

ELISA kits for blood coagulation – 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 name	abID	SimpleStep	Sensitivity (pg/ml)	Range (pg/ml)	Assay Time	Validated Sample Types							
						CS	S	HP	EP	CP	U	Sv	M
Alpha 2 Macroglobulin	ab108883		2	2-500	4h	●						●	●
Alpha 2 Macroglobulin	ab108888		1000	1250-40000	3h	●	●	●	●	●			
Thrombin	ab108909		0.3	0.313-20	4h	●							
Tissue Factor	ab220653	✓	0.0036	0.01563-1	1h 30m	●		●	●	●	●	●	
Factor V	ab137976		0.9	0.938-60	4h	●		●	●	●	●		
Factor VII	ab108829		6	6.25-200	4h	●	●	●	●	●		●	
Factor VII	ab190810	✓	0.292	0.315-20	1h 30m		●	●	●	●			
Factor IX	ab108831		0.4	1.56-100	4h	●	●	●	●	●			●
Factor IX	ab188393	✓	0.23	0.78-50	1h 30m		●	●	●	●			
Factor X	ab108832		1.5	1.563-100	4h	●	●	●	●	●	●	●	●
Factor XI	ab108834		1.5	1.56-50	4h	●	●	●	●	●			
Factor XII	ab108835		0.02	6.25-100	4h	●	●	●	●	●	●		●
Factor XII	ab192144	✓	0.7	0.94-60	1h 30m	●	●	●	●	●			
Factor XIII	ab108836		2.1	5-80	4h	●	●	●	●	●	●	●	●
Plasma kallikrein	ab171015		1.2	1.25-80	4h	●	●	●	●	●		●	●
Plasma kallikrein	ab202405	✓	0.154	0.47-30	1h 30m	●	●	●	●	●			
Kininogen	ab108875		100	125-4000	3h	●	●	●	●	●			
Kininogen	ab108876		0.7	1.563-100	4h	●					●	●	●
Kininogen	ab189574	✓	0.0087	0.02344-1.5	1h 30m	●	●	●	●	●			
tPA / Tissue Plasminogen Activator	ab108914		0.03	0.031-2	4h	●	●	●	●	●	●	●	●
tPA / Tissue Plasminogen Activator	ab190812	✓	0.0035	0.078-5	1h 30m	●	●	●	●	●			
PAI1 + tPA	ab192151	✓	0.016	0.039-2.5	1h 30m	●	●	●	●	●	●		
uPA	ab108917		0.2	0.25-8	4h	●	●	●	●	●	●		
uPA	ab119611		0.005	0.0625-4		●	●	●	●				
Angiostatin	ab99973		20	31.25-2000		●	●	●					
Plasminogen	ab196262	✓	0.165	0.468-30	1h 30m	●	●	●	●	●	●	●	
Protein C	ab137987		3	3.13-200	4h	●					●	●	●
Protein C	ab137988		60	94-6000	3h		●	●	●	●			
Protein S	ab125969		250	250-8000	4h	●	●	●	●	●			
Protein S	ab190808	✓	0.4	3.12-200	1h 30m		●	●	●	●			
Alpha 1 Antitrypsin	ab108798		18	39-40000	3h		●	●	●	●		●	●
Alpha 1 Antitrypsin	ab108799		0.35	0.391-100	4h	●	●	●			●	●	●
Alpha 1 Antitrypsin	ab189579	✓	13	31.25-2000	1h 30m	●		●		●			
Thrombin-Antithrombin Complex	ab108907		1	1.481-120	4h	●	●	●	●	●			●
Antithrombin III	ab220073		30	63-4000	2h 40m		●	●	●	●			
Antithrombin III	ab222507	✓	0.33	3.13-200	1h 30m	●	●	●	●	●	●		
PAI1	ab184863	✓	0.021	0.09375-6	1h 30m	●	●	●	●	●			
TFPI / Tissue Factor Pathway Inhibitor	ab108904		0.12	0.313-20	5h	●	●	●	●	●			●
TFPI / Tissue Factor Pathway Inhibitor	ab188392	✓	0.0149	0.0781-5	1h 30m	●	●	●	●	●	●		
Thrombomodulin	ab46508		0.31	0.625-20	1h 45m	●	●	●	●	●			
Thrombomodulin	ab214029	✓	0.003	0.02-1.3	1h 30m	●	●	●	●	●	●		
VWF / Von Willebrand Factor	ab223864	✓	0.079	0.469-30	1h 30m		●	●	●	●			

Key: CS=cell culture supernatant, S=serum, HP=heparin plasma, EP=EDTA plasma, CP=citrate plasma, U=urine, Sv=saliva, M=milk
 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 blood coagulation

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.

Mouse



Target name	abID	SimpleStep	Sensitivity (pg/ml)	Range (pg/ml)	Assay Time	Validated Sample Types					
						CS	S	HP	EP	CP	U
Alpha 2 Macroglobulin	ab108887		0.07	0.078-20	4h	●	●	●	●	●	●
Prothrombin	ab157526		0.386	1-500		●	●		●	●	
Prothrombin / Thrombin Total	ab157527		0.52	1-500	1h 50m	●	●		●	●	
Tissue Factor	ab214091	✓	0.0037	0.025-1.6	1h 30m	●					●
uPA Total	ab198512		0.01	0.01-10					●	●	
Plasminogen Total	ab198511		0.38	1-500		●	●	●	●	●	●
Alpha 1 Antitrypsin	ab205088		1.492	6.25-200	2h 10m		●	●	●	●	●
Antithrombin III	ab108800		3	3.13-100	4h	●	●	●	●	●	●
Thrombin-Antithrombin Complex	ab137994		0.01	0.031-8	5h	●		●	●	●	
PAI1 Total	ab157529		0.006	0.05-50	1h 50m		●			●	
PAI1	ab197752	✓	0.2	0.298-5	1h 30m	●	●			●	
PAI1	ab201283 *	✓	0.159	0.298-5	1h 30m	●	●			●	●
TFPI / Tissue Factor Pathway Inhibitor	ab217776	✓	0.0017	0.019-1.2	1h 30m	●	●	●	●	●	●
Thrombomodulin	ab209880	✓	0.0175	0.094-6	1h 30m	●	●	●	●	●	●
VWF / Von Willebrand Factor	ab208980	✓	0.037	0.11-7	1h 30m	●	●	●	●	●	

*also reacts with rat

Rat



Target name	abID	SimpleStep	Sensitivity (pg/ml)	Range (pg/ml)	Assay Time	Validated Sample Types					
						CS	S	HP	EP	CP	U
Alpha 2 Macroglobulin	ab157730		2.23	6.25-400		●	●	●	●		
Kininogen	ab157742		4.971	6.25-400		●	●	●	●		
tPA / Tissue Plasminogen Activator	ab198510		0.043	0.05-50					●	●	
Plasminogen	ab157740		3.24	3.13-200		●	●	●	●		
PAI1	ab198509		0.079	0.1-50			●			●	
PAI1	ab201283 †	✓	0.159	0.298-5	1h 30m	●	●			●	●

†also reacts with mouse

Key: CS=cell culture supernatant, S=serum, HP=heparin plasma, EP=EDTA plasma, CP=citrate plasma, U=urine, Sv=saliva, M=milk
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).