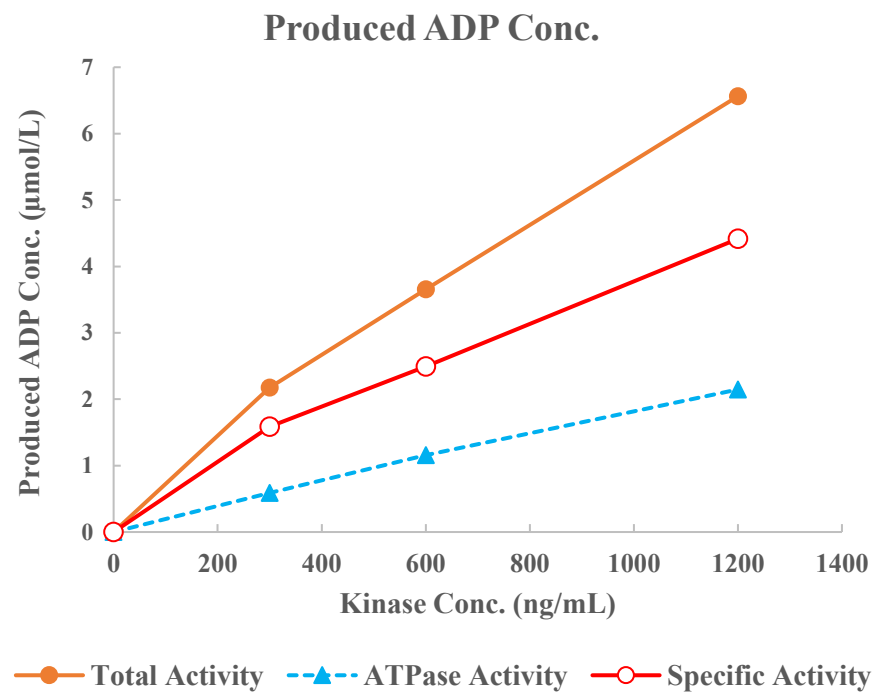
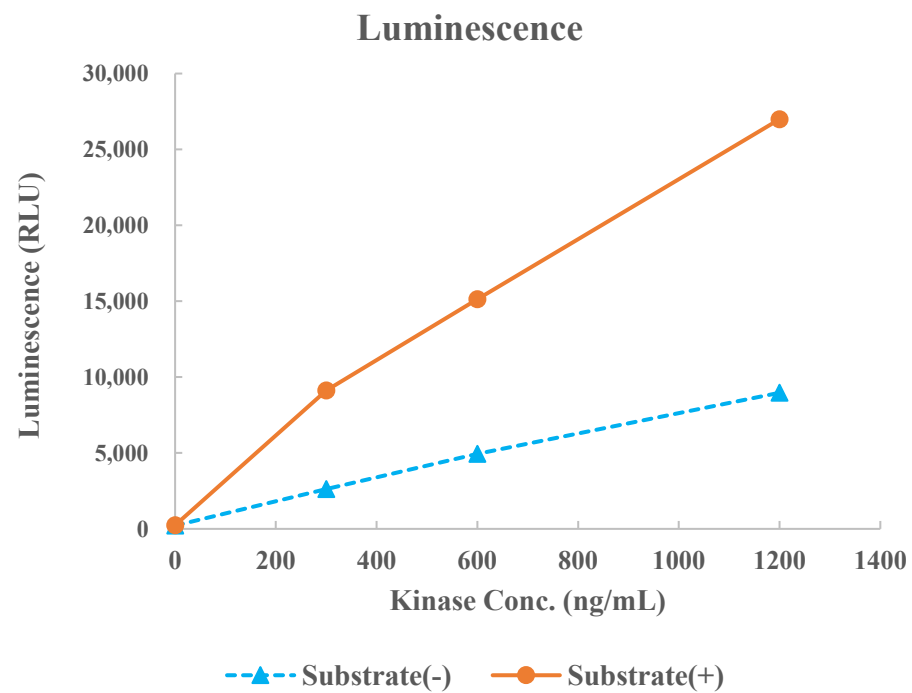


DAPK1



Kinase	DAPK1, 02-134
Substrate	Biotin-DAPK1tide, BTSU17
Substrate Conc.	10 μmol/L
ATP Conc.	10 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour