


List of Kinases for the WildType+ 2Point-Profiler

No.	Kinase	No.	Kinase	No.	Kinase	No.	Kinase	No.	Kinase	No.	Kinase	No.	Kinase	No.	Kinase
1	ABL1 wt	53	CDC42BPB	105	DYRK3	157	FYN wt	209	MAP3K11	261	NEK6	313	PKC-mu	365	SRMS
2	ABL1 E255K	54	CDC7/DBF4	106	DYRK4	158	GRK2	210	MAP3K7/MAP3K7IP1	262	NEK7	314	PKC-nu	366	SRPK1
3	ABL1 F317I	55	CDK1/CycA2	107	EEF2K	159	GRK3	211	MAP3K9	263	NEK9	315	PKC-theta	367	SRPK2
4	ABL1 G250E	56	CDK1/CycB1	108	EGF-R wt	160	GRK4	212	MAP4K2	264	NIK	316	PKC-zeta	368	STK17A
5	ABL1 H396P	57	CDK1/CycE1	109	EGF-R d746-750	161	GRK5	213	MAP4K4	265	NLK	317	PKC-zeta wt aa184-592 (PKM-zeta)	369	STK23
6	ABL1 M351T	58	CDK12 wt/CycK	110	EGF-R d747-749/A750P	162	GRK6	214	MAP4K5	266	NPM1 ALK	318	PKMYT1	370	STK25
7	ABL1 Q252H	59	CDK12 R722C/CycK	111	EGF-R d747-752/P753S	163	GRK7	215	MAPKAPK2	267	NPM1 ALK F1174L	319	PKN3	371	STK33
8	ABL1 T315I	60	CDK16/CycY	112	EGF-R d752-759	164	GSG2	216	MAPKAPK3	268	p38-alpha	320	PLK1	372	STK39
9	ABL1 Y253F	61	CDK19/CycC	113	EGF-R G719C	165	GSK3-alpha	217	MAPKAPK5	269	p38-beta	321	PLK3	373	SYK aa1-635
10	ABL2	62	CDK2/CycA2	114	EGF-R G719S	166	GSK3-beta	218	MARK1	270	p38-delta	322	PRK1	374	TAOK2
11	ACK1	63	CDK2/CycE1	115	EGF-R L858R	167	HCK	219	MARK2	271	p38-gamma	323	PRK2	375	TAOK3
12	ACV-R1	64	CDK3/CycC	116	EGF-R L861Q	168	HIPK1	220	MARK3	272	PAK1	324	PRKD2	376	TBK1
13	ACV-R1B	65	CDK3/CycE1	117	EGF-R T790M	169	HIPK2	221	MARK4	273	PAK2	325	PRKG1	377	TEC
14	ACV-R2A	66	CDK4/CycD1	118	EGF-R T790M/L858R	170	HIPK3	222	MATK	274	PAK3	326	PRKG2	378	TGFB-R1
15	ACV-R2B	67	CDK4/CycD3	119	E1F2AK2	171	HIPK4	223	MEK1 wt	275	PAK4	327	PRKX	379	TGFB-R2
16	ACV-RL1	68	CDK5/p25NCK	120	E1F2AK3	172	HRI	224	MEK2	276	PAK6	328	PYK2	380	TIE2 wt
17	AKT1 aa106-480	69	CDK5/p35NCK	121	EML4 ALK	173	IGF1-R	225	MEK5	277	PAK7	329	RAF1 Y340D/Y341D (untagged)*	381	TIE2 R849W
18	AKT2 aa107-481	70	CDK6/CycD1	122	EML4 ALK F1174L	174	IKK-alpha	226	MEKK2	278	PASK	330	RET wt	382	TIE2 Y1108F
19	AKT3 aa106-479	71	CDK6/CycD3	123	EPHA1	175	IKK-beta	227	MEKK3	279	PBK	331	RET E762Q	383	TIE2 Y897S
20	ALK wt (GST-HIS-tag)	72	CDK7/CycH/MAT1	124	EPHA2	176	IKK-epsilon	228	MELK	280	PDGFR-alpha wt	332	RET G691S	384	TLK1
21	ALK C1156Y (GST-HIS-tag)	73	CDK8/CycC	125	EPHA3	177	INS-R	229	MERTK	281	PDGFR-alpha D842V	333	RET M918T	385	TLK2
22	ALK F1174L (GST-HIS-tag)	74	CDK9/CycK	126	EPHA4	178	INSR-R	230	MET wt	282	PDGFR-alpha T674I	334	RET R749T	386	TNK1
23	ALK F1174S (GST-HIS-tag)	75	CDK9/CycT1	127	EPHA5	179	IRAK1	231	MET D1228H	283	PDGFR-alpha V561D	335	RET R813Q	387	TRK-A wt
24	ALK L1196M (GST-HIS-tag)	76	CHK1	128	EPHA6	180	IRAK4 (untagged)	232	MET D1228N	284	PDGFR-beta	336	RET S891A	388	TRK-B
25	ALK R1275Q (GST-HIS-tag)	77	CHK2	129	EPHA7	181	ITK	233	MET F1200I	285	PDK1	337	RET V804L	389	TRK-C
26	AMPK-alpha1 aa1-550	78	CK1-alpha1	130	EPHA8	182	JAK1	234	MET M1250T	286	PHKG1	338	RET V804M	390	TSF1
27	ARK5	79	CK1-delta	131	EPHB1	183	JAK2	235	MET Y1230A	287	PHKG2	339	RET Y791F	391	TSK2
28	ASK1	80	CK1-epsilon	132	EPHB2	184	JAK3	236	MET Y1230C	288	PI4K2A	340	RIPK2	392	TSSK1
29	Aurora-A	81	CK1-gamma1	133	EPHB3	185	JNK1	237	MET Y1230D	289	PI4K2B	341	RIPK5	393	TTBK1
30	Aurora-B	82	CK1-gamma2	134	EPHB4	186	JNK2	238	MET Y1230H	290	PI4KB	342	ROCK1	394	TTBK2
31	Aurora-C	83	CK1-gamma3	135	ERBB2 wt	187	JNK3	239	MET Y1235D	291	PIK3C2A	343	ROCK2	395	TTK
32	AXL	84	CK2-alpha1	136	ERBB4	188	KIT wt	240	MINK1	292	PIK3C2B	344	RON	396	TXK
33	BLK	85	CK2-alpha2	137	ERK1	189	KIT A829P	241	MKK4	293	PIK3C2G	345	ROS	397	TYK2
34	BMPR1A	86	CLK1	138	ERK2	190	KIT D816H	242	MKK6 S207D/T211D*	294	PIK3C3	346	RPS6KA1	398	TYRO3
35	BMX	87	CLK2	139	ERK5	191	KIT D816V	243	MKK7	295	PIK3CA/PIK3R1	347	RPS6KA2	399	VEGF-R1
36	B-RAF wt	88	CLK3	140	ERK7	192	KIT T670I	244	MKNK1	296	PIK3CB wt/PIK3R1	348	RPS6KA3	400	VEGF-R2
37	B-RAF V600E	89	CLK4	141	FAK aa2-1052	193	KIT V559D	245	MKNK2	297	PIK3CD/PIK3R1	349	RPS6KA4	401	VEGF-R3
38	BRK	90	COT	142	FER	194	KIT V559D/T670I	246	MLK2	298	PIK3CG	350	RPS6KA5	402	VRK1
39	BRSK1	91	CSF1-R	143	FES	195	KIT V559D/V654A	247	MST1	299	PIM1	351	RPS6KA6	403	VRK2
40	BRSK2	92	CSK	144	FGF-R1 wt	196	KIT V560G	248	MST2	300	PIM2	352	S6K	404	WEE1
41	BTK	93	DAPK1	145	FGF-R1 V561M	197	KIT V654A	249	MST3	301	PIM3	353	S6K-beta	405	WNK1
42	BUB1B	94	DAPK2	146	FGF-R2	198	LCK	250	MST4	302	PIP5K1A	354	SAK	406	WNK2
43	CAMK1D	95	DAPK3	147	FGF-R3 wt	199	LIMK1	251	mTOR	303	PIP5K1C	355	SGK1	407	WNK3
44	CAMK2A	96	DCAMK2	148	FGF-R3 G697C	200	LIMK2	252	MUSK	304	PKA	356	SGK2	408	YES
45	CAMK2B	97	DDR2 wt	149	FGF-R3 K650E	201	LRKK2 wt	253	MYLK	305	PKC-alpha	357	SGK3	409	ZAK
46	CAMK2D	98	DDR2 N456S	150	FGF-R3 K650M	202	LRKK2 G2019S	254	MYLK2	306	PKC-beta1	358	SIK1	410	ZAP70
47	CAMK2G	99	DDR2 T654M	151	FGF-R4	203	LRKK2 I2020T	255	MYLK3	307	PKC-beta2	359	SIK2		
48	CAMK4	100	DMPK	152	FGR	204	LRKK2 R1441C	256	NEK1	308	PKC-delta	360	SIK3		
49	CAMKK1	101	DNA-PK	153	FLT3 wt	205	LYN	257	NEK11	309	PKC-epsilon	361	SLK		
50	CAMKK2	102	DYRK1A	154	FLT3 D835Y	206	LYN	258	NEK2	310	PKC-eta	362	SNARK		
51	CCDC6-RET	103	DYRK1B	155	FLT3 ITD	207	MAP3K10	259	NEK3	311	PKC-gamma	363	SNK		
52	CDC42BPA	104	DYRK2	156	FRK	208	MAP3K10	260	NEK4	312	PKC-iota	364	SRC (GST-HIS-tag)		

 Protein Kinases

 Lipid Kinases