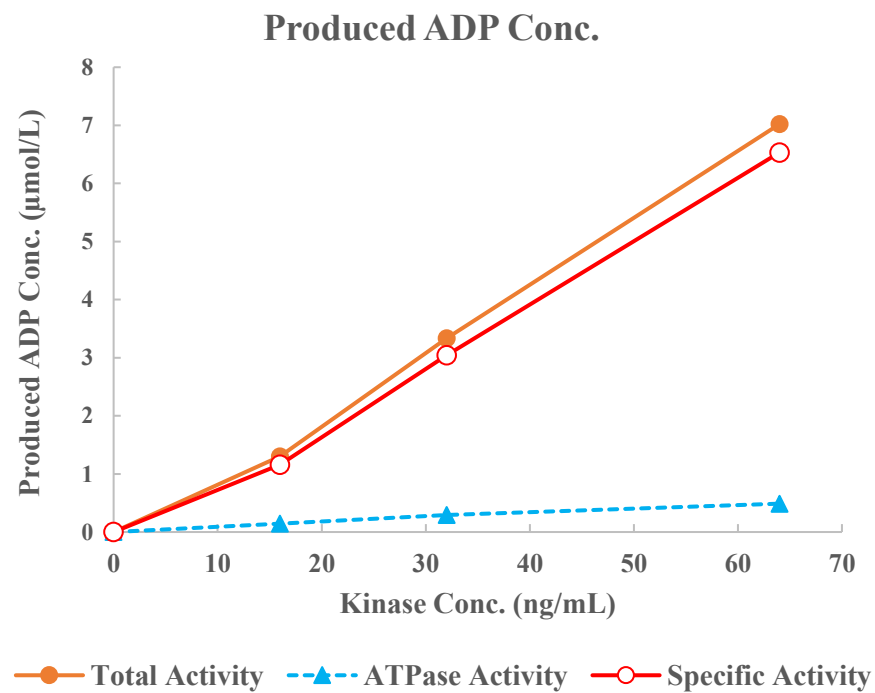
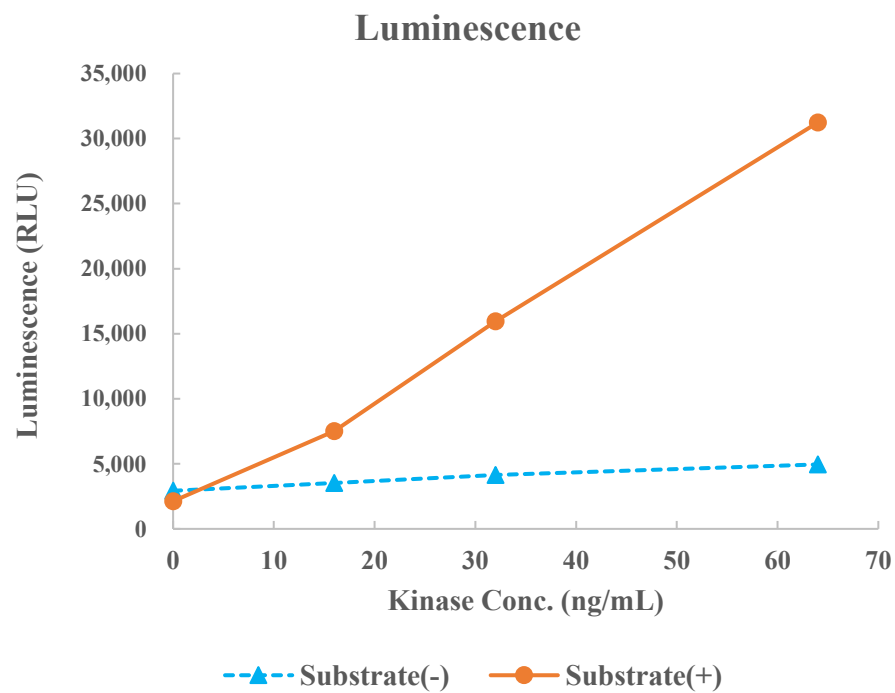


FGFR3 [V555M]



Kinase	FGFR3 [V555M], 08-543
Substrate	Biotin-CSKtide, BTSU02
Substrate Conc.	10 μmol/L
ATP Conc.	50 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour