAB415-500	500 µg
	Detection (Biotin)	BAF415	50 µg
	Protein	415-ML-010	10 µg
Rat GM-CSF	Capture	MAB518	500 µg
	Detection (Biotin)	BAF518	50 µg
	Protein	518-GM-005	5 µg
Human gp130	Capture	MAB628-500	500 µg
	Detection (Biotin)	BAF228	50 µg
	Protein	228-GP-010	10 µg
Mouse gp130	Capture	MAB4682-500	500 µg
	Detection (Biotin)	BAF468	50 µg
	Protein	468-MG-100	100 µg
Mouse Granzyme B	Capture	AF1865	100 µg
	Detection (Biotin)	BAF1865	50 µg
	Protein	1865-SE-010	10 µg
Mouse Gremlin	Capture	MAB956-500	500 µg
	Detection (Biotin)	BAF956	50 µg
	Protein	956-GR-050	50 µg
Mouse H60	Capture	MAB1155	500 µg
	Detection (Biotin)	BAF1155	50 µg
	Protein	1155-H6-050	50 µg
Human HAI-1	Capture	MAB10481	500 µg
	Detection (Biotin)	BAM10482	250 µg
	Protein	1048-PI-010	10 µg
Mouse HAI-1	Capture	MAB1141	500 µg
	Detection (Biotin)	BAF1141	50 µg
	Protein	1141-PI-010	10 µg
Human HAI-2	Capture	MAB1106	500 µg
	Detection (Biotin)	BAF1106	50 µg
	Protein	1106-PI-010	10 µg

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Human HB-EGF	Capture	AF-259-NA	250 µg
	Detection (Biotin)	BAF259	50 µg
	Protein	259-HE-050	50 µg
Human HGF	Capture	MAB694-500	500 µg
	Detection (Biotin)	BAF294	100 µg
	Protein	294-HG-005	5 µg
Mouse HGF	Capture	AF2207	100 µg
	Detection (Biotin)	BAF2207	50 µg
	Protein	2207-HG-025	25 µg
Canine HGF R/c-MET	Capture	AF4140	100 µg
	Detection (Biotin)	BAF4140	50 µg
	Protein	4140-ME-050	50 µg
Human HGF R/c-MET	Capture	MAB3581-500	500 µg
	Detection (Biotin)	BAF358	50 µg
	Protein	358-MT-100	100 µg
Mouse HGF R/c-MET	Capture	MAB5271	500 µg
	Detection (Biotin)	BAF527	50 µg
	Protein	527-ME-100	100 µg
Human HVEM/TNFRSF14	Capture	MAB356-500	500 µg
	Detection (Biotin)	BAF356	50 µg
	Protein	356-HV-100	100 µg
Human ICAM-1/CD54	Capture	MAB720-500	500 µg
	Detection (Biotin)	BAF720	50 µg
	Protein	ADP4-050	50 µg
Mouse ICAM-1/CD54	Capture	MAB796-500	500 µg
	Detection (Biotin)	BAF796	50 µg
	Protein	796-IC-050	50 µg
Rat ICAM-1/CD54	Capture	MAB5832-500	500 µg
	Detection (Biotin)	BAF583	50 µg
	Protein	583-IC-050	50 µg
Human ICAM-3/CD50	Capture	BBA15	200 µg
	Detection (Biotin)	BAF813	50 µg
	Protein	715-IC-050	50 µg
Mouse IFN-α	Capture	22100-1	250 µg
	Detection	32100-1	3000 UN
	Protein	12100-1	3000 UN
Bovine IFN-γ	Capture	MAB23001	500 µg
	Detection (Biotin)	BAF2300	50 µg
	Protein	2300-BG-025	25 µg
Canine IFN-γ	Capture	MAB781-500	500 µg
	Detection (Biotin)	BAM7811	250 µg
	Protein	781-CG-050	50 µg
Cotton Rat IFN-γ	Capture	AF565	100 µg
	Detection (Biotin)	BAF565	50 µg
	Protein	565-RG-025	25 µg
Equine IFN-γ	Capture	AF1586	100 µg
	Detection (Biotin)	BAF1586	50 µg
	Protein	1586-HG-025	25 µg
Human IFN-γ	Capture	MAB2852-500	500 µg
	Detection (Biotin)	BAF285	50 µg
	Protein	285-IF-100	100 µg

For research use only. Not for use in diagnostic procedures.



Analyte	Reagent	Catalog #	Size
Mouse IFN- $\gamma$	Capture	MAB785-500	500 $\mu$ g
	Detection (Biotin)	BAF485	50 $\mu$ g
	Protein	485-MI-100	100 $\mu$ g
Rat IFN- $\gamma$	Capture	MAB5851	500 $\mu$ g
	Detection (Biotin)	BAF585	50 $\mu$ g
	Protein	585-IF-100	100 $\mu$ g
Rhesus Macaque IFN- $\gamma$	Capture	AF961	100 $\mu$ g
	Detection (Biotin)	BAF961	50 $\mu$ g
	Protein	961-RM-025	25 $\mu$ g
Mouse IFN- $\gamma$ R1	Capture	MAB10262	500 $\mu$ g
	Detection (Biotin)	BAF1026	50 $\mu$ g
	Protein	1026-GR-100	100 $\mu$ g
Human IGFBP-1	Capture	MAB675-500	500 $\mu$ g
	Detection (Biotin)	BAF871	50 $\mu$ g
	Protein	871-B1-025	25 $\mu$ g
Human IGFBP-2	Capture	MAB6741-500	500 $\mu$ g
	Capture	MAB6741R-10	100 $\mu$ g
	Detection (Biotin)	BAF674	50 $\mu$ g
	Protein	674-B2-025	25 $\mu$ g
Mouse IGFBP-2	Capture	MAB7971	500 $\mu$ g
	Detection (Biotin)	BAF797	50 $\mu$ g
	Protein	797-B2-025	25 $\mu$ g
Human IGFBP-3	Capture	MAB305-500	500 $\mu$ g
	Detection (Biotin)	BAF675	50 $\mu$ g
	Protein	675-B3-025	25 $\mu$ g
Mouse IGFBP-3	Capture	MAB775	500 $\mu$ g
	Detection (Biotin)	BAF775	50 $\mu$ g
	Protein	775-B3-025	25 $\mu$ g
Human IGFBP-4	Capture	MAB8041	500 $\mu$ g
	Capture	MAB8041R-10	100 $\mu$ g
	Detection (Biotin)	BAF804	50 $\mu$ g
	Protein	804-GB-025	25 $\mu$ g
Human IGFBP-5	Capture	MAB8751-500	500 $\mu$ g
	Detection (Biotin)	BAF875	50 $\mu$ g
	Protein	875-B5-025	25 $\mu$ g
Mouse IGFBP-5	Capture	AF578	100 $\mu$ g
	Detection (Biotin)	BAF578	50 $\mu$ g
	Protein	578-B5-025	25 $\mu$ g
Human IGFBP-6	Capture	MAB8761	500 $\mu$ g
	Detection (Biotin)	BAF876	50 $\mu$ g
	Protein	876-B6-025	25 $\mu$ g
Mouse IGFBP-6	Capture	MAB7761	500 $\mu$ g
	Detection (Biotin)	BAF776	50 $\mu$ g
	Protein	776-B6-025	25 $\mu$ g
Human IGF-I	Capture	MAB291-500	500 $\mu$ g
	Detection (Biotin)	BAF291	50 $\mu$ g
Mouse IGF-I	Capture	MAB791-500	500 $\mu$ g
	Detection (Biotin)	BAF791	50 $\mu$ g
	Protein	791-MG-050	50 $\mu$ g
Human IGF-I R	Capture	MAB391-500	500 $\mu$ g
	Detection (Biotin)	BAF391	50 $\mu$ g
	Protein	391-GR-050	50 $\mu$ g

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Mouse IGF-II	Capture	MAB792-500	500 µg
	Detection (Biotin)	BAF792	50 µg
	Protein	792-MG-050	50 µg
Human IGF-II R	Capture	MAB24471	500 µg
	Detection (Biotin)	BAF2447	50 µg
	Protein	2447-GR-050	50 µg
Human IL-1α/IL-1F1	Capture	MAB200-500	500 µg
	Capture	MAB200R-100	100 µg
	Detection (Biotin)	BAF200	50 µg
	Protein	200-LA-002	2 µg
Mouse IL-1α/IL-1F1	Capture	MAB400-500	500 µg
	Detection (Biotin)	BAF400	50 µg
	Protein	400-ML-005	5 µg
Rat IL-1α/IL-1F1	Capture	MAB500	500 µg
	Detection (Biotin)	BAF500	50 µg
	Protein	500-RL-005	5 µg
Feline IL-1β/IL-1F2	Capture	AF1796	100 µg
	Detection (Biotin)	BAF1796	50 µg
	Protein	1796-FL-010	10 µg
Human IL-1β/IL-1F2	Capture	MAB601-500	500 µg
	Capture	MAB601R-100	100 µg
	Detection (Biotin)	BAF201	50 µg
	Protein	201-LB-005	5 µg
Mouse IL-1β/IL-1F2	Capture	MAB401-500	500 µg
	Detection (Biotin)	BAF401	50 µg
	Protein	401-ML-005	5 µg
Porcine IL-1β/IL-1F2	Capture	MAB6811-500	500 µg
	Detection (Biotin)	BAF681	50 µg
	Protein	681-PI-010	10 µg
Rat IL-1β/IL-1F2	Capture	AF-501-NA	100 µg
	Detection (Biotin)	BAF501	50 µg
	Protein	501-RL-010	10 µg
Equine IL-1ra/IL-1F3	Capture	AF2466	100 µg
	Detection (Biotin)	BAF2466	50 µg
	Protein	2466-RA-050	50 µg
Human IL-1ra/IL-1F3	Capture	MAB280	1 mg
	Detection (Biotin)	BAF280	50 µg
	Protein	280-RA-010	10 µg
Mouse IL-1ra/IL-1F3	Capture	AF-480-NA	100 µg
	Detection (Biotin)	BAF480	50 µg
	Protein	480-RM-010	10 µg
Human IL-1 RI	Capture	MAB269-500	500 µg
	Detection (Biotin)	BAF269	50 µg
	Protein	269-1R-100	100 µg
Human IL-1 RII	Capture	MAB663-500	500 µg
	Detection (Biotin)	BAF263	50 µg
	Protein	263-2R-050	50 µg
Bovine IL-2	Capture	AF2465	100 µg
	Detection (Biotin)	BAF2465	50 µg
	Protein	2465-BL-020	20 µg
Canine IL-2	Capture	AF1815	100 µg
	Detection (Biotin)	BAF1815	50 µg
	Protein	1815-CL-020	20 µg

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Equine IL-2	Capture	AF1613	100 µg
	Detection (Biotin)	BAF1613	50 µg
	Protein	1613-IL-020	20 µg
Feline IL-2	Capture	AF1890	100 µg
	Detection (Biotin)	BAF1890	50 µg
	Protein	1890-FL-020	20 µg
Human/Primate IL-2	Capture	MAB602-500	500 µg
	Detection (Biotin)	BAF202	50 µg
	Protein	202-IL-010	10 µg
Rat IL-2	Capture	AF-502-NA	100 µg
	Detection (Biotin)	BAF502	50 µg
	Protein	502-RL-010	10 µg
Human IL-3	Capture	MAB603-500	500 µg
	Capture	MAB603R-100	100 µg
	Detection (Biotin)	BAF203	50 µg
	Protein	203-IL-010	10 µg
Mouse IL-3	Capture	MAB403-500	500 µg
	Detection (Biotin)	BAF403	50 µg
	Protein	403-ML-010	10 µg
Mouse IL-3 Rβ	Capture	MAB5491	500 µg
	Detection (Biotin)	BAF549	50 µg
	Protein	549-R3-100	100 µg
Cotton Rat IL-4	Capture	MAB584	500 µg
	Detection (Biotin)	BAF584	50 µg
	Protein	584-R4-010	10 µg
Equine IL-4	Capture	AF1809	100 µg
	Detection (Biotin)	BAF1809	50 µg
	Protein	1809-EL-020	20 µg
Feline IL-4	Capture	MAB984	500 µg
	Detection (Biotin)	BAF984	50 µg
	Protein	984-FL-025	25 µg
Human IL-4	Capture	MAB604-500	500 µg
	Detection (Biotin)	BAF204	50 µg
	Protein	204-IL-010	10 µg
Mouse IL-4	Capture	MAB404-500	500 µg
	Capture	MAB404R-500	500 µg
	Detection (Biotin)	BAF404	50 µg
	Protein	404-ML-010	10 µg
Porcine IL-4	Capture	AF654	100 µg
	Detection (Biotin)	BAF654	50 µg
	Protein	654-P4-025	25 µg
Rat IL-4	Capture	MAB504-500	500 µg
	Detection (Biotin)	BAF504	50 µg
	Protein	504-RL-005	5 µg
Feline IL-5	Capture	MAB1795	500 µg
	Detection (Biotin)	BAM17951	250 µg
	Protein	1795-FL-010	10 µg
Human IL-5	Capture	MAB405-500	500 µg
	Capture	MAB405R-100	100 µg
	Detection (Biotin)	BAM6051	250 µg
	Protein	205-IL-005	5 µg

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Mouse IL-5	Capture	MAB405-500	500 µg
	Capture	MAB405R-100	100 µg
	Detection (Biotin)	BAM705	50 µg
	Protein	405-ML-005	5 µg
Canine IL-6	Capture	AF1609	100 µg
	Detection (Biotin)	BAF1609	50 µg
	Protein	1609-CL-025	25 µg
Human/Primate IL-6	Capture	MAB206-500	500 µg
	Detection (Biotin)	BAF206	50 µg
	Protein	206-IL-010	10 µg
Mouse IL-6	Capture	MAB406-500	500 µg
	Detection (Biotin)	BAF406	50 µg
	Protein	406-ML-005	5 µg
Porcine IL-6	Capture	AF686	100 µg
	Detection (Biotin)	BAF686	50 µg
	Protein	686-PI-025	25 µg
Rat IL-6	Capture	MAB506-500	500 µg
	Detection (Biotin)	BAF506	50 µg
	Protein	506-RL-010	10 µg
Human IL-6 Rα	Capture	MAB227-500	500 µg
	Capture	MAB227R-100	100 µg
	Detection (Biotin)	BAF227	50 µg
	Protein	227-SR-025	25 µg
Mouse IL-6 Rα	Capture	AF1830	100 µg
	Detection (Biotin)	BAF1830	50 µg
	Protein	1830-SR-025	25 µg
Human IL-7	Capture	MAB207-100	100 µg
	Detection (Biotin)	BAF207	50 µg
	Protein	207-IL-005	5 µg
Mouse IL-7	Capture	AF407	100 µg
	Detection (Biotin)	BAF407	50 µg
	Protein	407-ML-025	25 µg
Mouse IL-7 Rα/CD127	Capture	MAB4774	500 µg
	Detection (Biotin)	BAM47741	250 µg
	Protein	747-MR-050	50 µg
Canine IL-8/CXCL8	Capture	MAB16081	500 µg
	Detection (Biotin)	BAM16082	100 µg
	Protein	1608-CL-025	25 µg
Feline IL-8/CXCL8	Capture	MAB22772	500 µg
	Detection (Biotin)	BAM22771	250 µg
	Protein	2277-FL-025	25 µg
Human IL-8/CXCL8	Capture	MAB208-500	500 µg
	Detection (Biotin)	BAF208	50 µg
	Protein	208-IL-010	10 µg
Porcine IL-8/CXCL8	Capture	MAB5351-500	500 µg
	Detection (Biotin)	BAF535	50 µg
	Protein	535-IN-025	25 µg
Canine IL-10	Capture	MAB7351	500 µg
	Detection (Biotin)	BAF735	50 µg
	Protein	735-CL-010	10 µg
Equine IL-10	Capture	AF1605	100 µg
	Detection (Biotin)	BAF1605	50 µg
	Protein	1605-IL-010	10 µg

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Feline IL-10	Capture	MAB7361	500 µg
	Detection (Biotin)	BAF736	50 µg
	Protein	736-FL-010	10 µg
Human IL-10	Capture	MAB2172-100	100 µg
	Detection (Biotin)	BAF217	100 µg
	Protein	1064-IL-010	10 µg
Mouse IL-10	Capture	MAB417-500	500 µg
	Detection (Biotin)	BAF417	50 µg
	Protein	417-ML-005	5 µg
Rat IL-10	Capture	MAB519-500	500 µg
	Detection (Biotin)	BAF519	50 µg
	Protein	522-RL-005	5 µg
Human IL-11	Capture	MAB618-500	500 µg
	Detection (Biotin)	BAF218	50 µg
	Protein	218-IL-005	5 µg
Mouse IL-11	Capture	MAB4181	500 µg
	Detection (Biotin)	BAF418	50 µg
	Protein	418-ML-005	5 µg
Feline IL-12/IL-23 p40	Capture	AF1954	100 µg
	Detection (Biotin)	BAF1954	50 µg
	Protein	2117-FL-025	25 µg
Human IL-12 p70	Capture	MAB611-500	500 µg
	Capture	MAB611R-100	100 µg
	Detection (Biotin)	BAF219	100 µg
	Protein	219-IL-005	5 µg
Mouse IL-12 p70	Capture	MAB419-500	500 µg
	Detection (Biotin)	BAF419	50 µg
	Protein	419-ML-010	10 µg
Porcine IL-12	Capture	MAB9121	500 µg
	Detection (Biotin)	BAM9122	100 µg
	Protein	912-PL-025	25 µg
Human IL-12/IL-23 p40	Capture	MAB609-500	500 µg
	Detection (Biotin)	BAF219	100 µg
	Protein	309-IL-002	2 µg
Mouse IL-12/IL-23 p40 Allele-specific	Capture	MAB499-500	500 µg
	Detection (Biotin)	BAF499	50 µg
	Protein	499-ML-025	25 µg
Mouse IL-12/IL-23 p40 Nonallele-specific	Capture	MAB4991-500	500 µg
	Detection (Biotin)	BAF499	50 µg
	Protein	499-ML-025	25 µg
Human IL-13	Capture	MAB213-500	500 µg
	Detection (Biotin)	BAF213	100 µg
	Protein	213-ILB-005	5 µg
Mouse IL-13	Capture	MAB413-500	500 µg
	Detection (Biotin)	BAF413	50 µg
	Protein	413-ML-005	5 µg
Human IL-13 Rα1	Capture	MAB146	500 µg
	Detection (Biotin)	BAF152	50 µg
	Protein	146-IR-100	100 µg
Human IL-13 Rα2	Capture	MAB6141-500	500 µg
	Detection (Biotin)	BAF614	50 µg
	Protein	614-INS-100	100 µg

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Human/Primate IL-15	Capture	MAB647-500	500 µg
	Detection (Biotin)	BAM247	50 µg
	Protein	247-IL-005	5 µg
Mouse IL-15	Capture	MAB447-500	500 µg
	Detection (Biotin)	BAF447	50 µg
	Protein	447-ML-010	10 µg
Human IL-16	Capture	MAB316-500	500 µg
	Detection (Biotin)	BAF316	50 µg
	Protein	316-IL-002	2 µg
Human/Primate IL-17	Capture	MAB317-500	500 µg
	Detection (Biotin)	BAF317	50 µg
	Protein	317-ILB-050	50 µg
Mouse IL-17	Capture	MAB721-100	100 µg
	Detection (Biotin)	BAF421	50 µg
	Protein	421-ML-025	25 µg
Human IL-17B	Capture	MAB1248-500	500 µg
	Detection (Biotin)	BAM12481	250 µg
	Protein	1248-IB-025	25 µg
Mouse IL-17B	Capture	AF1709	100 µg
	Detection (Biotin)	BAF1709	50 µg
	Protein	1709-ML-025	25 µg
Human IL-17C	Capture	MAB1234	500 µg
	Detection (Biotin)	BAF1234	50 µg
	Protein	1234-IL-025	25 µg
Human IL-17D	Capture	MAB1504	500 µg
	Detection (Biotin)	BAF1504	50 µg
	Protein	1504-IL-025	25 µg
Mouse IL-17E/IL-25	Capture	MAB13991-50	500 µg
	Detection (Biotin)	BAF1399	50 µg
	Protein	1399-IL-025	25 µg
Human/Primate IL-17F	Capture	AF1335	100 µg
	Detection (Biotin)	BAF1335	50 µg
	Protein	1335-IL-025	25 µg
Mouse IL-17F	Capture	AF2057	100 µg
	Detection (Biotin)	BAF2057	50 µg
	Protein	2057-IL-025	25 µg
Human IL-17 RA/IL-17 R	Capture	MAB1771	500 µg
	Detection (Biotin)	BAF177	50 µg
	Protein	177-IR-100	100 µg
Human IL-17 RB	Capture	MAB12071	500 µg
	Detection (Biotin)	BAM1207	50 µg
	Protein	1207-BR-050	50 µg
Mouse IL-17 RB	Capture	MAB10401	500 µg
	Detection (Biotin)	BAF1040	50 µg
	Protein	1040-BR-050	50 µg
Human IL-18/IL-1F4	Capture	D044-3	100 µg
	Detection (Biotin)	D045-6	50 µL
	Protein	B001-5	25 µg
Mouse IL-18/IL-1F4	Capture	D047-3	100 µg
	Detection (Biotin)	D048-6	50 µL
	Protein	B002-5	25 µg

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Human IL-18 BPa	Capture	MAB1192	500 µg
	Detection (Biotin)	BAF119	50 µg
	Protein	119-BP-100	100 µg
Human IL-20	Capture	MAB11022	500 µg
	Detection (Biotin)	BAF1102	50 µg
	Protein	1102-IL-025	25 µg
Mouse IL-20	Capture	MAB12042	500 µg
	Detection (Biotin)	BAF1204	50 µg
	Protein	1204-ML-025	25 µg
Mouse IL-21	Capture	AF594	100 µg
	Detection (Biotin)	BAF594	50 µg
	Protein	594-ML-010	10 µg
Human IL-22	Capture	MAB7822-500	500 µg
	Detection (Biotin)	BAM7821	100 µg
	Protein	782-IL-010	10 µg
Mouse IL-22	Capture	AF582	100 µg
	Detection (Biotin)	BAF582	50 µg
	Protein	582-ML-010	10 µg
Mouse IL-2	Capture	MAB702-500	500 µg
	Detection (Biotin)	BAF402	50 µg
	Protein	402-ML-020	20 µg
Human IL-24	Capture	MAB19652	500 µg
	Detection (Biotin)	BAF1965	50 µg
	Protein	1965-IL-025	25 µg
Mouse IL-23	Capture	MAB1887	500 µg
	Detection (Biotin)	BAF499	50 µg
	Protein	1887-ML-010	10 µg
Human IL-27	Capture	AF2526	100 µg
	Detection (Biotin)	BAF2526	50 µg
	Protein	2526-IL-010	10 µg
Human IL-28A/IFN-λ2	Capture	MAB15871-50	500 µg
	Detection (Biotin)	BAF1587	50 µg
	Protein	1587-IL-025	25 µg
Mouse IL-28A/IFN-λ2	Capture	MAB17892-50	500 µg
	Detection (Biotin)	BAM17891	250 µg
	Protein	1789-ML-025	25 µg
IL-29/IFN-λ1	Capture	MAB15981-50	500 µg
	Detection (Biotin)	BAF1598	50 µg
	Protein	1598-IL-025	25 µg
Human IL-31	Capture	AF2824	100 µg
	Detection (Biotin)	BAF2824	50 µg
	Protein	2824-IL-010	10 µg
Human IL-33	Capture	AF3625	100 µg
	Detection (Biotin)	BAF3625	50 µg
	Protein	3625-IL-010	10 µg
Mouse IL-33	Capture	AF3626	100 µg
	Detection (Biotin)	BAF3626	50 µg
	Protein	3626-ML-010	10 µg
Human IL-34	Capture	MAB52651	500 µg
	Detection (Biotin)	BAM52652	250 µg
	Protein	5265-IL-010	10 µg

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Human Proinsulin	Capture	MAB13363	500 µg
	Detection (Biotin)	BAM13362	250 µg
	Protein	1336-PN-050	50 µg
Rat Jagged 1	Capture	AF599	100 µg
	Detection (Biotin)	BAF599	50 µg
	Protein	599-JG-100	100 µg
Mouse JAM-A	Capture	AF1077	100 µg
	Detection (Biotin)	BAF1077	50 µg
	Protein	1077-JM-050	50 µg
Human Kallikrein 3/PSA	Capture	MAB13442-50	500 µg
	Detection (Biotin)	BAF1344	50 µg
	Protein	1344-SE-010	10 µg
Human Kallikrein 14	Capture	AF2626	100 µg
	Detection (Biotin)	BAF2626	50 µg
	Protein	2626-SE-010	10 µg
Mouse Kremen-1	Capture	MAB16471	500 µg
	Detection (Biotin)	BAF1647	50 µg
	Protein	1647-KR-025	25 µg
Human LAG-3	Capture	MAB23192	500 µg
	Detection (Biotin)	BAM23191	250 µg
	Protein	2319-L3-050	50 µg
Human Layilin	Capture	MAB26371	500 µg
	Detection (Biotin)	BAF2637	50 µg
	Protein	2536-LA-025	25 µg
Human LDL R	Capture	MAB2148-500	500 µg
	Detection (Biotin)	BAF2148	50 µg
	Protein	2148-LD-025	25 µg
Human Legumain	Capture	MAB21992	500 µg
	Detection (Biotin)	BAF2199	50 µg
	Protein	2199-CY-010	10 µg
Human Leptin/OB	Capture	MAB398-500	500 µg
	Detection (Biotin)	BAM398	250 µg
	Protein	398-LP-01M	1 mg
Mouse Leptin/OB	Capture	AF498	100 µg
	Detection (Biotin)	BAF498	50 µg
	Protein	498-OB-01M	1 mg
Mouse Leptin R	Capture	MAB497	500 µg
	Detection (Biotin)	BAF497	50 µg
	Protein	497-LR-100	100 µg
Human LIF	Capture	MAB250-500	500 µg
	Detection (Biotin)	BAF250	50 µg
Human LIGHT/TNFSF14	Capture	MAB664	500 µg
	Detection (Biotin)	BAF664	50 µg
	Protein	664-LI-025	25 µg
Mouse Limitin	Capture	MAB5971	500 µg
	Detection (Biotin)	BAF597	50 µg
	Protein	1535-LM-010	10 µg
Human LIMPII/SR-B2	Capture	AF1966	100 µg
	Detection (Biotin)	BAF1966	50 µg
	Protein	1966-LM-050	50 µg

For research use only. Not for use in diagnostic procedures.



Analyte	Reagent	Catalog #	Size
Human Lipocalin-2/NGAL	Capture	MAB17571-50	500 µg
	Capture	MAB17571R-1	100 µg
	Detection (Biotin)	BAF1757	50 µg
	Protein	1757-LC-050	50 µg
Mouse Lipocalin-2/NGAL	Capture	MAB18571-50	500 µg
	Detection (Biotin)	BAM1857	250 µg
	Protein	1857-LC-050	50 µg
Rat Lipocalin-2	Capture	AF3508	100 µg
	Detection (Biotin)	BAF3508	50 µg
	Protein	3508-LC-050	50 µg
Mouse LIX	Capture	MAB433-500	500 µg
	Detection (Biotin)	BAF433	50 µg
	Protein	433-MC-025	25 µg
Rat LIX	Capture	MAB543	500 µg
	Detection (Biotin)	BAF543	50 µg
	Protein	543-RL-025	25 µg
Human LOX-1/OLR1	Capture	MAB17981-50	500 µg
	Detection (Biotin)	BAF1798	50 µg
	Protein	1798-LX-050	50 µg
Mouse LOX-1/OLR1	Capture	MAB15641	500 µg
	Detection (Biotin)	BAF1564	50 µg
	Protein	1564-LX-050	50 µg
Human Lymphotoxin- $\alpha$ /TNF- $\beta$	Capture	MAB621	1 mg
	Capture	MAB621R-100	100 µg
	Detection (Biotin)	BAF211	50 µg
	Protein	211-TBB-010	10 µg
Human LYVE-1	Capture	MAB20891	500 µg
	Detection (Biotin)	BAF2089	50 µg
	Protein	2089-LY-025	25 µg
Mouse MAdCAM-1	Capture	MAB9931	500 µg
	Detection (Biotin)	BAF993	50 µg
	Protein	993-MC-050	50 µg
Rat MAG/Siglec-4a	Capture	MAB5381	500 µg
	Detection (Biotin)	BAF538	50 µg
	Protein	538-MG-100	100 µg
Human Marapsin/Pancreasin	Capture	AF1988	100 µg
	Detection (Biotin)	BAF1988	50 µg
	Protein	1988-SE-010	10 µg
Mouse Marapsin/Pancreasin	Capture	MAB19891	500 µg
	Detection (Biotin)	BAF1989	50 µg
	Protein	1989-SE-010	10 µg
Human MBL	Capture	AF2307	100 µg
	Detection (Biotin)	BAF2307	50 µg
	Protein	2307-MB-050	50 µg
Mouse MBL-2	Capture	AF2208	100 µg
	Detection (Biotin)	BAF2208	50 µg
	Protein	2208-MB-050	50 µg
Human M-CSF	Capture	MAB616-500	500 µg
	Detection (Biotin)	BAF216	50 µg
	Protein	216-MC-005	5 µg
Mouse M-CSF	Capture	MAB416-500	500 µg
	Detection (Biotin)	BAF416	50 µg
	Protein	416-ML-010	10 µg

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Human M-CSF R/CD115	Capture	MAB3292-500	500 µg
	Detection (Biotin)	BAF329	50 µg
	Protein	329-MR-100	100 µg
Human Mer	Capture	MAB8912-500	500 µg
	Detection (Biotin)	BAF891	50 µg
	Protein	891-MR-100	100 µg
Mouse Meteorin	Capture	MAB34752	500 µg
	Detection (Biotin)	BAM34751	250 µg
	Protein	3475-MN-025	25 µg
Mouse MFG-E8	Capture	MAB2805-500	500 µg
	Detection (Biotin)	BAF2805	50 µg
	Protein	2805-MF-050	50 µg
Human MICA	Capture	MAB13002-50	500 µg
	Detection (Biotin)	BAF1300	50 µg
	Protein	1300-MA-050	50 µg
Human MICB	Capture	MAB1599-500	500 µg
	Detection (Biotin)	BAF1599	50 µg
	Protein	1599-MB-050	50 µg
Human MMP-1	Capture	AF901	100 µg
	Detection (Biotin)	BAF901	50 µg
	Protein	901-MP-010	10 µg
Human/Primate MMP-3	Capture	AF513	100 µg
	Detection (Biotin)	BAF513	50 µg
	Protein	513-MP-010	10 µg
Human/Primate MMP-7	Capture	MAB9072-500	500 µg
	Detection (Biotin)	BAF907	50 µg
	Protein	907-MP-010	10 µg
Human MMP-8	Capture	MAB908-500	500 µg
	Detection (Biotin)	BAF908	50 µg
	Protein	908-MP-010	10 µg
Human/Primate MMP-9	Capture	MAB936-500	500 µg
	Detection (Biotin)	BAF911	50 µg
	Protein	911-MP-010	10 µg
Mouse Pro-MMP-9	Capture	MAB9092	500 µg
	Detection (Biotin)	BAM909	250 µg
	Protein	909-MM-010	10 µg
Human MMP-10	Capture	MAB9103	500 µg
	Detection (Biotin)	BAF910	50 µg
	Protein	910-MP-010	10 µg
Human MSP	Capture	MAB352	500 µg
	Detection (Biotin)	BAM3521	250 µg
	Protein	352-MS-010	10 µg
Human/Primate NCAM-1/CD56	Capture	MAB2408	500 µg
	Detection (Biotin)	BAF2408	50 µg
	Protein	2408-NC-050	50 µg
Human Nephilysin/CD10	Capture	AF1182	100 µg
	Detection (Biotin)	BAF1182	50 µg
	Protein	1182-ZNC-010	10 µg
Mouse Nephilysin/CD10	Capture	AF1126	100 µg
	Detection (Biotin)	BAF1126	50 µg
	Protein	1126-ZN-010	10 µg

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Human NRG1- $\beta$ 1/HRG1- $\beta$ 1	Capture	MAB3771-500	500 $\mu$ g
	Detection (Biotin)	BAF377	50 $\mu$ g
	Protein	377-HB-050	50 $\mu$ g
Human $\beta$ -NGF	Capture	MAB256-500	500 $\mu$ g
	Detection (Biotin)	BAF256	50 $\mu$ g
	Protein	256-GF-100	100 $\mu$ g
Rat $\beta$ -NGF	Capture	AF-556-NA	100 $\mu$ g
	Detection (Biotin)	BAF556	50 $\mu$ g
	Protein	556-NG-100	100 $\mu$ g
Human Nidogen-1/Entactin	Capture	MAB2570-500	500 $\mu$ g
	Detection (Biotin)	BAF2570	50 $\mu$ g
	Protein	2570-ND-050	50 $\mu$ g
Mouse Nope	Capture	AF1394	100 $\mu$ g
	Detection (Biotin)	BAF1394	50 $\mu$ g
	Protein	1394-NP-050	50 $\mu$ g
Human Notch-1	Capture	MAB5317-500	500 $\mu$ g
	Detection (Biotin)	BAM53171	250 $\mu$ g
	Protein	3647-TK-050	50 $\mu$ g
Human NOV/CCN3	Capture	AF1640	100 $\mu$ g
	Detection (Biotin)	BAF1640	50 $\mu$ g
	Protein	1640-NV-050	50 $\mu$ g
Mouse NOV/CCN3	Capture	MAB19762	500 $\mu$ g
	Detection (Biotin)	BAM19761	250 $\mu$ g
	Protein	1976-NV-050	50 $\mu$ g
Human NrCAM	Capture	MAB20341	500 $\mu$ g
	Detection (Biotin)	BAF2034	50 $\mu$ g
	Protein	2034-NR-050	50 $\mu$ g
Human NT-3	Capture	MAB267-500	500 $\mu$ g
	Detection (Biotin)	BAF267	50 $\mu$ g
	Protein	267-N3-005	5 $\mu$ g
Human NT-4	Capture	MAB268	500 $\mu$ g
	Detection (Biotin)	BAF268	50 $\mu$ g
	Protein	268-N4-005	5 $\mu$ g
Human Oncostatin M/OSM	Capture	MAB295-500	500 $\mu$ g
	Detection (Biotin)	BAF295	50 $\mu$ g
	Protein	295-OM-010	10 $\mu$ g
Human Osteoactivin/GPNMB	Capture	AF2550	100 $\mu$ g
	Detection (Biotin)	BAF2550	50 $\mu$ g
	Protein	2550-AC-050	50 $\mu$ g
Mouse Osteoactivin/GPNMB	Capture	AF2330	100 $\mu$ g
	Detection (Biotin)	BAF2330	50 $\mu$ g
	Protein	2330-AC-050	50 $\mu$ g
Human Osteopontin/OPN	Capture	MAB14332-50	500 $\mu$ g
	Detection (Biotin)	BAF1433	50 $\mu$ g
	Protein	1433-OP-050	50 $\mu$ g
Mouse Osteopontin/OPN	Capture	AF808	100 $\mu$ g
	Detection (Biotin)	BAF808	50 $\mu$ g
	Protein	441-OP-050	50 $\mu$ g
Human Osteoprotegerin/TNFRSF11B	Capture	MAB8051-500	500 $\mu$ g
	Detection (Biotin)	BAF805	50 $\mu$ g
	Protein	185-OS-025	25 $\mu$ g

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Mouse Osteoprotegerin/TNFRSF11B	Capture	MAB4591	500 µg
	Detection (Biotin)	BAF459	50 µg
	Protein	459-MO-100	100 µg
Mouse OX40 Ligand/TNFSF4	Capture	MAB12362	500 µg
	Detection (Biotin)	BAM12361	250 µg
	Protein	1236-OX-025	25 µg
Human PD-1	Capture	AF1086	100 µg
	Detection (Biotin)	BAF1086	50 µg
	Protein	1086-PD-050	50 µg
Human PDGF-AB	Capture	MAB1739-500	500 µg
	Detection (Biotin)	BAF221	50 µg
	Protein	222-AB-010	10 µg
Human PDGF-BB	Capture	385-PR-100/C	100 µg
	Detection (Biotin)	BAF220	50 µg
	Protein	220-BB-010	10 µg
Human Pentraxin 3/TSG-14	Capture	MAB1826-500	500 µg
	Detection (Biotin)	BAF1826	50 µg
	Protein	1826-TS-025	25 µg
Mouse Pentraxin 3/TSG-14	Capture	MAB21661	500 µg
	Detection (Biotin)	BAF2166	50 µg
	Protein	2166-TS-025	25 µg
Mouse Periostin/OSF-2	Capture	MAB2955	500 µg
	Detection (Biotin)	BAF2955	50 µg
	Protein	2955-F2-050	50 µg
Human PGLYRP1/PGRP-S	Capture	AF2590	100 µg
	Detection (Biotin)	BAF2590	50 µg
	Protein	2590-PG-050	50 µg
Human PIGF	Capture	MAB264-500	500 µg
	Detection (Biotin)	BAF264	50 µg
	Protein	264-PG-010	10 µg
Mouse PIGF-2	Capture	MAB4651	500 µg
	Detection (Biotin)	BAF465	50 µg
	Protein	465-PL-010	10 µg
Mouse PRDC	Capture	AF2069	100 µg
	Detection (Biotin)	BAF2069	50 µg
	Protein	2069-PR-050	50 µg
Mouse Progranulin	Capture	MAB2557	500 µg
	Detection (Biotin)	BAM25572	50 µg
	Protein	2557-PG-050	50 µg
Human Prolactin	Capture	AF682	100 µg
	Detection (Biotin)	BAF682	50 µg
	Protein	682-PL-050	50 µg
Mouse Prolactin	Capture	AF1445	100 µg
	Detection (Biotin)	BAF1445	50 µg
	Protein	1445-PL-050	50 µg
Mouse Prostatin/Prss8	Capture	AF2968	100 µg
	Detection (Biotin)	BAF2968	50 µg
	Protein	2968-SE-010	10 µg
Canine RAGE	Capture	100-B-001	1 µg
	Capture	AF4750	100 µg
	Detection (Biotin)	BAF4750	50 µg
	Protein	4750-RG-050	50 µg

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Human RAGE	Capture	MAB11451-50	500 µg
	Detection (Biotin)	BAF1145	50 µg
	Protein	1145-RG-050	50 µg
Mouse RAGE	Capture	MAB11792	500 µg
	Detection (Biotin)	BAM11791	250 µg
	Protein	1179-RG-050	50 µg
Rat RAGE	Capture	AF1616	100 µg
	Detection (Biotin)	BAF1616	50 µg
	Protein	1616-RG-050	50 µg
Human RANK/TNFRSF11A	Capture	MAB683	500 µg
	Detection (Biotin)	BAF683	50 µg
	Protein	683-RK-100	100 µg
Human RBP4/Retinol Binding Protein 4	Capture	MAB33781-50	500 µg
	Detection (Biotin)	BAM33782	250 µg
	Protein	3378-LC-050	50 µg
Mouse Renin 1	Capture	AF4277	100 µg
	Detection (Biotin)	BAF4277	50 µg
	Protein	4277-AS-020	20 µg
Human Resistin	Capture	MAB13591-50	500 µg
	Detection (Biotin)	BAM1359	100 µg
	Protein	1359-RN-005	5 µg
Mouse Resistin	Capture	MAB1069	500 µg
	Detection (Biotin)	BAF1069	50 µg
	Protein	1069-RN-050	50 µg
Rat ROBO1	Capture	MAB17491	500 µg
	Detection (Biotin)	BAF1749	50 µg
	Protein	1749-RB-050	50 µg
Canine SCF/c-kit Ligand	Capture	AF2278	100 µg
	Detection (Biotin)	BAF2278	50 µg
	Protein	2278-SC-025	25 µg
Human SCF/c-kit Ligand	Capture	MAB655-500	500 µg
	Detection (Biotin)	BAF255	50 µg
	Protein	255-SC-010	10 µg
Mouse SCF/c-kit Ligand	Capture	AF-455-NA	100 µg
	Detection (Biotin)	BAF455	50 µg
	Protein	455-MC-010	10 µg
Human E-Selectin/CD62E	Capture	BBA16	200 µg
	Capture	BBA2	1 mg
	Detection (Biotin)	BBA8	100 µg
	Protein	ADP1-050	50 µg
Mouse E-Selectin/CD62E	Capture	MAB5751-500	500 µg
	Detection (Biotin)	BAM5752	100 µg
	Protein	575-ES-100	100 µg
Human L-Selectin/CD62L	Capture	BBA24	200 µg
	Detection (Biotin)	BAF728	50 µg
	Protein	ADP2-050	50 µg
Mouse L-Selectin/CD62L	Capture	MAB5762	500 µg
	Detection (Biotin)	BAM5761	250 µg
	Protein	576-LS-100	100 µg
Rat L-Selectin/CD62L	Capture	MAB1534	500 µg
	Detection (Biotin)	BAF1534	50 µg
	Protein	1534-LS-050	50 µg

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Mouse P-Selectin/CD62P	Capture	MAB737	500 µg
	Detection (Biotin)	BAF737	50 µg
	Protein	737-PS-050	50 µg
Human Serpin A4/Kallistatin	Capture	MAB16691	500 µg
	Detection (Biotin)	BAF1669	50 µg
Human Serpin E1/PAI-1	Capture	MAB1786-500	500 µg
	Detection (Biotin)	BAF1786	50 µg
	Protein	1786-PI-010	10 µg
Human/Mouse sFRP-3	Capture	MAB1921	500 µg
	Detection (Biotin)	BAM192	250 µg
	Protein	192-SF-010	10 µg
Human Siglec-5	Capture	MAB10721	500 µg
	Detection (Biotin)	BAM10722	250 µg
	Protein	1072-SL-050	50 µg
Human Siglec-9	Capture	MAB1139-500	500 µg
	Detection (Biotin)	BAF1139	50 µg
	Protein	1139-SL-050	50 µg
Mouse SLAM/CD150	Capture	MAB4330	100 µg
	Detection (Biotin)	BAF4330	50 µg
	Protein	4330-SL-050	50 µg
Mouse Sonic Hedgehog	Capture	MAB4641-500	500 µg
	Detection (Biotin)	BAM4642	50 µg
	Protein	461-SH-025	25 µg
Human SOST	Capture	MAB1406-500	500 µg
	Detection (Biotin)	BAM14061	250 µg
	Protein	1406-ST-025	25 µg
Human/Primate ST2/IL-33 R	Capture	MAB523-500	500 µg
	Detection (Biotin)	BAF523	50 µg
	Protein	523-ST-100	100 µg
Mouse ST2/IL-33 R	Capture	MAB10041-50	500 µg
	Detection (Biotin)	BAF1004	50 µg
	Protein	1004-MR-050	50 µg
Human Syndecan-1/CD138	Capture	AF2780	100 µg
	Detection (Biotin)	BAF2780	50 µg
	Protein	2780-SD-050	50 µg
Human TACE/ADAM17	Capture	MAB9304	500 µg
	Detection (Biotin)	BAM9303	250 µg
	Protein	930-ADB-010	10 µg
Human TACI/TNFRSF13B	Capture	MAB174-500	500 µg
	Detection (Biotin)	BAF174	50 µg
	Protein	174-TC-050	50 µg
Mouse TACI/TNFRSF13B	Capture	MAB1041	500 µg
	Detection (Biotin)	BAF1041	50 µg
	Protein	1041-TC-050	50 µg
Mouse Testican 3/SPOCK	Capture	MAB2346	500 µg
	Detection (Biotin)	BAF2346	50 µg
	Protein	2346-PI-050	50 µg
Human TFPI	Capture	MAB29741	500 µg
	Detection (Biotin)	BAF2974	50 µg
	Protein	2974-PI-010	10 µg
Human TGF-α	Capture	AF-239-NA	100 µg
	Detection (Biotin)	BAF239	50 µg
	Protein	239-A-100	100 µg

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Human LAP TGF- $\beta$ 1	Capture	MAB2461-500	500 $\mu$ g
	Detection (Biotin)	BAM2462	50 $\mu$ g
	Protein	246-LP-025	25 $\mu$ g
Human TGF- $\beta$ 1	Capture	341-BR-050	50 $\mu$ g
	Capture	MAB1835-500	500 $\mu$ g
	Capture	MAB240-500	500 $\mu$ g
	Detection (Biotin)	BAF240	50 $\mu$ g
	Protein	240-B-002	2 $\mu$ g
Human TGF- $\beta$ 2	Capture	MAB612-500	500 $\mu$ g
	Detection (Biotin)	BAF302	50 $\mu$ g
	Protein	302-B2-002	2 $\mu$ g
Human TGF- $\beta$ 3	Capture	MAB643-500	500 $\mu$ g
	Detection (Biotin)	BAF243	50 $\mu$ g
	Protein	243-B3-002	2 $\mu$ g
Human TGF- $\beta$ RII	Capture	AF-241-NA	250 $\mu$ g
	Detection (Biotin)	BAF241	50 $\mu$ g
	Protein	241-R2-025	25 $\mu$ g
Human TGF- $\beta$ RIII	Capture	MAB2421	500 $\mu$ g
	Detection (Biotin)	BAF242	50 $\mu$ g
	Protein	242-R3-100	100 $\mu$ g
Human Thrombomodulin/BDCA-3	Capture	MAB3947-100	100 $\mu$ g
	Detection (Biotin)	BAM39471	50 $\mu$ g
	Protein	3947-PA-010	10 $\mu$ g
Human Thrombopoietin/Tpo	Capture	MAB288-500	500 $\mu$ g
	Detection (Biotin)	BAF288	50 $\mu$ g
	Detection (Biotin)	BAM288	250 $\mu$ g
	Protein	288-TPN-025	25 $\mu$ g
Mouse Thrombopoietin/Tpo	Capture	MAB488	1 mg
	Detection (Biotin)	BAF488	50 $\mu$ g
	Protein	488-TO-005	5 $\mu$ g
Human Thrombospondin-1	Capture	MAB3074-500	500 $\mu$ g
	Detection (Biotin)	BAF3074	50 $\mu$ g
	Protein	3074-TH-050	50 $\mu$ g
Human Thrombospondin-2	Capture	MAB16351-50	500 $\mu$ g
	Detection (Biotin)	BAF1635	50 $\mu$ g
	Protein	1635-T2-050	50 $\mu$ g
Human Tie-2	Capture	MAB3132-500	500 $\mu$ g
	Detection (Biotin)	BAM3313	250 $\mu$ g
	Protein	313-TI-100	100 $\mu$ g
Human TIM-1/KIM-1/HAVCR	Capture	AF1750	100 $\mu$ g
	Detection (Biotin)	BAF1750	50 $\mu$ g
	Protein	1750-TM-050	50 $\mu$ g
Mouse TIM-1/KIM-1/HAVCR	Capture	MAB1817-500	500 $\mu$ g
	Detection (Biotin)	BAF1817	50 $\mu$ g
	Protein	1817-TM-050	50 $\mu$ g
Rat TIM-1/KIM-1/HAVCR	Capture	MAB3689-500	500 $\mu$ g
	Detection (Biotin)	BAF3689	50 $\mu$ g
	Protein	3689-TM-050	50 $\mu$ g
Human TIMP-1	Capture	MAB970-500	500 $\mu$ g
	Detection (Biotin)	BAF970	50 $\mu$ g
	Protein	970-TM-010	10 $\mu$ g

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Mouse TIMP-1	Capture	MAB980-500	500 µg
	Detection (Biotin)	BAF980	50 µg
	Protein	980-MT-010	10 µg
Rat TIMP-1	Capture	MAB5802	500 µg
	Detection (Biotin)	BAF580	50 µg
	Protein	580-RT-010	10 µg
Human TIMP-2	Capture	MAB9711-500	500 µg
	Detection (Biotin)	BAF971	50 µg
	Protein	971-TM-010	10 µg
Human TIMP-3	Capture	MAB973-500	500 µg
	Detection (Biotin)	BAM9731	250 µg
	Protein	973-TM-010	10 µg
Human TIMP-4	Capture	MAB974	500 µg
	Detection (Biotin)	BAF974	50 µg
	Protein	974-TSF-010	10 µg
Human TLR2	Capture	AF2616	100 µg
	Detection (Biotin)	BAF2616	50 µg
	Protein	2616-TR-050	50 µg
Bovine TNF-α	Capture	AF2279	100 µg
	Detection (Biotin)	BAF2279	50 µg
	Protein	2279-BT-025	25 µg
Canine TNF-α	Capture	MAB1507	500 µg
	Detection (Biotin)	BAF1507	50 µg
	Protein	1507-CT-025	25 µg
Guinea Pig TNF-α	Capture	MAB5035	500 µg
	Detection (Biotin)	BAM50351	250 µg
	Protein	5035-TG-025	25 µg
Human TNF-α	Capture	MAB610-500	500 µg
	Capture	MAB610R-100	100 µg
	Detection (Biotin)	BAF210	50 µg
Mouse TNF-α	Capture	AF-410-NA	100 µg
	Detection (Biotin)	BAF410	50 µg
	Protein	410-MT-010	10 µg
Porcine TNF-α	Capture	MAB6902	500 µg
	Detection (Biotin)	BAM6903	250 µg
	Protein	690-PT-025	25 µg
Rat TNF-α	Capture	MAB510-500	500 µg
	Capture	MAB510R-100	100 µg
	Detection (Biotin)	BAF510	50 µg
	Protein	510-RT-010	10 µg
Rhesus Macaque TNF-α	Capture	MAB10702	500 µg
	Detection (Biotin)	BAF1070	50 µg
	Protein	1070-RM-025	25 µg
Human TNF RI/TNFRSF1A	Capture	MAB625-500	500 µg
	Detection (Biotin)	BAF225	50 µg
	Protein	636-R1-025	25 µg
Mouse TNF RI/TNFRSF1A	Capture	MAB425	500 µg
	Detection (Biotin)	BAF425	50 µg
	Protein	425-R1-050	50 µg
Human TNF RII/TNFRSF1B	Capture	MAB726-500	500 µg
	Detection (Biotin)	BAF726	50 µg
	Protein	1089-R2-025	25 µg

For research use only. Not for use in diagnostic procedures.



Analyte	Reagent	Catalog #	Size
Mouse TNF RII/TNFRSF1B	Capture	MAB426	500 µg
	Detection (Biotin)	BAF426	50 µg
	Protein	426-R2-050	50 µg
Human/Primate TRAIL/TNFSF10	Capture	MAB3751-500	500 µg
	Detection (Biotin)	BAF375	50 µg
	Protein	375-TL-010	10 µg
Mouse TRAIL/TNFSF10	Capture	MAB1121	500 µg
	Detection (Biotin)	BAF1121	50 µg
	Protein	1121-TL-010	10 µg
Human TRAIL R3/TNFRSF10C	Capture	MAB6301-500	500 µg
	Detection (Biotin)	BAF630	50 µg
	Protein	630-TR-100	100 µg
Human TRAIL R4/TNFRSF10D	Capture	MAB633	500 µg
	Detection (Biotin)	BAF633	50 µg
	Protein	633-TR-100	100 µg
Mouse TRANCE/TNFSF11/RANK L	Capture	AF462	100 µg
	Detection (Biotin)	BAF462	50 µg
	Protein	462-TR-010	10 µg
Human Trappin-2/Elafin	Capture	AF1747	100 µg
	Detection (Biotin)	BAF1747	50 µg
	Protein	1747-PI-010	10 µg
Human TREM-1	Capture	AF1278	100 µg
	Detection (Biotin)	BAF1278	50 µg
	Protein	1278-TR-050	50 µg
Mouse TREM-1	Capture	AF1187	100 µg
	Detection (Biotin)	BAF1187	50 µg
	Protein	1187-TR-025	25 µg
Mouse TREML1/TLT-1	Capture	MAB24241	500 µg
	Detection (Biotin)	BAF2424	50 µg
	Protein	2424-TL-050	50 µg
Mouse TROY/TNFRSF19	Capture	AF723	100 µg
	Detection (Biotin)	BAF723	50 µg
	Protein	723-TR-100	100 µg
Mouse Tryptase ε/BSSP-4	Capture	MAB20592	500 µg
	Detection (Biotin)	BAM20591	250 µg
	Protein	2059-SE-010	10 µg
Human TSLP	Capture	AF1398	100 µg
	Detection (Biotin)	BAF1398	50 µg
	Protein	1398-TS-010	10 µg
Mouse TSLP	Capture	MAB5551-500	500 µg
	Detection (Biotin)	BAF555	50 µg
	Protein	555-TS-010	10 µg
Human TWEAK/TNFSF12	Capture	MAB1090-500	500 µg
	Detection (Biotin)	BAF1090	50 µg
	Protein	1090-TW-025	25 µg
Mouse TWEAK/TNFSF12	Capture	MAB1237	500 µg
	Detection (Biotin)	BAF1237	50 µg
	Protein	1237-TW-025	25 µg
Mouse TWEAK R/TNFRSF12	Capture	AF1610	100 µg
	Detection (Biotin)	BAF1610	50 µg
	Protein	1610-TW-050	50 µg

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Human Ubiquitin+1	Capture	MAB7031	500 µg
	Detection (Biotin)	BAM7032	250 µg
	Protein	703-UB-025	25 µg
Human u-Plasminogen Activator/Urokinase	Capture	MAB1310-500	500 µg
	Detection (Biotin)	BAF1310	50 µg
	Protein	1310-SE-010	10 µg
Human uPAR	Capture	MAB807-500	500 µg
	Detection (Biotin)	BAF807	50 µg
	Protein	807-UK-100	100 µg
Human VCAM-1/CD106	Capture	MAB809-500	500 µg
	Detection (Biotin)	BAF809	50 µg
	Protein	809-VR-050	50 µg
Mouse VCAM-1/CD106	Capture	MAB6432-500	500 µg
	Detection (Biotin)	BAM6433	100 µg
	Protein	643-VM-050	50 µg
Canine VEGF	Capture	MAB1603	500 µg
	Detection (Biotin)	BAF1603	50 µg
	Protein	1603-CV-010	10 µg
Human/Primate VEGF	Capture	MAB293-500	500 µg
	Capture	MAB293R-100	100 µg
	Detection (Biotin)	BAF293	50 µg
	Protein	293-VE-010	10 µg
Mouse VEGF	Capture	AF-493-NA	100 µg
	Detection (Biotin)	BAF493	50 µg
	Protein	493-MV-005	5 µg
Rat VEGF	Capture	MAB564-500	500 µg
	Detection (Biotin)	BAF564	50 µg
	Protein	564-RV-010	10 µg
Human VEGF-D	Capture	MAB2861-500	500 µg
	Detection (Biotin)	BAM286	250 µg
	Protein	622-VD-005	5 µg
Mouse VEGF-D	Capture	MAB469	500 µg
	Detection (Biotin)	BAF469	50 µg
	Protein	469-VD-025	25 µg
Mouse VEGF R1/Flt-1	Capture	AF471	100 µg
	Detection (Biotin)	BAF471	50 µg
	Protein	471-F1-100	100 µg
Human VEGF R2/KDR	Capture	MAB3573-500	500 µg
	Detection (Biotin)	BAF357	50 µg
	Protein	357-KD-050	50 µg
Mouse VEGF R2/Flk-1	Capture	MAB4431-500	500 µg
	Detection (Biotin)	BAF644	50 µg
	Protein	443-KD-050	50 µg
Human VEGF R3/Flt-4	Capture	MAB349-500	500 µg
	Detection (Biotin)	BAM3492	250 µg
	Protein	349-F4-050	50 µg
Mouse VEGF R3/Flt-4	Capture	MAB7431	500 µg
	Detection (Biotin)	BAF743	50 µg
	Protein	743-R3-100	100 µg
Human WIF-1	Capture	MAB1342	500 µg
	Detection (Biotin)	BAM1341	250 µg
	Protein	1341-WF-050	50 µg

For research use only. Not for use in diagnostic procedures.

Analyte	Reagent	Catalog #	Size
Human WISP-1/CCN4	Capture	MAB16271	500 µg
	Detection (Biotin)	BAF1627	50 µg
	Protein	1627-WS-050	50 µg
Human XCL1/Lymphotactin	Capture	MAB6951	500 µg
	Detection (Biotin)	BAF695	50 µg
	Protein	695-LT-025	25 µg
Mouse XCL1/Lymphotactin	Capture	MAB486	500 µg
	Detection (Biotin)	BAF486	50 µg
	Protein	486-LT-025	25 µg

For research use only. Not for use in diagnostic procedures.