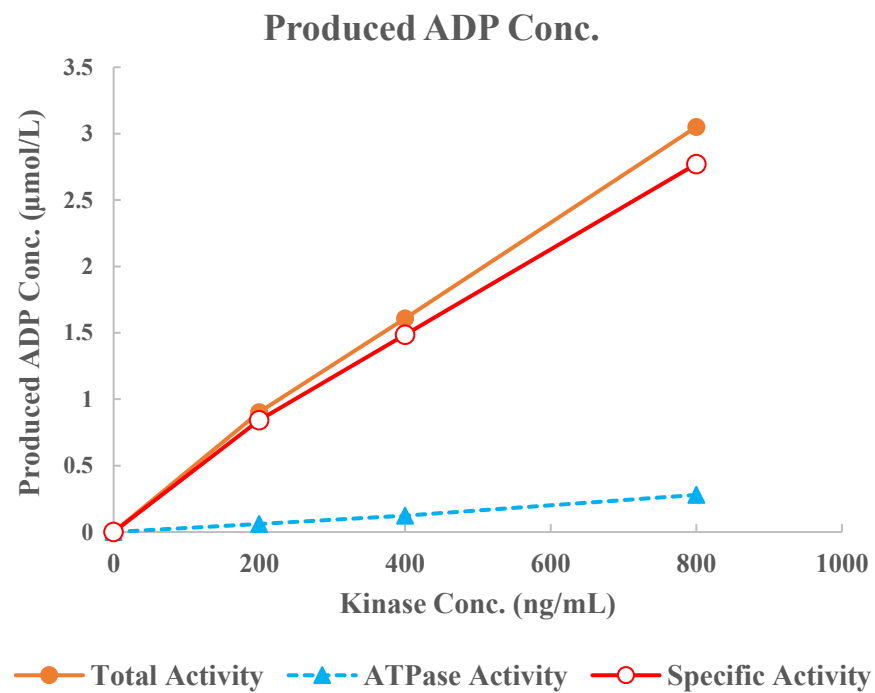
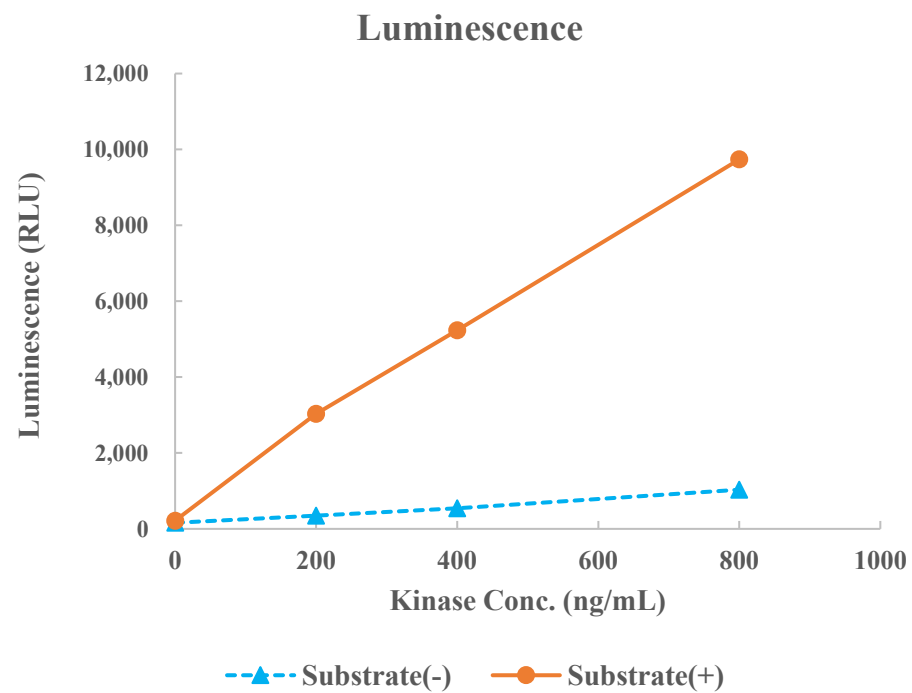


ZAP70



Kinase	ZAP70, 08-177
Substrate	Biotin-Blk/Lyntide, BTSU04
Substrate Conc.	10 μmol/L
ATP Conc.	5 μmol/L
Metal	Mg, Mn
Metal Conc.	5, 1 mmol/L
Additives	-
Reaction Time	1 hour