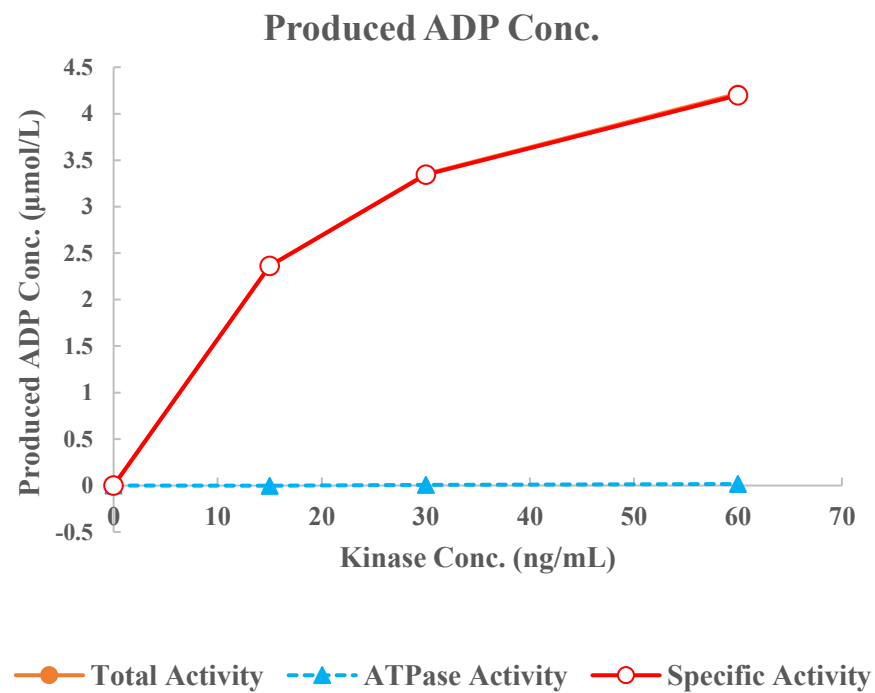
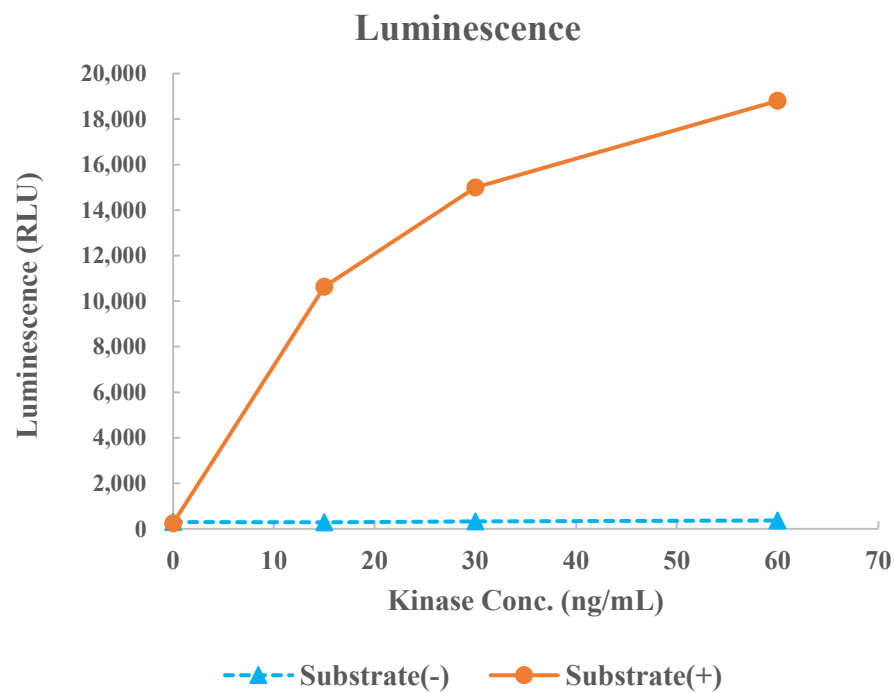


CaMK2δ(CAMK2D)



Kinase	CaMK2δ(CAMK2D), 02-111
Substrate	Biotin-GS peptide, BTSU05
Substrate Conc.	10 μmol/L
ATP Conc.	5 μmol/L
Metal	Mg, Ca
Metal Conc.	5, 1 mmol/L
Additives	10 μg/mL Calmodulin
Reaction Time	1 hour