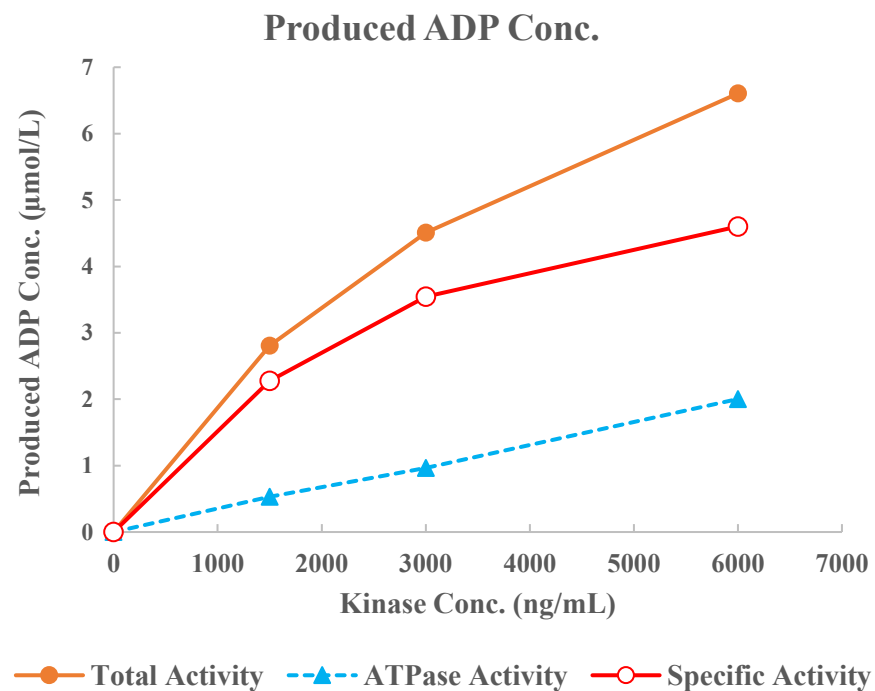
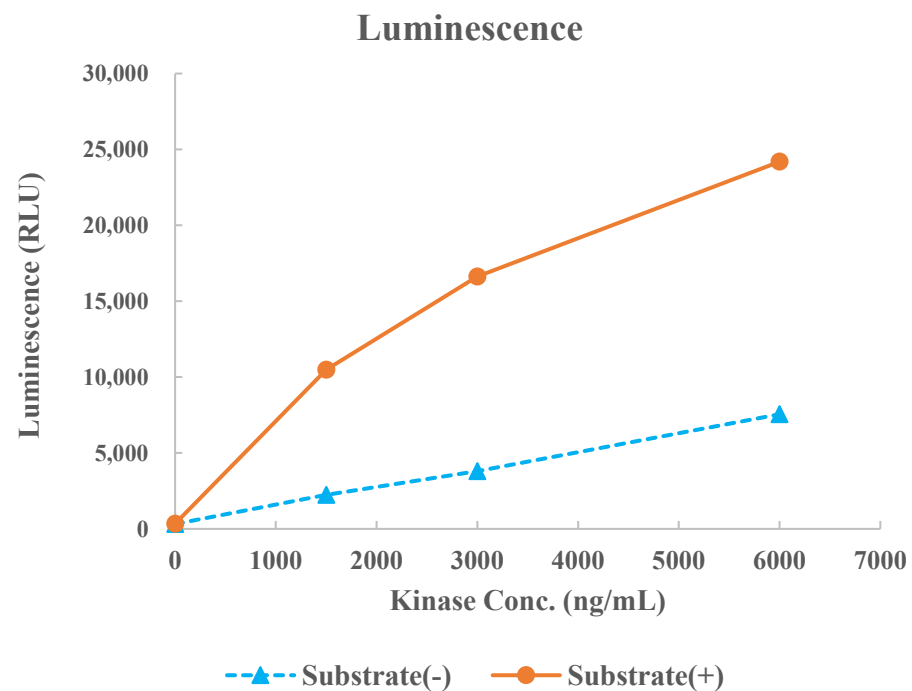


BUB1/BUB3



Kinase	BUB1/BUB3, 05-187
Substrate	Biotin-H2A peptide, BTSU47
Substrate Conc.	10 μmol/L
ATP Conc.	10 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour