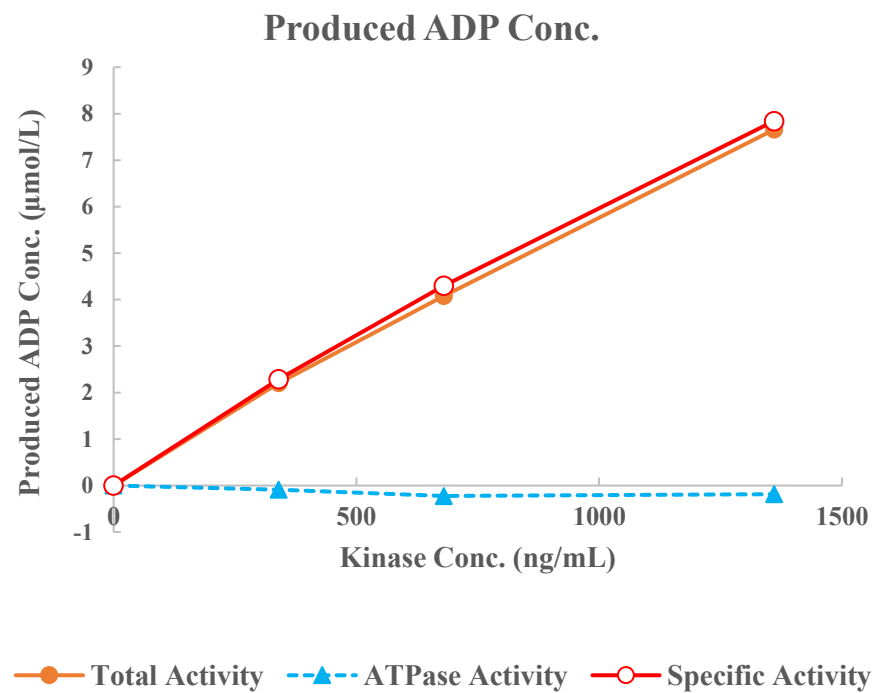
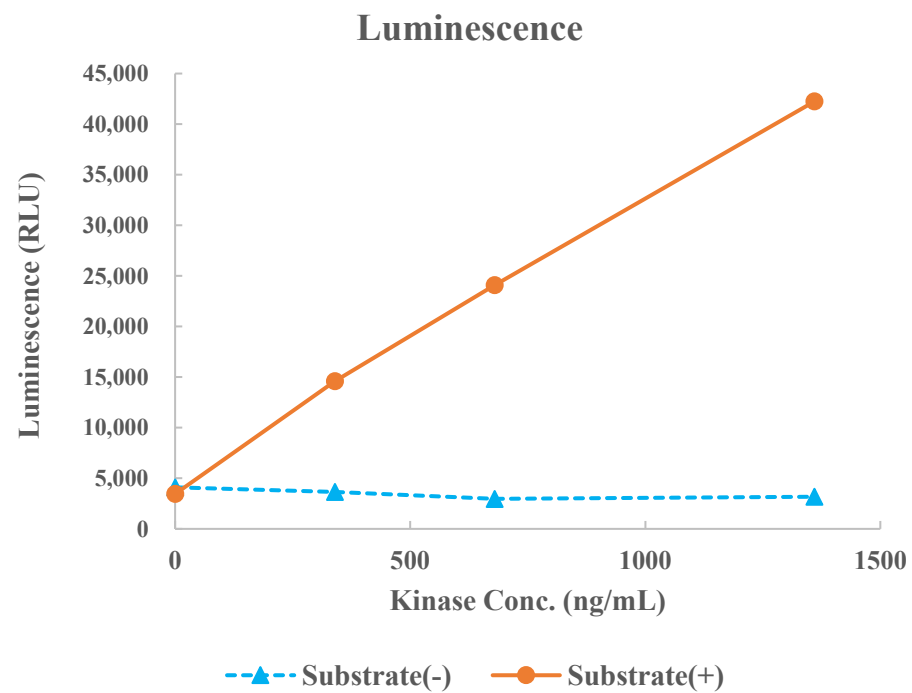


EML4-ALK



Kinase	EML4-ALK, 08-516
Substrate	Biotin-Srcptide, BTSU01
Substrate Conc.	50 μmol/L
ATP Conc.	50 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour