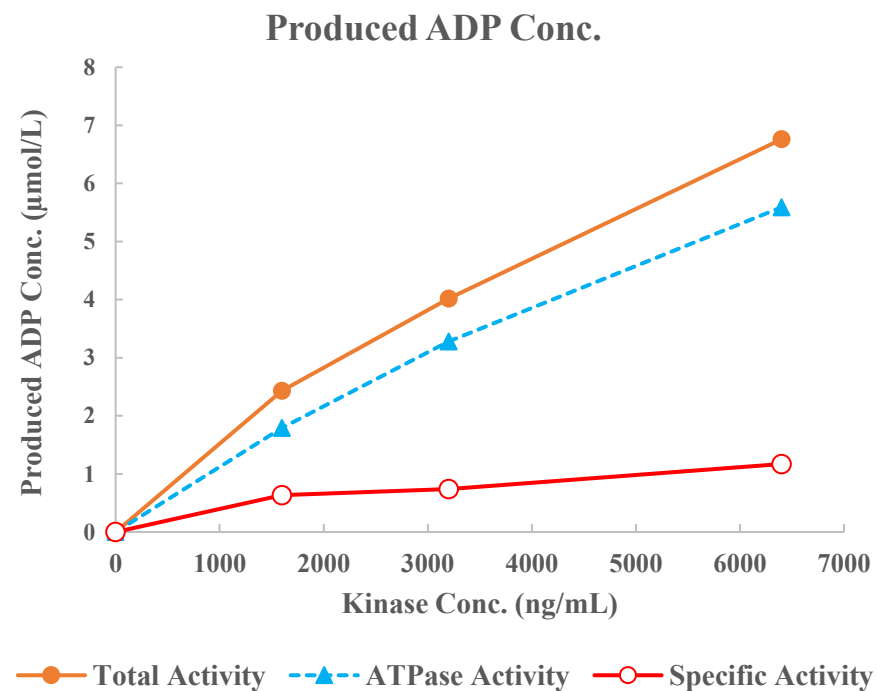
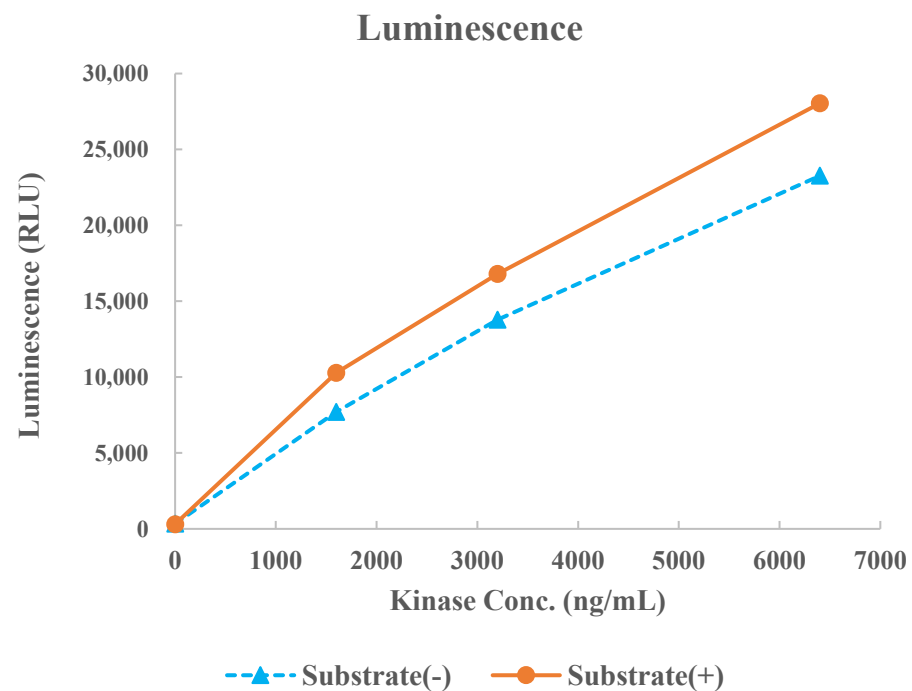


CDC7/ASK



Kinase	CDC7/ASK, 05-109
Substrate	Biotin-MCM2 peptide, BTSU33
Substrate Conc.	10 μmol/L
ATP Conc.	10 μmol/L
Metal	Mg
Metal Conc.	10 mmol/L
Additives	-
Reaction Time	1 hour