

List of Kinases for Cellular Phosphorylation Assays



No.	Kinase	Cell Line	No.	Kinase	Cell Line	No.	Kinase	Cell Line	No.	Kinase	Cell Line
1	AKT1 (h)	Rat1-myrAKT1 (r)	29	FYN (h)	Rat1 (r)						
2	ALK (h)	Karpas-299 (h)	30	IGF1-R (h)	MEF-IGF1-R (m)						
3	Aurora-B (h)	HT-29 (h)	31	KIT (h)	M07e (h)						
4	AXL (h)	MEF-AXL (m)	32	MET (h)	MKN45 (h)						
5	BCR-ABL	K562	33	MET D1288H Mutant (h)	Rat1 (r)						
6	B-RAF-VE (h)	Rat1-B-RAF-VE (r)	34	MET D1228N Mutant (h)	Rat1 (r)						
7	EGF-R (h)	A431 (h)	35	MET F1200I Mutant (h)	Rat1 (r)						
8	EGF-R D746-750/T790M Mutant (h)	Rat1 (r)	36	MET M1250T Mutant (h)	Rat1 (r)						
9	EGF-R D746-750/C797S Mutant (h)	Rat1 (r)	37	MET Y1230H Mutant (h)	Rat1 (r)						
10	EGF-R D746-750/T790M/C797S Mutant (h)	Rat1 (r)	38	MET Y1230A Mutant (h)	Rat1 (r)						
11	EGF-R G719S Mutant (h)	Rat1 (r)	39	MET Y1230C Mutant (h)	Rat1 (r)						
12	EGF-R L858R Mutant (h)	Rat1 (r)	40	MET Y1230D Mutant (h)	Rat1 (r)						
13	EGF-R L861Q Mutant (h)	Rat1 (r)	41	MET control wt (h)	Rat1 (r)						
14	EGF-R T790M Mutant (h)	Rat1 (r)	42	MNK1 (h)	Karpas-299 (h)						
15	EGF-R T790M/C797S/L858R (h)	Rat1 (r)	43	PDGFR-beta (m)	NIH3T3 (m)						
16	EGF-R T790/L858R Mutant (h)	Rat1 (r)	44	PIM1	HEK293T (h)						
17	EGF-R Δ752-759 Mutant (h)	Rat1 (r)	45	PIM2	HEK293T (h)						
18	EGF-R Δ747-749/A750P Mutant (h)	Rat1 (r)	46	PIM3	HEK293T (h)						
19	EGF-R control wt (h)	Rat1 (r)	47	ROCK (r)	A7r5 (r)						
20	EPHB4 (h)	MEF-EPHB4 (m)	48	RON (h)	T47D (h)						
21	ERBB2 (h)	NIH3T3-ERBB2 (m)	49	S6K (h)	Karpas-299 (h)						
22	ERBB4 (h)	T47D (h)	50	SRC (h)	MEF-SRC (m)						
23	FAK (h)	MEF-FAK (m)	51	TIE2 (h)	CHO-TIE2 (ha)						
24	FGF-R2	Kato III	52	VEGF-R2 (h)	HUE (h)						
25	FLT3-wt (h)	MEF-FLT3-wt (m)	53	VEGF-R3 (h)	MEF-VEGF-R3 (m)						
26	FLT3-DY (h)	MEF-FLT3-DY (m)									
27	FLT3-ITD (h)	MEF-FLT3-ITD (m)									
28	Haspin (GSG2)	HT-29 (h)									

In brackets: origin of kinase and cell line (h: human; ha: hamster; m: mouse; r: rat)

 Cell Lines