

## List of Recombinant Kinase Substrates for Sale



No.	Substrate	No.	Substrate	No.	Substrate	No.	Substrate
1	ATF2						
2	ELK1*						
3	ERK2 K54R						
4	JNK1 K55R/K56R*						
5	JUN						
6	MEK1 K97M						
7	p38-alpha K53A						
8	RB-CTF						
9	RB-S6P						
10	RB-ER-CDC25tide*						
11	RB-ER-CHKtide						
12	RB-ER-GSK3(14-27)						
13	RB-ER-IRStide						
14	RB-ER-NTRK3tide						

\*not recommended for use in ATP consumption assays