

PhosphoSens[®] Technology - Sox-based Homogeneous, Kinetic & Fluorescence-based Kinase Assays
List of Assays by Target Kinase

Revised May 16, 2017

Currently 167 assays with 100 Tyrosine Kinases & 67 Serine/Threonine Kinases (actively being expanded)

Summary of PhosphoSens [®] Assays in Terms of Assay Strength based on Reaction Rates				
Categories	Very Strong (VS)	Strong (S)	Medium (M)	Weak (W) ¹
Initial Reaction Rate (RFU/pmole kinase/min)	>1000	500-999	250-499	<250
# of Assays by Category	134	18	15	0
% of Assays by Category	80%	11%	9%	0%

¹Improvements in progress

Pricing: \$295 for a 96 assay kit, including all required reagents besides kinase. Bulk pricing starts at 1 mg (net) - please inquire.

Special Offer: Mention this listing and receive 10% off your first order (kit or bulk sensor).

#	Recombinant Kinase	Source	Kinase Type	Best Sox-based Substrate	Substrate Mr (Da)	Working [Substrate], μ M	Strength of Assay	Status
Tyrosine Kinases (Wild-type or mutant forms)								
1	ABL1	Carna BS	Tyr	AQT0032	1673.8	10	VS	Available
2	ABL1 (E255K)	Carna BS	Tyr	AQT0032	1673.8	10	VS	Available
3	ABL1 (T315I)	Carna BS	Tyr	AQT0032	1673.8	10	VS	Available
4	ABL2	Carna BS	Tyr	AQT0032	1673.8	10	VS	Available
5	ALK	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
6	ALK (C1156Y)	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
7	ALK (F1174L)	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
8	ALK (G1269A)	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
9	ALK (I202R)	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
10	ALK (R1275Q)	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
11	ALK (T1151-L1152 INS T)	Carna BS	Tyr	AQT0026	1367.3	10	S	Available
12	AXL	Carna BS	Tyr	AQT0101	1384.6	10	VS	Available
13	BLK	Carna BS	Tyr	AQT0032	1673.8	10	VS	Available
14	BMX	Carna BS	Tyr	AQT0102	1268.2	10	VS	Available
15	BTK	Carna BS	Tyr	AQT0101	1384.6	10	VS	Available
16	EGFR	BPS BS	Tyr	AQT0099	1383.5	10	VS	Available
17	EGFR	Carna BS	Tyr	AQT0099	1383.5	10	VS	Available
18	EGFR (d746-750)	Carna BS	Tyr	AQT0099	1383.5	10	VS	Available
19	EGFR (d746-750/T790M)	Carna BS	Tyr	AQT0099	1383.5	10	VS	Available
20	EGFR (L851Q)	Carna BS	Tyr	AQT0099	1383.5	10	VS	Available
21	EGFR (L858R)	Carna BS	Tyr	AQT0099	1383.5	10	VS	Available
22	EGFR (T790M)	Carna BS	Tyr	AQT0099	1383.5	10	VS	Available
23	EGFR (T790M/C797S/L858R)	BPS BS	Tyr	AQT0099	1383.5	10	VS	Available
24	EGFR (T790M/L858R)	BPS BS	Tyr	AQT0099	1383.5	10	VS	Available
25	EPHA1	Carna BS	Tyr	AQT0025	2008.1	10	VS	Available
26	EPHA2	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
27	EPHA3	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
28	EPHA4	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
29	EPHA5	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
30	EPHA6	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
31	EPHA7	Carna BS	Tyr	AQT0025	2008.1	10	M	Available
32	EPHA8	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
33	EPHB1	Carna BS	Tyr	AQT0025	2008.1	10	VS	Available

34	EPHB2	Carna BS	Tyr	AQT0025	2008.1	10	VS	Available
35	EPHB3	Carna BS	Tyr	AQT0025	2008.1	10	VS	Available
36	EPHB4	Carna BS	Tyr	AQT0025	2008.1	10	VS	Available
37	FAK (+ 0.5 nM Src)	Carna BS	Tyr	AQT0061	2122.9	20	VS	Available
38	FER	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
39	FES	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
40	FGFR1	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
41	FGFR2	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
42	FGFR3	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
43	FGFR3 (G697C)	ProKinase	Tyr	AQT0026	1367.3	10	VS	Available
44	FGFR3 (K650E)	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
45	FGFR3 (K650M)	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
46	FGFR3 (V555L)	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
47	FGFR3 (V555M)	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
48	FGFR4	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
49	FGR	Carna BS	Tyr	AQT0032	1673.8	10	VS	Available
50	FLT1	Carna BS	Tyr	AQT0001	1431.3	10	VS	Available
51	FLT3	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
52	FLT4	Carna BS	Tyr	AQT0001	1431.3	10	VS	Available
53	FMS	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
54	FRK	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
55	FYNa	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
56	FYNb	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
57	HCK	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
58	HER2	BPS BS	Tyr	AQT0001	1431.3	10	S	Available
59	HER2 (725YVMA726)	ProKinase	Tyr	AQT0094	2042.1	10	S	Available
60	HER4	Carna BS	Tyr	AQT0001	1431.3	10	VS	Available
61	IGF1R	Carna BS	Tyr	AQT0025	2008.1	10	VS	Available
62	INSR	Carna BS	Tyr	AQT0101	1384.6	10	VS	Available
63	IRR	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
64	ITK	Carna BS	Tyr	AQT0101	1384.6	10	S	Available
65	JAK1	Thermo Fisher	Tyr	AQT0001	1431.3	10	S	Available
66	JAK2	Carna BS	Tyr	AQT0001	1431.3	10	VS	Available
67	JAK3	Carna BS	Tyr	AQT0001	1431.3	10	VS	Available
68	KDR	Carna BS	Tyr	AQT0007	1332.2	10	VS	Available
69	KIT	Carna BS	Tyr	AQT0102	1268.2	10	VS	Available
70	LCK	Carna BS	Tyr	AQT0025	2008.1	10	VS	Available
71	LTK	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
72	LYNa	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
73	LYNb	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
74	MER	Carna BS	Tyr	AQT0101	1384.6	10	VS	Available
75	MET	Carna BS	Tyr	AQT0104	1361.5	10	VS	Available
76	NPM1-ALK	Carna BS	Tyr	AQT0025	2008.1	10	S	Available
77	PDGFR α	Carna BS	Tyr	AQT0001	1431.3	10	S	Available
78	PDGFR β	Carna BS	Tyr	AQT0025	2008.1	10	VS	Available
79	PYK2	Carna BS	Tyr	AQT0101	1384.6	10	VS	Available
80	RET	Carna BS	Tyr	AQT0009/102	1268.2	10	VS	Available
81	RET (G691S)	Carna BS	Tyr	AQT0009/102	1268.2	10	VS	Available
82	RET (M918T)	Carna BS	Tyr	AQT0009/102	1268.2	10	VS	Available
83	RET (S891A)	Carna BS	Tyr	AQT0009/102	1268.2	10	VS	Available
84	RET (Y791F)	Carna BS	Tyr	AQT0009/102	1268.2	10	VS	Available
85	RET-CCDC6	SignalChem	Tyr	AQT0009/102	1268.2	10	VS	Available
86	RON	Carna BS	Tyr	AQT0001	1431.3	10	S	Available
87	ROS	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
88	SRC	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
89	SRM	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
90	SYK	Carna BS	Tyr	Y7/AQT0016	1375.2	10	VS	Available
91	TEC	Carna BS	Tyr	AQT0101	1384.6	10	VS	Available
92	TIE2	Carna BS	Tyr	AQT0025	2008.1	10	VS	Available

93	TRKA	BPS BS	Tyr	AQT0096	1353.3	10	VS	Available
94	TRKB	BPS BS	Tyr	AQT0096	1353.3	10	VS	Available
95	TRKC	BPS BS	Tyr	AQT0096	1353.3	10	VS	Available
96	TXK	Carna BS	Tyr	AQT0104	1361.5	10	VS	Available
97	Tyk2	Thermo Fisher	Tyr	AQT0001	1431.3	10	M	Available
98	TYRO3	Carna BS	Tyr	AQT0104	1361.5	10	VS	Available
99	YES	Carna BS	Tyr	AQT0026	1367.3	10	VS	Available
100	ZAP70	Carna BS	Tyr	AQT0001	1431.3	10	VS	Available

Serine/Threonine Kinases (rapidly being expanded)

#	Recombinant Kinase	Source	Kinase Type	Best Sox-based Substrate	Substrate Mr (Da)	Working [Substrate], μ M	Strength of Assay	Status
1	AKT1	Carna BS	Ser/Thr	AQT0076	1469.4	10	VS	Available
2	AKT2	Carna BS	Ser/Thr	AQT0076	1469.4	10	VS	Available
3	AKT3	Carna BS	Ser/Thr	AQT0076	1469.4	10	VS	Available
4	AMPK α 1 β 1 γ 1	SignalChem	Ser/Thr	L2-1B12	2463.6	10	M	Production
5	AURORA A	Carna BS	Ser/Thr	AUR sensor	1398.4	10	VS	Available
	CDC2 /CycB1	Carna BS	Ser/Thr	Assay development in progress, please inquire				
	CDC2L6 /CycC	Carna BS	Ser/Thr					
	CDC7/ASK	Carna BS	Ser/Thr					
	CDK1/CycA2	ProQinase	Ser/Thr					
	CDK1/CycB1	ProQinase	Ser/Thr					
	CDK1/CycE1	ProQinase	Ser/Thr					
	CDK2/cyclin A1	Thermo Fisher	Ser/Thr					
	CDK2/CycA2	Carna BS	Ser/Thr					
	CDK2/CycE1	Carna BS	Ser/Thr					
	CDK2/cyclin O	Thermo Fisher	Ser/Thr					
	CDK3 /CycE1	Carna BS	Ser/Thr					
	CDK4/cyclin D1	Thermo Fisher	Ser/Thr					
	CDK4 /CycD3	Carna BS	Ser/Thr					
	CDK5 /p25	Carna BS	Ser/Thr					
	CDK5 /p35	ProQinase	Ser/Thr					
	CDK6/CycD1	Carna BS	Ser/Thr					
	CDK6/CycD3	Carna BS	Ser/Thr					
	CDK7 /CycH/MAT1	Carna BS	Ser/Thr					
	CDK8 /CycC	Carna BS	Ser/Thr					
	CDK9/CycK	Carna BS	Ser/Thr					
	CDK9/CycT1	Carna BS	Ser/Thr					
	CDK11/cyclin C	Thermo Fisher	Ser/Thr					
	CDK12 wt/CycK	ProQinase	Ser/Thr					
	CDK13/CycK	ProQinase	Ser/Thr					
	CDK14 (PFTK1)/cyclin Y	Thermo Fisher	Ser/Thr					
	CDK16/CycY	ProQinase	Ser/Thr					
	CDK17/cyclin Y	Thermo Fisher	Ser/Thr					
	CDK17/p35NCK	ProQinase	Ser/Thr					
	CDK18/cyclin Y	Thermo Fisher	Ser/Thr					
	CDK19/CycC	ProQinase	Ser/Thr					
	CDKL5	Thermo Fisher	Ser/Thr					
6	CDPK1 (toxoplasma gondii)	Academic	Ser/Thr	AQT0124	1748.0	10	VS	Available
7	CDPK3 (toxoplasma gondii)	Academic	Ser/Thr	AQT0124	1748.0	10	VS	Available
8	CHEK1	Carna BS	Ser/Thr	CHEK2-S2	2234.3	10	S	Available
9	CHEK2	Carna BS	Ser/Thr	CHEK2-S2	2234.3	10	VS	Available
10	DAPK1	Carna BS	Ser/Thr	L2-1B12	2463.9	10	VS	Production
11	DAPK2	Carna BS	Ser/Thr	L2-2A05	2476.9	10	S	Production
12	DAPK3	SignalChem	Ser/Thr	AQT0175	2367.5	10	VS	Available
13	DRAK1	Carna BS	Ser/Thr	AQT0173	2383.8	10	S	Available
14	DRAK2	SignalChem	Ser/Thr	AQT0174	2383.5	10	VS	Available
15	ERK1	Carna BS	Ser/Thr	AQT0082	1025.0	10	VS	Available

16	ERK2	Carna BS	Ser/Thr	AQT0082	1025.0	10	VS	Available
17	GCK (MAP4K2)	SignalChem	Ser/Thr	L1-A05	1403.4	10	VS	Production
18	GLK (MAP4K3)	SignalChem	Ser/Thr	AQT0135	2564.1	10	M	Available
19	HGK (MAP4K4)	SignalChem	Ser/Thr	AQT0135	2564.1	10	VS	Available
20	HIPK2	SignalChem	Ser/Thr	AQT0134	2357.5	10	M	Available
21	HPK1 (MAP4K1)	Thermo Fisher	Ser/Thr	AQT0135	2564.1	10	VS	Available
22	IKK β	Thermo Fisher	Ser/Thr	AQT-ST LIB10 2F12	2472.4	10	VS	Production
23	IKK ϵ	Thermo Fisher	Ser/Thr	AQT-ST LIB10 2F10	2067.9	10	VS	Production
24	KHS1 (MAP4K5)	SignalChem	Ser/Thr	L1-D07	1354.4	10	VS	Production
25	MAPKAPK2	SignalChem	Ser/Thr	AQT0150	2413.5	10	VS	Available
26	MARK2	SignalChem	Ser/Thr	L2-1B12	2463.6	10	M	Production
27	MEK1 (Cascade)	Carna BS	Ser/Thr	AQT0082	1025.0	10	VS	Available
28	MEK2 (Cascade)	Carna BS	Ser/Thr	AQT0082	1025.0	10	VS	Available
29	MINK (MAP4K6)	SignalChem	Ser/Thr	L1-D07	1354.4	10	VS	Production
30	MLCK	SignalChem	Ser/Thr	AQT0151	2342.5	10	M	Available
31	MLK1	SignalChem	Ser/Thr	AQT0136	1413.5	10	S	Available
32	MST1	SignalChem	Ser/Thr	AQT0135	2564.1	10	S	Available
33	MST2	SignalChem	Ser/Thr	AQT0135	2563.7	10	VS	Available
34	MST3	SignalChem	Ser/Thr	L2-1B12	2463.6	10	M	Production
35	MST4	SignalChem	Ser/Thr	L2-1B12	2463.6	10	M	Production
36	MYLK2	SignalChem	Ser/Thr	AQT0134	2357.5	10	VS	Available
37	MYLK4	SignalChem	Ser/Thr	AQT0151	2342.5	10	M	Available
38	NEK2	SignalChem	Ser/Thr	AQT0136	1413.5	10	M	Available
39	p38 γ	Carna BS	Ser/Thr	AQT0082	1025.0	10	S	Available
40	p38 δ	Carna BS	Ser/Thr	AQT0082	1025.0	10	M	Available
41	p70S6K	Carna BS	Ser/Thr	AQT0074	2037.1	10	VS	Available
42	PHKG2	Carna BS	Ser/Thr	AQT0074	2037.1	10	M	Available
43	PIM1	Carna BS	Ser/Thr	PIM sensor	1897.9	10	VS	Production
44	PKA α	Carna BS	Ser/Thr	AQT0128	1336.3	10	VS	Available
45	PKA β	Carna BS	Ser/Thr	AQT0128	1336.3	10	VS	Available
46	PKC α (alpha)	Carna BS	Ser/Thr	ST8	1237.4	10	VS	Available
47	PKC β 1 (beta1)	Carna BS	Ser/Thr	ST8	1237.4	10	S	Available
48	PKC β 2 (beta2)	Carna BS	Ser/Thr	ST8	1237.4	10	VS	Available
49	PKC δ (delta)	Carna BS	Ser/Thr	AQT0127	1603.71	20	M	Available
50	PKC γ (gamma)	Carna BS	Ser/Thr	ST8	1237.4	10	VS	Available
51	PKC ϵ (epsilon)	Carna BS	Ser/Thr	ST8	1237.4	10	VS	Available
52	PKC η (eta)	Carna BS	Ser/Thr	ST8	1237.4	20	S	Available
53	PKC θ (theta)	Carna BS	Ser/Thr	ST8	1237.4	10	VS	Available
54	PKC ι (iota)	Carna BS	Ser/Thr	AQT0127	1603.7	20	VS	Production
55	PKC ζ (zeta)	Carna BS	Ser/Thr	AQT0127	1603.71	20	S	Available
56	PLK1	SignalChem	Ser/Thr	L2-1B12	2463.6	10	M	Production
57	Raf, b- (Cascade)	Carna BS	Ser/Thr	AQT0082	1025.0	10	VS	Available
58	Raf, c- (Cascade)	Carna BS	Ser/Thr	AQT0082	1025.0	10	VS	Available
59	ROCK1	Carna BS	Ser/Thr	AQT0134	2357.5	10	S	Available
60	RSK1	Carna BS	Ser/Thr	PKA-S4	1336.3	10	VS	Production
61	RSK2	Carna BS	Ser/Thr	PKA-S4	1336.3	10	VS	Production
62	RSK3	Carna BS	Ser/Thr	PKA-S4	1336.3	10	VS	Production
63	SGK1	Carna BS	Ser/Thr	AQT0080	1571.5	10	VS	Available
64	SGK2	Carna BS	Ser/Thr	AQT0080	1571.5	10	VS	Available
65	SGK3	Carna BS	Ser/Thr	AQT0080	1571.5	10	VS	Available
66	TBK1	Carna BS	Ser/Thr	AQT-ST LIB10 2F10	2067.9	10	VS	Production
67	TNIK (MAP4K)	SignalChem	Ser/Thr	AQT0135	2564.1	10	VS	Available