

ELISA kits for T cell-secreted cytokines (human)

Our range of kits includes SimpleStep ELISA® kits (SSE) that contain recombinant monoclonal antibodies for specific and sensitive results, and a single-wash protocol to reduce assay time to 90 minutes.



Target	abID	Product Name	SimpleStep	Sensitivity (pg/ml)	Range (pg/ml)	Assay Time	Validated Sample Types								Does not react with	
							CS	S	HP	EP	CP	U	CL	TL		
Amphiregulin	ab222504	Human Amphiregulin ELISA Kit	✓	2.4	18 - 600	1h 30m	●	●		●	●					Mouse
MIP1a	ab214569	Human MIP1a ELISA Kit	✓	0.4	1.56 - 100	1h 30m	●	●	●	●	●					Rat, Others
MIP1b	ab216173	Human MIP1b ELISA Kit	✓	0.96	4.69 - 300	1h 30m	●	●	●	●	●					Mouse, Rat
RANTES	ab174446	Human RANTES ELISA Kit (CCL5)	✓	2	3.125 - 200	1h 30m	●	●	●	●	●					
MIP3a	ab178015	Human MIP3a ELISA Kit (CCL20)	✓	13	39.06 - 2500	1h 30m	●	●	●	●	●					
GM-CSF	ab174448	Human GM-CSF ELISA Kit (CSF2)	✓	1.6	1.4 - 1000	1h 30m	●	●	●	●	●					
IFN γ	ab46048	Human IFN gamma High Sensitivity ELISA Kit		< 0.69	0.78 - 25	2h 20m	●	●	●	●	●					
IFN γ	ab174443	Human IFN gamma ELISA Kit	✓	470	468 - 30000	1h 30m	●	●	●	●	●					
IL-1α	ab178008	Human IL-1α ELISA Kit	✓	0.5	3.1 - 200	1h 30m	●	●	●	●	●					
IL-2	ab46054	Human IL-2 ELISA Kit (Interleukin-2) High Sensitivity		< 0.97	1.87 - 60	2h	●	●	●	●	●					
IL-2	ab174444	Human IL-2 ELISA Kit	✓	9	15.63 - 2000	1h 30m	●	●	●	●	●					Mouse
IL-4	ab46063	Human IL-4 ELISA Kit (Interleukin-4) High Sensitivity		0.31	0.31 - 10	4h	●	●	●	●	●					
IL-4	ab215089	Human IL-4 ELISA Kit	✓	1.08	6.25 - 400	1h 30m	●	●	●	●	●					
IL-5	ab215536	Human IL-5 ELISA Kit	✓	3.3	4.7 - 300	1h 30m	●	●	●	●	●	●				Mouse, Rat, Others
IL-6	ab46042	Human IL-6 ELISA Kit (Interleukin-6) High Sensitivity		< 0.8	1.56 - 50	3h 45m	●	●	●	●	●					
IL-6	ab178013	Human IL-6 ELISA Kit	✓	12	78.12 - 5000	1h 30m	●	●	●	●	●					
IL-9	ab100577	Human IL-9 ELISA Kit (Interleukin-9)		< 100	40 - 30000	4h 45m	●		●	●	●					
IL-10	ab46059	Human IL-10 ELISA Kit (Interleukin-10) High Sensitivity		< 1.3	1.56 - 50	3h	●	●	●	●	●					
IL-10	ab185986	Human IL-10 ELISA Kit (Interleukin-10)	✓	6.6	6.86 - 5000	1h 30m	●	●	●	●	●					
IL-13	ab178014	Human IL-13 ELISA Kit	✓	39	78 - 5000	1h 30m	●	●	●	●	●					
IL-17A	ab216167	Human IL-17A ELISA Kit	✓	1.1	1.56 - 100	1h 30m	●	●	●	●	●					Mouse, Rat
IL-17F	ab100557	Human IL-17F ELISA Kit (Interleukin-17F)		< 20	13.72 - 10000	4h 45m	●	●	●	●	●					
IL-17F	ab119537	Human IL-17F ELISA Kit (Interleukin-17F)		15.5	31.3 - 2000	4h 10m	●	●	●		●					
IL-21	ab119542	Human IL-21 ELISA Kit (Interleukin-21)		20	78 - 5000	4h 10m	●	●		●	●					
IL-22	ab119543	Human IL-22 ELISA Kit (Interleukin-22)		5	31.3 - 2000	4h 10m	●	●	●	●	●					
IL-22	ab216170	Human IL-22 ELISA Kit	✓	2.61	13.12 - 840	1h 30m	●	●	●	●	●	●				Mouse, Rat, Others
IL-25	ab213793	Human IL-17E ELISA Kit (IL-25)		< 10	62.5 - 4000	3h 30m	●	●	●	●	●		●			
TNF β	ab119561	Human TNF beta ELISA Kit		4.6	15.6 - 1000	4h 10m	●	●	●	●	●					
TNF β	ab119576	Human TNF beta ELISA Kit		< 5	62.5 - 4000	3h 25m	●	●	●	●	●			●		
TGF β1	ab100647	Human TGF beta 1 ELISA Kit		< 80	82 - 60000	4h 45m	●	●	●	●	●					
TGF β1	ab108912	Human TGF beta 1 ELISA Kit		30	31 - 2000	5h	●									
TNF α	ab181421	Human TNF alpha ELISA Kit	✓	14	31.25 - 2000	1h 30m	●	●	●	●	●					Mouse
Lymphotactin	ab100584	Human Lymphotactin ELISA Kit		< 150	69 - 50000	4h 45m	●	●	●	●	●					

Key: CS=cell culture supernatant, S=serum, HP=heparin plasma, EP=EDTA plasma, CP=citrate plasma, U=urine, CL=cell lysate, TL=tissue lysate
 All products in the list are coated ELISA kits, and are available as 1x96 tests. Overall, listed ELISA kits have intra- and inter-assay precision lower than 10% (CV).

ELISA kits for T cell-secreted cytokines (mouse)

Our range of kits includes SimpleStep ELISA® kits (SSE) that contain recombinant monoclonal antibodies for specific and sensitive results, and a single-wash protocol to reduce assay time to 90 minutes.



Target	abID	Product Name	SimpleStep	Sensitivity (pg/ml)	Range (pg/ml)	Assay Time	Validated Sample Types								Does not react with	
							CS	S	HP	EP	CP	U	CL	TL		
Amphiregulin	ab100668	Mouse Amphiregulin ELISA Kit		< 2	1.1 - 800	4h 45m	●	●	●	●	●					
MIP1a	ab200017	Mouse MIP1a ELISA Kit (CCL3)	✓	1.95	15.6 - 1000	1h 30m	●	●	●	●	●	●				Human, Others
MIP1b	ab215115	Mouse MIP1b ELISA Kit	✓	2.1	5 - 350	1h 30m	●	●	●	●	●					Rat, Human, Others
RANTES	ab215537	Mouse RANTES ELISA Kit*	✓	5.04	7.81 - 500	1h 30m	●	●								Human, Others
MIP3a	ab100728	Mouse MIP3 alpha ELISA Kit (CCL20)		< 3	2.06 - 1500	4h 45m	●	●	●	●	●					
GM-CSF	ab201276	Mouse GM-CSF ELISA Kit (CSF2)	✓	3.1	12.5 - 800	1h 30m	●									
IFN γ	ab46081	Mouse Interferon gamma ELISA Kit		< 15	31.25 - 1000	3h 45m		●								
IFN γ	ab100689	Mouse Interferon gamma ELISA Kit (IFNG)		< 5	2.74 - 2000	4h 45m	●	●	●	●	●					
IFN γ	ab100690	Mouse Interferon gamma ELISA Kit (IFNG)		< 10	2.74 - 2000	4h 45m							●	●		
IL-1a	ab199076	Mouse IL-1a ELISA Kit	✓	0.62	4.69 - 300	1h 30m	●	●				●				
IL-2	ab100706	Mouse IL-2 ELISA Kit (Interleukin-2)		< 1	0.625 - 40	4h 45m	●	●	●	●	●					
IL-2	ab223588	Mouse IL-2 ELISA Kit	✓	3.7	15.625 - 1000	1h 30m	●	●		●	●					Rat, Human
IL-4	ab100710	Mouse IL-4 ELISA Kit (Interleukin-4)		< 1	0.819 - 200	4h 45m							●	●		
IL-4	ab221833	Mouse IL-4 ELISA Kit	✓	6.3	18.75 - 1200	1h 30m	●	●					●			
IL-5	ab204523	Mouse IL-5 ELISA Kit	✓	1.2	3.9 - 250	1h 30m	●	●	●	●	●					Human, Others
IL-6	ab100712	Mouse IL-6 ELISA Kit (Interleukin-6)		< 2	0.82 - 600	4h 45m	●	●	●	●	●				●	
IL-6	ab100713	Mouse IL-6 ELISA Kit (Interleukin-6)		< 2	0.82 - 600	4h 45m							●	●		
IL-6	ab222503	Mouse IL-6 ELISA Kit*	✓	11.3	15.6 - 1000	1h 30m	●	●	●	●	●					Human
IL-9	ab222505	Mouse IL-9 ELISA Kit	✓	1.67	3.1 - 200	1h 30m	●	●	●		●	●				Rat, Human, Others
IL-10	ab46103	Mouse IL-10 ELISA kit (Interleukin-10)		< 20	31.25 - 1000	3h 45m	●	●			●					
IL-10	ab100697	Mouse IL-10 ELISA kit (Interleukin-10)		< 45	78.13 - 5000	4h 45m	●	●	●	●	●					
IL-10	ab108870	Mouse IL-10 ELISA kit (Interleukin-10)		30	31 - 2000	5h	●	●			●				●	
IL-13	ab219634	Mouse IL-13 ELISA Kit	✓	1.2	1.95 - 125	1h 30m	●	●	●	●	●					Rat, Human
IL-17A	ab199081	Mouse IL-17A ELISA Kit	✓	0.5	6.25 - 400	1h 30m	●	●			●					
IL-17F	ab204522	Mouse IL-17F ELISA Kit	✓	0.1	9.375 - 600	1h 30m	●	●	●	●	●					Rat, Human, Others
IL-21	ab100707	Mouse IL-21 ELISA Kit (Interleukin-21)		< 6	8.19 - 2000	4h 45m	●		●	●	●					
IL-21	ab171180	Mouse IL-21 ELISA Kit (Interleukin-21)		< 10	62.5 - 4000	3h 55m	●	●								
TGF β	ab119557	Mouse TGF beta 1 ELISA Kit		7.8	31.3 - 2000	5h	●	●		●	●					
TNF α	ab208348	Mouse TNF alpha ELISA Kit	✓	9.1	46.88 - 3000	1h 30m	●	●	●	●	●					
Lymphotoxin	ab100720	Mouse Lymphotoxin ELISA Kit (SCM-1)		< 40	46.9 - 3000	4h 45m	●	●	●	●	●					

Key: CS=cell culture supernatant, S=serum, HP=heparin plasma, EP=EDTA plasma, CP=citrate plasma, U=urine, CL=cell lysate, TL=tissue lysate
 All products in the list are coated ELISA kits, and are available as 1x96 tests. Overall, listed ELISA kits have intra- and inter-assay precision lower than 10% (CV).

ELISA kits for T cell-secreted cytokines (rat)

Our range of kits includes SimpleStep ELISA® kits (SSE) that contain recombinant monoclonal antibodies for specific and sensitive results, and a single-wash protocol to reduce assay time to 90 minutes.



Target	abID	Product Name	SimpleStep	Sensitivity (pg/ml)	Range (pg/ml)	Assay Time	Validated Sample Types								Does not react with	
							CS	S	HP	EP	CP	U	CL	TL		
MIP1a	ab213916	Rat MIP-1 alpha ELISA Kit (CCL3)		< 1	7.8 - 500	3h 30m	●	●	●	●	●					
MIP1b	ab171336	Rat MIP-1 beta ELISA Kit (CCL4)		< 10	15.6 - 1000	3h 25m	●		●	●	●					
RANTES	ab213907	Rat Rantes ELISA Kit		< 1	15.6 - 1000	3h 30m	●	●	●	●	●					
MIP3a	ab213904	Rat MIP-3 alpha ELISA Kit (CCL20)		< 1	7.8 - 500	3h 30m	●	●	●	●	●					
GM-CSF	ab100762	Rat GM-CSF ELISA Kit		< 10	12.35 - 3000	4h 45m	●	●	●	●	●					
IFN γ	ab46107	Rat IFN gamma ELISA Kit		< 10	31.25 - 1000	3h 45m	●	●								
IFN γ	ab113349	Rat IFN gamma ELISA Kit		< 100	123 - 30000	4h 45m	●		●	●	●					
IL-1A	ab113350	Rat IL-1 alpha ELISA Kit (Interleukin-1 alpha)		< 15	11 - 8000	4h 45m	●	●	●	●	●					
IL-2	ab221834	Rat IL-2 ELISA Kit	✓	4.9	31.25 - 2000	1h 30m	●	●								Mouse
IL-3	ab46073	Rat IL-4 ELISA Kit (Interleukin-4)		< 15	31.25 - 1000	4h 45m	●	●								
IL-4	ab100770	Rat IL-4 ELISA Kit (Interleukin-4)		< 1.5	0.66 - 160	4h 45m	●	●	●	●	●					
IL-5	ab100771	Rat IL-4 ELISA Kit (Interleukin-4)		< 1.5	0.66 - 160	4h 45m							●	●		
IL-6	ab100772	Rat IL-6 ELISA Kit (Interleukin-6)		< 30	40.96 - 10000	4h 45m	●		●	●	●					
IL-7	ab119548	Rat IL-6 ELISA Kit (Interleukin-6)		12	31 - 2000	3h 10m	●	●								
IL-10	ab100765	Rat IL-10 ELISA Kit (Interleukin-10)		< 10	8.23 - 6000	4h 45m							●	●		
IL-10	ab214566	Rat IL-10 ELISA Kit	✓	4.89	31.3 - 2000	1h 30m	●	●			●					Mouse, Human, Others
IL-13	ab100766	Rat IL-13 ELISA Kit (Interleukin-13)		< 20	27.43 - 20000	4h 45m							●	●		
IL-17A	ab214028	Rat IL-17A ELISA Kit	✓	1.1	6.25 - 400	1h 30m	●	●			●					Human
IL-17F	ab213913	Rat IL-17F ELISA Kit		< 10	31.2 - 2000	3h 30m	●	●	●	●	●		●	●		
TGF β1	ab119558	Rat TGF beta 1 ELISA Kit		8	31.3 - 2000	5h	●	●	●	●	●				●	
TNF α	ab46070	Rat TNF alpha ELISA Kit		< 15	31.25 - 1000	3h 45m	●	●								
TNF α	ab100784	Rat TNF alpha ELISA Kit		< 25	82.3 - 20000	4h 45m	●	●	●	●	●					
TNF α	ab100785	Rat TNF alpha ELISA Kit		< 25	82.3 - 20000	4h 45m	●						●	●		
TNF α	ab108913	Rat TNF alpha ELISA Kit		15	16 - 1000	4h 30m	●									

Key: CS=cell culture supernatant, S=serum, HP=heparin plasma, EP=EDTA plasma, CP=citrate plasma, U=urine, CL=cell lysate, TL=tissue lysate
 All products in the list are coated ELISA kits, and are available as 1x96 tests. Overall, listed ELISA kits have intra- and inter-assay precision lower than 10% (CV).