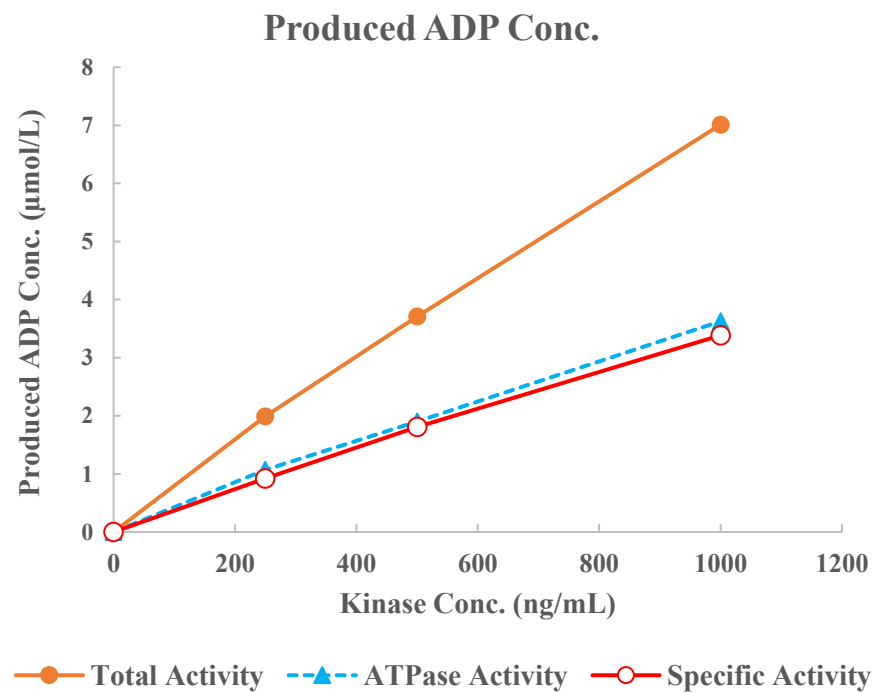
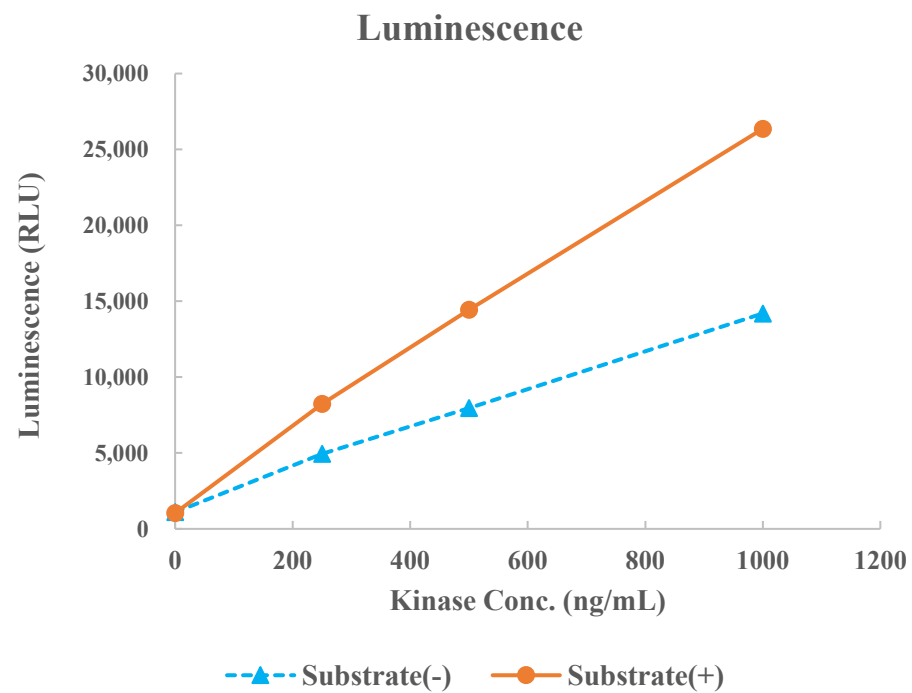


MINK(MINK1)



Kinase	MINK(MINK1), 07-139
Substrate	Biotin-Modified Erktide, BTSU03
Substrate Conc.	10 μmol/L
ATP Conc.	50 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour