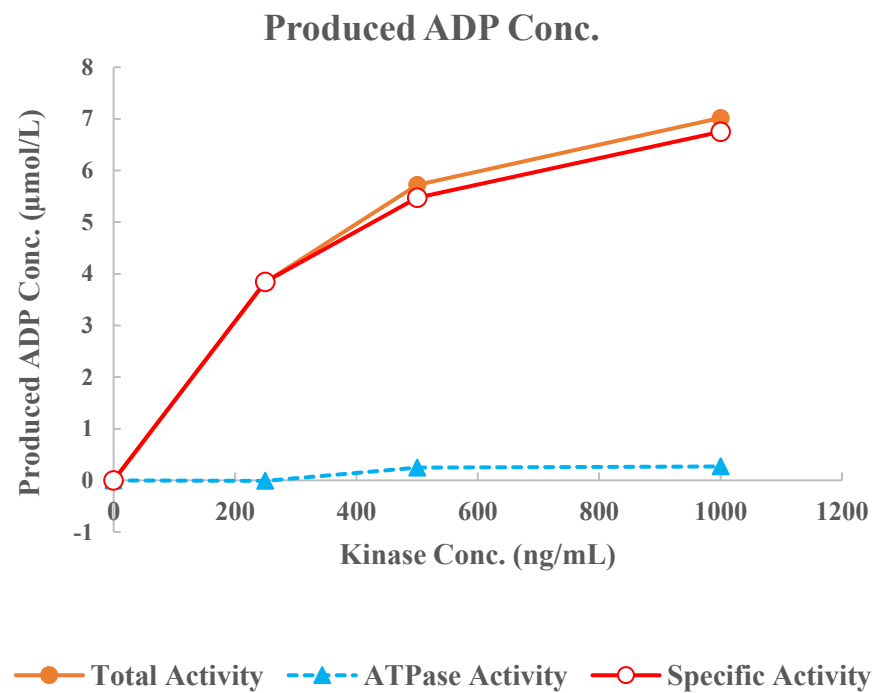
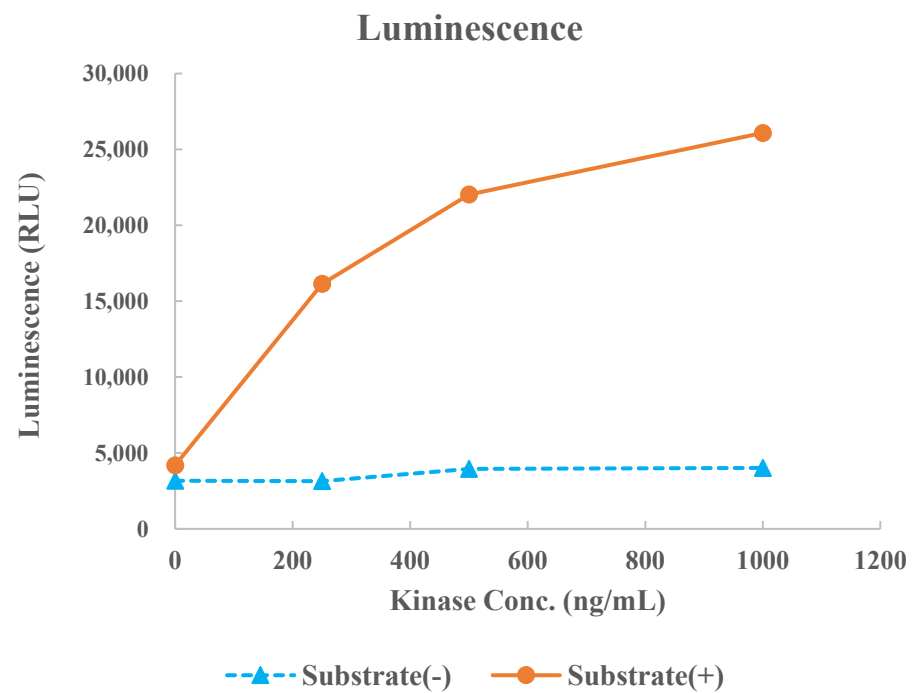


EPHB2



Kinase	EPHB2, 08-129
Substrate	Biotin-Blk/Lyntide, BTSU04
Substrate Conc.	10 μmol/L
ATP Conc.	100 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour