

List of Recombinant Additional Kinase Variants for Sale



No.	Kinase	No.	Kinase	No.	Kinase	No.	Kinase
1	AKT1 aa1-480	29	p38-gamma				
2	CDC7	30	p38-delta				
3	CDK1	31	PKC-zeta aa184-592 K281M (PKM-zeta)				
4	CDK2						
5	CDK3						
6	CDK4						
7	CDK5						
8	CDK6						
9	CDK6/CycD1						
10	CDK6/CycD3						
11	CDK7						
12	CDK8						
13	CDK9						
14	CDK12 Y901C/CycK						
15	CDK16						
16	ERK2						
17	ERK7						
18	FAK						
19	JNK1						
20	JNK2						
21	JNK3						
22	MAPKAPK3						
23	MAPKAPK5						
24	MEK1						
25	MKK4						
26	MKK7						
27	p38-alpha						
28	p38-beta						