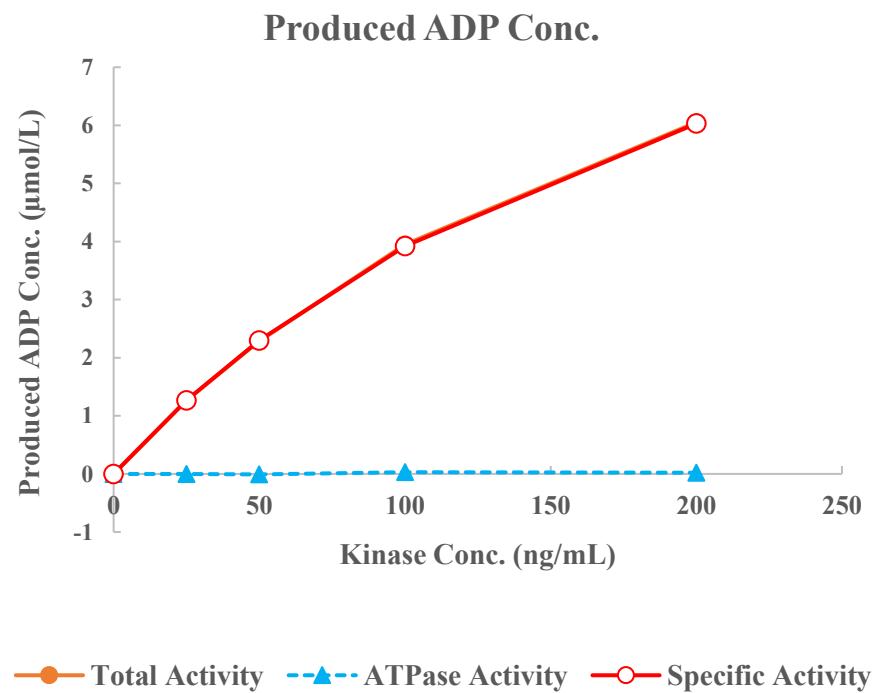
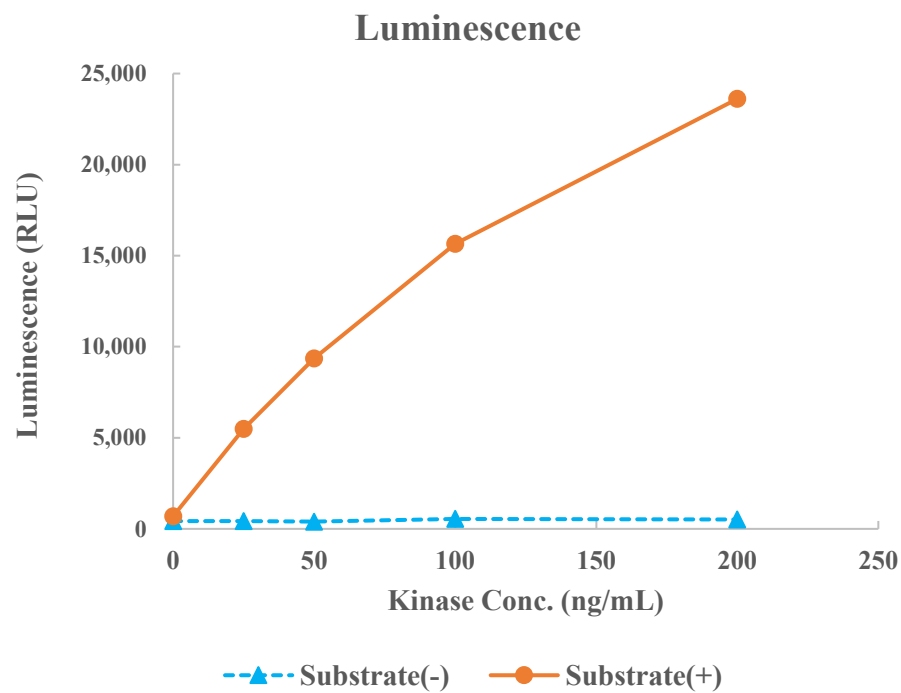


CaMK1δ(CAMK1D)



Kinase	CaMK1δ(CAMK1D), 02-106
Substrate	Biotin-Synapsin peptide, BTSU45
Substrate Conc.	10 μmol/L
ATP Conc.	10 μmol/L
Metal	Mg, Ca
Metal Conc.	5, 1 mmol/L
Additives	10 μg/mL Calmodulin
Reaction Time	1 hour