

# List of Kinases for the Kinase WildType 2Point-Profiler and the WildType IC50-Profiler

No.	Kinase	No.	Kinase	No.	Kinase	No.	Kinase	No.	Kinase	No.	Kinase	No.	Kinase
1	ABL1 wt	51	CDK5/p25NCK	101	EPHB2	151	KIT wt	201	NEK3	251	RAF1 Y340D/Y341D*	301	TSK2
2	ABL2	52	CDK5/p35NCK	102	EPHB3	152	LCK	202	NEK4	252	RET wt	302	TSSK1
3	ACK1	53	CDK6/CycD1	103	EPHB4	153	LIMK1	203	NEK6	253	RIPK2	303	TTBK1
4	ACV-R1	54	CDK6/CycD3	104	ERBB2 wt	154	LIMK2	204	NEK7	254	RIPK5	304	TTBK2
5	ACV-R1B	55	CDK7/CycH/MAT1	105	ERBB4	155	LRRK2 wt	205	NEK9	255	ROCK1	305	TTK
6	ACV-R2A	56	CDK8/CycC	106	ERK1	156	LTK	206	NIK	256	ROCK2	306	TXK
7	ACV-R2B	57	CDK9/CycK	107	ERK2	157	LYN	207	NLK	257	RON	307	TYK2
8	ACV-RL1	58	CDK9/CycT1	108	ERK5	158	MAP3K1	208	p38-alpha	258	ROS	308	TYRO3
9	AKT1 aa106-480	59	CHK1	109	ERK7	159	MAP3K10	209	p38-beta	259	RPS6KA1	309	VEGF-R1
10	AKT2 aa107-481	60	CHK2	110	FAK aa2-1052	160	MAP3K11	210	p38-delta	260	RPS6KA2	310	VEGF-R2
11	AKT3 aa106-479	61	CK1-alpha1	111	FER	161	MAP3K7/MAP3K7IP1	211	p38-gamma	261	RPS6KA3	311	VEGF-R3
12	ALK wt (GST-HIS-tag)	62	CK1-delta	112	FES	162	MAP3K9	212	PAK1	262	RPS6KA4	312	VRK1
13	AMPK-alpha1 aa1-550	63	CK1-epsilon	113	FGF-R1 wt	163	MAP4K2	213	PAK2	263	RPS6KA5	313	VRK2
14	ARK5	64	CK1-gamma1	114	FGF-R2	164	MAP4K4	214	PAK3	264	RPS6KA6	314	WEE1
15	ASK1	65	CK1-gamma2	115	FGF-R3 wt	165	MAP4K5	215	PAK4	265	S6K	315	WNK1
16	Aurora-A	66	CK1-gamma3	116	FGF-R4	166	MAPKAPK2	216	PAK6	266	S6K-beta	316	WNK2
17	Aurora-B	67	CK2-alpha1	117	FGR	167	MAPKAPK3	217	PAK7	267	SAK	317	WNK3
18	Aurora-C	68	CK2-alpha2	118	FLT3 wt	168	MAPKAPK5	218	PASK	268	SGK1	318	YES
19	AXL	69	CLK1	119	FRK	169	MARK1	219	PBK	269	SGK2	319	ZAK
20	BLK	70	CLK2	120	FYN wt	170	MARK2	220	PDGFR-alpha wt	270	SGK3	320	ZAP70
21	BMPR1A	71	CLK3	121	GRK2	171	MARK3	221	PDGFR-beta	271	SIK1		
22	BMX	72	CLK4	122	GRK3	172	MARK4	222	PK1	272	SIK2		
23	B-RAF wt	73	COT	123	GRK4	173	MATK	223	PHKG1	273	SIK3		
24	BRK	74	CSF1-R	124	GRK5	174	MEK1 wt	224	PHKG2	274	SLK		
25	BRSK1	75	CSK	125	GRK6	175	MEK2	225	PIM1	275	SNARK		
26	BRSK2	76	DAPK1	126	GRK7	176	MEK5	226	PIM2	276	SNK		
27	BTK	77	DAPK2	127	GS2	177	MEKK2	227	PIM3	277	SRC (GST-HIS-tag)		
28	BUB1B	78	DAPK3	128	GSK3-alpha	178	MEKK3	228	PKA	278	SRMS		
29	CAMK1D	79	DCAMK2	129	GSK3-beta	179	MELK	229	PKC-alpha	279	SRPK1		
30	CAMK2A	80	DDR2 wt	130	HCK	180	MERTK	230	PKC-beta1	280	SRPK2		
31	CAMK2B	81	DMPK	131	HIPK1	181	MET wt	231	PKC-beta2	281	STK17A		
32	CAMK2D	82	DNA-PK	132	HIPK2	182	MINK1	232	PKC-delta	282	STK23		
33	CAMK2G	83	DYRK1A	133	HIPK3	183	MKK4	233	PKC-epsilon	283	STK25		
34	CAMK4	84	DYRK1B	134	HIPK4	184	MKK6 S207D/T211D*	234	PKC-eta	284	STK33		
35	CAMKK1	85	DYRK2	135	HRI	185	MKK7	235	PKC-gamma	285	STK39		
36	CAMKK2	86	DYRK3	136	IGF1-R	186	MKNK1	236	PKC-iota	286	SYK aa1-635		
37	CDC42BPA	87	DYRK4	137	IKK-alpha	187	MKNK2	237	PKC-mu	287	TAOK2		
38	CDC42BPB	88	EEF2K	138	IKK-beta	188	MLK4	238	PKC-nu	288	TAOK3		
39	CDC7/DBF4	89	EGF-R wt	139	IKK-epsilon	189	MST1	239	PKC-theta	289	TBK1		
40	CDK1/CycA2	90	EIF2AK2	140	INS-R	190	MST2	240	PKC-zeta	290	TEC		
41	CDK1/CycB1	91	EIF2AK3	141	INSR-R	191	MST3	241	PKMYT1	291	TGFB-R1		
42	CDK1/CycE1	92	EPHA1	142	IRAK1	192	MST4	242	PLK1	292	TGFB-R2		
43	CDK16/CycC	93	EPHA2	143	IRAK4 (untagged)	193	mTOR	243	PLK3	293	TIE2 wt		
44	CDK19/CycC	94	EPHA3	144	ITK	194	MUSK	244	PRK1	294	TLK1		
45	CDK2/CycA2	95	EPHA4	145	JAK1	195	MYLK	245	PRK2	295	TLK2		
46	CDK2/CycE1	96	EPHA5	146	JAK2	196	MYLK2	246	PRKD2	296	TNK1		
47	CDK3/CycC	97	EPHA6	147	JAK3	197	MYLK3	247	PRKG1	297	TRK-A wt		
48	CDK3/CycE1	98	EPHA7	148	JNK1	198	NEK1	248	PRKG2	298	TRK-B		
49	CDK4/CycD1	99	EPHA8	149	JNK2	199	NEK11	249	PRKX	299	TRK-C		
50	CDK4/CycD3	100	EPHB1	150	JNK3	200	NEK2	250	PYK2	300	TSF1		

\* constitutively active

 Protein Kinases