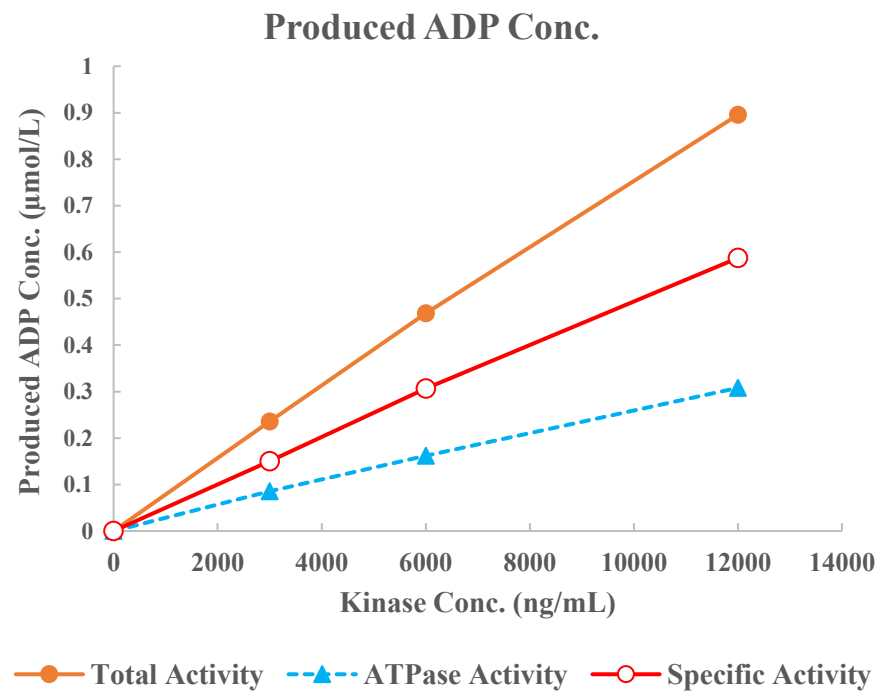
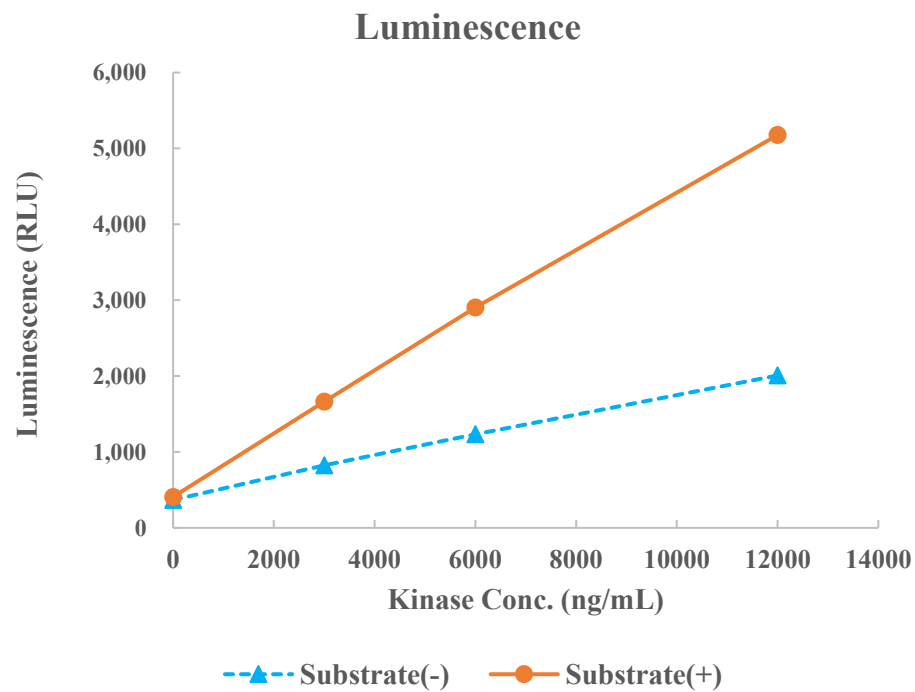


MRCK α (CDC42BPA)



Kinase	MRCK α (CDC42BPA), 01-107
Substrate	Biotin-DAPK1tide, BTSU17
Substrate Conc.	50 μ mol/L
ATP Conc.	10 μ mol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour