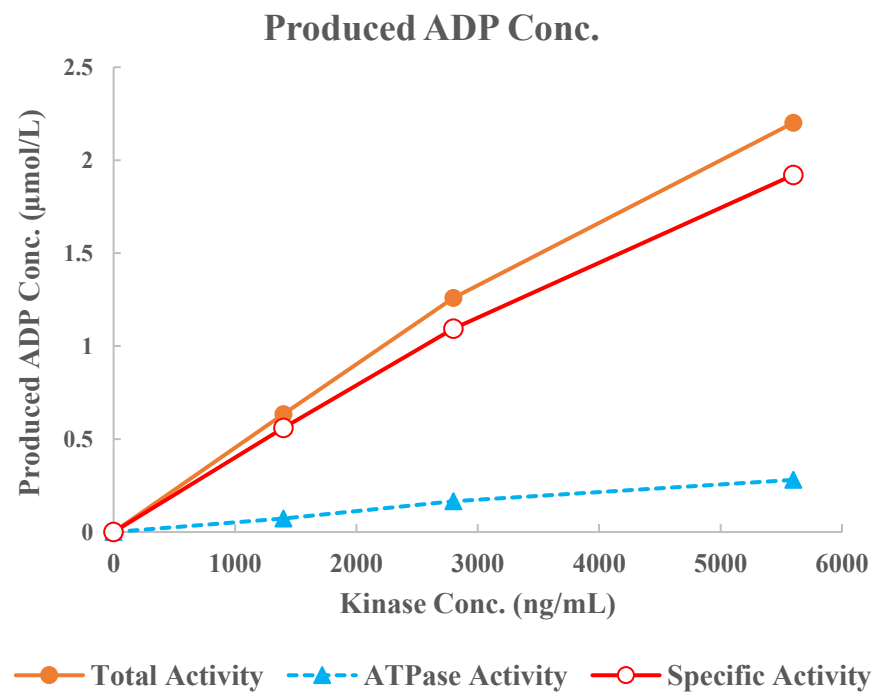
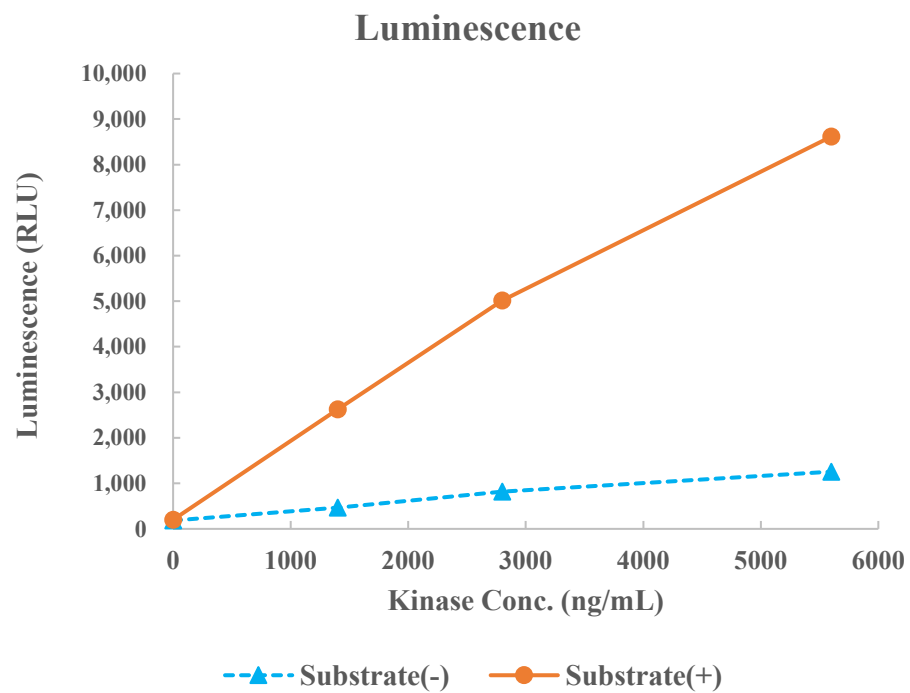


PDGFR α (PDGFRA) [T674I]



Kinase	PDGFR α (PDGFRA) [T674I], 08-503
Substrate	Biotin-CSKtide, BTSU02
Substrate Conc.	10 μ mol/L
ATP Conc.	10 μ mol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour