

# List of Active Recombinant Protein Kinases for Sale



No.	Kinase	No.	Kinase	No.	Kinase	No.	Kinase	No.	Kinase
1	ABL1 wt	54	CDK4/CycD3	107	EPHB1	160	KIT T670I	213	NPM1 ALK
2	ABL1 E255K	55	CDK5/p25NCK	108	EPHB2	161	KIT V559D	214	NPM1 ALK F1174L
3	ABL1 F317I	56	CDK5/p35NCK	109	EPHB3	162	KIT V559D/T670I	215	p38-alpha
4	ABL1 G250E	57	CDK6/CycD1	110	EPHB4	163	KIT V559D/V654A	216	p38-beta
5	ABL1 H396P	58	CDK6/CycD2	111	ERBB2 wt	164	KIT V560G	217	p38-delta
6	ABL1 M351T	59	CDK6/CycD3	112	ERBB2 775YVMA776	165	KIT V654A	218	p38-gamma
7	ABL1 Q252H	60	CDK7/CycH/MAT1	113	ERBB4	166	LCK	219	PAK1
8	ABL1 T315I	61	CDK8/CycC	114	ERK1	167	LYN	220	PAK2
9	ABL1 Y253F	62	CDK9/CycK	115	ERK2	168	MAP3K1	221	PAK3
10	ABL2	63	CDK9/CycT1	116	ERK5	169	MAP4K4	222	PAK4
11	ACK1	64	CDK12 wt/CycK	117	ERK7	170	MAPKAPK2	223	PAK6
12	ACV-R1	65	CDK12 R722C/CycK	118	FAK aa2-1052	171	MAPKAPK3	224	PAK7
13	AKT1 aa106-480	66	CDK13/CycK	119	FAK aa411-686	172	MAPKAPK5	225	PBK
14	AKT1 aa1-480	67	CDK16/CycY	120	FER	173	MARK1	226	PDGFR-alpha wt
15	AKT2 aa1-481	68	CDK17/p35NCK	121	FGF-R1 wt	174	MARK3	227	PDGFR-alpha D842V
16	AKT2 aa107-481	69	CDK18/CycY	122	FGF-R1 V561M	175	MASTL	228	PDGFR-alpha T674I
17	AKT3 aa1-479	70	CDK19/CycC	123	FGF-R2	176	MEK1 wt	229	PDGFR-alpha V561D
18	AKT3 aa106-479	71	CDK20/CycH	124	FGF-R3 wt	177	MEK1 F53L	230	PDGFR-beta
19	ALK wt (GST-HIS-tag)	72	CDK20/CycT1	125	FGF-R3 G697C	178	MEK1 S218E/S222E	231	PDK1
20	ALK wt (HIS-tag)	73	CHK1	126	FGF-R3 K650E	179	MEK5	232	PIM1
21	ALK C1156Y (GST-HIS-tag)	74	CK1-alpha1	127	FGF-R3 K650M	180	MEKK2	233	PIM2
22	ALK F1174L (GST-HIS-tag)	75	CK2-alpha1	128	FGF-R4	181	MEKK3	234	PIM3
23	ALK F1174S (GST-HIS-tag)	76	CK2-alpha2	129	FGR	182	MELK	235	PKC-alpha
24	ALK G1202R (GST-HIS-tag)	77	CLK1	130	FLT3 wt	183	MERTK	236	PKC-beta1
25	ALK L1196M (GST-HIS-tag)	78	CLK3	131	FLT3 D835Y	184	MET wt	237	PKC-beta2
26	ALK R1275Q (GST-HIS-tag)	79	COT	132	FLT3 ITD	185	MET D1228H	238	PKC-delta
27	AMPK-alpha1 aa1-312	80	CSF1-R	133	FRK	186	MET D1228N	239	PKC-epsilon
28	AMPK-alpha1 aa1-550	81	CSK	134	FYN wt	187	MET F1200I	240	PKC-eta
29	ARK5	82	DAPK1	135	FYN Y531F	188	MET G1163R	241	PKC-gamma
30	Aurora-A	83	DDR2 wt	136	GSG2	189	MET L1195V	242	PKC-iota
31	Aurora-B	84	DDR2 N456S	137	GSK3-beta	190	MET M1250T	243	PKC-mu
32	Aurora-C	85	DDR2 T654M	138	HCK	191	MET Y1230A	244	PKC-nu
33	AXL	86	EGF-R wt	139	HIPK3	192	MET Y1230C	245	PKC-theta
34	BLK	87	EGF-R C797S	140	HRI	193	MET Y1230D	246	PKC-zeta
35	B-RAF wt	88	EGF-R d746-750	141	IGF1-R	194	MET Y1230H	247	PKC-zeta wt aa184-592 (PKM-zeta)
36	B-RAF V600E	89	EGF-R d747-749/A750P	142	IKK-beta	195	MET Y1235D	248	PKMYT1
37	BMPR1B	90	EGF-R d747-752/P753S	143	IKK-epsilon	196	MINK1	249	PLK1
38	BRK	91	EGF-R d752-759	144	INS-R	197	MKK4	250	PLK3
39	BRSK2	92	EGF-R G719C	145	IRAK4 (GST-HIS-tag)	198	MKK6 S207D/T211D	251	PRK1
40	BTK	93	EGF-R G719S	146	IRAK4 (untagged)	199	MKK7	252	PYK2
41	CCDC6-RET	94	EGF-R L858R	147	ITK	200	MLK4	253	RAF1 Y340D/Y341D (GST-HIS-tag)
42	CDC42BPB	95	EGF-R L861Q	148	JAK1 aa583-1154 wt	201	MST1	254	RAF1 Y340D/Y341D (untagged)
43	CDC7/DBF4	96	EGF-R T790M	149	JAK1 aa583-1154 S729C	202	MST2	255	RET wt
44	CDK1/CycA2	97	EGF-R T790M/L858R	150	JAK1 aa850-1154 wt	203	MST4	256	RET E762Q
45	CDK1/CycB1	98	EML4 ALK	151	JAK2	204	MUSK	257	RET G691S
46	CDK1/CycE1	99	EML4 ALK F1174L	152	JAK3	205	MYLK2	258	RET M918T
47	CDK2/CycA2	100	EPHA1	153	JNK1	206	NEK1	259	RET R749T
48	CDK2/CycD1	101	EPHA2	154	JNK2	207	NEK2	260	RET R813Q
49	CDK2/CycE1	102	EPHA3	155	JNK3	208	NEK3	261	RET S891A
50	CDK3/CycC	103	EPHA4	156	KIT wt	209	NEK6	262	RET V804E
51	CDK3/CycE1	104	EPHA5	157	KIT A829P	210	NEK7	263	RET V804L
52	CDK4/CycD1	105	EPHA6	158	KIT D816H	211	NIK	264	RET V804M
53	CDK4/CycD2	106	EPHA7	159	KIT D816V	212	NLK	265	RET Y791F