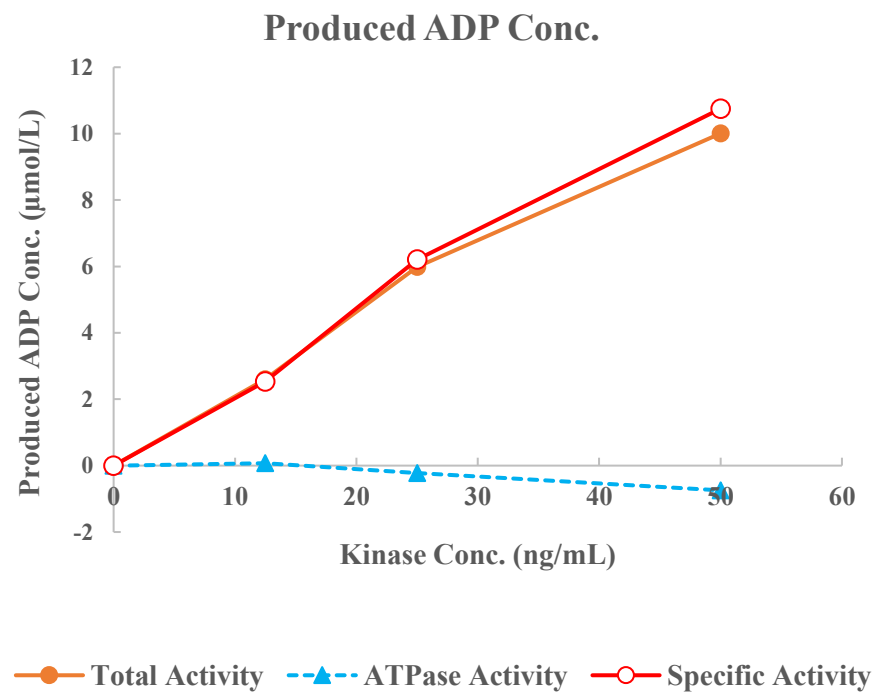
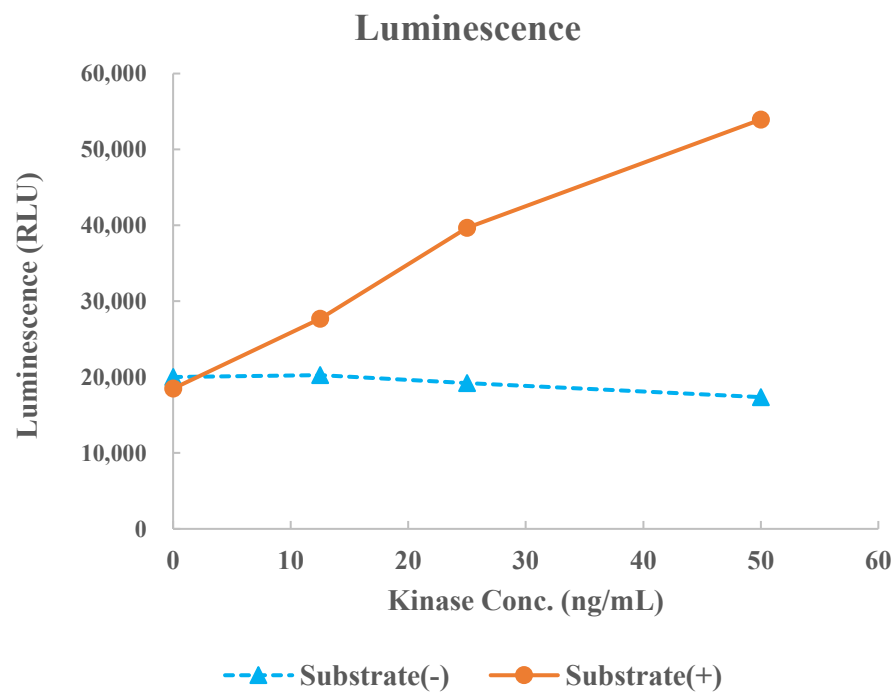


PIM1



Kinase	PIM1, 02-054
Substrate	Biotin-S6K2 peptide, BTSU14
Substrate Conc.	10 μmol/L
ATP Conc.	500 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour