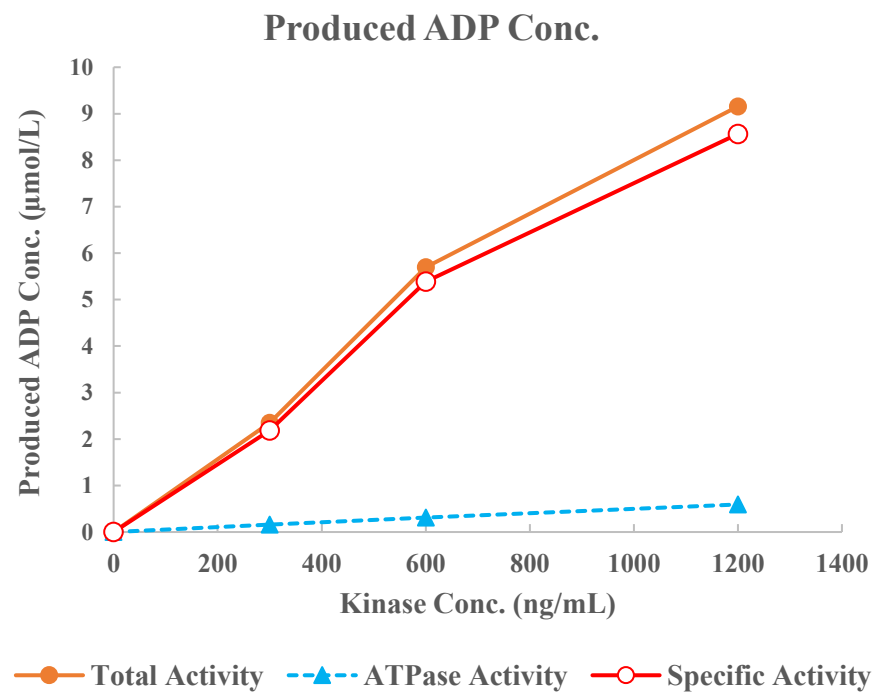
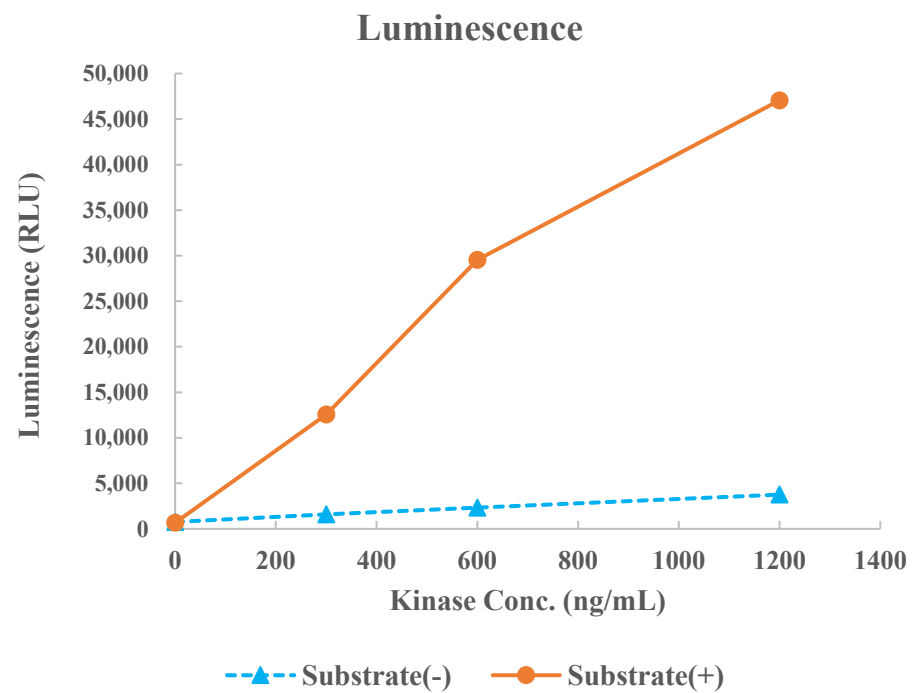


MUSK



Kinase	MUSK, 08-153
Substrate	Biotin-CSKtide, BTSU02
Substrate Conc.	50 μmol/L
ATP Conc.	10 μmol/L
Metal	Mg, Mn
Metal Conc.	5, 1 mmol/L
Additives	-
Reaction Time	1 hour