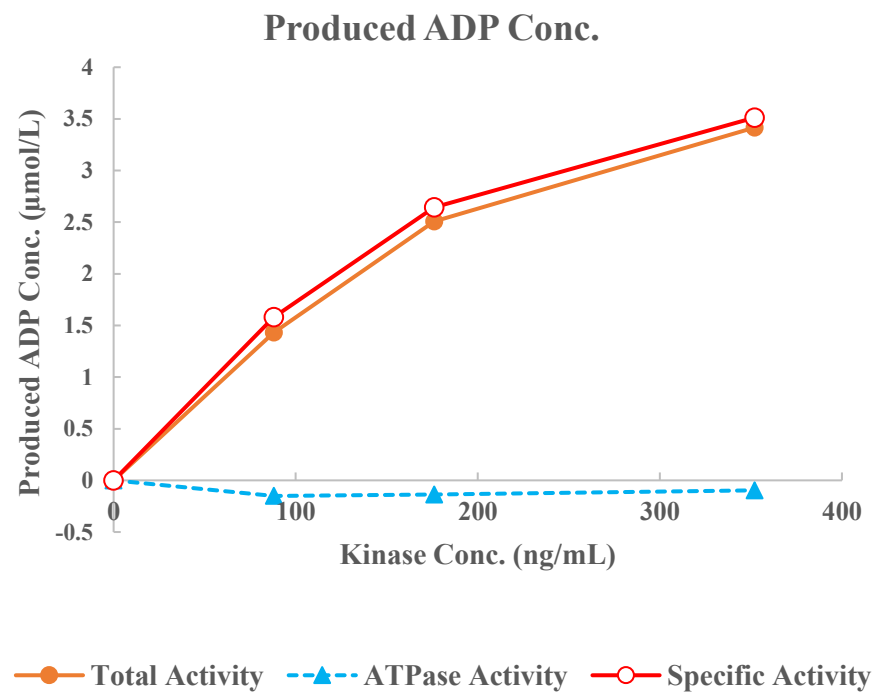
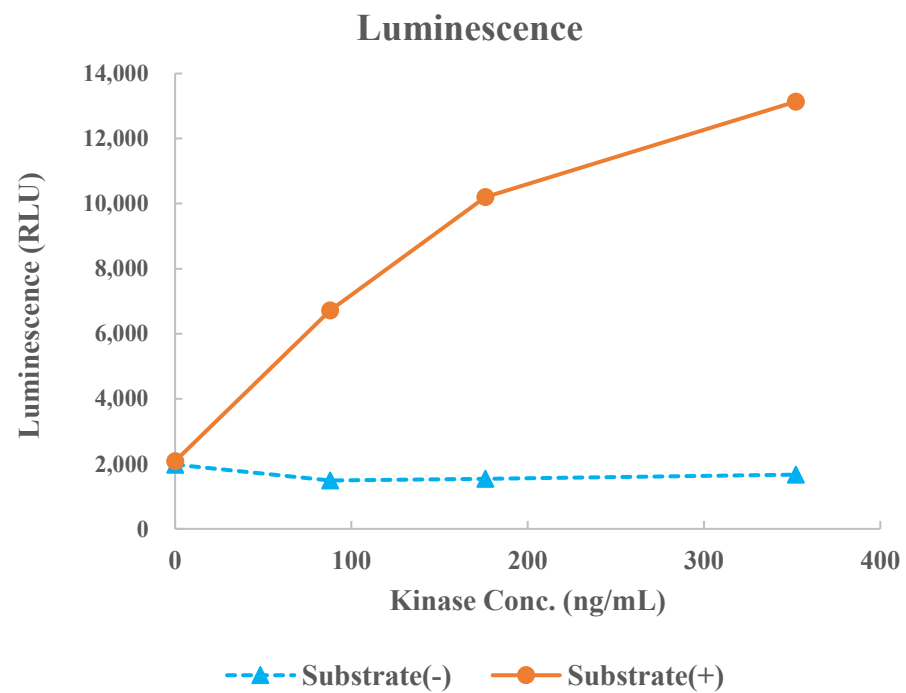


EPHB4



Kinase	EPHB4, 08-131
Substrate	Biotin-Blk/Lyntide, BTSU04
Substrate Conc.	10 μmol/L
ATP Conc.	50 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour