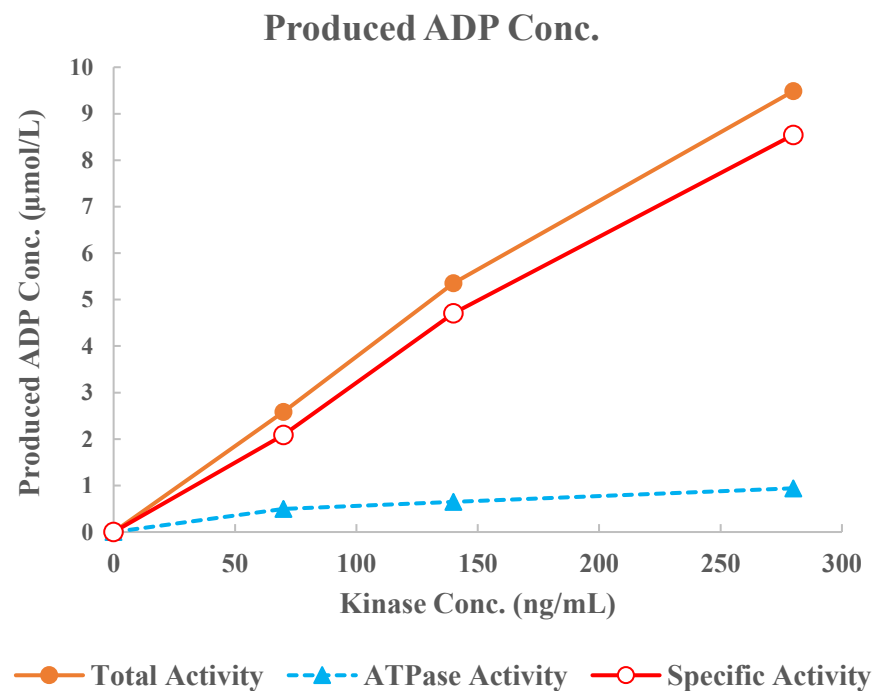
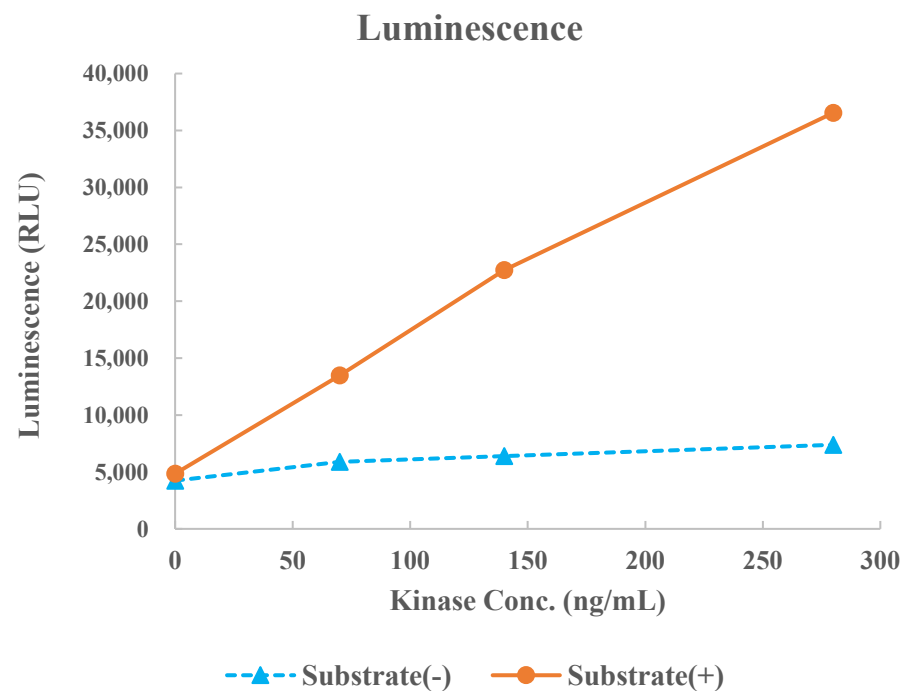


PIM3



Kinase	PIM3, 02-156
Substrate	Biotin-S6K2 peptide, BTSU14
Substrate Conc.	10 μmol/L
ATP Conc.	150 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour