

# ELISA kits for Ras signalling pathway (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.

Kits validated on cell culture supernatant (CS), serum (S), heparin plasma (HP), EDTA plasma (EP), citrate plasma (CP), urine (U), saliva (Sv) and milk (M)

Target Name	abID	SimpleStep	Sensitivity (ng/mL)	Range (ng/mL)	Assay Time	validated samples							
						CS	S	HP	EP	CP	U	Sv	M
Angiopoietin 1	<a href="#">ab99972</a>		< 0.03	0.074 - 18	5h	•	•	•	•	•			
Angiopoietin 2	<a href="#">ab99971</a>		< 0.01	0.004 - 3	5h	•	•	•	•	•			
Angiopoietin 4	<a href="#">ab193693</a>		0.25	0.25 - 200	5h	•	•	•	•	•			
C-Peptide	<a href="#">ab178641</a>		0.01	0.7 - 1.9	2h 30m		•	•	•	•			
c-kit	<a href="#">ab45924</a>		< 0.095	0.312 - 10	1h 45	•	•	•	•	•			
EGF	<a href="#">ab100504</a>		< 0.001	0.0008 - 0.2	5h	•	•	•	•	•			
EGF	<a href="#">ab217772</a>	✓	0.001	0.0015 - 0.1	1h 30m	•		•	•	•	•	•	•
EGFR	<a href="#">ab100505</a>		< 0.004	0.0027 - 2	5h	•	•	•	•	•			
soluble EGFR	<a href="#">ab193764</a>	✓	0.001	0.078 - 5	1h 30m	•	•	•	•	•			
Fas Ligand	<a href="#">ab100515</a>		< 0.002	0.0013 - 1	5h	•	•	•	•	•			
Fas Ligand	<a href="#">ab45892</a>		0.012	0.062 - 2	3h 45	•	•	•	•	•			
FGF1*	<a href="#">ab219636</a>	✓	0.00197	0.019 - 1.2	1h 30m	•	•				•	•	
FGF4	<a href="#">ab100516</a>		< 0.05	0.025 - 18	5h	•		•	•	•			
FGF6	<a href="#">ab100518</a>		< 0.006	0.008 - 6	5h	•		•	•	•			
FGF7	<a href="#">ab100519</a>		< 0.005	0.055 - 0.4	5h	•	•	•	•	•			
FGF7	<a href="#">ab183362</a>	✓	0.0072	0.031 - 2	1h 30m		•	•	•	•			
FGF9	<a href="#">ab100520</a>		< 0.2	0.358 - 35	5h	•		•	•	•			
FGF19	<a href="#">ab193703</a>		0.03	0.03 - 8	5h		•	•	•	•			
FGF21	<a href="#">ab125966</a>		0.03	0.03 - 2	4h	•	•	•	•	•			
FGF21	<a href="#">ab222506</a>	✓	0.00033	0.031 - 2	1h 30m		•	•	•	•			•
FGF basic	<a href="#">ab99979</a>		< 0.05	0.102 - 10	5h	•	•	•	•	•			
acid FGF	<a href="#">ab193691</a>		0.012	0.012 - 4	5h	•	•	•	•	•			
HGF	<a href="#">ab100534</a>		< 0.003	0.027 - 2	5h	•	•	•	•	•			
IGF1	<a href="#">ab211651</a>	✓	0.051	0.008 - 5	1h 30m	•	•	•	•	•		•	
IGF1	<a href="#">ab100545</a>		< 0.2	0.1 - 30	5h	•	•	•	•	•			
IGF 1	<a href="#">ab108873</a>		> 0.3	0.375 - 24	5h	•	•	•	•	•			
IGF1 R	<a href="#">ab100546</a>		< 0.006	0.008 - 6	5h	•	•	•	•	•			
Insulin	<a href="#">ab200011</a>	✓	<b>1.1 pmol/l</b>	<b>2.1 – 133.6 pmol/l</b>			•	•					
Insulin	<a href="#">ab100578</a>		<b>&lt; 28 pmol/l</b>	<b>32 - 2083 pmol/l</b>	5h	•	•	•	•	•			
Proinsulin	<a href="#">ab213977</a>		<b>0.17 pmol/l</b>	<b>1.6 - 50 pmol/l</b>	3h		•		•				
M-CSF	<a href="#">ab100590</a>		< 0.005	0.0028 - 2	5h	•	•	•	•	•			
PDGF BB	<a href="#">ab184860</a>	✓	0.002	0.002 - 1.5	1h 30m		•	•	•	•			
PDGF BB	<a href="#">ab100624</a>		< 0.001	0.0005 - 0.4	5h	•	•	•	•	•			

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). \* Reacts with mouse - \*\* Reacts with mouse and rat – SQ: semi quantitative

# ELISA kits for Ras signalling pathway (human) (continued)

Kits validated on cell culture supernatant (CS), serum (S), heparin plasma (HP), EDTA plasma (EP), citrate plasma (CP), urine (U), saliva (S) and milk (M)

Target Name	abID	SimpleStep	Sensitivity (ng/mL)	Range (ng/mL)	Assay Time	validated samples							
						CS	S	HP	EP	CP	U	Sv	M
PDGF AA	<a href="#">ab100622</a>		<0.04	0.04 - 30	5h	•	•	•	•	•			
PDGF AB	<a href="#">ab100623</a>		< 0.01	0.01 - 2.5	5h	•	•	•	•	•			
PDGF receptor beta	<a href="#">ab100626</a>		< 0.025	0.025 - 18	5h	•	•	•	•	•			
PDGF receptor alpha	<a href="#">ab155438</a>		< 0.4	0.49 - 120	5h	•	•	•	•	•			
PLGF	<a href="#">ab100629</a>		< 0.002	0.0013 - 1	5h	•	•	•	•	•			
SCF	<a href="#">ab100636</a>		< 0.002	0.0027 - 2	5h	•	•	•	•	•			
SCF	<a href="#">ab176109</a>		0.0005	0.016 - 1	5h		•	•	•	•			
SCF	<a href="#">ab108901</a>		0.1	0.156 - 40	5h	•	•	•	•	•			
TIE2	<a href="#">ab100650</a>		< 0.020	0.021 - 15	5h	•	•	•	•	•			
VEGF	<a href="#">ab100662</a>		< 0.010	0.008 - 6	5h	•	•	•	•	•			
VEGF	<a href="#">ab222510</a>		0.0027	0.012 - 0.8				•	•	•	•	•	•
VEGFA	<a href="#">ab119566</a>		0.0079	0.016 - 1		•	•	•	•	•			
VEGFC	<a href="#">ab100664</a>		< 0.015	0.014 - 10	3h 30m	•	•	•	•	•			
VEGF R1	<a href="#">ab119567</a>		0.03	0.16 - 10	3h 30m	•	•	•	•	•			
VEGF R2	<a href="#">ab195210</a>	✓	0.146	0.391 - 25	1h 30m	•	•	•	•	•			
VEGF R2	<a href="#">ab119613</a>		< 0.03	0.156 - 10	3h 30m	•	•	•	•	•			
VEGF R3	<a href="#">ab213476</a>	✓	0.01	0.117 - 7.5	1h 30m	•	•	•	•	•			
VEGF R3	<a href="#">ab100665</a>		< 0.07	0.034 - 25	5h	•	•	•	•	•			
VEGF R3	<a href="#">ab100666</a>		< 0.015	0.012 - 3	5h	•	•	•	•	•			

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). \* Reacts with mouse - \*\* Reacts with mouse and rat – SQ: semi quantitative

# ELISA kits for Ras signalling pathway (human) (continued)

Kits validated on cell lysate (CL), tissue lysate (TL), cell culture extract (CE), tissue extract (TE), cell culture media (CM)

Target Name	abID	SimpleStep	Sensitivity (ng/mL)	Range (ng/mL)	Assay Time	Validated samples				
						CL	TL	CE	TE	CM
AKT 1/2/3 pS473 + AKT1** (SQ)	ab176657	✓	0.03		1h 30m	•	•			
AKT 1/2/3 pT308 + AKT1* (SQ)	ab176658	✓	0.03		1h 30m	•	•			
akt (ps473) + total akt** (SQ)	ab126433				5h	•				
AKT 1/2/3 (pS473)*	ab176635	✓	0.03	1 - 100	1h 30m	•	•			
AKT 1/2/3 (pT308)*	ab176636	✓	0.03	1 - 100	1h 30m	•	•			
Akt (pS473)** (SQ)	ab126432				5h	•				
AKT1*	ab176637	✓	0.03	0.5 - 5	1h 30m	•	•			
AKT1	ab214023	✓	0.027	0.185 - 12	1h 30m			•	•	
AKT2**	ab208986	✓	0.005	0.078 - 5	1h 30m			•	•	
AKT3	ab216946	✓	0.041	0.625 - 40	1h 30m	•		•	•	
Soluble EGFR	ab193764	✓	0.001	0.078 - 5	1h 30m					•
EGFR (pY992) (SQ)	ab126442				5h		•			
EGFR (pY1045) (SQ)	ab126436				5h		•			
EGFR (pY1068) (SQ)	ab126438				5h		•			
EGFR (pS1070) (SQ)	ab126434				5h		•			
EGFR (pY1086) (SQ)	ab126440				5h		•			
EGFR (pY992) + total EGFR (SQ)	ab126443				5h		•			
EGFR (pY1068) + total EGF (SQ)	ab126439				5h		•			
EGFR (pY1045) + total EGFR (SQ)	ab126437				5h		•			
EGFR (pS1070) + total EGFR (SQ)	ab126435				5h		•			
EGFR (pY1086) + total EGFR (SQ)	ab126441				5h		•			
EGFR (pY992) + total EGFR (SQ)	ab126443				5h		•			
EphA2	ab219051	✓	0.0009	0.009 - 0.6	1h 30m					•
ERK 1/2 (pT202/Y204 + Total)* (SQ)	ab176660	✓		ERK1/2 (pT202/Y204): 100 ERK 1/2 total: 600	1h 30m	•	•			
ERK 1/2 (pT202/Y204)**	ab176640	✓	0.1	0.2 - 20	1h 30m	•	•			
ERK1/2**	ab176641	✓	0.6	0.6 - 60	1h 30m	•	•			
Erk1 (pT202/pY204) + Erk2 (pT185/pY187) + total Erk1/2** (SQ)	ab126445				5h	•				

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). \* Reacts with mouse - \*\* Reacts with mouse and rat – SQ: semi quantitative

# ELISA kits for Ras signalling pathway (human) (continued)

Kits validated on cell cell lysate (CL), tissue lysate (TL), cell culture extract (CE), tissue extract (TE), cell culture media (CM)

Target Name	abID	SimpleStep	Sensitivity (ng/mL)	Range (ng/mL)	Assay Time	validated sample types				
						CL	TL	CE	TE	CM
Erk1 (pT202/pY204) + Erk2 (pT185/pY187)**	ab126444				5h	•				
ERK1	ab215545	✓	0.0129	0.094 - 6	1h 30m			•	•	
FGF1**	ab219636	✓	0.00197	0.019 - 1.2	1h 30m				•	•
FGF4	ab100517		< 0.05	0.025 - 18	5h			•	•	
FGFR3	ab214027	✓	0.061	0.312 - 20	1h 30m			•	•	
FGF basic	ab99980		< 0.05	0.102 - 10	5h				•	•
HGF	ab100535		< 0.008	0.008 - 6	5h			•	•	
JNK 1/2 (pT183/Y185 + Total)* (SQ)	ab176662	✓	JNK1/2 (pT183/Y185): 0.5 JNK1/2 (Total): 0.1		1h 30m		•	•		
JNK 1/2 (pT183/Y185)*	ab176645	✓	0.5	1 - 100	1h 30m		•	•		
JNK 1/2*	ab176646	✓	0.1	0.2 - 20	1h 30m		•	•		
JNK2	ab218265	✓	0.081	0.234 - 15					•	•
Met (pY1234/pY1235) + total Met** (SQ)	ab126451				5h			•	•	
Met (pY1234/pY1235)* (SQ)	ab126450				5h			•		
Mek1 (pS217/221) (SQ)	ab126446				5h			•		
Mek1 (pS217/221) + total Mek1 (SQ)	ab126448				5h			•		
NF kappaB p65 (pS536 + Total)* (SQ)	ab176663	✓	0.5		1h 30m		•	•		
NF kappaB p65*	ab176648	✓	5000	5000 - 600000	1h 30m		•	•		
NF kappaB p65 (pS536)*	ab176647	✓	5000	5000 - 600000	1h 30m		•	•		
PDGF BB	ab184860	✓	0.002	0.002 - 1.5	1h 30m					•
PDGF BB	ab100625		< 0.002	0.0005 - 0.5	5h			•	•	
SCF	ab176109		0.0005	0.016 - 1	5h					
VEGF	ab100663		< 0.010	0.008 - 6	5h			•	•	
ZAP70	ab173192	✓	< 0.5	1.6 - 100	1h 30m		•	•	•	

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). \* Reacts with mouse - \*\* Reacts with mouse and rat – SQ: semi quantitative

# ELISA kits for Ras signalling pathway (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.

Kits validated on cell culture supernatant (CS), serum (S), heparin plasma (HP), EDTA plasma (EP), citrate plasma (CP), urine (U), saliva (S) and milk (M)

Target Name	abID	SimpleStep	Sensitivity (ng/mL)	Range (ng/mL)	Assay Time	validated samples							
						CS	S	HP	EP	CP	U	Sv	M
AKT2**	<a href="#">ab208986</a>	✓	0.005	0.078 - 5	1h 30m	•							
Angiopoietin 2***	<a href="#">ab209883</a>	✓	0.017	0.07 - 4.4	1h 30m	•	•	•	•	•			
Angiopoietin 2	<a href="#">ab171335</a>		<0.01	0.015 - 10	3h30m	•	•	•	•		•		
EGF	<a href="#">ab100679</a>		<0.003	0.003 - 0.8	5h	•	•	•	•	•			
EGFR	<a href="#">ab201275</a>	✓	0.06	1.875 - 120	1h 30m	•	•	•	•	•			
EphA2	<a href="#">ab207622</a>	✓	0.012	0.023 - 1.5	1h 30m	•	•	•	•	•			
FGF basic	<a href="#">ab100670</a>		<0.008	0.019 - 0.6	5h	•	•	•	•	•	•		
FGF1***	<a href="#">ab223587</a>	✓		0.016 - 1	1h 30m	•	•	•	•	•	•		
FGF1**	<a href="#">ab219636</a>	✓	0.00197	0.019 - 1.2	1h 30m	•	•				•	•	
FGF-21	<a href="#">ab212160</a>	✓	0.001	0.009 - 0.6	1h 30m	•	•	•	•				
FGF23	<a href="#">ab213863</a>		<0.01	0.015 - 1	3h30m	•	•	•	•				
HGF	<a href="#">ab100686</a>		<0.4	0.614 - 60	5h	•	•	•	•	•			
HGF***	<a href="#">ab223862</a>	✓	0.008	0.062 - 4	1h 30m		•	•	•	•			
IGF1	<a href="#">ab100695</a>		<0.004	0.003 - 2	4h45m	•	•	•	•				
IGF1	<a href="#">ab108874</a>		0.1	0.156 - 10	5h	•	•	•	•				
M-CSF	<a href="#">ab199084</a>	✓	0.004	0.047 - 3	1h 30m	•	•				•		
M-CSF	<a href="#">ab155457</a>		<0.001	0.001 - 0.35	5h	•	•	•	•	•			
M-CSFR	<a href="#">ab213858</a>		<0.01	0.062 - 4	3h 30m	•	•	•	•				
NGF R***	<a href="#">ab205569</a>	✓	0.002	0.01 - 0.7	1h 30m	•	•	•	•	•			
PDGF BB	<a href="#">ab224879</a>	✓			1h 30m	•	•	•	•	•			
PLGF	<a href="#">ab197748</a>	✓	0.0005	0.005 - 0.3	1h 30m	•	•				•		
PLGF-2	<a href="#">ab155459</a>		<0.001	0.002 - 0.4	5h	•	•	•	•	•			
SCF***	<a href="#">ab100470</a>		<0.005	0.003 - 2	5h	•	•	•	•	•			
SCF	<a href="#">ab197750</a>	✓	0.001	0.015 - 1	1h 30m	•	•	•	•	•			
VEGF	<a href="#">ab100751</a>		<0.002	0.004 - 1	5h	•	•	•	•				
VEGF	<a href="#">ab209882</a>	✓	0.0003	0.006 - 0.37	1h 30m	•							
VEGF R3	<a href="#">ab203367</a>	✓	0.018	0.047 - 3	1h 30m	•	•	•	•	•			
VEGF A	<a href="#">ab119565</a>		0.02	0.062 - 4	4h 30m	•	•	•	•	•			
VEGF-B	<a href="#">ab213897</a>		<0.002	0.031 - 2	3h 30m	•	•	•					
VEGF	<a href="#">ab209882</a>	✓	0.0003	0.006 - 0.37	1h 30m	•							

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). \* Reacts with mouse - \*\* Reacts with mouse and rat – SQ: semi quantitative

# ELISA kits for Ras signalling pathway (mouse) (continued)

Kits validated on cell lysate (CL), tissue lysate (TL), cell culture extract (CE), tissue extract (TE), cell culture media (CM)

Target Name	abID	SimpleStep	Sensitivity (ng/mL)	Range (ng/mL)	Assay Time	Validated samples				
						CL	TL	CE	TE	CM
Akt (pS473)** (SQ)	ab126432				5h	•				
Akt (pS473) + total Akt** (SQ)	ab126433				5h	•				
AKT 1/2/3 (pS473)* (SQ)	ab176635	✓	0.03	1 - 100	1h 30m	•	•			
AKT 1/2/3 (pT308)* (SQ)	ab176636	✓	0.03	1 - 100	1h 30m	•	•			
AKT 1/2/3 pS473 + AKT1** (SQ)	ab176657	✓	0.03		1h 30m	•	•			
AKT 1/2/3 pT308 + AKT1* (SQ)	ab176658	✓	0.03		1h 30m	•	•			
AKT1* (SQ)	ab176637	✓	0.03	0.5 - 5	1h 30m	•	•			
AKT2**	ab208986	✓	0.005	0.078 - 5	1h 30m			•	•	
Bad	ab214089	✓	0.018	0.078 - 5	1h 30m			•	•	
c-Kit	ab203361	✓	0.03	0.23 - 15	1h 30m			•	•	
ERK 1/2 (pT202/Y204 + Total)* (SQ)	ab176660	✓	ERK1/2 (pT202/Y204): 100 ERK 1/2 total: 600		1h 30m	•	•			
Erk1 (pT202/pY204) + Erk2 (pT185/pY187)** (SQ)	ab126444				5h	•				
Erk1 (pT202/pY204) + Erk2 (pT185/pY187) + total Erk1/2** (SQ)	ab126445				5h	•				
ERK1/2* (SQ)	ab176641	✓	0.6	0.6 - 60	1h 30m					
FGF basic	ab100670		<0.008	0.019 - 0.6	5h					
FGF1**	ab219636	✓	0.00197	0.019 - 1.2	1h 30m				•	
FGF1***	ab223587	✓		0.016 - 1	1h 30m		•			
FGF23	ab213863		<0.01	0.015 - 1	3h30m		•			
HGF	ab100687		<0.1	0.25 - 60	5h			•	•	
HGF***	ab223862	✓	0.008	0.062 - 4	1h 30m			•	•	
IGF1	ab100695		<0.004	0.003 - 2	5h		•			
JNK 1/2 (pT183/Y185 + Total)* (SQ)	ab176662	✓	JNK1/2 (pT183/Y185): 0.5 JNK1/2 (Total): 0.1		1h 30m	•	•			
JNK 1/2 (pT183/Y185)* (SQ)	ab176645	✓	0.5	1 - 100	1h 30m	•	•			
M-CSF	ab199084	✓	0.004	0.047 - 3	1h 30m			•		
M-CSFR	ab213858		<0.01	0.062 - 4	3h 30m	•	•			
Mek1 (pS217/221)** (SQ)	ab126446				5h	•				
Mek1 (pS217/221) + total Mek1** (SQ)	ab126448				5h	•				
Met (pY1234/pY1235)* (SQ)	ab126450				5h	•				

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). \* Reacts with mouse - \*\* Reacts with mouse and rat – SQ: semi quantitative

## ELISA kits for Ras signalling pathway (mouse) (continued)

Kits validated on cell lysate (CL), tissue lysate (TL), cell culture extract (CE), tissue extract (TE), cell culture media (CM)

Target Name	abID	SimpleStep	Sensitivity (ng/mL)	Range (ng/mL)	Assay Time	validated sample types				
						CL	TL	CE	TE	CM
Met (pY1234/pY1235) + total Met** (SQ)	<a href="#">ab126451</a>				5h	•	•			
NF kappaB p65* (SQ)	<a href="#">ab176648</a>	✓	5000	5000 - 600000	1h 30m	•	•			
NF kappaB p65 (pS536 + Total)* (SQ)	<a href="#">ab176663</a>	✓	0.5		1h 30m	•	•			
NF kappaB p65 (pS536)* (SQ)	<a href="#">ab176647</a>	✓	5000	5000 - 600000	1h 30m	•	•			
NGF R***	<a href="#">ab205569</a>	✓	0.002	0.01 - 0.7	1h 30m			•	•	
SHP-2**	<a href="#">ab206309</a>	✓	0.24	0.63 - 40	1h 30m			•	•	
VEGF	<a href="#">ab209882</a>	✓	0.0003	0.006 - 0.37	1h 30m			•	•	
VEGF	<a href="#">ab100752</a>		<0.002	0.004 - 1	5h			•	•	

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). \* Reacts with mouse - \*\* Reacts with mouse and rat – SQ: semi quantitative

## ELISA kits for Ras signalling pathway (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.

Kits validated on cell culture supernatant (CS), serum (S), heparin plasma (HP), EDTA plasma (EP), citrate plasma (CP), urine (U), saliva (Sv) and milk (M)

Target Name	abID	SimpleStep	Sensitivity (ng/mL)	Range (ng/mL)	Assay Time	validated samples							
						CS	S	HP	EP	CP	U	Sv	M
AKT2**	<a href="#">ab208986</a>	✓	0.005	0.078 - 5	1h 30m	•							
Angiopoietin-1	<a href="#">ab213920</a>		<0.01	0.16 - 10	3h 30m	•	•	•	•				
EGF	<a href="#">ab193734</a>		0.008	0.008 - 0.2	5h	•	•	•	•	•			
Fas Ligand	<a href="#">ab100759</a>		<0.09	0.082 - 60		•		•	•	•			
FGF1	<a href="#">ab213914</a>		<0.01	0.032 - 2		•	•	•	•				
FGF1*	<a href="#">ab223587</a>	✓		0.016 - 1	1h 30m	•	•	•	•	•	•		
FGF1**	<a href="#">ab219636</a>	✓	0.00197	0.019 - 1.2	1h 30m	•	•				•	•	
FGF-21	<a href="#">ab223589</a>	✓	0.007	0.031 - 2	1h 30m	•	•	•	•	•			
HGF*	<a href="#">ab223862</a>	✓	0.008	0.062 - 4	1h 30m		•	•	•	•			
IGF1	<a href="#">ab213902</a>		<0.005	0.063 - 4	3h 30m	•	•	•	•				
PDGF-AA	<a href="#">ab155464</a>		<0.09	0.094 - 6	5h		•	•	•	•	•		
PDGF-AB	<a href="#">ab213906</a>		<0.003	0.031 - 2	3h 30m	•	•	•	•				
SCF*	<a href="#">ab100740</a>		<0.005	0.003 - 2	5h	•	•	•	•	•			
SCF	<a href="#">ab213908</a>		<0.01	0.031 - 2	3h 30m	•	•		•				
VEGF	<a href="#">ab100786</a>		<0.002	0.082 - 0.2	5h	•	•	•	•	•			
VEGF R3	<a href="#">ab213926</a>		0.01	0.16 - 10	3h 30m	•	•						

Kits validated on cell cell lysate (CL), tissue lysate (TL), cell culture extract (CE), tissue extract (TE), cell culture media (CM)

Target Name	abID	SimpleStep	Sensitivity (ng/mL)	Range (ng/mL)	Assay Time	Validated samples				
						CL	TL	CE	TE	CM
Akt (pS473)** (SQ)	<a href="#">ab126432</a>				5h	•				
Akt (pS473) + total Akt ** (SQ)	<a href="#">ab126433</a>				5h	•				
AKT 1/2/3 pS473 + AKT1** (SQ)	<a href="#">ab176657</a>	✓	0.03		1h 30m	•	•			
AKT2**	<a href="#">ab208986</a>	✓	0.005	0.078 - 5	1h 30m				•	•
Angiopoietin-1	<a href="#">ab213920</a>		<0.01	0.16 - 10	3h 30m	•	•			
Erk1 (pT202/pY204) + Erk2 (pT185/pY187)** (SQ)	<a href="#">ab126444</a>				5h	•				
Erk1 (pT202/pY204) + Erk2 (pT185/pY187) + total Erk1/2** (SQ)	<a href="#">ab126445</a>				5h	•				
FGF1**	<a href="#">ab219636</a>	✓	0.00197	0.019 - 1.2	1h 30m					•
FGF1*	<a href="#">ab223587</a>	✓		0.016 - 1	1h 30m				•	
FGF1	<a href="#">ab213914</a>		<0.01	0.032 - 2	3h 30	•	•			

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). \* Reacts with mouse - \*\* Reacts with mouse and rat – SQ: semi quantitative



## ELISA kits for Ras signalling pathway (rat) (continued)

Kits validated on cell cell lysate (CL), tissue lysate (TL), cell culture extract (CE), tissue extract (TE), cell culture media (CM)

Target Name	abID	SimpleStep	Sensitivity (ng/mL)	Range (ng/mL)	Assay Time	validated sample types				
						CL	TL	CE	TE	CM
IGF-1	ab213902		<0.005	0.063 - 4	3h 30m	•	•			
JNK2**	ab218265	✓	0.081	0.234 - 15				•	•	
Mek1 (pS217/221)** (SQ)	ab126446				5h	•				
Mek1 (pS217/221) + total Mek1** (SQ)	ab126448				5h	•				
Met (pY1234/pY1235) + total Met** (SQ)	ab126451				5h	•	•			
PDGF-AA	ab155464		<0.09	0.094- 6	5h					•
PDGF-AB	ab213906		<0.003	0.031 - 2	3h 30m	•	•			
VEGF	ab100787		<0.002	0.082 - 0.2	5h			•	•	

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). \* Reacts with mouse - \*\* Reacts with mouse and rat – SQ: semi quantitative