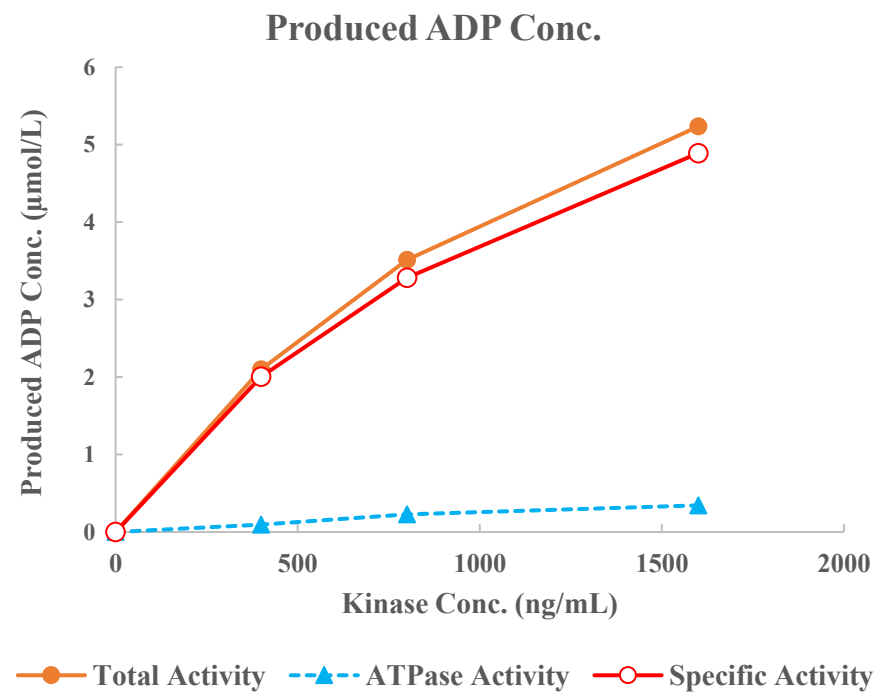
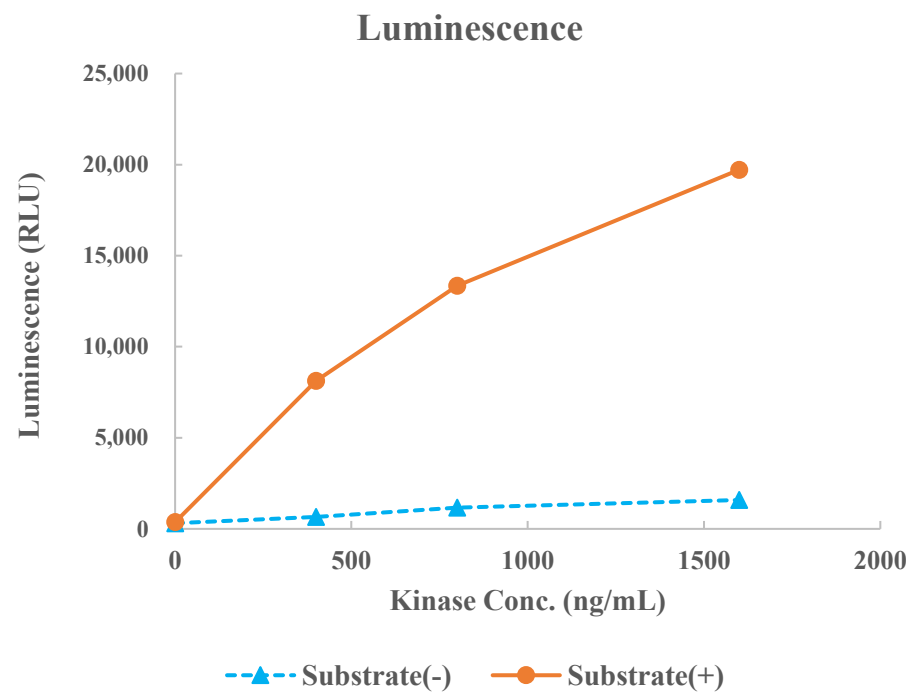


CK1δ(CSNK1D)



Kinase	CK1δ(CSNK1D), 03-103
Substrate	Biotin-CKtide, BTSU12
Substrate Conc.	10 μmol/L
ATP Conc.	10 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour